

Subcloning Notebook

Vital tools
and techniques
for transferring
your inserts
between vectors



Promega

Need Help? Get Answers!

If you need help with any product or protocol featured in this notebook, just give a call. All Promega Technical Service scientists have life science degrees and extensive lab experience. Most have post-grad or post-doc lab experience. They understand you and what you face each day. Even if your question involves a product that isn't ours, feel free to call.



Scientists Helping Scientists Succeed!

Visit the Promega Web site for 24/7 support. Find online protocols, product FAQs, application information, references, publications, calculators and many other valuable support tools. It's all easy to access and understand.

- ***Troubleshooting experiments***
- ***Automation installation and support***
- ***Global availability***
- ***Multi-language capabilities***



www.promega.com

Toll Free in USA: 800-356-5926

Contact your local Promega Branch Office or Distributor

Subcloning Notebook

Table of Contents

Chapter 1: Classic Subcloning

Basic Steps for Subcloning	3
Subcloning Strategy	4–9
Restriction Digestion	10–12
Double Enzyme Digests	13
Partial Restriction Digestion	14
Creating Blunt Ends	15–17
Dephosphorylating Vectors	18–22
Ligation	23–26
Purifying Vector and Insert	27–29
Gel Electrophoresis	30
DNA Markers	31
Ordering Information	32–34

Chapter 2: PCR Subcloning

Introduction	35
T-Vector Systems	36–37
Giving Blunt-Ended DNA an A-tail for T-Vector Subcloning	38
Subcloning with RE Sites	39
Subcloning using PCR Primers Containing Restriction Sites	40
Ordering Information	41–42

Chapter 3: Transforming Bacteria

Properties of <i>E. coli</i> Strains for Subcloning	43
Ready-to-Use Competent Cells	44–45
Making Your Own Competent Cells	46
Determining Transformation Efficiency of Competent Cells	47
Transforming Ligation Reactions	47
Media and Solutions	48

Chapter 4: Screening for Recombinants

Introduction	49
Colony PCR	50
Go Directly to Gel	51
Screening by Plasmid Minipreps and RE Digests	52
Plasmid Minipreps	53
Troubleshooting Subcloning Experiments	54–55
Ordering Information	56

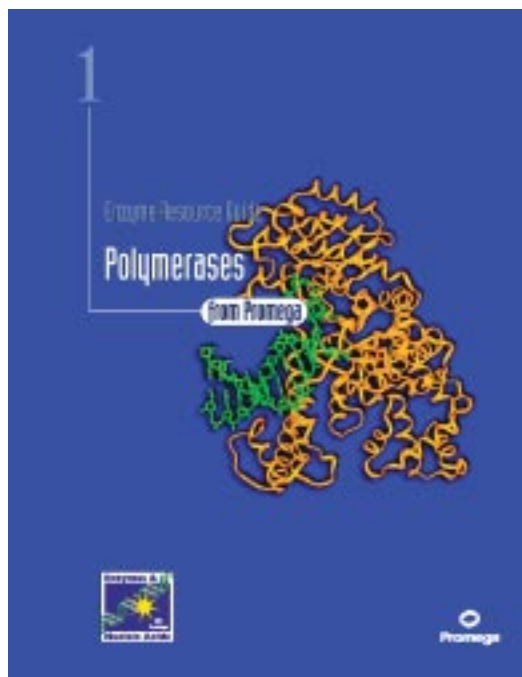
Chapter 5: Technical Appendix

Restriction Enzyme Activity In Promega 10X Buffers, Reaction Temperature and Heat Inactivation	57–58
Isoschizomers	59–60
Compatible Ends	61
Site-Specific Methylation Sensitivity of Promega Restriction Enzymes	62
Restriction Enzyme Buffer Composition	63
Copy Number of Commonly Used Plasmids	63
Star Activity	63
Genotypes of Frequently Used Bacterial Strains	64
Genetic Markers in <i>E. coli</i>	65–66
Nucleic Acid Calculations	67
Formulas for DNA Molar Conversions	67

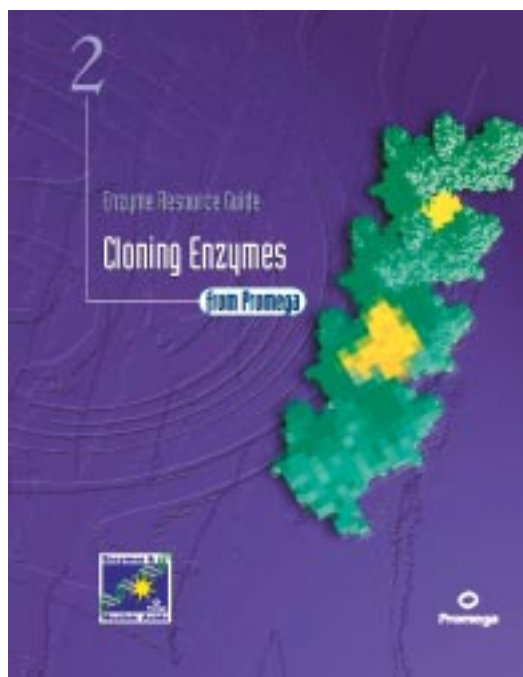
Classic Subcloning

Starving for more in-depth knowledge of the enzymes mentioned in this notebook?

Savor these tasty offerings at the Promega Web site:



The **Polymerases Guide** features mesophilic and thermophilic DNA polymerases, RNA polymerases, reverse transcriptases and terminal transferase. The guide includes background information, a list of applications for each polymerase, enzyme properties and other product information such as the quality control assays performed to ensure high-quality enzymes. Figures and tables enhance the text reference information. The guide is available in pdf format online at: www.promega.com/guides/polym_guide/

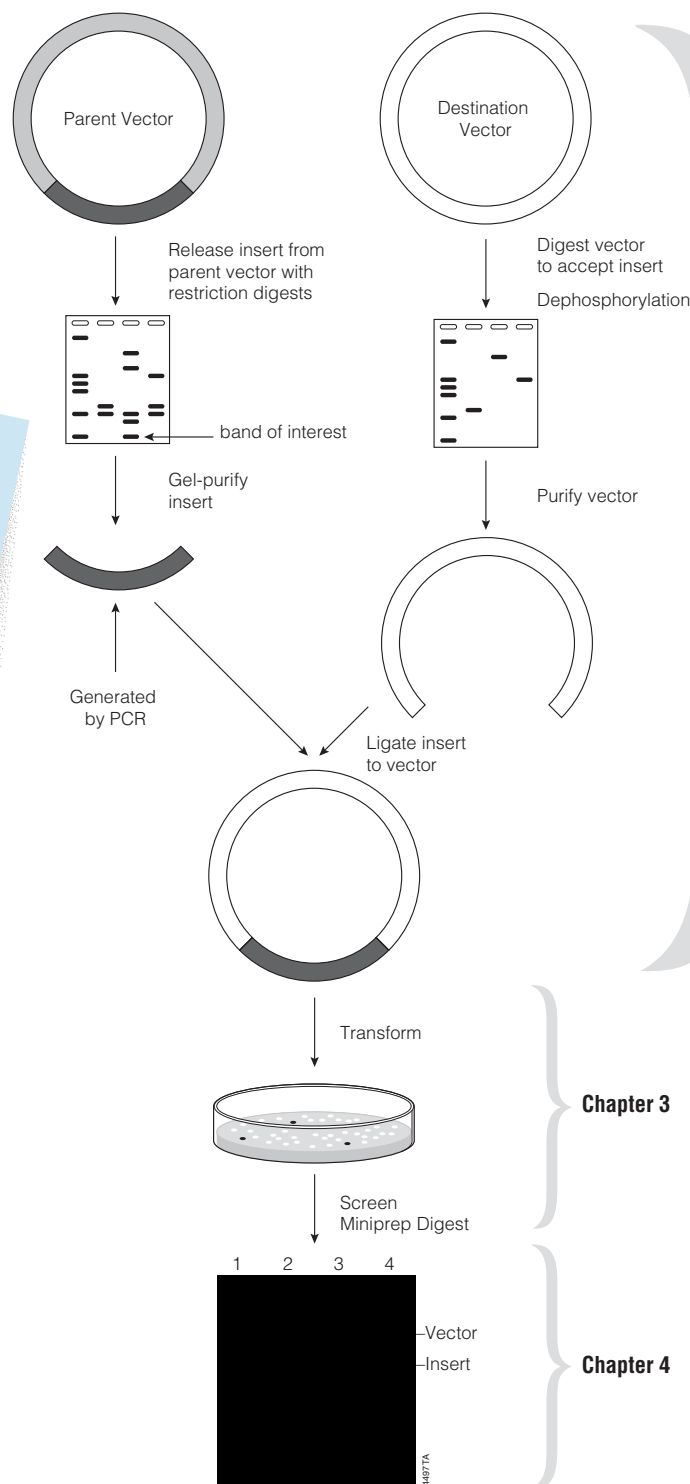


Cloning Enzymes, in the Enzyme Resource Guide series, highlights those enzymes important in nucleic acid cloning procedures. Enzymes that modify nucleic acids provide the foundation for many molecular biology techniques. Specific features of the in vivo functions of these enzymes have been exploited in vitro to provide many of the protocols currently used in nucleic acid manipulations. The guide is available in pdf format online at: www.promega.com/guides/cloning_guide/

Classic Subcloning

Basic Steps for Subcloning

Subcloning is a basic procedure in molecular biology required to move inserts from one vector to another to gain the desired functionality to study your insert. Essentially all subcloning reactions proceed the same way as illustrated in the figure below. You release and purify your insert from the parent vector, ligate this insert into a prepared destination vector, transform this ligation reaction into competent bacterial cells. Then you screen the transformed cells for the insert. This *Subcloning Notebook* will guide you through every step in the process.



Dephosphorylation reduces the chance of vector self-ligation to virtually zero.

Chapters 1 and 2

Gel isolation of vector reduces background by eliminating uncut vector from the transformation.

Chapter 3

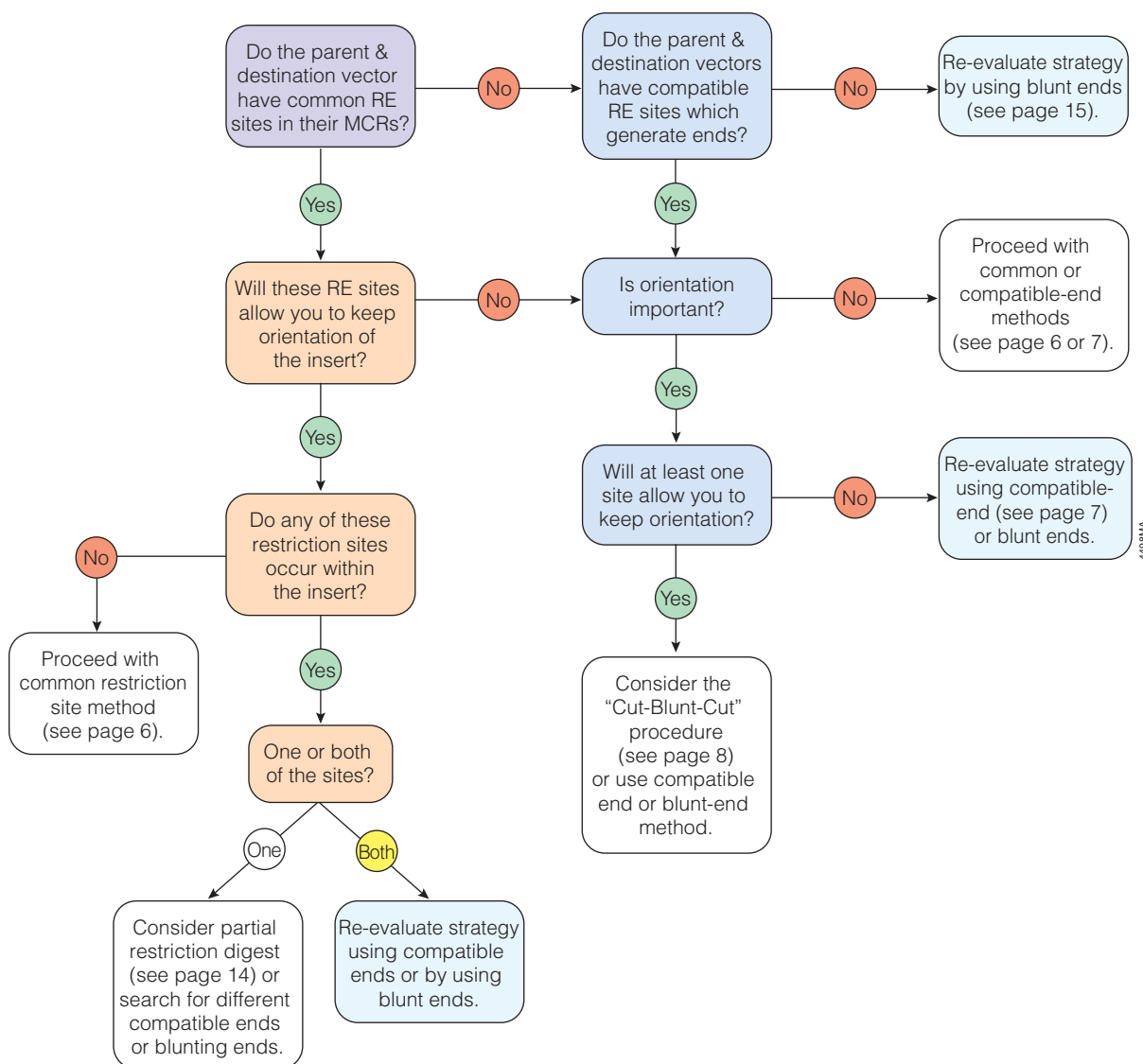
Chapter 4

Classic Subcloning

Subcloning Strategy

Before you begin your subcloning, you need to know: The restriction enzyme (RE) sites available for subcloning in your parent vector multiple cloning region (or in the insert if you need to digest the insert); the RE sites available in the destination vector multiple cloning region (MCR); and if these same sites also occur in your insert. Once you know this information, you can begin to ask questions about which subcloning strategy to use.

See the Compatible Ends Table on page 61 for a listing of overhangs compatible with Promega enzymes.

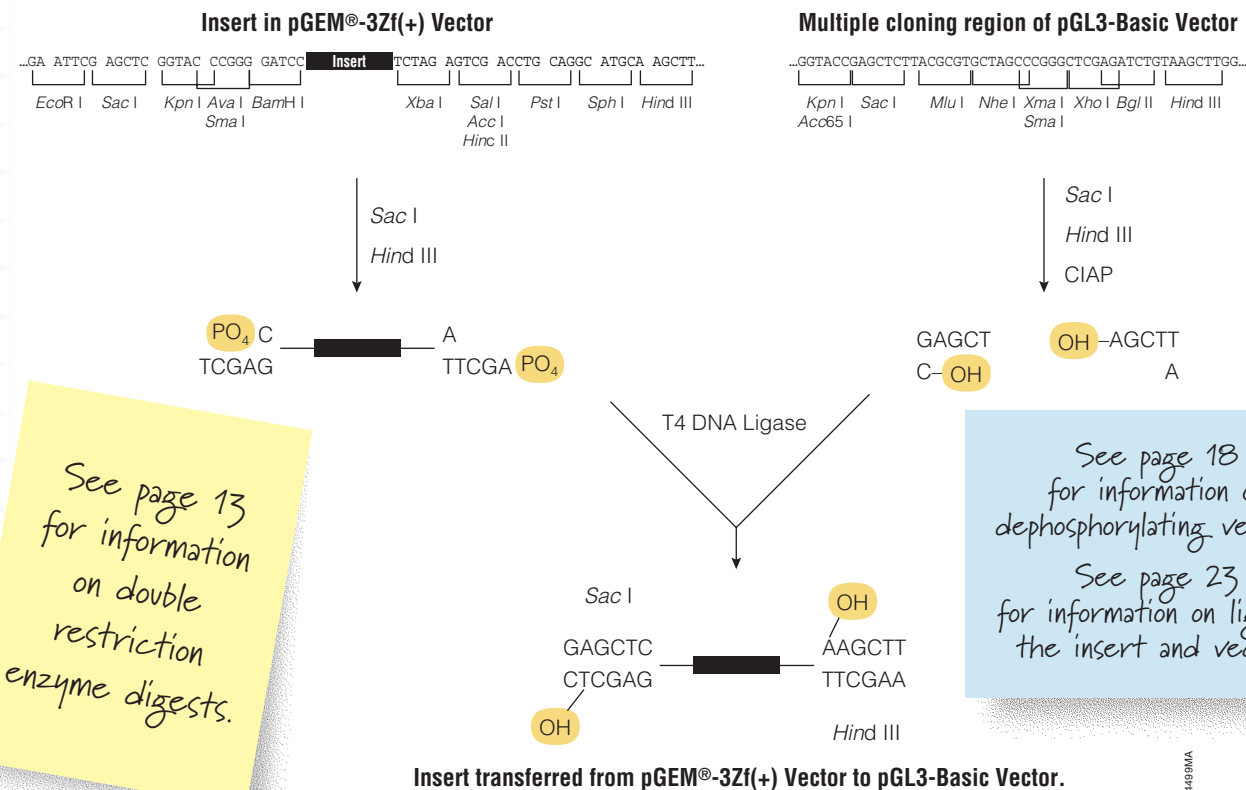


Classic Subcloning

Subcloning Strategy: Common Restriction Sites

If your parent and destination vector multiple cloning regions contain common restriction sites and neither of these restriction sites occur within your insert, you have a very straightforward subcloning process. You digest your parent and destination vectors with the same two enzymes followed by dephosphorylation of the destination vector. The insert and the dephosphorylated vector are then separated on an agarose gel and purified using a system such as the Wizard® SV Gel and PCR Clean-Up System (see page 28) and ligated.

The T4 DNA Ligase will join the DNA through reforming the bond between the 5'-PO₄ coming from the insert and the 3'-OH of the vector. The vector has been dephosphorylated so the second bond will not be formed in vitro (indicated by the OH). These nicks will be repaired in the bacteria upon transformation.

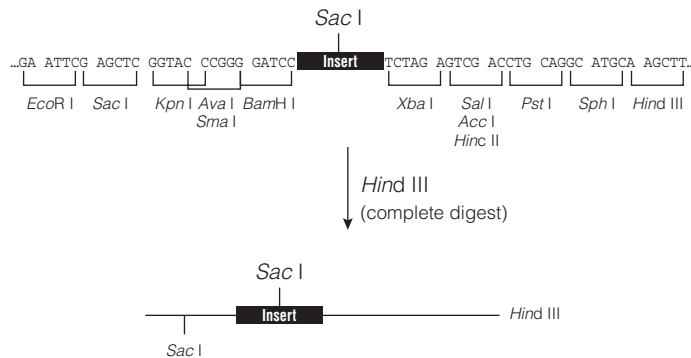


Classic Subcloning

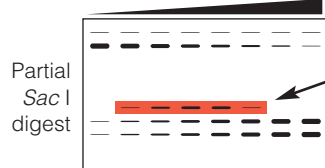
Subcloning Strategy: Common Restriction Sites with Partial Digests

Having a restriction site in both the multiple cloning region and the insert does not exclude the use of this site for subcloning. A partial restriction digest strategy can be employed.

Insert in pGEM®3Zf(+) Vector

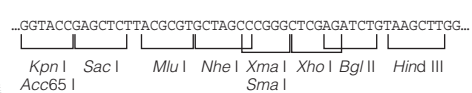


Sac I Concentration



Gel isolate the band you want.

Multiple cloning region of pGL3-Basic Vector



*Sac*I
*Hind*III
(complete digest with both enzymes)
CIAP



T4 DNA Ligase



Insert transferred from pGEM®-3Zf(+) Vector to pGL3-Basic Vector.

See page 14 for information on partial restriction digests.

45201MA

Classic Subcloning

Subcloning Strategy: Moving Inserts with Compatible Restriction Sites

If you don't have common restriction sites in the parent and destination vector multiple cloning regions, you may have compatible restriction sites. Compatible restriction sites have the same overhang sequence and can be ligated together. In this example, *Xba* I and *Nhe* I both produce the same 5' overhang sequence. Cut sites from these two are exactly matching and ligate well. However, neither the *Xba* I or *Nhe* I sites are regenerated in the ligation. A table of compatible ends is present on page 61 of this Notebook. Compatible end ligation is straightforward after the enzymes are identified.

Xba I or *Spe* I is compatible with the *Nhe* I site of the pGL3-Basic Vector.

Xba I and *Sal* I have better buffer compatibility for the double digest than *Spe* I and *Sal* I.

Sal I is compatible with the *Xho* I site in the pGL3-Basic Vector.

Insert in pGEM®-9Zf(-) Vector

...TAT GCATCACTAG TAAGC TTTGC TCTAG A **Insert** GAATT CGTCG ACGAG CTC...

Nsi I Spe I Hind III Xba I EcoR I Sal I Sac I

Xba I

Sal I

Gel Isolation of insert

PO₄ CTAGA
T

G
CAGCT PO₄

Multiple cloning region of pGL3-Basic Vector

...GGTACCGAGCTCTTACGCGTCTAGCCCGGCTCGAGATCTGTAAGCTTGG...

Kpn I Sac I Mlu I *Nhe* I *Xma* I *Xho* I Bgl II Hind III
Acc65 I Sma I

Nhe I

Xho I

Dephosphorylation
Gel Isolation

G
CGATC OH

OH TCGAG
C

T4 DNA Ligase

GCTAGA
CGATCT

GTCGAG
CAGCTC

~~*Xba* I~~

~~*Nhe* I~~

OH

~~*Xho* I~~

~~*Sal* I~~

Insert transferred from pGEM-9Zf(-) Vector to pGL3-Basic Vector.

See the Compatible End Table on page 61 of the Technical Appendix for a listing of compatible ends to Promega enzymes.

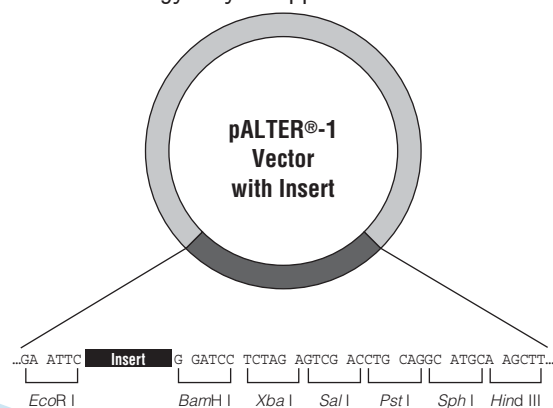
In this example, none of the restriction sites used for the compatible-end subcloning are regenerated in the final ligation product.

4501MA

Classic Subcloning

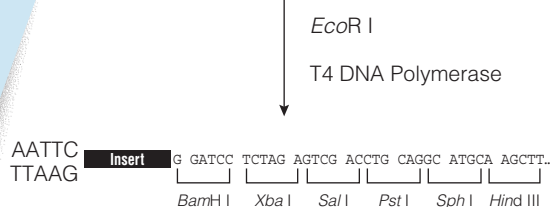
Subcloning Strategy: Moving Inserts with Only One Common Site

You've looked for common sites or compatible sites and you can find only one match on one side of your insert. What do you do about the other side of the insert? You can use a method commonly referred to as "cut-blunt-cut". Any restriction site can be made blunt through the action of T4 DNA Polymerase. Simply digest the parent vector and blunt that site with T4 DNA Polymerase (protocols on page 16), run the products on a gel, purify and proceed with the common or compatible end restriction enzyme digestion. In this example, the destination vector has *Sma* I site, which leaves a blunt end. Most vectors have at least one blunt-ended restriction site that can accept the newly created blunt end from the insert. If you don't have such a site or the site would not be in the correct orientation, the same "cut-blunt-cut" strategy may be applied to the destination vector as well.



The pALTER®-1 Vector is used with the Altered Sites® in vitro Site-Directed Mutagenesis System.

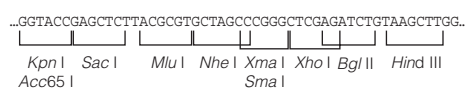
See page 16 for T4 DNA Polymerase procedure.



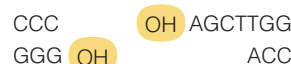
Hind III



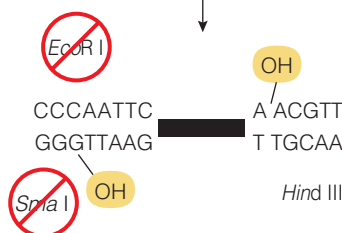
Multiple cloning region of pGL3-Basic Vector



Sma I
Hind III
CIAP
Gel Isolation



T4 DNA Ligase



This may commonly be referred to as the "Cut-Blunt-Cut" strategy.

The cut-blunt-cut strategy can also be used on destination vectors as well. If you don't have a ready-to-use blunt site, make one!

Insert transferred from pALTER®-1 Vector to pGL3-Basic Vector.

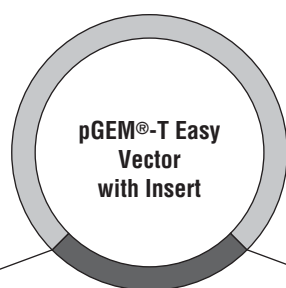
4502MA

Classic Subcloning

Subcloning Strategy: Blunt-End Method

You can't find a single common site or compatible site in the parent or destination vector. What do you do? Many people resort to amplifying the insert with restriction sites in the primers to provide the compatibility, but this strategy may cause some problems (i.e., introduction of mutations, difficulty digesting PCR products [see page 40]). Another method involves straight blunt-end cloning. You cut out your insert with whichever enzymes you desire. Treat with T4 DNA Polymerase to blunt either 5' or 3' overhangs and ligate into the destination vector opened with a blunt-end cutter or made blunt by T4 DNA Polymerase. Remember though, this method will not retain orientation of your insert so you will have to screen for orientation by methods like those outlined on page 50.

The pGEM-T Easy Vector is designed for direct cloning of PCR products. See page 37.



Multiple cloning site of pGL3 Basic Vector

...GGTACCGAGCTCTTACGCGTGCTAGCCGGGCTCGAGATCTGTAAGCTTGG...
Kpn I Sac I Mlu I Nhe I Xma I Xho I Bgl II Hind III
Acc65 I Sma I

...TCCC GGCCG CCATG GCGGC CGCGG GAATT CGAT
...AGGG CCGGC GGTAC CGCCG GCGCC CTTAA GCTA
BstZ I Nco I Not I BstZ I Sac II EcoR I
ATCAC TAGTG AATTC GCGGC CGCCT GCAG...
TAGTG ATCAC TTAAG CGCCG GCGGA CGTC...

See page 16 for T4 DNA Polymerase procedure.

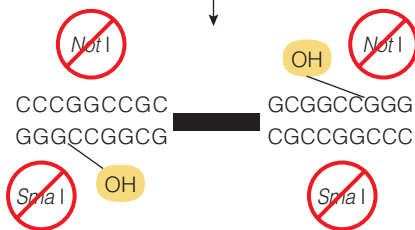
Not I
T4 DNA Polymerase
dNTPs

PO₄ GGCCGC CCGGCG Insert GCGGCC CGCCGG PO₄

Sma I
CIAP
Gel Isolation

CCC GGG OH
GGG OH

T4 DNA Ligase



Insert transferred from pGEM-T Easy Vector to pGL3-Basic Vector.

The blunt-end method will not maintain the orientation of your insert.

If there is no blunt-ended RE site in your destination vector, you can use T4 DNA Polymerase to make the cut vector blunt-ended.

Classic Subcloning

Restriction Digestion

Restriction endonucleases (RE), also referred to as restriction enzymes, are proteins that recognize short, specific (often palindromic) DNA sequences. Type II REs cleave double-stranded DNA (dsDNA) at specific sites within or adjacent to their recognition sequences. Many restriction enzymes will not cut DNA that is methylated on one or both strands of the recognition site, although some require substrate methylation (see page 62).

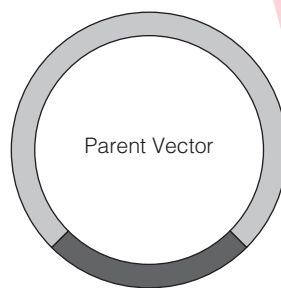
Restriction digestion is one of the most common reactions performed in molecular biology. For a digestion with a single RE the reaction is very simple:

Nuclease-Free Water	14µl
10X Restriction Buffer	2µl
Acetylated BSA (1mg/ml)	2µl
DNA (~1µg)	1µl
Restriction Enzyme (10u)	1µl
Final Volume	20µl

Mix by pipetting and collect the contents at the bottom of the tube. Incubate at the appropriate temperature for the enzyme for 1–4 hours. Add 4µl of 6X Blue/Orange Loading Dye and analyze digested DNA by gel electrophoresis.

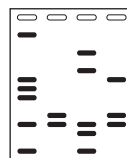
Preparing an insert for transfer from one vector to another usually requires a **double digest** (digest with two different REs). If both restriction enzymes work in the same restriction enzyme buffer, the reaction is straightforward. **Simply add 1µl of the second restriction enzyme and adjust the amount of water used.**

Remember, restriction enzymes are commonly stabilized in 50% glycerol solution. Do not exceed 5% glycerol in final digest with the two enzymes. Glycerol concentrations >5% may lead to star activity (see page 63).



Parent Vector

Release insert from parent vector with restriction digests.



Gel-purify insert.



Gel isolation is a necessity in subcloning. You get the insert you need.

Learn more about the history and enzymology of restriction enzymes with the Promega Restriction Enzyme Resource located at: www.promega.com/guides



Look at these search tools to help you plan your experiments. www.promega.com/guides/re_guide/default.htm

Classic Subcloning

Restriction Digestion

What is supplied with Promega Restriction Enzymes?

Each RE has specific requirements for optimal activity. Ideal storage and assay conditions favor the highest activity and highest fidelity in a particular enzyme's function. Conditions such as temperature, pH, enzyme cofactors, salt composition and ionic strength affect enzyme activity and stability.

Each Promega Restriction Enzyme is supplied with:

- The optimal reaction buffer
This may be from the 4-CORE® System (Reaction Buffers A, B, C, D) or one of the other optimal buffers (Reaction Buffers E–L). This buffer always yields 100% activity for the enzyme that it accompanies, and serves as the specific reaction buffer for single digests.
- MULTI-CORE™ Buffer
This is designed for broad compatibility and is provided with enzymes that have 25% or greater activity in this buffer. The MULTI-CORE™ Buffer is useful for multiple digests because it generally yields more activity for more enzyme combinations than any of the other buffers, but sometimes using the MULTI-CORE™ Buffer can compromise enzyme activity. Multiple digests using REs with significantly different buffer requirements may require a sequential reaction with the addition of RE buffer or salt before the second enzyme is used.
- 100X Acetylated BSA
We recommend adding 0.1mg/ml acetylated BSA to every reaction. The acetylated BSA improves the stability of the enzyme in the reaction.

For more information on the use of acetylated BSA in restriction digests, see "BSA and Restriction Enzyme Digestions" in *Promega Notes* 60 at:

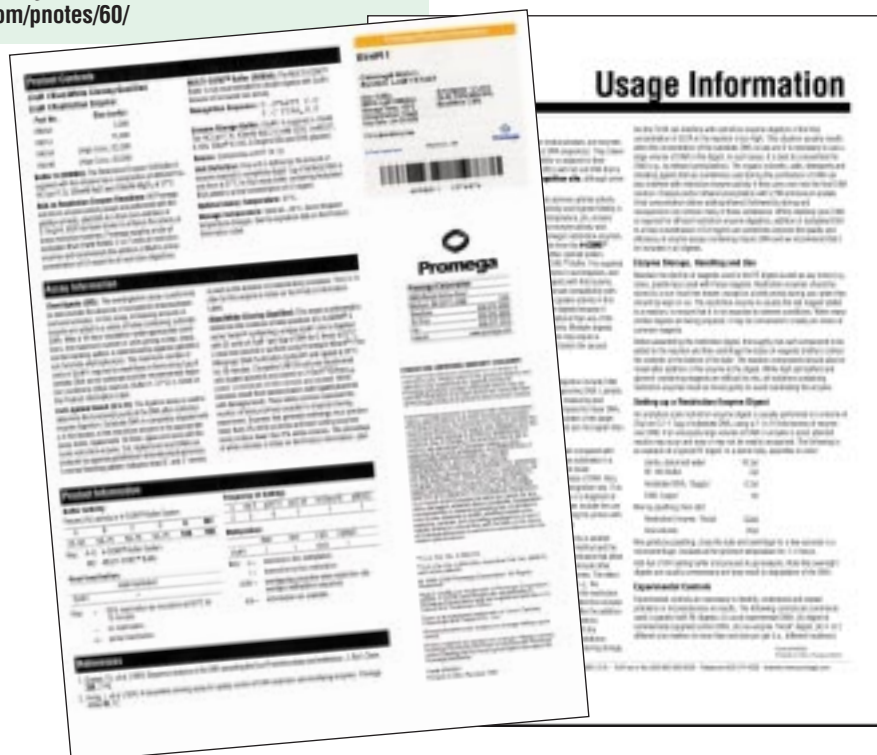
www.promega.com/pnotes/60/

For a recent review on restriction enzymes see:

Williams, R.J. (2003)
Restriction Endonucleases: Classification, properties and applications. Mol. Biotechnol. 23, 225–43.

Easily locate usage and lot information

Each enzyme comes with a Promega Product Information Sheet (PPI) that contains details of quality control assays performed, lot-specific information and usage information. The sheet also has protocol information and references. The lot-specific information is printed on a removable sticker that can be pasted into a notebook or logbook, making your record keeping easier.



Classic Subcloning

Restriction Digestion: Other Considerations

Do both enzymes work at the same temperature?

The majority of restriction enzymes work best at 37°C, but those isolated from thermophilic bacteria require higher temperatures for maximal activity (e.g., *Bst*X I and *Bst*Z I work best at 50°C). Some work below 37°C like *Sma* I (25°C) and *Csp* I (30°C). If you must work with two enzymes with different optimum temperatures, you can use the sequential digest method (assemble all components, perform for the lower-temperature digest first, then digest at the higher temperature second). Usually an hour at each temperature will work fine.

When working with an enzyme that requires a temperature above 37°C, evaporation of the reaction can lead to increased glycerol concentration, which can in turn lead to star activity. Evaporation can be avoided in such reactions by applying a few drops of molecular biology grade mineral oil above the reaction. Clean up with the Wizard® SV Gel and PCR Clean-Up System to remove the mineral oil and recover the pure DNA.

Do my enzymes exhibit methylation sensitivity?

An often overlooked reason for a restriction enzyme failure is sensitivity to *dam* and *dcm* methylation. Many common bacterial strains like JM109, XL1-Blue, and DH5α™ are positive for these two genes. The *dam* gene encodes a DNA adenosine methylase that methylates the N6 position of the adenine residue in the sequence: 5'...GATC...3', a common sequence within many restriction sites. The *dcm* gene encodes a DNA cytosine methylase that methylates the C5 position of the internal cytosine residue in the sequence: 5'...CCAGG...3'. Some restriction enzymes are sensitive to these methylations and will not cut their recognition sequence if the methylation occurs within the recognition site (e.g., *Bcl* I and *dam* methylation) or overlaps the recognition site (e.g., the ATCGAT recognition site falling within the context of ...GATCGAT... or ...ATCGATC... for *dam* methylation).

See the tables on pages 57–58 for optimal reaction temperatures of Promega Restriction Enzymes.

See the table on page 62 for methylation sensitivities of Promega Restriction Enzymes.

- Need to digest a piece of DNA with a *dam* or *dcm* sensitive enzyme?
- Check to see if the enzyme has an isoschizomer or neoschizomer. The isoschizomer or neoschizomer may not be sensitive to the methylation.
 - Transform the plasmid into a *dam/dcm* minus bacterial strain like JM110.

See the tables on pages 59–60 for listings of isoschizomers and neoschizomers.

Classic Subcloning

Double Enzyme Digests

Double Digests with a Common Buffer

In many cases, the enzymes are not supplied with the same reaction buffer, and another buffer may be appropriate. In these cases, activities in other buffers must be assessed by consulting buffer activity charts like those on pages 57–58. In this chart, all Promega REs are tested in Buffers A, B, C, D and MULTI-CORE™ Buffers. Promega Blue/White Cloning-Qualified REs are also assayed in Buffers E and H. Ideally you want to choose a buffer in which each enzyme retains at least 75%. For instance, if you were to perform a double digest with *EcoR* I (optimal in Buffer H) and *Bam*H I (optimal in Buffer E) you would choose in Buffer E because the *Bam*H I has 100% activity and *EcoR* I has 75–100% activity. Both enzymes will maintain acceptable levels of activity in this buffer. Promega has developed an online restriction enzyme compatible buffer search engine available at: www.promega.com/guides/re_guide/ to assist you in finding the right buffer for double digests with all Promega Restriction Enzymes.

A table that describes activity of Promega REs in Promega Restriction Buffers is located on pages 57–58 of this notebook and in the Promega Catalog Appendix.



Double Digests without a Common Buffer

Some enzymes just do not partner well [e.g., double digest with *Pst* I (optimal in Buffer H) and *Spe* I (optimal in Buffer B)]. A review of the tables on pages 57–58 shows that the best-case scenario is provided by Buffer B. *Spe* I of course is optimal in B (100%) but *Pst* I has only 50–75% activity. Three choices are available.

Sequential Method: Perform sequential digests: First digest with *Spe* I in Buffer B, purify DNA, and then perform the *Pst* I digest in Buffer H.

Incubate Longer: Assemble the reaction as usual in Buffer B and incubate 2–4 hours.

Add More Enzyme: Add 1.5–2.0μl of *Pst* I and incubate 1–2 hours.

All three methods work. The first scenario seems intensive, but systems like the Wizard® SV Gel and PCR Clean-Up System make the process very easy (see page 28). The entire reaction can be cleaned and eluted in 15μl of water, and the buffer, enzyme and BSA can be added to bring the reaction to 20μl for the second optimal digest. This is really your only option if both enzymes have no compatibility (i.e., activity in buffer less than 25%).

The second and third methods may provide alternatives to performing sequential digests, depending on the enzymes involved. The second method simply takes more time. The activities in the tables on pages 57–58 are based on a 1-hour incubation. Longer incubation can improve the percent cleavage of the template. This is useful if the two enzymes have a buffer capable of at least 50% activity for both enzymes. The third method is tricky, especially if one of the enzymes is prone to star activity in higher glycerol concentrations. Remember, restriction enzymes are usually stabilized by 50% glycerol so they do not freeze in –20°C storage. Star activity (see page 63) may occur when the digestion glycerol concentration in the reaction rises above 5%. This method is usually only acceptable for two enzymes that have more than 50% activity in the same buffer.

Compare conditions for two Promega REs quickly online. See the RE resource tools at: www.promega.com/techserv/tools

Classic Subcloning

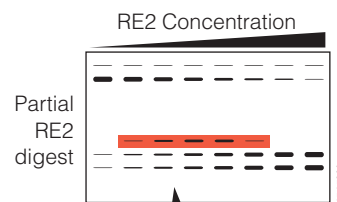
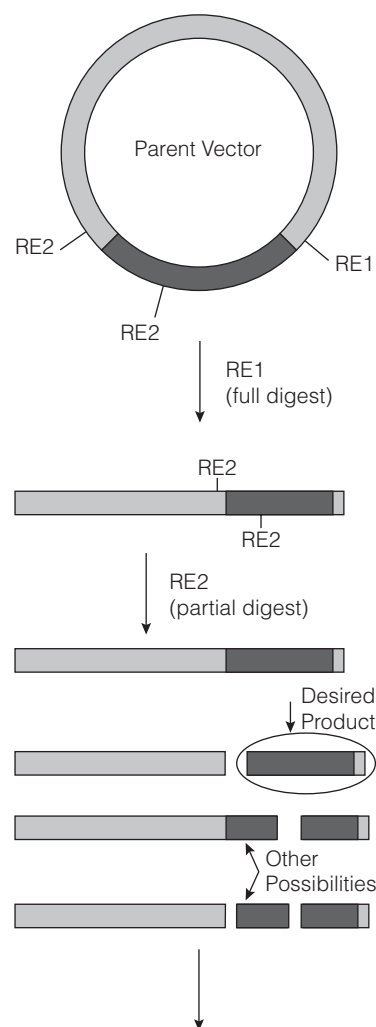
Partial Restriction Digestion

Controlling Cut Frequency in Restriction Digestion

The presence of a restriction recognition site in the insert and the multiple cloning region does not necessarily preclude use of that restriction site in a subcloning strategy. Under normal restriction digest conditions, the enzyme is in excess so that all recognition sites in the plasmid can be cleaved. You can manipulate the restriction digest conditions such that you will digest only a subset of sites. Many strategies have been employed to do partial digests: Decreasing reaction temperature, using a non-optimal buffer, and decreasing units of enzyme. The method presented here uses dilutions of enzyme in the optimal buffer.

A key to doing partial digests is to have a way in which you can differentiate partial digests from complete digests. In other words, you must have a discernable base pair-size difference on the agarose gel so you can cut out the band and perform gel isolation to purify the fragment for ligation into the destination vector. In the following example, the parent vector is first linearized and a partial digest performed on the linearized vector.

1. Digest 10µg of parent vector to completion to linearize (i.e., RE1; 50µl reaction).
2. Purify vector with the Wizard® SV Gel and PCR Clean-Up System directly from the reaction. Elute in 20µl nuclease-free water.
3. On ice, create serial dilutions of RE2 in 1X RE Buffer containing 0.1mg/ml Acetylated BSA (e.g. to yield 5, 2.5, 1.25, 0.625, 0.313, 0.156, 0.078, 0.039u of RE per 18µl of solution).
4. Add 2µl of the purified vector to each tube.
5. Incubate all reactions at 37°C for 30–45 minutes.
6. Add loading dye to each reaction and analyze digests by agarose gel electrophoresis.
7. Identify and cut bands from the gel containing the DNA fragment of interest.
8. Purify insert using the Wizard® SV Gel and PCR Clean-Up System. Elute in 15–20µl nuclease-free water.
9. Proceed to ligation reaction.



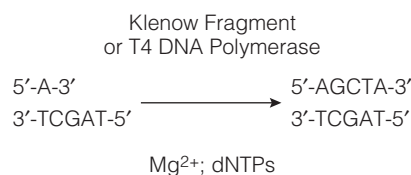
Gel isolate the band you want!

Classic Subcloning

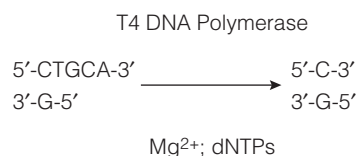
Creating Blunt Ends

Turning an Overhang into a Blunt End

Occasionally you encounter a subcloning application where the choice of restriction sites you can use is limited or where no restriction sites exist in common between vectors and insert. Blunt-ended ligation is an option in these situations. Most vectors contain a blunt cutter like *EcoR* V or *Sma* I in the multiple cloning region, but the parent vector containing your insert may not contain a blunt-cutter site. A blunting reaction can come in handy. Two enzymes are commonly used to generate blunt ends: T4 DNA Polymerase (see page 16) and the Klenow Fragment of DNA Polymerase I (see page 17). The T4 DNA Polymerase is useful for blunting both 5' and 3' overhangs. Klenow works best with 5' overhangs.



5' Overhang Fill-In Reaction



3' Overhang Blunting Reaction

4508MA

Classic Subcloning

Creating Blunt Ends

Blunting a 5' Overhang

T4 DNA Polymerase Method

T4 DNA Polymerase has excellent activity in Promega Restriction Enzyme Buffers B, C, E, and MULTI-CORE™, displaying more than 70% activity. The protocol below is for an integrated blunting reaction following the restriction digestion, and has been tested with the buffers listed above. The following protocol works from a 50µl digestion. The 50µl digestion is recommended to reduce the concentration of glycerol coming from both the restriction enzymes and the T4 DNA Polymerase. Reducing the glycerol concentration prevents potential star activity that may be associated with some restriction enzymes.

1. Digest DNA (0.5–2.0µg) in a 50µl volume.*
2. Add 5u of T4 DNA Polymerase/µg DNA.
3. Add dNTPs to a final concentration of 100µM (e.g., 0.5µl of dNTP Mix [Cat.# U1511]).
4. Incubate at 37°C for 10 minutes.
5. Purify DNA with the Wizard® SV Gel and PCR Clean-Up System direct purification protocol. If both ends of the DNA are being blunted in this reaction, use gel electrophoresis followed by the gel purification protocol to purify the DNA from the enzymes.

*Restriction digest should contain 0.1µg/µl acetylated BSA.

Blunting a 3' Overhang

T4 DNA Polymerase Method

T4 DNA Polymerase has excellent activity in Promega Restriction Enzyme Buffers B, C, E, and MULTI-CORE™, displaying more than 70% activity. The protocol below is for an integrated blunting reaction following the restriction digestion and has been tested with the buffers listed above. The following protocol works from a 50µl digestion. The 50µl digestion is recommended to reduce the concentration of glycerol coming from both the restriction enzymes and the T4 DNA Polymerase. Reducing the glycerol concentration prevents potential star activity that may be associated with some restriction enzymes.

1. Digest DNA (0.5–2.0µg) in a 50µl volume.*
2. Add 5u of T4 DNA Polymerase/µg DNA.
3. Add dNTPs to a final concentration of 100µM (e.g., 0.5µl of dNTP Mix [Cat.# U1511]).
4. Incubate at 37°C for 5 minutes.
5. Purify DNA with the Wizard® SV Gel and PCR Clean-Up System direct purification protocol. If both ends of the DNA are being blunted in this reaction, use gel electrophoresis followed by the gel purification protocol to purify the DNA from the enzymes.

*Restriction digest should contain 0.1µg/µl acetylated BSA.

T4 DNA Polymerase
Cat.# M4211 100u
5–10u/µl
Cat.# M4215 500u
5–10u/µl

See the Product
Information Sheet at:
www.promega.com/tbs

Note: With high concentrations of dNTPs (i.e., 100µM), degradation of the DNA will stop at the duplex DNA. However, if the dNTPs are exhausted, the highly active exonuclease activity (200 times more active than that of DNA polymerase I) of T4 DNA Polymerase will degrade the dsDNA.

Classic Subcloning

Creating Blunt Ends

Blunting a 5' Overhang

Klenow Polymerase Method

Following the restriction enzyme digestion that generated the 5'-protruding ends, purify the DNA from the reaction with a system like the Wizard® SV Gel and PCR Clean-Up System (see page 28 for more information).

1. Assemble the following reaction:

DNA template	1–4µg
10X Klenow Buffer	2µl
Acetylated BSA (10µg/µl)	0.2µl
dNTPs (1mM each)*	0.8µl
Klenow Polymerase	1µl
Nuclease-Free Water	to 20µl

* A 1:10 dilution of the dNTP Mix (Cat.# U1511) in water.

2. Incubate at ambient room temperature for 10 minutes.
3. Purify the DNA from the reaction using the Wizard® SV Gel and PCR Clean-Up System with the direct purification protocol. If both ends of the DNA are being blunted in this reaction, use gel electrophoresis followed by the gel purification protocol.

Note: Promega Restriction Enzyme Buffers A, B, C, D, E, and H may be substituted for the 10X Klenow Buffer, but polymerase activity is 27–43% of the 10X Klenow Buffer.

Note: This method will not work for 3' overhangs.

DNA Polymerase I Large
(Klenow) Fragment

Cat.# M2201 150v
5–10v/µl

Cat.# M2206 500v
5–10v/µl

See the Product
Information Sheet at:
www.promega.com/tbs

Classic Subcloning

Dephosphorylating Vectors to Limit Self-Ligation

Preventing vector self-ligation is critical for reducing subcloning background. The efficiency of ligating the plasmid to itself is far better than ligating a separate piece of DNA into the vector and is the favored reaction. Removing the 5' phosphates of the linearized vector will prevent T4 DNA Ligase from recircularizing the vector. Calf Intestinal Alkaline Phosphatase is the classic enzyme for vector dephosphorylation. The enzyme can be used on 5' recessed ends (i.e., results from an enzyme leaving a 3' overhang), 5' overhangs and blunt-ends. After dephosphorylation, the enzyme must be removed either by direct purification or gel electrophoresis and gel isolation with DNA purification systems like the Wizard® SV Gel and PCR Clean-Up System. Shrimp Alkaline Phosphatase can be used in place of Calf Intestinal Alkaline Phosphatase and offers the advantage of simple heat denaturation to inactivate the enzyme without the need for further purification.

Multiple cloning region of pGL3-Basic Vector

...GGTACCGAGCTCTTACGCGTGTAGCCCGGCTCGAGATCTGTAAGCTTGG...
Kpn I Sac I Mlu I Nhe I Xma I Xho I Bgl II Hind III
Acc65 I Sma I

Hind III
Dephosphorylation

...A OH-AGCTT...
...TTCGA-OH A...

4507MA

Dephosphorylation can reduce the chance of vector self-ligation to virtually zero.

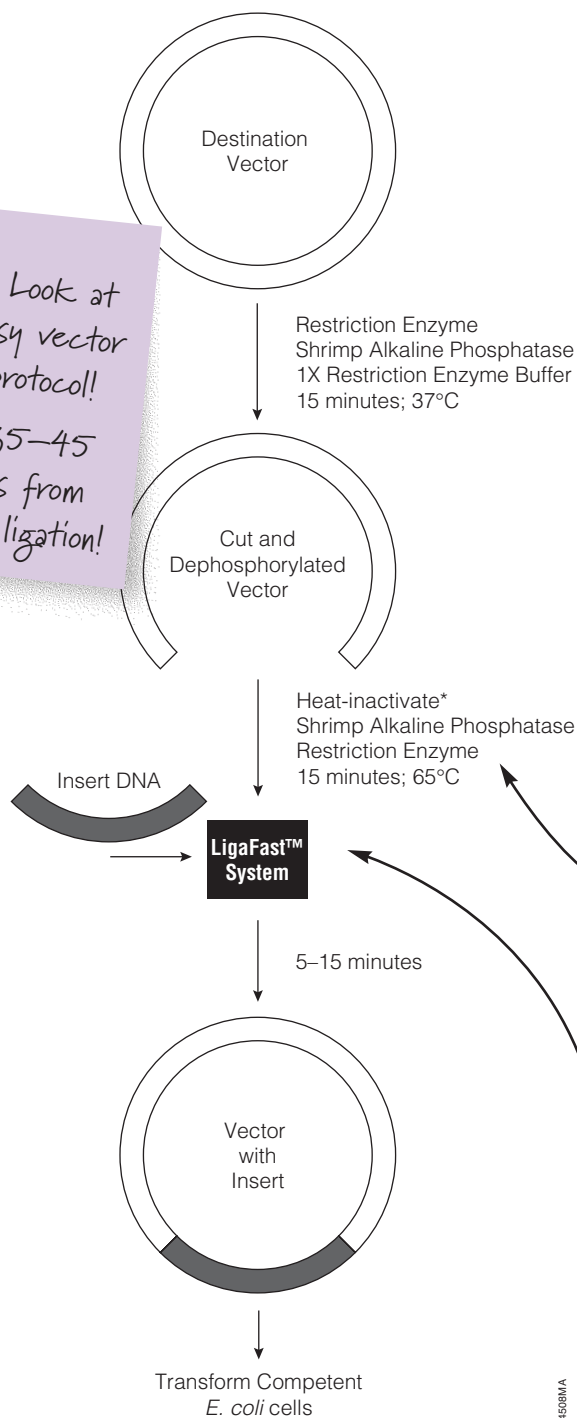
Is it necessary to dephosphorylate linearized vectors before performing the insert ligation?

If the plasmid vector being used was linearized with a single restriction enzyme (generating either a blunt or overhanging end), then dephosphorylation of the vector is a prerequisite to reduce religated vector background. However, if the vector was cut with two different restriction enzymes that leave incompatible ends (this does not include two different enzymes that each leave blunt ends), then dephosphorylation may be omitted. One exception to this is when the selected restriction sites lie close to one another in the vector. In this case, it is still advisable to dephosphorylate the vector, because you cannot be certain from looking at the digested plasmids on the gel if both enzymes cut the plasmid to completion. The presence of a small amount of singly cut plasmid vector in the subsequent ligation reaction can dramatically increase background, which could make it difficult to identify your desired recombinant.

Gel purification of the processed destination vector before ligation ensures that uncut and partially cut vectors are removed from the subcloning reaction.

Classic Subcloning

Dephosphorylating Vectors: Shrimp Alkaline Phosphatase



* Not all restriction enzymes can be heat-inactivated.

Streamlined Restriction Digestion, Dephosphorylation and Ligation Procedure

1. Combine restriction digestion and dephosphorylation of DNA vector in 1X restriction enzyme buffer. Use 15 units of restriction enzyme/ μg vector and 10 units Shrimp Alkaline Phosphatase (SAP)/ μg vector in a final volume of 30–50 μl . Incubate at 37°C for 15 minutes. This is a sufficient amount of SAP to completely dephosphorylate the vector regardless of overhang type (5', 3', or blunt) in any Promega RE buffer.
2. Heat-inactivate both restriction enzyme and SAP for 15 minutes at 65°C.
Note: Not all restriction enzymes can be heat inactivated (see pages 57–58).
3. Centrifuge and remove 1–2 μl of vector for ligation with appropriate DNA insert using T4 DNA Ligase and 2X Rapid Ligation Buffer from LigaFast™ Rapid DNA Ligation System at 15°C for 5 minutes (3' or 5' ends) or 15 minutes for blunt ends in a final reaction volume of 10–50 μl . We recommend starting with a 1:2 molar ratio of vector:insert DNA.
4. Transform the ligated material directly into competent *E. coli* cells.

If your restriction enzyme cannot be heat-inactivated, use the Wizard® SV Gel and PCR Clean-Up System for direct purification. Full purification in just 15 minutes, and you can elute the DNA in as little as 15 μl of water.

Contains the Promega Blue/White Cloning-Qualified T4 DNA Ligase and 2X Rapid Ligation Buffer.

Five-minute ligations for sticky ends; 15-minute ligations for blunt ends.

Wow! Look at this easy vector prep protocol! Only 35–45 minutes from start to ligation!

Classic Subcloning

Dephosphorylating Vectors: Shrimp Alkaline Phosphatase

Dephosphorylation of Purified DNA

1. Purify vector from restriction digest using the Wizard® SV Gel and PCR Clean-Up System.
2. Combine the following:

DNA (1–2µg)	Xµl
10X SAP Buffer	3–5µl
SAP (1u/µl)	1µl/µg DNA
Nuclease-Free Water	to 30–50µl
3. Incubate at 37°C for 15 minutes (works for both 5' and 3' overhangs or blunt ends).
4. Inactivate SAP by heating to 65°C for 15 minutes or purify with the Wizard® SV Gel and PCR Clean-Up System. Proceed to ligation.

SAP Activity in Promega RE Buffers

Buffer	% Activity of SAP
A	20%
B	20%
C	25%
D	35%
E	20%
F	60%
G	30%
H	30%
J	30%
K	20%
L	30%
MULTI-CORE™ Buffer	10%

Using the protocol above with MULTI-CORE™ Buffer in place of SAP Buffer and blunt-ended ligation, greater than 90% of the transformants contained inserts.

Shrimp Alkaline Phosphatase
Cat.# M8201 500u
1u/µl
See the Product
Information Sheet at:
www.promega.com/tbs

This protocol is designed to handle most situations with 5', 3' and blunt ends on the DNA.

Below are the minimal unit requirements for the various ends in 1X SAP Buffer:

5' Overhang: 0.015u SAP/pmol ends

Blunt Overhang: 0.03u SAP/pmol ends

3' Overhang: 0.4u SAP/pmol ends

Classic Subcloning

Dephosphorylating Vectors: Calf Intestinal Alkaline Phosphatase

Dephosphorylation Immediately After Restriction Digestion

1. Add the following components directly to the digested DNA. The CIAP may be diluted on ice in 1X CIAP Buffer immediately before use. Discard any unused, diluted enzyme.

CIAP 10X Reaction Buffer	10µl
CIAP (0.01u/pmol of ends*)	1–2µl
Nuclease-Free Water	to 100µl

*For pmol of ends, simply multiply the pmol of DNA by 2. For example, 1µg of a 1kb DNA fragment will convert to 1.52pmol of DNA and converts to 3pmol of ends.

Note: Dilution of the standard CIAP (1u/µl) is not absolutely necessary, but these are the conditions under which we test the enzyme.

2. Incubate using one of the following conditions, depending on the type of ends present:

5' Overhangs: Incubate for 30 minutes at 37°C. Add another 0.01u CIAP/pmol ends and incubate an additional 30 minutes at 37°C.

3' Overhangs or Blunt Ends: Incubate for 15 minutes at 37°C, then for 15 minutes at 56°C. Add another 0.01u CIAP/pmol ends and repeat incubations at both temperatures.

3. Purify DNA using the Wizard® SV Gel and PCR Clean-Up System and proceed to ligation.

The CIAP Buffer must be added to the reaction for efficient dephosphorylation. The diluted CIAP needs the Zn^{2+} from the buffer to work effectively.

Calculating pmol of DNA from micrograms of DNA.

$$\mu\text{g DNA} \times \frac{\text{pmol}}{660\text{pg}} \times \frac{10^6\text{pg}}{1\mu\text{g}} \times \frac{1}{N} = \text{pmol DNA}$$

N is the number of nucleotides and 660pg/pmol is the average molecular weight of a nucleotide pair.

Online calculators for this equation and many other useful equations are available on the Promega BioMath page: www.promega.com/biomath

Classic Subcloning

Dephosphorylating Vectors: Calf Intestinal Alkaline Phosphatase

Dephosphorylation of Purified DNA

1. Dilute sufficient CIAP for immediate use in CIAP 1X Reaction Buffer to a final concentration of 0.01u/μl. Each pmol of DNA ends will require 0.01u CIAP.

2. Assemble the following reaction:

DNA (up to 10pmol of ends)	40μl
CIAP 10X Reaction Buffer	5μl
diluted CIAP (0.01u/μl)	up to 5μl
Nuclease-Free Water to	50μl

See previous page for calculation of pmol of ends.

Note: Diluting the standard CIAP (1u/μl) is not absolutely necessary, but these are the conditions under which we test the enzyme.

3. Incubate using one of the following conditions, depending on the type of ends present:

5' Overhangs: Incubate for 30 minutes at 37°C, add another 0.01u/pmol of ends of CIAP and repeat incubation.

3' Overhangs or Blunt Ends: Incubate for 15 minutes at 37°C then for 15 minutes at 56°C. Add another 0.01u CIAP/pmol ends and repeat incubations at both temperatures.

4. Purify DNA using the Wizard® SV Gel and PCR Clean-Up System and proceed to ligation.

Alkaline Phosphatase, Calf Intestinal
Cat.# M1821 1,000u
1u/μl
Cat.# M2825 1,000u
20u/μl

See the Product
Information Sheet at:
www.promega.com/tbs

Calf Intestinal Alkaline Phosphatase must be removed prior to the ligation reaction. The Wizard® SV Gel and PCR Clean-Up System can do the purification in 15 minutes, and the dephosphorylated vector can be eluted from the membrane in as little as 15μl of water.

Classic Subcloning

Ligation: Ligating Vector and Insert

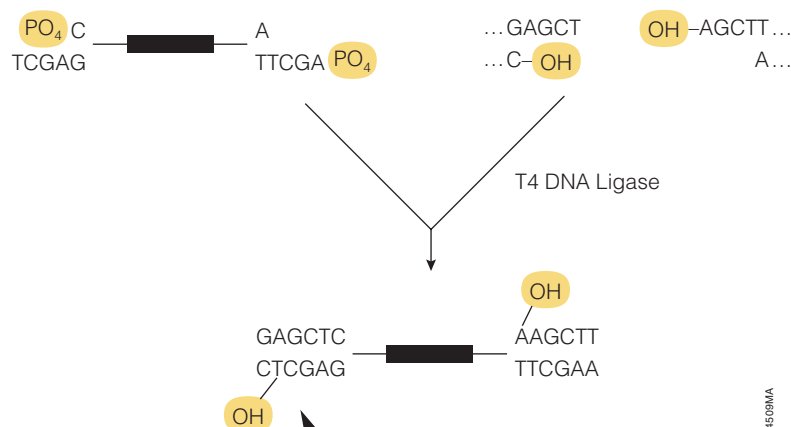
Molecular biologists have exploited DNA ligases to insert pieces of DNA into vectors for decades. The enzyme most commonly used is derived from bacteriophage T4. T4 DNA Ligase is about 400-fold more active than *E. coli* DNA ligase for ligating blunt ends, and thus is the enzyme of choice for all molecular biology requirements. Promega offers **T4 DNA Ligase** in standard or high-concentrate form (see page 25), with the standard Ligase Buffer or with the 2X Rapid Ligation Buffer offered in the **LigaFast™ Rapid DNA Ligation System** (see page 24). The LigaFast™ System allows rapid, 5-minute ligations for 5' or 3' overhang cohesive ends or 15-minute ligations for blunt ends.

How Does DNA Ligase Work?

DNA ligases are responsible for joining gaps that form in DNA during replication, DNA repair and recombination (1). DNA ligases catalyze the formation of a phosphodiester bond between adjacent nucleotides with the concomitant hydrolysis of ATP to AMP and inorganic phosphate. DNA ligases will only form this covalent linkage in a duplex molecule (e.g., at a nick in dsDNA or when joining cohesive- or blunt-ended dsDNAs; 2). The ligation mechanism occurs in three stages. First is the formation of an enzyme-nucleotide intermediate through transfer of an adenylyl group (AMP) from ATP to the ϵ -amine group of a lysine residue in the enzyme. This results in the release of pyrophosphate from ATP. Second, the adenylyl group is transferred from the enzyme to the 5'-phosphate of the DNA, thereby activating it. Third, a phosphodiester bond is formed by nucleophilic attack of the 3'-hydroxyl group of the DNA with concomitant release of AMP.

1. Okazaki, R. *et al.* (1968) *Proc. Natl. Acad. Sci. USA* **59**, 598.

2. Higgins, N.P. and Cozzarelli, R. (1989) In: *Recombinant DNA Methodology* Wu, R., Grossman, L. and Moldave, K., eds. Academic Press, Inc., San Diego, California.



These nicks will be repaired within the host bacteria upon transformation.

Classic Subcloning

Ligation

LigaFast™ Rapid DNA Ligation System

We recommend starting with a 1:2 molar ratio of vector:insert DNA when cloning a fragment into a plasmid vector. The following example illustrates the conversion of molar ratios to mass ratios for a 3.0kb plasmid and a 0.5kb insert DNA fragment

$$\frac{\text{ng of vector} \times \text{kb size of insert}}{\text{kb size of vector}} \times \text{molar ratio of} \frac{\text{insert}}{\text{vector}} = \text{ng of insert}$$

Example:

How much 0.5kb insert DNA should be added to a ligation in which 100ng of 3kb vector will be used? The desired vector:insert ratio will be 1:2.

$$\frac{100\text{ng vector} \times 0.5\text{kb insert}}{3\text{kb vector}} \times \frac{2}{1} = 33.3\text{ng insert}$$

The following ligation reaction of a 3kb vector and a 0.5kb insert DNA uses the 1:2 vector:insert ratio. Typical ligation reactions use 100–200ng of vector DNA.

1. Assemble the following reaction in a sterile microcentrifuge tube:

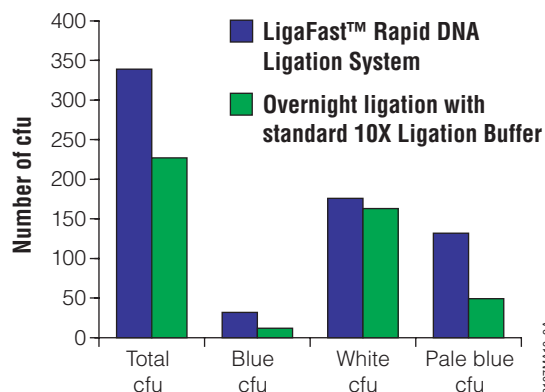
vector DNA	100ng
insert DNA	33ng
2X Rapid Ligation Buffer	5μl
T4 DNA Ligase (3u/μl)	1μl
nuclease-free water to	10μl

2. Incubate the reaction at room temperature for 5 minutes for cohesive-ended ligations, or 15 minutes for blunt-ended ligations.

LigaFast™ Rapid DNA Ligation System
Cat.# M8221 30 reactions
Cat.# M8225 150 reactions

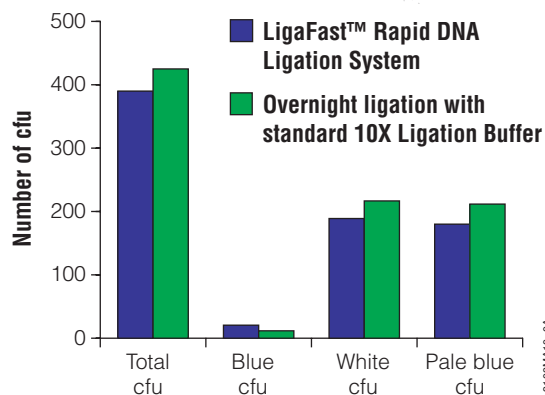
See the Product
Information Sheet at:
www.promega.com/tbs

Blunt-ended
Ligation = 15 minutes



Comparison of overnight ligations and the LigaFast™ Rapid DNA Ligation System using blunt-ended DNA inserts. Experiment performed with blunt-end insert ligated into an *EcoR* V-cut, dephosphorylated pGEM® Vector. Ligations were performed under standard conditions (see pages 24 and 25) using 4°C overnight for the T4 DNA Ligase (3u with standard 10X Ligation Buffer) or 15 minutes at room temperature for the LigaFast™ System. Ligated DNA was transformed into High Competency JM109 cells and plated on indicator media. White and pale blue colonies were confirmed to contain recombinant vector by restriction enzyme analysis.

Sticky-ended
Ligation = 5 minutes



Comparison of overnight ligations and the LigaFast™ Rapid DNA Ligation System using a DNA insert with 5' overhangs. Experiment performed with blunt-end insert ligated into an *Sa*I-cut, dephosphorylated pGEM® Vector. Ligations were performed under standard conditions using 4°C overnight for the T4 DNA Ligase (3u in with standard 10X Buffer) or 5 minutes at room temperature for the LigaFast™ System. Ligated DNA was transformed into High Competency JM109 cells and plated on indicator media. White and pale blue colonies were confirmed to contain recombinant vector by restriction enzyme analysis.

Classic Subcloning

Ligation

T4 DNA Ligase

We recommend using a 1:1, 1:3 or 3:1 molar ratio of vector:insert DNA when cloning a fragment into a plasmid vector.

The following ligation reaction of a 3.0kb vector and a 0.5kb insert DNA uses the 1:3 vector:insert ratio. Typical ligation reactions use 100–200ng of vector DNA.

1. Assemble the following reaction in a sterile microcentrifuge tube:

vector DNA	100ng
insert DNA	50ng
Ligase 10X Buffer	1µl
T4 DNA Ligase (3u/µl)	1µl
Nuclease-Free Water to	10µl

2. Incubate the reaction:

22–25°C	3 hours	Cohesive ends
4°C	Overnight	Cohesive ends
15°C	4–18 hours	Blunt ends

Ligation temperature and duration vary widely in the scientific literature. These are the conditions we use when testing the enzyme.

Ligase Buffers contain ATP to drive the reaction.

Try to avoid multiple freeze-thaw cycles of the buffer.

Dispense the buffer into smaller volumes to minimize the freeze-thaw cycles on each aliquot.

**T4 DNA Ligase
Blue/White Cloning Qualified**

Cat.# M1801	100v 1–3v/µl
Cat.# M1804	500v 1–3v/µl
Cat.# M1794	500v 10–20v/µl

See the Product
Information Sheet at:
www.promega.com/tbs

Standard T4 DNA Ligase methods are more forgiving toward dilute DNA concentrations. Vector and insert can make up 80% of the final volume.

Classic Subcloning

Ligation: Control Reaction

Controls help ensure that everything is functioning normally in your subcloning reaction. If something does go wrong, you can use your controls to figure out where a problem might have occurred.

When ligating insert and vector, you can do a control ligation of vector with no insert. Carry this reaction through transformation and plating. The number of colonies you see can be a good indicator of how a ligation reaction performed and how many background colonies you will have on your plate.

Quick Checks of T4 DNA Ligase

You can always do a quick test of your ligase by simply taking 1 µg of a DNA digest marker (e.g., Lambda DNA *Hind* III Markers [Cat.# G1711]) and performing a 15- to 30-minute ligation reaction under normal conditions. Run the ligation reaction on a gel in comparison to the standard marker. You should see DNA of much higher molecular weight on the gel in comparison to the marker.

Another quick test is to cut a plasmid with a single restriction enzyme. Add this vector to a ligation reaction and transform.

Questions on Subcloning? Call Promega Technical Services

The Promega Worldwide Technical Service Group, Field Applications Specialists, and Distributors are committed to providing you with the highest quality products available to ensure your success. Each of these individuals has an extensive background in molecular biology research, hands-on bench experience with Promega products, and training in problem solving and troubleshooting. Additionally, the full resources of our R&D, Quality Assurance and Production Scientists are available to help increase your laboratory's productivity.

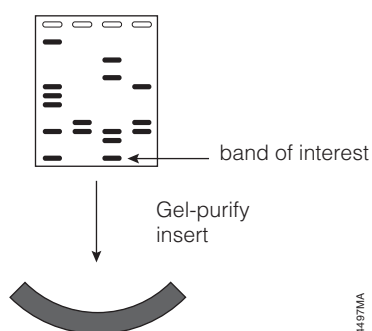
Contact Promega Technical Services directly or through your Branch Office or by email at: **techserv@promega.com**



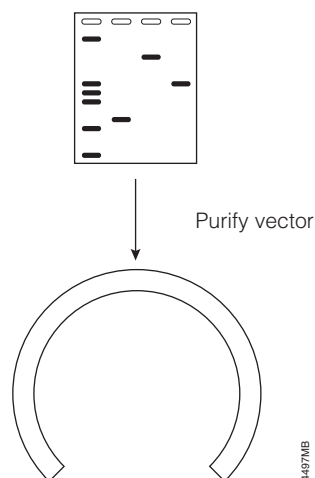
Classic Subcloning

Purifying Vector and Insert

Purification of the insert and destination vector are absolutely critical for success in subcloning applications. Years ago, each step called for phenol:chloroform extractions followed by ethanol precipitation to remove enzymes such as calf intestinal alkaline phosphatase from enzymatic vector manipulations. Guanidine-based nucleic acid clean-up systems greatly simplified the removal of enzymes. Gel isolation methods further improved the efficiency of subcloning by segregating the wanted reactants from the unwanted reactants.



Gel isolation is a practical necessity in subcloning. You get the insert you need.



Gel isolation of vector reduces background by eliminating uncut vector from the ligation.

Classic Subcloning

Purifying Vector and Insert

Wizard® SV Gel and PCR Clean-Up System

The Wizard SV Gel and PCR Clean-Up System is designed to extract and purify DNA fragments directly from PCR^(a) or from agarose gels. Fragments of 100bp to 10kb can be recovered from standard or low-melt agarose gels in either Tris acetate (TAE) buffer or Tris borate buffer (TBE). Up to 95% recovery is achieved, depending upon the DNA fragment size. This membrane-based system, which can bind up to 40µg of DNA, allows recovery of isolated DNA fragments or PCR products in as little as 15 minutes, depending on the number of samples processed and the protocol used. Samples can be eluted in as little as 15µl of nuclease-free water. The purified DNA can be used for automated fluorescent sequencing, cloning, labeling, restriction enzyme digestion or in vitro transcription/ translation without further manipulation.

Process up to 10 gel slices (3.5g total) on a single column with sequential loading.

Capture up to 40µg of DNA on a single column!

From start to purified DNA in 15 minutes!

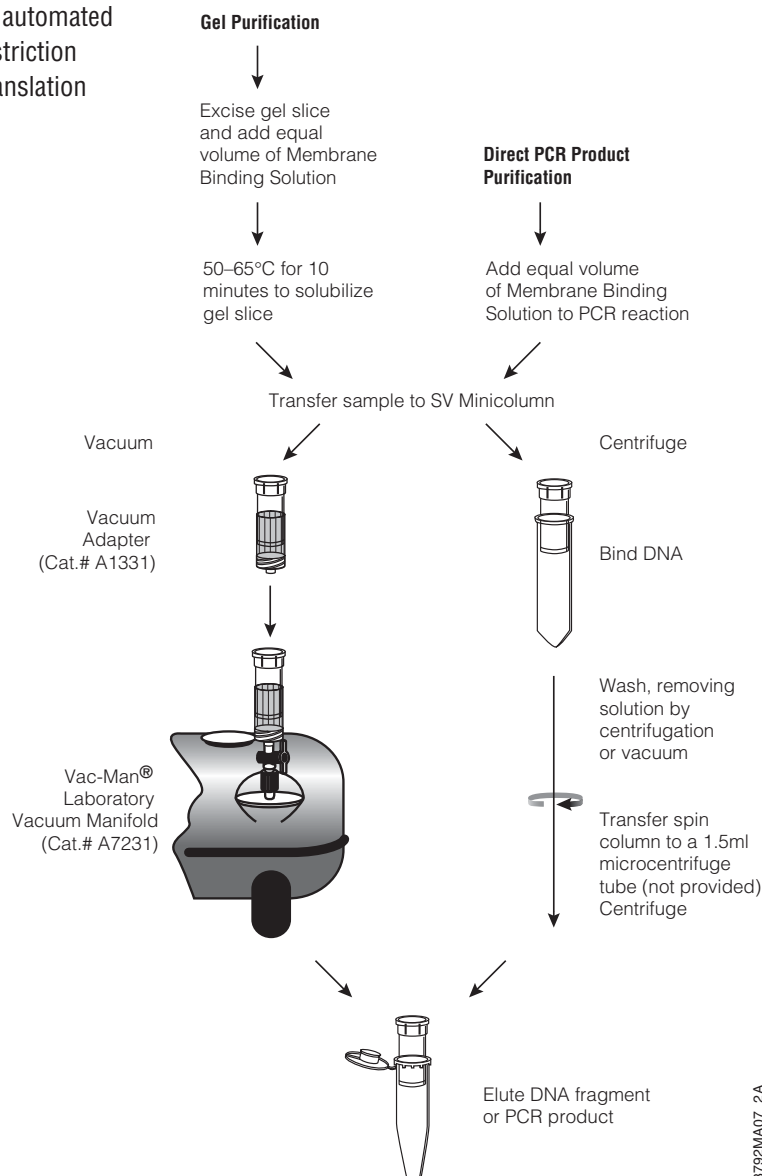
Wizard® SV Gel and PCR Clean-Up System

Cat.# A9281 50 preps

Cat.# A9282 250 preps

Protocol available at:

www.promega.com/tbs/tb308/tb308.html

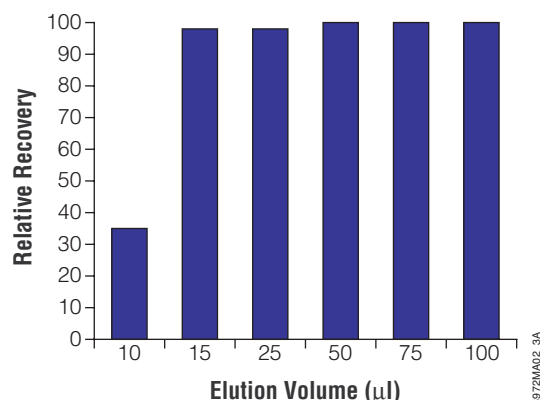


Flow chart of DNA fragment gel purification or direct PCR product purification using the Wizard SV Gel and PCR Clean-Up System.

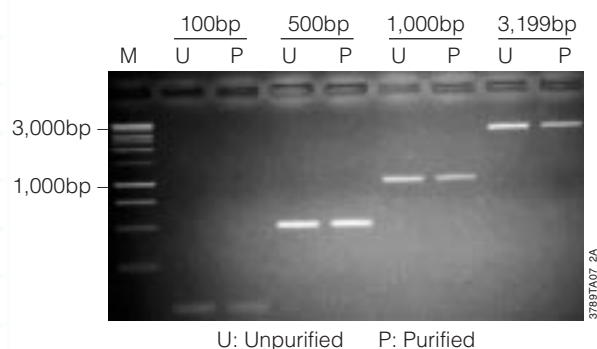
3792MA07_2A

Classic Subcloning

Purifying Vector and Insert



Elution volume versus recovery for a 700bp PCR product. One hundred percent is based on recovery with 50µl elution. Adapted from Table 4 in Betz, N. and Strader, T. (2002) Clean Up with Wizard® SV for Gel and PCR. *Promega Notes* **82**, 2–5.

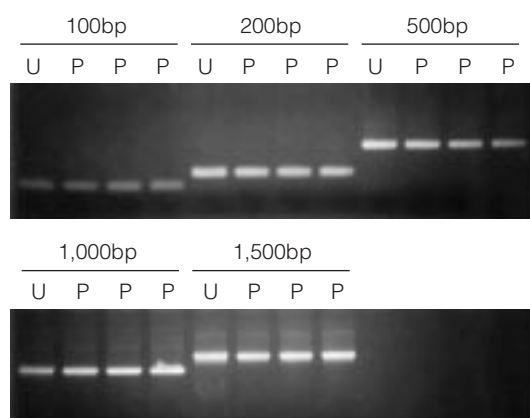


Recovery of various sized unpurified (U) and purified (P) PCR products. Purified lanes were extracted from a 1% agarose gel run with TAE buffer.

Wizard SV Gel and PCR Clean-Up System can remove ethidium bromide and tough enzymes like calf intestinal alkaline phosphatase. See Buross, M. and Betz, N. (2002) Removal of ethidium bromide and calf intestinal alkaline phosphatase using the Wizard SV Gel and PCR Clean-Up System. This can be viewed online at: www.promega.com/enotes/applications/ap0045_tabs.htm

Concentrate DNA by eluting in as little as 15µl.

Linear DNA as big as 10kb can be purified with the system with up to 95% recovery.



Recovery comparison of various sized unpurified (U) and purified (P) PCR products directly purified from PCR amplifications.

Wizard SV Gel and PCR Clean-Up System is tested for purification from up to 3% agarose gels.

Gel Percentages and Resolution of Linear DNA on Agarose Gels.

% Agarose	Resolution
0.8	800bp–10kb+
1.0	400bp–8kb+
1.2	300bp–7kb
1.5	200bp–4kb
2.0	100bp–3kb
3.0	100bp–1kb

Adapted from Brown, T.A. (1998) In: *Molecular Biology LABFAX II: Gene Analysis*. 2nd ed. Academic Press, **90**.

Classic Subcloning

Gel Electrophoresis

Agarose Gel Electrophoresis of DNA

Running double-stranded, linear DNA (like plasmid DNA from restriction enzyme digests) on an agarose gel is a routine activity in molecular biology laboratories. The basic method is very straightforward:

1. Set up the minigel apparatus as recommended by the manufacturer.
2. Weigh the required amount of agarose and add it to the appropriate amount of TAE or TBE 1X Buffer in a flask or bottle. For example, to prepare a 1% agarose gel, add 1.0g of agarose to 100ml of buffer. **Note:** The volume of buffer and agarose should not exceed half the volume of the container.
3. Heat the mixture in a microwave oven or on a hot plate for the minimum time required to allow all the agarose to dissolve. Interrupt the heating at regular intervals and swirl the container to mix the contents. Do not allow the solution to boil over.

CAUTION: The container and contents will be hot! Swirling may also cause solution to boil vigorously. Use adequate precautions.

4. Cool the solution to 50–60°C and pour the gel. Allow the gel to form completely (typically, 30 minutes at room temperature is sufficient). Remove the comb from the gel, place it in the electrophoresis chamber and add a sufficient volume of TAE or TBE 1X buffer to just cover the surface of the gel.
5. Load samples with 1X Blue/Orange Loading Dye into the wells.
6. Connect the gel apparatus to an electrical power supply and apply an appropriate voltage to the gel. For minigels, typical gradients used are between 1–5 volts/cm. Higher voltages and shorter runs will decrease the resolution of the gel and may also cause overheating that may melt the agarose.
7. After electrophoresis is complete, remove the gel and stain it by soaking it in a solution of 0.5µg/ml ethidium bromide for 30 minutes at room temperature. **Note:** Ethidium bromide may also be incorporated in the gel and electrophoresis buffer, at a concentration of 0.5µg/ml, during gel preparation. This eliminates the need for post-electrophoretic staining but may interfere with accurate size determination of DNA fragments. **CAUTION:** Always wear gloves when working with ethidium bromide.

8. Place the gel on a UV lightbox and photograph the gel according to the specification recommended for your camera and film type. **CAUTION:** Use protective eyewear when working with a UV light source.

Note: You may wish to destain or rinse the gel in fresh 1X running buffer prior to viewing it on the UV lightbox.

Recipes

Nearly all of these reagents can be purchased premade including the agarose gels. Here are the directions if you wish to prepare your own reagents.

Blue/Orange Loading Dye, 6X

(available from Promega [Cat.# G1881])

10mM	Tris-HCl, pH 7.5
50mM	EDTA
15%	Ficoll® 400
0.03%	bromophenol blue
0.03%	xylene cyanol FF
0.4%	orange G

One or more dyes can be left out of the recipe to create a custom loading dye.

TAE 50X Buffer (1L)

(Available in a 10X or 40X solution from Promega [Cat.# V4271 and V4281, respectively])

Dissolve 242g Tris base and 37.2g disodium EDTA, dihydrate in 900ml of deionized water. Add 57.1ml glacial acetic acid and adjust the final volume with water to 1 liter. Store at room temperature or 4°C.

TBE 10X Buffer (1L)

(Available in a 10X solution from Promega [Cat.# V4251])

Dissolve 108g of Tris base and 55g boric acid in 900ml deionized water. Add 40ml 0.5M EDTA (pH 8.0) and increase the final volume to 1L. Store at room temperature or 4°C.

Ethidium bromide can detect as little as 1ng of dsDNA in a band.
Brown, T.A. (1998)
In: Molecular Biology LABFAX
II: Gene Analysis. 2nd ed.
Academic Press, 101.

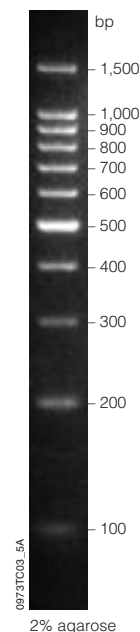
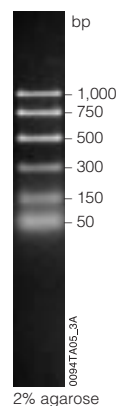
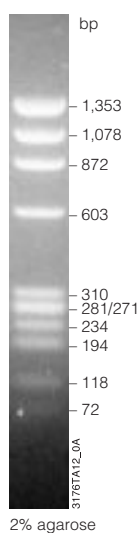
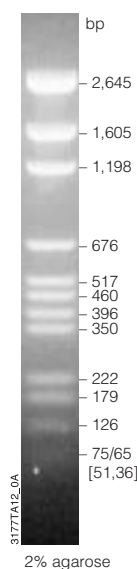
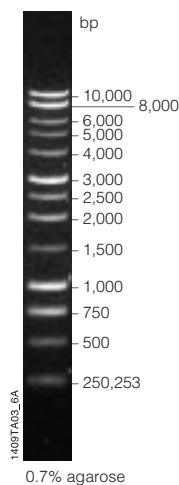
Classic Subcloning

DNA Markers

DNA markers should always be run on agarose gels to aid in identifying bands of interest. This is especially true if you are performing applications such as partial restriction digestion. Promega offers a wide variety of DNA markers to fit your needs. Below is a sampling of marker options available from Promega. BenchTop Markers come premixed with Blue/Orange Loading Dye ready to load onto the gel. As the name implies, you can store them on your benchtop, no need to freeze and thaw every time you need it. Conventional markers are pure DNA solutions and come with a tube of 6X Blue/Orange Loading Dye for use with the marker and your samples.

Each of these markers is available in a ready-to-use BenchTop version or in a conventional version.

1kb DNA Ladder BenchTop Cat.# G7541 Conventional Cat.# G5711	pGEM® DNA Markers BenchTop Cat.# G7521 Conventional Cat.# G1741	ϕ X174 DNA/ Hae III Markers BenchTop Cat.# G7511 Conventional Cat.# G1761	PCR Markers BenchTop Cat.# G7531 Conventional Cat.# G3161	100bp DNA Ladder BenchTop Cat.# G8291 Conventional Cat.# G2101
--	---	---	---	--



Classic Subcloning: Ordering Information

	Enzyme	Heat Inactivated	Buffer	Recognition Site	Size (u)	Conc. (u/μl)	Cat.#
●	<i>Aat</i> II	+	J	GACGT▼C	50	3–5	R6541
					250	3–5	R6545
●	<i>Acc</i> I	–	G	GT▼(A/C)(T/G)AC	100	3–10	R6411
					500	3–10	R6415
	<i>Acc</i> III	–	F	T▼CCGGA	200	10	R6581
●	<i>Acc</i> 65 I (<i>Kpn</i> I)	+	D	G▼GTACC	1,500	10	R6921
	<i>Acc</i> B7 I	+	E	CCANNNN▼NTGG	200	10	R7081
	<i>Age</i> I	+	K	A▼CCGGT	100	3–10	R7251
	<i>Alu</i> I	+	B	AG▼CT	500	10	R6281
	<i>Alw</i> 26 I	+	C	GTCTC(N) ₅ ▼	100	8–12	R6761
				GTCTC(N) ₅	500	8–12	R6765
	<i>Alw</i> 44 I	+	C	G▼TGAC	1,000	10	R6771
●	<i>Apa</i> I	+	A	GGGCC▼C	5,000	10	R6361
					25,000	40–80	R4364
●	<i>Ava</i> I	+/-	B	C▼(T/C)CG(A/G)G	200	8–12	R6091
					1,000	8–12	R6095
	<i>Ava</i> II (<i>Sin</i> I)	+	C	G▼G(A/T)CC	100	1–10	R6131
					1,000	1–10	R6135
	<i>Bal</i> I	+	G	TGG▼CCA	50	2–10	R6691
					250	2–10	R6695
●	<i>Bam</i> HI	+	E	G▼GATCC	2,500	10	R6021
					12,500	10	R6025
					12,500	40–80	R4024
					50,000	40–80	R4027
	<i>Ban</i> I	–	G	G▼G(T/C)(A/G)CC	200	8–12	R6891
	<i>Ban</i> II	+	E	G(A/G)GC(T/C)▼C	1,000	8–12	R6561
	<i>Bbu</i> I (<i>Sph</i> I)	+	A	GCA TG▼C	200	10	R6621
					1,000	40–80	R4624
●	<i>Bcl</i> I	–	C	T▼GATCA	1,000	10	R6651
					5,000	40–80	R4654
●	<i>Bgl</i> I	+	D	GCCNNNN▼NGGC	1,000	10	R6071
					5,000	10	R6077
					5,000	40–80	R4074
	<i>Bgl</i> II	–	D	A▼GATCT	500	10	R6081
					2,500	10	R6085
					10,000	10	R6087
					2,500	40–80	R4084
	<i>Bsa</i> MI	–	D	GAATGCN▼	500	10	R6991
	<i>Bsp</i> 1286 I	+	A	G(G/A/T)GC(C/A/T)▼C	500	10	R6741
	<i>Bsr</i> S I	–	D	ACTGGN	500	10	R7241
●	<i>Bss</i> HI	–	H	G▼CGCGC	100	10	R6831
					500	10	R6835
	<i>Bst</i> 98 I	–	D	C▼TTAAG	500	8–12	R7141
	<i>Bst</i> E II	–	D	G▼GTNACC	2,000	10	R6641
	<i>Bst</i> O I	–	C	CC▼(A/T)GG	2,000	10	R6931
●	<i>Bst</i> X I	+/-	D	CCANNNNN▼NTGG	250	8–12	R6471
					1,000	8–12	R6475
	<i>Bst</i> Z I	–	D	C▼GGCCG	500	10	R6881
●	<i>Bsu</i> 36 I	–	E	CC▼TNAGG	500	10	R6821
	<i>Cfo</i> I (<i>Hha</i> I)	+/-	B	GCG▼C	3,000	10	R6241
●	<i>Cla</i> I	+	C	AT▼CGAT	500	10	R6551
					2,500	10	R6555
●	<i>Csp</i> I	+	K	CG▼G(A/T)CCG	100	10	R6671
					500	10	R6675




















	Enzyme	Heat Inactivated	Buffer	Recognition Site	Size (u)	Conc. (u/μl)	Cat.#
●	<i>Csp</i> 45 I	+	B	TT▼CGAA	2,500	10	R6571
	<i>Dde</i> I	+/-	D	C▼TNAG	200	10	R6291
					1,000	10	R6295
	<i>Dpn</i> I (<i>Sau</i> 3A I)	+	B	G ^{me} A▼TC	200	10	R6231
	<i>Dra</i> I	+	B	TTT▼AAA	2,000	10	R6271
	<i>Ecl</i> HK I	+	E	GACNNN▼NNGTC	100	10	R7111
●	<i>Eco</i> 47 III	+	D	AGC▼GCT	50	2–5	R6731
●	<i>Eco</i> 52 I (<i>Bst</i> Z I)	+	L	C▼GGCCG	50	1–5	R6751
●	<i>Eco</i> CR I (<i>Sac</i> I)	+	B	GAG▼CTC	1,000	10	R6951
					5,000	40–80	R4954
●	<i>Eco</i> R I	+	H	G▼AATTC	5,000	12	R6011
					15,000	12	R6017
					25,000	40–80	R4014
					50,000	40–80	R4017
●	<i>Eco</i> R V	+	D	GAT▼ATC	2,000	10	R6351
					10,000	10	R6355
					10,000	40–80	R4354
	<i>Fok</i> I	+	B	GGATG(N) ₉ GGATG(N) ₍₁₃₎ ▼	100	2–10	R6781
	<i>Hae</i> II	–	B	(A/G)GCGC▼(T/C)	1,000	10	R6661
	<i>Hae</i> III	–	C	GG▼CC	2,500	10	R6171
					10,000	10	R6175
					12,500	40–80	R4174
	<i>Hha</i> I (<i>Cfo</i> I)	+	C	GCG▼C	1,000	10	R6441
●	<i>Hinc</i> II	+	B	GT(T/C)▼(A/G)AC	200	10	R6031
					1,000	10	R6035
					5,000	10	R6037
					1,000	40–80	R4034
●	<i>Hind</i> III	+	E	A▼AGCTT	5,000	10	R6041
					15,000	10	R6045
					25,000	40–80	R4044
					50,000	40–80	R4047
	<i>Hinf</i> I	–	B	G▼ANTC	1,000	10	R6201
					5,000	10	R6205
					5,000	40–80	R4204
	<i>Hpa</i> I	–	J	GTT▼AAC	100	3–10	R6301
					500	3–10	R6305
	<i>Hpa</i> II (<i>Msp</i> I)	–	A	C▼CGG	1,000	10	R6311
					5,000	10	R6315
	<i>Hsp</i> 92 I	+	F	G(A/G)▼CG(T/C)C	500	10	R7151
	<i>Hsp</i> 92 II	+	K	CATG▼	1,000	10	R7161
	I- <i>Ppo</i> I (Intron-Encoded Endonuclease)	+	10X	CTCTCTTAA▼GGTAGC I- <i>Ppo</i> I	10,000	100–200	R7031
●	<i>Kpn</i> I ^(b) (<i>Acc</i> 65 I)	+/-	J	GGTAC▼C	2,500	8–12	R6341
					10,000	8–12	R6345
					12,500	40–80	R4344
	<i>Mbo</i> II	+	B	GAAGA(N) ₈ GAAGA(N) ₇ ▲	100	2–10	R6723
	<i>Mlu</i> I	+/-	D	A▼CGCGT	1,000	10	R6381
●	<i>Msp</i> I (<i>Hpa</i> II)	+	B	C▼CGG	2,000	10	R6401
					10,000	10	R6405
					10,000	40–80	R4404
	<i>Msp</i> A1 I	+	C	C(A/C)G▼C(G/T)G	1,000	10	R7021
	Turbo™ <i>Nae</i> I ^(c)	+	Turbo™	GCC▼GGC	250	4	R7231

● Indicates Genome Qualified.

● Indicates Blue/White Cloning Qualified.







Promega Subcloning Notebook

Classic Subcloning: Ordering Information

Enzyme	Heat Inactivated	Buffer	Recognition Site	Size (u)	Conc. (u/μl)	Cat.#
<i>Nae</i> I (<i>Ngo</i> M IV)	+	A	GCC▼GGC	250	4	R7131
				1,000	4	R7135
 Turbo™ <i>Nar</i> I(c)	+	Turbo™	GG▼CGCC	200	10	R7261
 <i>Nar</i> I	+	G	GG▼CGCC	200	10	R6861
<i>Nci</i> I	+	B	CC▼(C/G)GG	1,000	10	R7061
 <i>Nco</i> I	+	D	C▼CATGG	200	10	R6513
				1,000	10	R6515
<i>Nde</i> I	+	D	CA▼TATG	500	10	R6801
<i>Nde</i> II	+	D	▼GATC	200	10	R7291
(<i>Dpn</i> I, <i>Sau</i> 3A I)				1,000	10	R7295
<i>Ngo</i> M IV (<i>Nae</i> I)	+	MULTI-CORE™	G▼CCGGC	500	10	R7171
 <i>Nhe</i> I	+	B	G▼CTAGC	250	10	R6501
				1,250	10	R6505
  <i>Not</i> I	+	D	GC▼GGCCGC	200	10	R6431
				1,000	10	R6435
				1,000	40–80	R4434
 <i>Nru</i> I	+	K	TCG▼CGA	200	10	R7091
 <i>Nsi</i> I	+/-	D	ATGCA▼T	250	10	R6531
 <i>Pst</i> I	+/-	H	CTGCA▼G	3,000	10	R6111
				15,000	10	R6115
				15,000	40–80	R4114
				50,000	40–80	R4117
<i>Pvu</i> I	–	D	CGAT▼CG	100	2–10	R6321
				500	2–10	R6325
<i>Pvu</i> II	+	B	CAG▼CTG	1,000	8–12	R6331
				5,000	8–12	R6335
<i>Rsa</i> I	+	C	GT▼AC	1,000	10	R6371
				5,000	40–80	R4374
 <i>Sac</i> I (<i>Eco</i> CR I)	+	J	GAGCT▼C	1,000	10	R6061
				5,000	10	R6065
				5,000	40–80	R4064
 <i>Sac</i> II	+	C	CCGC▼GG	500	10	R6221
  <i>Sal</i> I	+	D	G▼TCGAC	2,000	10	R6051
				10,000	10	R6055
				10,000	40–80	R4054
<i>Sau</i> 3A I (<i>Dpn</i> I, <i>Nde</i> II)	+	B	▼GATC	100	3–10	R6191
				500	3–10	R6195
<i>Sca</i> I	+	K	AGT▼ACT	1,000	8–12	R6211
				5,000	40–80	R4214
  <i>Sfi</i> I(d)	+	B	GGCCNNNN▼NGGCC	250	10	R6391
				1,250	40–80	R4394
 <i>Sgf</i> I	+/-	C	GCGAT▼CGC	250	8–12	R7103
				1,250	40–80	R5104
<i>Sin</i> I (<i>Ava</i> II)	+	A	G▼G(A/T)CC	200	8–12	R6141
				1,000	40–80	R4144
  <i>Sma</i> I (<i>Xma</i> I)	+	J	CCC▼GGG	1,000	8–12	R6121
				5,000	8–12	R6125
				5,000	40–80	R4124
<i>Sna</i> B I	–	B	TAC▼GTA	100	2–10	R6791
				500	2–10	R6795
 <i>Spe</i> I	+	B	A▼CTAGT	200	10	R6591
				1,000	10	R6595

 Indicates Genome Qualified.

 Indicates Blue/White Cloning Qualified.

Enzyme	Heat Inactivated	Buffer	Recognition Site	Size (u)	Conc. (u/μl)	Cat.#
 <i>Sph</i> I (<i>Bbu</i> I)	+	K	GCATG▼C	200	10	R6261
				1,000	10	R6265
<i>Ssp</i> I	+	E	AAT▼ATT	500	10	R6601
				2,500	40–80	R4604
<i>Stu</i> I	+	B	AGG▼CCT	400	10	R6421
<i>Sty</i> I	+	F	C▼C(A/T)(T/A)GG	2,000	10	R6481
<i>Taq</i> I	–s	E	T▼CGA	1,000	10	R6151
				10,000	10	R6155
				5,000	40–80	R4154
<i>Tru</i> 9 I	–	F	T▼TAA	200	8–12	R7011
 <i>Tth</i> 111 I	–	B	GACN▼NNGTC	500	8–12	R6841
 <i>Vsp</i> I	+	D	AT▼TAAT	500	8–12	R6851
  <i>Xba</i> I	–	D	T▼CTAGA	2,000	8–12	R6181
				10,000	8–12	R6185
				10,000	40–80	R4184
<i>Xho</i> I	+	D	C▼TCGAG	3,000	10	R6161
				10,000	10	R6165
				15,000	40–80	R4164
 <i>Xho</i> II	+	C	(A/G)▼GATC(T/C)	100	5–10	R6811
				500	5–10	R6815
<i>Xma</i> I (<i>Sma</i> I)	+	B	C▼CCGGG	50	1–5	R6491
				250	1–5	R6495
<i>Xmn</i> I	+	B	GAANN▼NNTTC	500	10	R7271
				2,500	10	R7273

Product	Size	Conc. (mg/ml)	Cat.#
BSA, (Bovine Serum Albumin) Acetylated	400μl	1	R9461
	1ml	10	R3961
MULTI-CORE™ Buffer Pack	3 × 1ml	—	R9991
4-CORE® Buffer Pack (1 each A–D)	4 × 1ml	—	R9921

For Laboratory Use.

Turbo™ Enzymes are provided with a reaction buffer containing a noncleavable affector sequence that facilitates efficient digestion of slow and resistant sites.

Restriction enzymes are shown to be heat inactivated (+) if they show >95% loss of activity after a 15 minute incubation at 65°C.

Enzymes followed by another enzyme name in parentheses indicate that the enzyme is an isoschizomer or neoschizomer of the enzyme in parentheses.

Restriction Enzyme Buffer Composition (1X).

Buffer	pH (at 37°C)	Tris-HCl (mM)	MgCl ₂ (mM)	NaCl (mM)	KCl (mM)	DTT (mM)
A	7.5	6	6	6	–	1
B	7.5	6	6	50	–	1
C	7.9	10	10	50	–	1
D	7.9	6	6	150	–	1
E	7.5	6	6	100	–	1
F	8.5	10	10	100	–	1
G	8.2	50	5	–	–	–
H	7.5	90	10	50	–	–
J	7.5	10	7	–	50	1
K	7.4	10	10	–	150	–
L	9.0	10	3	100	–	–

MULTI-CORE™ Buffer (1X): 25mM Tris-Acetate (pH 7.8 @ 25°C), 100mM potassium acetate, 10mM magnesium acetate, 1mM DTT.

1. For each 10°C rise in temperature between 0°C and 25°C, the pH of Tris buffers decreases 0.31 pH units.

2. For each 10°C rise in temperature between 25°C and 37°C, the pH of Tris buffers decreases 0.25 pH units.

Classic Subcloning: Ordering Information

Enzymes	Size	Conc.	Cat. #
T4 DNA Polymerase ^(d)	100u	5–10u/μl	M4211
	500u	5–10u/μl	M4215
DNA Polymerase I Large (Klenow) Fragments	150u	5–10u/μl	M2201
	500u	5–10u/μl	M2206
Shrimp Alkaline Phosphatase	500u	1u/μl	M8201
Alkaline Phosphatase, Calf Intestinal	1,000u	1u/μl	M1821
	1,000u	20u/μl	M2825
LigaFast™ Rapid DNA Ligation System	30 reactions	—	M8221
	150 reactions	—	M8225
T4 DNA Ligase	100u	1–3u/μl	M1801
	500u	1–3u/μl	M1804
	500u	10–20u/μl	M1794

For Laboratory Use.

Purification Systems	Size	Cat. #
Wizard® SV Gel and PCR Clean-Up System* (ready for spin protocol)	50 preps	A9281
	250 preps	A9282
Vac-Man® Laboratory Vacuum Manifold, 20-sample capacity (required for vacuum protocol)	1	A7231
Vacuum Adapters (required for vacuum protocol)	20	A1331

*For Laboratory Use.

Ready-to-Load BenchTop DNA Markers	Size	Cat. #
BenchTop 100bp DNA Ladder	50 lanes	G8291
BenchTop 1kb DNA Ladder	100 lanes	G7541
BenchTop PCR Markers	50 lanes	G7531
BenchTop pGEM® DNA Markers	50 lanes	G7521
BenchTop φX174 DNA/Hae III Markers	50 lanes	G7511

For Laboratory Use.

Conventional DNA Markers (supplied with 6X Blue/Orange Loading Dye)	Size	Cat. #
100bp DNA Ladder	50 lanes	G2101
1kb DNA Ladder	100 lanes	G5711
PCR Markers	50 lanes	G3161
pGEM® DNA Markers	50 lanes	G1741
φX174 DNA/Hae III Markers	50 lanes	G1761

For Laboratory Use.

Accessory Items	Size	Conc.	Cat. #
4-CORE® Buffer Pack*	4 × 1ml	—	R9921
MULTI-CORE™ Buffer Pack*	3 × 1ml	—	R9991
Bovine Serum Albumin, Acetylated*	400μl	1μg/μl	R9461
	1,000μl	10mg/ml	R3961
T4 DNA Ligase Buffer Pack*	3 × 500μl	—	C1263
CIAP Buffer Pack*	3 × 500μl	—	M1833
dNTP Mix*	200μl	10mM	U1511
	1,000μl	10mM	U1515
Agarose, LE, Analytical Grade	100g	—	V3121
	500g	—	V3125
Blue/Orange Loading Dye, 6X*	3 × 1ml	—	G1881
TAE Buffer, 10X	1,000ml	10X	V4271
TAE Buffer, 40X	1,000ml	40X	V4281
TBE Buffer, 10X	1,000ml	10X	V4251
Ethidium Bromide Solution, Molecular Grade	10ml	10mg/ml	H5041
Mineral Oil*	12ml	—	DY1151

*For Laboratory Use.

PCR Subcloning

Introduction

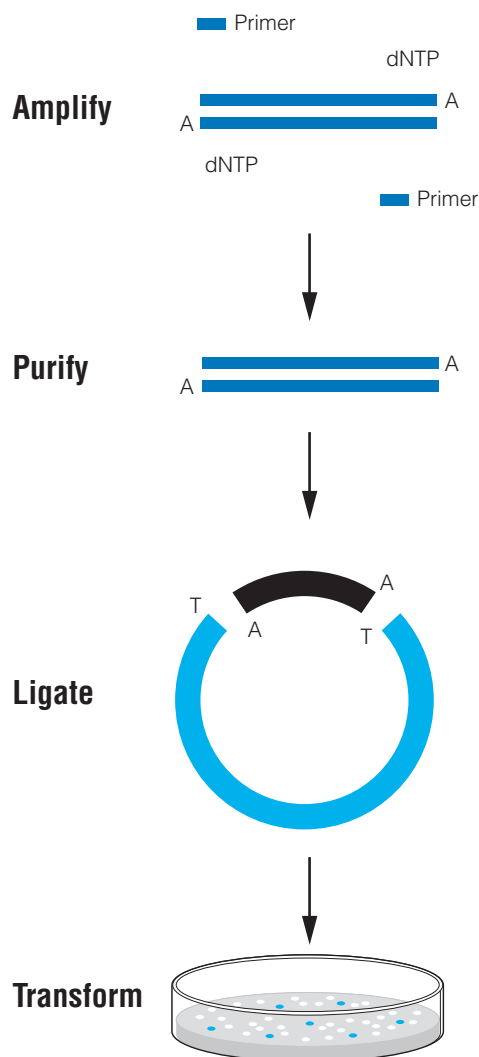
You may wish to subclone your PCR product into a plasmid cloning vector. When PCR was in its infancy, researchers found that subcloning PCR products by simple blunt-ended ligation into blunt-ended plasmid cloning vectors was not easy. Thermostable DNA polymerases, like *Taq* DNA polymerase, add a single nucleotide base extension to the 3' end of blunt DNA in a template-independent fashion (1,2). These polymerases usually add an adenine, leaving an "A overhang."

Historically, researchers have used several approaches to overcome the cloning difficulties presented by the presence of A overhangs on PCR products. One method involves treating the product with the Klenow fragment of *E. coli* DNA Polymerase I to create a blunt-ended fragment for subcloning. However this technique is not particularly efficient.

Another method commonly used by researchers is to add restriction enzyme recognition sites to the ends of the PCR primers (3). The PCR product is then digested and subcloned into the desired plasmid cloning vector in a desired orientation. Care must be exercised in primer design when using this method, as not all REs cleave efficiently at the ends of DNA, and you may not be able to use every RE you desire (4). Some REs require extra bases outside the recognition site (see page 40), adding further expense to the PCR primers as well as risk of priming to unrelated sequences in the genome.

A method of choice for cloning PCR products is T-Vector cloning. In essence, the plasmid cloning vector is treated to contain a 3' T overhang to match the 3' A overhang of the amplicon (5). The A-tailed amplicon is directly ligated to the T-tailed plasmid vector with no need for further enzymatic treatment of the amplicon other than the action of T4 DNA ligase. Promega has systems based on this technology for routine subcloning, and direct mammalian expression.

4 Simple Steps to Success



References

1. Clark, J.M. (1988) Novel non-template nucleotide addition reactions catalyzed by procaryotic and eucaryotic DNA polymerases. *Nucl. Acids Res.* **16**, 9677–86.
2. Mole, S.E., Iggo, R.D. and Lane, D.P. (1989) Using the polymerase chain reaction to modify expression plasmids for epitope mapping. *Nucl. Acids Res.* **17**, 3319.
3. Scharf, S.J., Horn, G.T. and Erlich, H.A. (1986) Direct cloning and sequence analysis of enzymatically amplified genomic sequences. *Science* **233**, 1076–8.
4. Kaufman, D.L. and Evans, G.A. (1990) Restriction endonuclease cleavage at the termini of PCR products. *BioTechniques* **9**, 304–6.
5. Mezei, L.M. and Storts, D.R. (1994) Cloning PCR Products. In: *PCR Technology Current Innovations*. Griffin, H.G. and Griffin, A.M. (eds). CRC Press, 21–7.

PCR Subcloning

T-Vector Systems

pGEM®-T and pGEM®-T Easy Vector Systems

The most basic need in PCR subcloning is a simple, general cloning vector. The pGEM-T and pGEM-T Easy Vector Systems^(e,f,g) are designed for just that purpose. The vectors are based on the pGEM-5Zf(+) Vector^(g) backbone. Each provide convenient T7 and SP6 promoters to serve as sequencing primer binding sites or for in vitro transcription of either strand of the insert with the appropriate RNA polymerase. The vectors have the *lacZ* α , allowing easy blue/white screening of the inserts with an appropriate bacterial strain (e.g., JM109, DH5 α [™], XL1 Blue, etc). To speed your research, these vectors are provided with 2X Rapid Ligation Buffer, allowing efficient ligation in just 1 hour with the supplied T4 DNA Ligase. You can either supply your own favorite *E. coli* cells or purchase the system with Promega JM109 Competent Cells. The choice is yours.

pGEM®-T Vector System I
(you supply competent cells)

Cat.# A3600

20 reactions

pGEM®-T Vector System II

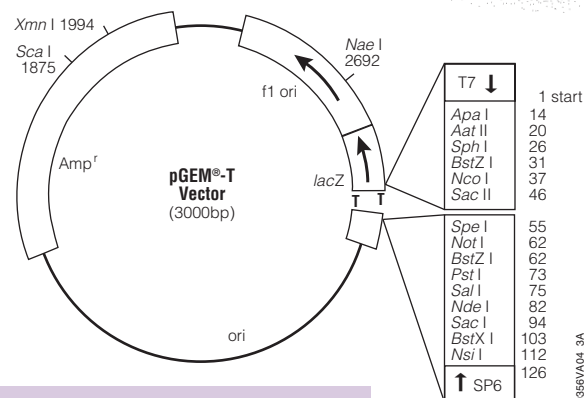
(supplied with High Efficiency JM109 Competent Cells)

Cat.# A3610

20 reactions

Protocol available at:

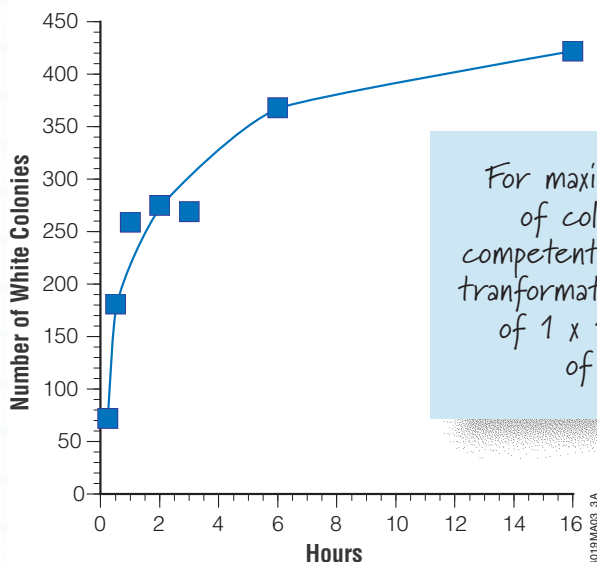
www.promega.com/tbs/tm042/tm042.html



Select recombinants by blue/white selection.

Sequence inserts with the following:

- SP6 Promoter Primer
- T7 Promoter Primer
- M13 Forward Primer
- M13 Reverse Primer

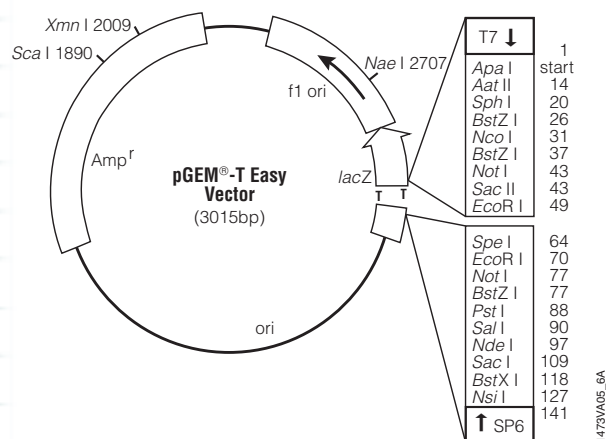


Number of white colonies (transformants) versus time of ligation. Control pGEM-T Easy ligation reactions were set up at room temperature (24°C) and allowed to proceed from 0.25 to 16 hours using the methods described in Technical Manual #TM042. Graph was adapted from Table 2 in Frackman, S. and Kephart, D. (1999) Rapid ligation for the pGEM-T and pGEM-T Easy Vector Systems. *Promega Notes* 71, 8–10.

Drop out insert with a single Bst Z I digest.

PCR Subcloning

T-Vector Systems



pGEM®-T Easy Vector System I
(you supply competent cells)
Cat.# A1360 20 reactions

pGEM®-T Easy Vector System II
(supplied with High Efficiency JM109 Competent Cells)
Cat.# A1380 20 reactions

Protocol available at:
www.promega.com/tbs/tm042/tm042.html

Sequence inserts with the following:

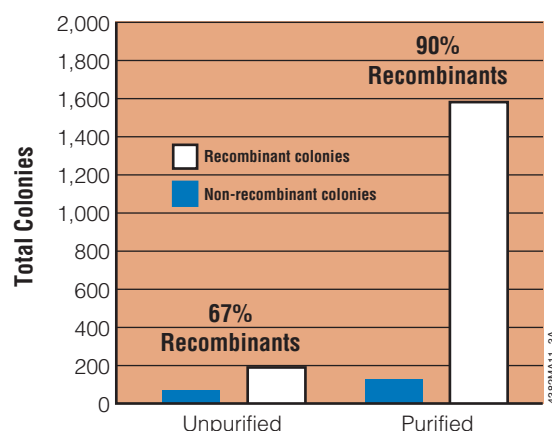
- SP6 Promoter Primer
- T7 Promoter Primer
- M13 Forward Primer
- M13 Reverse Primer

Drop out insert with a single BstZ I, EcoR I or Not I digest.

For maximum subcloning efficiency, purify the PCR product before subcloning. The presence of PCR primers and primer dimers can reduce subcloning efficiency.

Purification achieved with Wizard® SV Gel and PCR Clean-Up System.

For more information, see page 28.

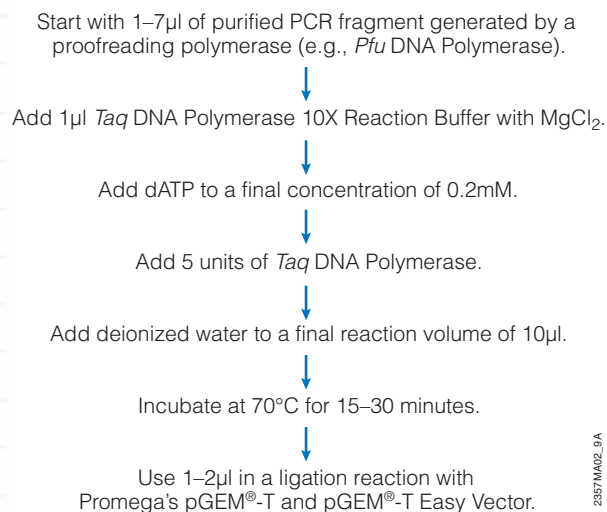


Cloning efficiency of a purified PCR product and an unpurified PCR product.

PCR Subcloning

Giving Blunt-Ended DNA an A-tail for T-Vector Subcloning

PCR amplicons generated with proofreading polymerases like *Pfu* or *Tli* DNA Polymerase are blunt-ended. Promega has developed an easy method to add an A-Tail to the DNA so that it can be used for T-Vector cloning.



Full details of the protocol are available in the *pGEM®-T* and *pGEM®-T Easy Vector Systems Technical Manual*, TM042. The proofreading enzyme must be removed using a system like the Wizard® SV Gel and PCR Clean-Up System prior to the A-tailing procedure. Any remaining proofreading enzyme in the PCR will remove the A-overhangs created by the *Taq* DNA polymerase.

Ends Left on PCR Products by Thermostable Enzymes.

Polymerase	Type of End*
<i>Taq</i> DNA Polymerase	3' A overhang
GoTaq® DNA Polymerase	3' A overhang
<i>Tfi</i> DNA Polymerase	3' A overhang
<i>Tth</i> DNA Polymerase	3' A overhang
<i>Pfu</i> DNA Polymerase	Blunt end
<i>Tli</i> DNA Polymerase	Blunt end
Long PCR mixes	Blunt end
Other Proofreading Polymerases	Blunt end

*All bases may be found at 3' overhang; adenine tends to be encountered most often.

For more information and techniques for cloning PCR DNA, check out Promega Frequently Asked Questions for the T-Vector cloning systems at: www.promega.com/faq

Promega recommends a guanidine-based purification method, like the Wizard SV Gel and PCR Clean-Up System, to remove proofreading polymerases.

PCR Subcloning

Subcloning with RE Sites

What PCR Cloning Controls Can Do for You

Each Promega PCR cloning system is provided with a control insert. The ligation and subsequent transformation of this positive control can give you a lot of information with regard to the ligation and transformation of your insert.

Typical Results

	Efficiency*	% White
Control insert	1110	92%
Control insert	1125	92%
No insert	92	—
No insert	109	—

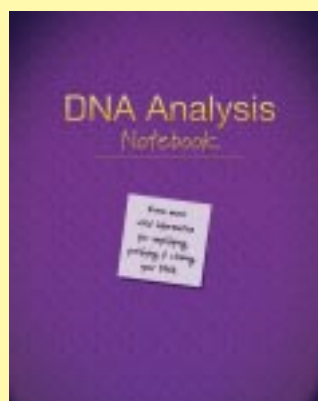
*cfu/ng control insert DNA; JM109 cells at 1.5×10^8 cfu/ μ g; pGEM®-T Easy Vector System II using room temperature ligation for 1 hour.

The total number of blue colonies obtained with positive control insert and no-insert controls should be approximately equal. The negative control may have some white colonies as well.

Need more information about PCR template preparation, PCR, PCR clean-up and PCR cloning? Request the DNA Analysis Notebook

Literature# BR129

www.promega.com/guides/dna_guide/default.htm



Interpreting Results from T-Cloning

Experimental insert looks like control insert in efficiency and percent white colonies.

Successful experiment. Greater than 80% of the white colonies should contain inserts

Experimental insert and control insert look like negative control.

Ligation has failed. Avoid multiple freeze/thaws of the ligation buffer. Ligase buffer contains ATP and could be damaged by freeze/thaw cycles. You may need to dispense the ligase buffer into smaller aliquots for your experimental needs.

No colonies with experimental insert, control insert or negative control.

Transformation has failed. Reassess the competent cells with an intact, supercoiled plasmid and determine the transformation efficiency. Use cells $>1 \times 10^8$ cfu/ μ g to insure >100 colonies from the control insert ligation.

Experimental insert has more blue colonies than control insert or negative control and fewer white colonies than control insert.

In-frame insertion, no interruption of α -fragment. Although the pGEM®-T Vector Control DNA will produce recombinants that generate white colonies, the insertion of other DNA fragments into the *lacZ* coding sequence may not result in white colonies unless the fragments disrupt the *lacZ* reading frame. Although this tends to occur most frequently with PCR products of 500bp or less, inserts of up to 2kb can result in blue colonies. Moreover, some insert DNAs can also result in pale blue colonies or “bull’s-eye” colonies with a blue center and a white perimeter. In one case in particular, we found that a 1.8kb insert when oriented in one direction produced white colonies and in the other produced bull’s-eye colonies [Knoche, K. and Kephart, D. (1999) Cloning blunt-end *Pfu* DNA polymerase-generated PCR fragments into pGEM®-T Vector Systems. *Promega Notes* 71, 10–13.].

PCR Subcloning

Subcloning Using PCR Primers Containing Restriction Sites

Frequently, the ends of insert DNA do not contain a suitable restriction enzyme site. The problem can be solved by using PCR to generate a site at the desired location. For this technique, the restriction enzyme site is designed into the 5'-end of the PCR primer. Because certain restriction enzymes inefficiently cleave recognition sequences located at the end of a DNA fragment, it is advisable to include at least four additional bases in front of the restriction recognition site. For the majority of restriction enzymes this will result in efficient cleavage.

Success in digesting PCR products can depend on the purity of the PCR product. Primers and primer dimers are present in overwhelming quantities when compared to the actual PCR product. Your PCR product will be competing with primers and primer dimers for the attention of the restriction enzyme, resulting in conditions favoring partial restriction digest. A simple clean-up of the reaction with the Wizard® SV Gel and PCR Clean-Up System can improve RE cleavage.

If you encounter a situation where the PCR product will not subclone, the digest may be adversely affected by proximity to the end of the PCR product. To improve the "placement" of the restriction site, the PCR product can be subcloned into the pGEM®-T Easy Vector. The restriction site should be readily cleavable in the context of the vector.

Ability of Restriction Enzymes to Cut PCR Products With RE Sites Near the End of the Fragment.

Enzyme	Distance (in bp) from the end of the PCR Fragment			
	0	1	2	3
<i>Apa</i> I	–	–	+/-	+
<i>Bam</i> H I	–	+/-	+	+
<i>Bst</i> X I	–	+/-	+	+
<i>Cla</i> I	–	+/-	+	+
<i>Eco</i> R I	–	+/-	+	+
<i>Eco</i> R V	–	+	+	+
<i>Hind</i> III	–	–	+	+
<i>Not</i> I	–	–	+	+
<i>Pst</i> I	–	–	+/-	+
<i>Sac</i> I	–	+/-	+	+
<i>Sal</i> I	+	+	+	+
<i>Sma</i> I	–	+/-	+	+
<i>Spe</i> I	+	+	+	+
<i>Xba</i> I	–	+/-	+	+
<i>Xho</i> I	–	–	+/-	+

PCR products in which the end of the restriction enzyme recognition sequence was flush with the end of the product or 1, 2, or 3 base pairs away from the end of the product were digested with a variety of enzymes. Purified PCR fragments (10–50ng) were digested with 0.5units of RE in 10µl of the appropriate reaction buffer for 45 minutes. Digestion is indicated as follows: Cleavable (+), not cleavable (–) and not reproducible (+/-). Data are the result of at least duplicate experiments and are reproduced by permission of Eaton Publishing. Taken from Simmermann, K. *et al.* (1998) Digestion of terminal restriction endonuclease recognition sites on PCR products. *BioTechniques* **24**, 582–4.

PCR Subcloning: Ordering Information

Basic PCR Cloning Systems

Product	Size	Cat. #
pGEM®-T Vector System I ^(e,f,g) Supplied with linearized, ligation-ready pGEM®-T Vector, 2X Rapid Ligation Buffer, T4 DNA Ligase and Positive Control Insert.	20 reactions	A3600
pGEM®-T Vector System II ^(e,f,g) Same contents as System I with 6 × 200µl JM109 High Efficiency Competent Cells.	20 reactions	A3610
pGEM®-T Easy Vector System I ^(e,f,g) Supplied with linearized, ligation-ready pGEM®-T Easy Vector, 2X Rapid Ligation Buffer, T4 DNA Ligase and Positive Control Insert.	20 reactions	A1360
pGEM®-T Easy Vector System II ^(e,f,g) Same contents as System I with 6 × 200µl JM109 High Efficiency Competent Cells. For Laboratory Use.	20 reactions	A1380

Sequencing Primers

Product	Conc.	Size	Cat. #
T7 Promoter Primer [5'-d(TAATACGACTCACTATAGGG)-3']	10µg/ml	2µg	Q5021
SP6 Promoter Primer [5'-d(TATTAGGTGACTATAG)-3']	10µg/ml	2µg	Q5011
pUC/M13 Primer, Forward (24 mer) [5'-d(CGCCAGGGTTTCCAGTCACGAC)-3']	10µg/ml	2µg	Q5601
pUC/M13 Primer, Reverse (22 mer) [5'-d(TCACACAGGAAACAGCTATGAC)-3']	10µg/ml	2µg	Q5421

Thermostable DNA Polymerases

Product	Conc.	Size	Cat. #
PCR Master Mix ^(h)	2X	100 reactions	M7502
	2X	1,000 reactions	M7505
PCR Master Mix contains dNTPs, buffer, Mg ²⁺ and <i>Taq</i> DNA Polymerase. A standard reaction contains 25µl of PCR Master Mix giving 1.5mM Mg ²⁺ , 200µM each dNTP and 1.25u of <i>Taq</i> DNA Polymerase in the final 50µl reaction.			
GoTaq® DNA Polymerase ⁽ⁱ⁾	100u	5u/µl	M3001
	500u	5u/µl	M3005
	2,500u	5u/µl	M3008
Supplied with 5X Green and 5X Colorless GoTaq® Reaction Buffer. Each contain 1.5mM MgCl ₂ in the final 1X concentration. Use the Green Buffer for direct gel analysis of amplification reactions. Use the Colorless Buffer for any reaction requiring absorbance or fluorescence measurements without prior PCR clean-up.			
<i>Taq</i> DNA Polymerase in Storage Buffer B ^(j) (Supplied with 10X Thermophillic Reaction Buffer and 25mM MgCl ₂ Solution.)	5u/µl	100u	M1661
	5u/µl	500u	M1665
	5u/µl	2,500u	M1668
<i>Taq</i> DNA Polymerase in Storage Buffer B ^(j) (Supplied with 10X Thermophillic Reaction Buffer containing 15mM MgCl ₂ .)	5u/µl	100u	M2661
	5u/µl	500u	M2665
	5u/µl	2,500u	M2668
<i>Taq</i> DNA Polymerase in Storage Buffer A ^(j) (Supplied with 10X Thermophillic Reaction Buffer and 25mM MgCl ₂ Solution.)	5u/µl	100u	M1861
	5u/µl	500u	M1865
	5u/µl	2,500u	M1868
<i>Taq</i> DNA Polymerase in Storage Buffer A ^(j) (Supplied with 10X Thermophillic Reaction Buffer containing 15mM MgCl ₂ .)	5u/µl	100u	M2861
	5u/µl	500u	M2865
	5u/µl	2,500u	M2868
For Laboratory Use.			

PCR Subcloning: Ordering Information

Thermostable DNA Polymerases

Product	Size	Conc.	Cat. #
<i>Taq</i> Bead™ Hot Start Polymerase ⁽¹⁾ *	100 reactions	1.25u/bead	M5661
(Supplied with 10X Thermophilic Reaction Buffer and 25mM MgCl ₂ Solution.)			
<i>T7</i> DNA Polymerase ⁽¹⁾ *	100u	5u/μl	M1941
(Supplied with <i>T7</i> 10X Reaction Buffer and 25mM MgSO ₄ Solution.)	1,000u	5u/μl	M1945
<i>T7h</i> DNA Polymerase ⁽¹⁾ *	100u	5u/μl	M2101
(Supplied with 10X Reverse Transcription Buffer, 10X Chelate Buffer, 10X Thermophilic Reaction Buffer, 25mM MgCl ₂ and 25mM MnCl ₂ .)	500u	5u/μl	M2105
<i>Pfu</i> DNA Polymerase ⁽¹⁾ *	100u	2–3u/μl	M7741
(Supplied with <i>Pfu</i> 10X Reaction Buffer containing MgSO ₄ . Not available in North America.)	500u	2–3u/μl	M7745
<i>T7i</i> DNA Polymerase ⁽¹⁾ *			
(Supplied with 10X Thermophilic Reaction Buffer and 25mM MgCl ₂ .)	50u	3u/μl	M7101

*For Laboratory Use.

PCR-qualified Nucleotides

Product	Conc.	Size	Cat. #
Set of dATP, dCTP, dGTP, dTTP ⁽¹⁾	100mM	10μmol each	U1330
	100mM	25μmol each	U1420
	100mM	40μmol each	U1240
	100mM	200μmol each	U1410
PCR Nucleotide Mix ⁽¹⁾	10mM	200μl	C1141
	10mM	1,000μl	C1145

Equal mixture of dATP, dCTP, dGTP and dTTP. Use 1μl per 50μl reaction for a final dNTP concentration of 200μM each.

For Laboratory Use.

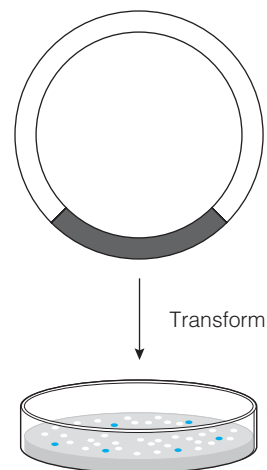
Transforming Bacteria

Properties of *E. coli* Strains for Subcloning

Common laboratory strains of *E. coli*, like JM109, DH5 α TM, and XL-1 Blue, are different from their wildtype counterparts. These strains carry some mutations designed to help you propagate plasmids. Typically laboratory strains have a mutation in the *recA* gene (*recA1*), a gene involved in recombination. The mutant gene limits recombination of the plasmid with the *E. coli* genome so that the plasmid inserts are more stable (the *recA1* mutation is more effective than the *recA13* mutation). Each of these strains also carries the *endA1* mutation that inactivates a nuclease that might copurify with plasmids during purification. This mutation helps you to purify higher quality plasmids. Special treatments must be performed on plasmids from strains that do not have this mutation (e.g., RR1, HB101, etc.) to eliminate the nuclease from the plasmid prep (e.g., the Alkaline Protease digestion in the Wizard[®] Plus SV Miniprep protocol).

Common laboratory strains of *E. coli* are typically defined as K strains or B strains based on the presence of the restriction and modification system that functions around *EcoK* I or *EcoB* I, respectively. In a wildtype K strain, the *E. coli* will have both the *EcoK* I restriction enzyme to cleave foreign DNA and *EcoK* I methylase to protect and mask host DNA recognition sequences. In B strains, the *EcoB* I restriction enzyme and methylase serve the same purpose. Strains like JM109, DH5 α TM and XL-1 Blue are K strains but carry the *hsdR17* (*r_K⁻*, *m_K⁺*) mutation. This mutation knocks out the *EcoK* I restriction enzyme but leaves the methylase intact. Therefore, these strains will not degrade plasmid DNA isolated from a B or K strain but will methylate it. This is useful if the DNA must be transferred to a K strain with an intact K restriction and methylation system.

If you wish to incorporate blue/white selection into your subcloning scheme, you need to transform *E. coli* carrying a *lacZ* Δ . This mutation deletes a portion of the β -galactosidase gene leaving what is termed the ω -fragment. The plasmid vector supplies this deleted portion, or α -fragment. Once inside the bacterium, the plasmid produces the α -fragment and the *E. coli* produces the ω -fragment, which combine to make a



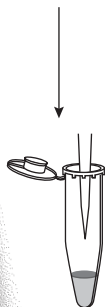
functional β -galactosidase. If grown on plate containing 5-bromo-4-chloro-3-indoyl- β -D-galactopyranoside (X-gal), the colony will turn blue as a result of β -galactosidase activity and indicate full complementation of the bacterium by the plasmid. This is termed α -complementation. Blue/White cloning methods use plasmids with a multiple cloning region within the coding sequence of the α -fragment. Disruption of the reading frame due to the presence of the insert will produce a non-functional α -fragment incapable of α -complementation. These disrupted plasmids are differentiated from the plasmids without insert by the color of the colony (white versus blue), hence the term blue/white selection. Strains like JM109, DH5 α TM and XL-1 Blue have the necessary deletion. One difference between these strains lies in how you get the bacterium to produce the ω -fragment. Both JM109 and XL-1 Blue have a second mutation call *lacI^q*. This mutation leads to increased production of the *lacI* repressor that stops transcription from the *lac* operon until substrate is present. To relieve this repression, these strains are grown on media containing the non-cleavable lactose analog, isopropyl- β -D-thiogalactopyranoside (IPTG). DH5 α TM does not have the *lacI^q* mutation and constantly produces a low level of the ω -fragment through leaky transcription of the *lac* operon and therefore does not require IPTG for blue/white selection.

Transforming Bacteria

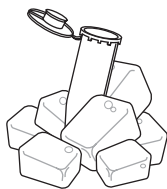
Ready-to-Use Competent Cells



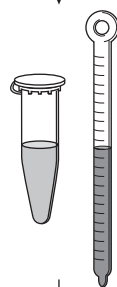
Thaw on crushed ice.



Add 1–50ng of DNA from ligation reaction.



Incubate on crushed ice for 30 minutes.

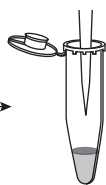


42°C

Heat shock at 42°C for 30 seconds.



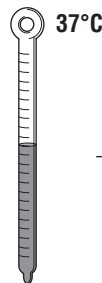
Place on crushed ice for 2 minutes.



Add 250µl room temperature SOC Media.



Incubate at 37°C; with shaking (225–250rpm) for 60 minutes.



37°C



Spread 100µl per plate. Incubate at 37°C overnight.

Use in blue/white selection cloning procedures!

Select96™ Competent Cells
Cat.# L3300 1 x 96 reactions

Full protocol available at:
www.promega.com/tbs/tb301/tb301.html

Single-use aliquots! Use one-at-a-time or all 96 at one time.



3722CA05_2A

4591MA

Transforming Bacteria

Making Your Own Competent Cells

Preparation of Competent Cells: Modified RbCl Method

This rubidium chloride protocol gives better transformation efficiencies than the CaCl_2 procedure for most strains. The procedure is an adaptation of one described in Hanahan, D. (1985) In: *DNA Cloning*, Volume 1, D. Glover, ed., IRL Press, Ltd., London, 109.

Materials to Be Supplied by the User

(Solution compositions are provided on page 48.)

- LB medium and plates
 - LB + 20mM MgSO_4
 - TFB1, ice-cold
 - TFB2, ice-cold
 - dry ice/isopropanol bath
1. Inoculate a single colony from an LB plate (for JM109, use M9 + B1 plate so that F' episome is maintained) into 2.5ml of LB medium in a plating tube. Incubate overnight at 37°C with shaking (approximately 225rpm).
 2. Subculture the overnight culture at a 1:100 dilution by inoculating 2.5ml into 250ml of LB supplemented with 20mM MgSO_4 . Grow the cells in a 1L flask until the OD_{600} reaches 0.4–0.6 (usually 5–6 hours, but the time may vary).
 3. Pellet the cells by centrifugation at $4,500 \times g$ for 5 minutes at 4°C. For a 250ml culture, use two 250ml centrifuge bottles in a large rotor.
 4. Gently resuspend the cell pellet in 0.4 original volume of ice-cold TFB1. For a 250ml subculture, use 100ml of TFB1 (50ml/bottle). Combine the resuspended cells in one bottle. For the remaining steps, keep the cells on ice and chill all pipettes, tubes and flasks.
 5. Incubate the resuspended cells on ice for 5 minutes at 4°C.
 6. Pellet the cells by centrifugation at $4,500 \times g$ for 5 minutes at 4°C.
 7. Gently resuspend the cells in 1/25 original volume of ice-cold TFB2. For a 250ml subculture, use 10ml of TFB2.
 8. Incubate the cells on ice for 15–60 minutes, then dispense 100µl/tube for storage at –70°C. Quick-freeze the tubes in a dry ice/isopropanol bath. JM109 competent cells prepared by this method are stable for 1 year.

Many *E. coli* strains carry episomes (e.g., F' and P2) expanding the capabilities of the bacterium for use in subcloning applications. For example, the XL1-Blue and JM109 strains carry the *lacI^qΔM15* mutation on the F' episome. The episomes are extrachromosomal, replicating plasmids with a selectable marker (page 64). When making competent cells from strains with episomes, the bacteria must first be plated on selective media. For XL1-Blue, colonies are selected on tetracycline plate since the episome contains the Tet^R gene. Due to this, however, the strain cannot be used with subcloning plasmids containing the Tet^R gene for selection. JM109 cells should be selected first on M9 minimal media containing thiamine (vitamin B1). The bacterial chromosome lacks the biosynthetic genes for proline synthesis (*proAB*) but the episome carries those genes. Colonies grown on the M9 + B1 plates (recipe on page 48) can then be processed into competent cells ready for blue/white selection.

Note: Competent cells may be conveniently quick-frozen using ice bath racks, which have an ice compartment bottom and a removable rack (American Scientific Products, Cat.# S9233-1). Set up an ice bath in one rack and an ethanol bath in another. Place the top-labeled tubes in the rack with ice, dispense 100µl cells per tube, then close the tubes. Add the dry ice to the ethanol bath, wait for it to stop bubbling, then transfer the rack and tubes to the dry ice bath for about 15 seconds. Drain the ethanol, wipe with a tissue, and transfer to an empty bottom compartment and place in a –70°C freezer. Do not get alcohol on the lips of the tubes. Liquid nitrogen also can be used for quick-freezing, but not with these racks. Use only plasticware designed for liquid nitrogen.

Note: Be careful not to get alcohol on the labels because it will remove them.

Transforming Bacteria

Determining Transformation Efficiency of Competent Cells

This is a general protocol for use with the procedure for producing competent cells that is provided on page 46. Please follow manufacturers' instructions when using purchased competent cells.

1. Thaw a 100µl aliquot of competent cells on ice.
2. Transfer 100µl of the cells to a 17 × 100mm polypropylene tube prechilled on ice.
3. Add 0.1ng of a supercoiled plasmid [e.g., pGEM®-3Zf(+) Vector] in a 10µl volume to the competent cells and gently mix by swirling the pipet tip (do not mix by pipetting).
4. Transfer the tubes from ice to a 42°C water bath and heat shock for 45–60 seconds. Place on ice immediately to cool for 2 minutes.
5. Add 890µl of SOC medium (giving a concentration of 0.1ng DNA/ml) and incubate for 45 minutes at 37°C with shaking (~150rpm).
6. Transfer 100µl of cells to 900µl of SOC medium (0.01ng DNA/ml) and plate 100µl of this (0.001ng/100µl) onto LB plates with the appropriate antibiotic. You may wish to plate 100µl of undiluted cells for determining efficiency as well. The 100µl aliquot will contain 0.01ng DNA.
7. Incubate the plates overnight in a 37°C incubator and count the number of colonies obtained. For example, if 200 colonies were obtained:

$$\frac{200\text{cfu}}{0.001\text{ng}} = 2 \times 10^5\text{cfu/ng} = 2 \times 10^8\text{cfu/}\mu\text{g DNA}$$

Competencies below 10^6 may not be useful for subcloning applications.

Transforming Ligation Reactions

This is a general protocol for use with the procedure for producing competent cells that is provided on page 46. Please follow manufacturers' instructions when using purchased competent cells.

1. Thaw a 100µl aliquot of competent cells on ice.
2. Transfer 100µl aliquot of the competent cells to a 17 × 100mm polypropylene tube prechilled on ice.
3. Add no more than 10ng of DNA in a maximum of 10µl from a ligation reaction to the cells and gently swirling the pipet tip (do not mix by pipetting). Incubate on ice for 30 minutes.
4. Transfer the tubes from ice to a 42°C water bath and heat-shock for 45–60 seconds. Place on ice immediately to cool for 2 minutes.
5. Add 1ml of LB or SOC medium and incubate for 45 minutes at 37°C with shaking (~150rpm).
6. Plate 100–200µl of the transformation mix, or an appropriate dilution onto selection plates. If you suspect low ligation efficiency, take the remaining cells and pellet by a quick 10–20 second spin in a microcentrifuge. Pour off media and resuspend pellet in about 200µl of SOC and plate.

Transformation Controls

Controls help you figure out where things may have gone wrong with the subcloning procedure. When transforming bacteria with your subcloning reaction DNA, also determine transformation and efficiency.

Transforming a ligation control (see page 26) of cut, dephosphorylated vector without insert can tell you how many background colonies you can expect in your actual vector + insert ligation.

Transforming more than 10ng of DNA from a ligation reaction may actually decrease transformation efficiency.

Transforming Bacteria

Media and Solutions

LB (Luria-Bertani) medium (1 liter)

- 10g tryptone
- 5g yeast extract
- 5g NaCl

Adjust pH to 7.5 with NaOH and autoclave.

LB plates (1 liter)

Add 15g agar to 1 liter of LB medium. Adjust to pH 7.5 with NaOH. Autoclave. Pour 30–35ml of medium into 85mm petri dishes. If necessary, flame the surface of the medium with a Bunsen burner to remove bubbles.

LB Medium

Allow the medium to cool to 55°C before adding antibiotic (either ampicillin, 125µg/ml final concentration; tetracycline, 12.5µg/ml final concentration; or chloramphenicol, 20µg/ml final concentration).

LB plates plus antibiotic (1 liter)

Add 15g agar to 1 liter of LB medium. Adjust to pH 7.5 with NaOH. Autoclave. Pour 30–35ml of medium into 85mm petri dishes. If necessary, flame the surface of the medium with a Bunsen burner to remove bubbles.

SOC medium

- 2.0g tryptone
- 0.5g yeast extract
- 1ml 1M NaCl
- 0.25ml 1M KCl
- 1ml Mg²⁺ stock
(1M MgCl₂ • 6H₂O, 1M MgSO₄ • 7H₂O) filter-sterilized
- 1ml 2M glucose, filter-sterilized

Add tryptone, yeast extract, NaCl and KCl to 97ml distilled water. Stir to dissolve. Autoclave and cool to room temperature. Add 2M Mg²⁺ stock and 2M glucose stock, each to a final concentration 20mM. Adjust to pH 7.0. Filter the complete medium through a 0.2µm filter unit.

M9 + B1 plates

- 6g Na₂HPO₄
- 3g KH₂PO₄
- 0.5g NaCl
- 1g NH₄Cl
- 15g Agar

Add deionized water to approximately 1L. Adjust to pH 7.4 with 10N NaOH. Autoclave and cool to 50°C. Then add the following sterile solutions:

- 2.0ml 1M MgSO₄
- 0.1ml 1M CaCl₂
- 10.0ml 20% Glucose (Filter Sterilized)
- 1.0ml 1M Thiamine-KCl (Filter Sterilized)

Pour 30–35ml of medium into 85mm petri dishes.

Bacterial Plates for Blue/White Selection.

Ampicillin Stock Solution

Dissolve at 50mg/ml in water, filter sterilize, store in aliquots at –20°C.

IPTG stock solution (0.1M)

1.2g IPTG (Cat.# V3951)

Add water to 50ml final volume. Filter-sterilize and store at 4°C.

X-Gal (2ml)

100mg X-gal (Cat.# V3941)

Dissolved at 50mg/ml in N,N'-dimethyl-formamide. Cover with aluminum foil and store at –20°C.

LB plates with ampicillin/IPTG/X-Gal

Add 15g agar to 1 liter of LB medium. Autoclave. Allow the medium to cool to 50°C before adding ampicillin to a final concentration of 100µg/ml then supplement with 0.5mM IPTG and 80µg/ml X-Gal and pour the plates. Pour 30–35ml of medium into 85mm petri dishes. Let the agar harden. Store at 4°C for up to 1 month or at room temperature for up to 1 week.

TFB1

- 30mM potassium acetate
- 10mM CaCl₂
- 50mM MnCl₂
- 100mM RbCl
- 15% glycerol

Adjust pH to 5.8 with 1M acetic acid. Be very careful as you approach 5.8; if the pH drops lower than 5.8, a black precipitate may form. Filter sterilize (0.2µm) and store at room temperature.

TFB2

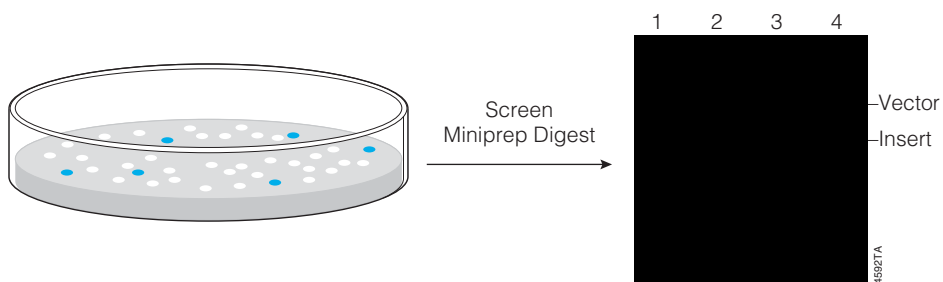
- 100mM MOPS or PIPES (pH 6.5)
- 75mM CaCl₂
- 10mM RbCl
- 15% glycerol

Adjust the pH to 6.5 with 1M KOH. Filter sterilize (0.2µm) and store at room temperature.

Screening for Recombinants

Introduction

Now that you've transformed your DNA and allowed the colonies to grow overnight, you need to determine if they contain the insert of interest. You can either screen them by colony PCR or the more traditional plasmid miniprep followed by restriction digestion. Colony PCR is the most rapid initial screen. A plasmid miniprep will take an extra day to grow up the culture but will provide a lot of material for further analysis. Some people do both and simply do not put the negative colonies that were identified through the colony PCR through the full miniprep procedure. The choice is yours.



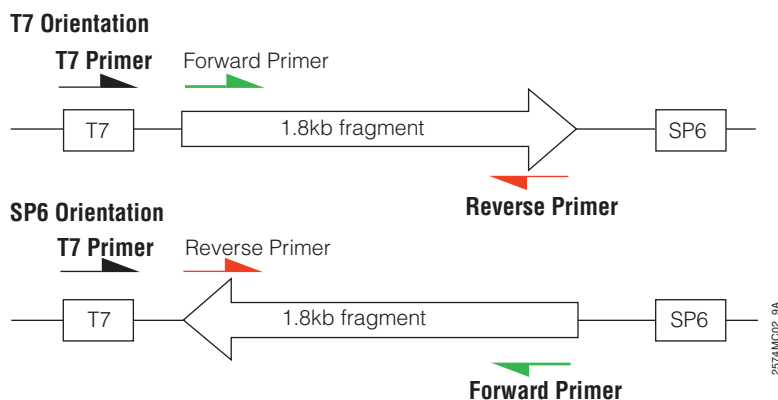
Screening for Recombinants

Colony PCR

Colony PCR involves lysing the bacteria and amplifying a portion of the plasmid. You can use either insert-specific primers or vector-specific primers to screen for recombinant plasmids. If your subcloning scheme will not maintain the orientation of the insert, you can use colony PCR to screen for orientation. Simply combine a vector-specific primer with an insert-specific primer.

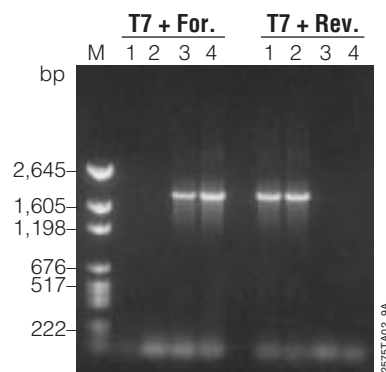
PCR cloning using the A-overhangs left by *Taq* DNA Polymerase^(†) and an appropriately T-tailed vector (e.g., pGEM[®]-T Easy Vector) is not a technique that will retain orientation. The orientation can be rapidly assessed with colony PCR using vector-specific primers and insert-specific primers as detailed below.

This technique was used when screening for orientation of a 1.8kb insert into the pGEM-T Easy Vector. Colony PCR was performed with the T7 Promoter Primer and either the insert-specific forward or reverse PCR primer. Eight white colonies were chosen from the cloning experiment for analysis. Clones with the T7 orientation will produce the fragment only with the T7 primer and reverse PCR primer, and clones in the opposite (SP6) orientation will only produce a product with the forward PCR primer as illustrated below.



Colony Prep for Colony PCR

1. Pick a well isolated colony and transfer to 50µl of sterile water. Part of the colony may be transferred to LB media containing the appropriate antibiotic for over night culture and miniprep if desired.
2. Boil for 10 minutes.
3. Centrifuge at $16,000 \times g$ for 5 minutes.
4. Use 5µl of the supernatant in a 50µl PCR.



Colony PCR. Colonies were suspended in 50µl sterile water, boiled for 10 minutes, centrifuged at $16,000 \times g$ for 5 minutes, and 5µl of the supernatant was used in each amplification. The DNA was amplified by PCR in 50µl volumes with 50pmol of each primer and 1.25 units of *Taq* DNA Polymerase (Cat.# M1661). After an initial denaturation of 2 minutes at 94°C, the amplification profile was 35 cycles of denaturation (94°C for 30 seconds), annealing (55°C for 1 minute) and extension (72°C for 2.5 minutes); PCR was concluded with 1 cycle of 72°C for 10 minutes. Amplification products (8µl) were analyzed on a 1% agarose gel containing ethidium bromide.

Screening for Recombinants

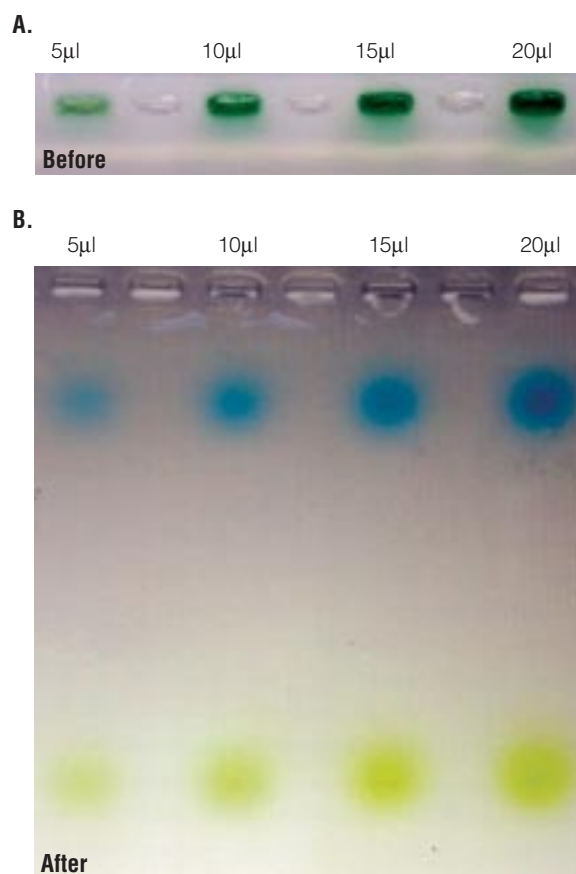
Go Directly to Gel

GoTaq® DNA Polymerase is the ideal choice for colony PCR applications. The enhanced buffer can handle the “dirty” template better than a conventional reaction buffer, and the Green GoTaq Reaction Buffer allows you to load the PCR products directly onto a gel after amplification. No need to add loading dye.

Colony PCR with GoTaq® DNA Polymerase Typical Reaction.

Nuclease-Free Water	to 50µl
5X Green GoTaq® Reaction Buffer	10µl
PCR Nucleotide Mix (Cat.# C1141)	1µl
GoTaq® DNA Polymerase (5u/µl)	1.25u
Downstream Primer	50pmol
Upstream Primer	50pmol
Colony Lysate	5µl

Assemble the reaction on ice in the order listed. A master mix containing everything but the colony lysate can be prepared and dispensed into reaction tubes ready to accept the colony lysate.



Amplification reactions using GoTaq DNA Polymerase with Green GoTaq Reaction Buffer. Panel A shows loaded wells of an agarose gel. Panel B shows the blue and yellow dyes after electrophoresis. Volumes of 5, 10, 15 and 20µl of the amplification reactions were loaded into a 1% agarose gel with TBE buffer and subjected to electrophoresis.

Cycling Conditions For GoTaq® Reactions.

Step	Temp	Time	Cycles
Initial			
Denaturation	94°C	2	1
Denaturation	94°C	0.5–1.0	
Annealing	42–65°C*	0.5–1.0	25–35
Extension	72°C	1 minute/kb	
Final Extension	72°C	5	1
Soak	4°C	Indefinite	1

Reactions are placed in a thermal cycler that has been preheated to 94°C.

*Annealing temperature should be optimized for each primer set based on the primer melting temperature (T_m). An online calculator for melting temperatures of primers in GoTaq® Reaction Buffer is available at:

www.promega.com/biomath

The extension time should be at least 1 minute/kilobase of target. Typically, anything smaller than 1kb uses a 1-minute extension.

GoTaq® DNA Polymerase
 Cat.# M3001 100v; 80 reactions
 Cat.# M3005 500v; 400 reactions
 Cat.# M3008 2,500v; 2,000 reactions
 Supplied with enzyme (5u/µl), 5X Green GoTaq® Reaction Buffer and 5X Colorless GoTaq® Reaction Buffer. Sufficient to give the indicated number of 50µl reactions using 1.25u of enzyme per reaction.
 Protocol available at:
www.promega.com/tbs/9pim300/9pim300.html

Screening for Recombinants

Screening by Plasmid Miniprep and RE Digests

The classic method for screening colonies involves performing a plasmid miniprep followed by restriction digestion. Well-isolated colonies are picked from a plate and transferred to culture medium containing the appropriate antibiotic for selection. Proper sterile technique is important. Many different culture media formulations are commonly used for minipreps. Promega recommends LB media supplemented with antibiotics (see page 48) for miniprep cultures to insure that the bacteria do not outgrow the ability of the antibiotic to select for the plasmid. If a rich medium like Terrific Broth is used, the bacteria can grow to very high cell densities and deplete the antibiotic. Once the antibiotic is depleted, the selection pressure to keep the plasmid is removed, and the plasmid may be lost.

You can inoculate the colony into 1–10ml of culture medium. If using a high-copy plasmid, 1–5ml (more typically, 1–2ml) is plenty. If you are using a low-copy plasmid, inoculate 10ml. Aerating the culture is very important for maximum cell density. A 17 × 100mm culture tube is fine for 1–2ml. If growing a larger volume, a 50ml sterile, disposable culture tube is better. Incubate the culture overnight (12–16 hours) with shaking (~250rpm). Remember, the greater the surface area, the greater the aeration. You can even grow miniprep cultures in sterile 25–50ml Erlenmeyer flasks.

Once the DNA is purified, a portion of the plasmid is screened by restriction digestion. For high-copy plasmids, you can obtain 4–10µg plasmid DNA per purification (1–5ml). For low-copy plasmids, you will obtain 1–3µg plasmid DNA per purification (10ml). Use 0.5–1µg of plasmid in your digest. Design the digest so that you can easily determine if your plasmid contains insert.

Note: Be sure to run uncut plasmid on the same gel for comparison.

Antibiotics: Mode of Action and Mechanism of Resistance.

Antibiotic	Mode of Action	Mechanism of Resistance	Working Concentration	Stock Solution
Ampicillin (Amp)	A derivative of penicillin that kills growing cells by interfering with bacterial cell wall synthesis.	The resistance gene (<i>bla</i>) specifies a periplasmic enzyme, β -lactamase, which cleaves the β -lactam ring of the antibiotic.	50–125µg/ml in water	50mg/ml
Chloramphenicol (Cm)	A bacteriostatic agent that interferes with bacterial protein synthesis by binding to the 50S subunit of ribosomes and preventing peptide bond formation.	The resistance gene (<i>cat</i>) specifies an acetyltransferase that acetylates, and thereby inactivates, the antibiotic.	20–170µg/ml in ethanol	34mg/ml
Kanamycin (Kan)	A bactericidal agent that binds to 70S ribosomes and causes misreading of messenger RNA.	The resistance gene (<i>kan</i>) specifies an enzyme (aminoglycoside phosphotransferase) that modifies the antibiotic and prevents its interaction with ribosomes.	30µg/ml in water	50mg/ml
Streptomycin (Sm)	A bactericidal agent that binds to the 30S subunit of ribosomes and causes misreading of the messenger RNA.	The resistance gene (<i>str</i>) specifies an enzyme that modifies the antibiotic and inhibits its binding to the ribosome.	30µg/ml in water	50mg/ml
Tetracycline (Tet)	A light-sensitive bacteriostatic agent that prevents bacterial protein synthesis by binding to the 30S subunit of ribosomes.	The resistance gene (<i>tet</i>) specifies a protein that modifies the bacterial membrane and prevents transport of the antibiotic into the cell.	10µg/ml in liquid culture; 12.5µg/ml in plates	12.5mg/ml in ethanol

Screening for Recombinants

Plasmid Minipreps

Wizard® Plus SV Minipreps DNA Purification System

The Wizard Plus SV Minipreps DNA Purification System^(h,i), a simple membrane-based system, provides a reliable method for rapidly isolating plasmid DNA. The entire procedure can be completed in 45 minutes or less. Work with up to 5ml of an overnight culture of a high-copy plasmid or up to 10ml of low-copy plasmid. Use the isolated DNA directly for applications such as automated fluorescent sequencing and restriction digests.

Vacuum protocol—
do 20 preps
at once.

Overnight culture

Centrifuge.

Remove culture media.
Resuspend cells.
Lyse cells.
Neutralize.

Clear lysate.

Transfer lysate.

Bind DNA.

Wash, removing solution by centrifugation or vacuum.

Elute plasmid DNA.

Transfer Spin Column to a Collection Tube. Centrifuge.

Spin protocol—
do as many preps
as your rotor
can hold.

Wizard® Plus SV Minipreps
Ready for spin protocols:

A1330

A1460

Ready for spin or vacuum protocols:

A1340

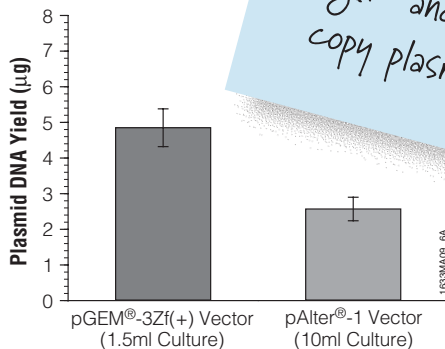
A1470

Protocol available at:
www.promega.com/tbs/tb225/tb225.html

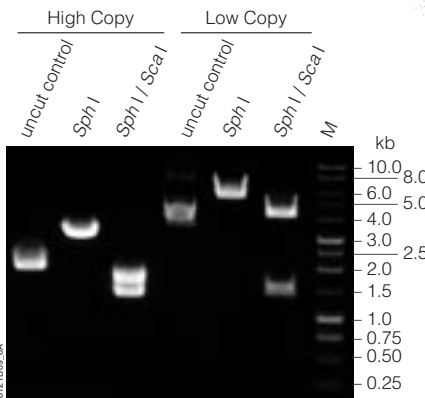
50 preps
250 preps

50 preps
250 preps

Works with
high- and low-
copy plasmids!



Plasmid DNA yield from high and low copy plasmids using the Wizard Plus SV Minipreps DNA Purification System. *E. coli* DH5α™ cells were transformed with either the pGEM®-3Zf(+) Vector (high-copy number plasmid) or the pALTER®-1 Vector (low-copy number plasmid) and grown in LB medium containing 50µg/ml of ampicillin (16 hours at 37°C, 200rpm). Plasmid DNA was isolated in sets of 42 on three consecutive days (126 total samples each) from 1.5ml (pGEM-3Zf(+) Vector) and 10ml (pALTER-1 Amp^r Vector) cultures using the Wizard® Plus SV Minipreps DNA Purification System.



Restriction enzyme digestion of high- and low-copy number plasmids. The plasmid vectors pGEM-3Zf(+) (high-copy) and pALTER-1 (low copy), isolated from *E. coli* DH5α using the Wizard Plus SV Minipreps DNA Purification System, were digested with 10 units of the indicated enzymes for 1 hour at 37°C. The digested samples were resolved on a 1% agarose gel and stained with ethidium bromide. The marker is the 1kb DNA Ladder (Cat.# G5711).

Screening for Recombinants

Troubleshooting Subcloning Experiments

Symptoms	Possible Causes	Comments
Few or no colonies obtained after transformation	Cells not competent	Competent cells may exhibit lower transformation efficiencies 5–6 weeks after preparation. To verify that bacteria are competent, perform a test transformation using a known amount of a standard supercoiled plasmid (see page 47).
	Unsuccessful ligation	Analyze samples of a linearized vector and the vector + insert ligation on an 0.8% agarose gel. If ligation was successful, the banding pattern of the ligation products should be different from that of the unligated sample.
	Inactive T4 DNA Ligase	Verify that the T4 DNA Ligase is active; perform a control ligation reaction with linear plasmid DNA.
	Inactive T4 DNA Ligase Buffer	Store T4 DNA Ligase 10X Reaction Buffer in small aliquots at -20°C to minimize freeze-thaw cycles of the buffer. Multiple freeze-thaw cycles may degrade the ATP in the buffer.
	Digested vector ends are not compatible with the fragment	Restriction enzyme sites that are adjacent within the multiple cloning region or near the ends can prove difficult to digest completely. See page 40 for a method to create PCR products with restriction sites in the primers.
	Excess ligation products added to competent cells	The added ligation products should not exceed 0.5% of the transformation reaction volume. Excess DNA ($>10\text{ng}$) may also inhibit the transformation.
High Background	Unsuccessful dephosphorylation of vector DNA	Attempt to religate the dephosphorylated vector. It should religate with low efficiency.
	Plates lack the correct antibiotic; the antibiotic is inactive	Perform a mock transformation with no DNA added. If colonies grow, discard the plates.
	Ratio of linearized, phosphorylated vector to insert DNA is too high	Reduce the amount of linearized vector in the reaction. Religation of the vector is favored when the vector:insert ratio is too high.

Screening for Recombinants

Troubleshooting Subcloning Experiments—Deletions of the Insert

During screening of your recombinant plasmids, you may encounter a situation where part of your insert is deleted. Perhaps careful examination of the subcloning strategy has not identified steps that may have led to this deletion. You attempt to sequence these clones and find that a portion of the insert has been deleted along with part of the vector. How did this happen? Two possibilities come to mind that could lead to such an event.

Symptoms	Possible Causes	Comments
Recombinants are isolated but contain deletions (usually unidirectional deletions of insert and part of vector)	Insert is unstable in the host strain	The insert may have been a substrate for recombination by recombinases in the host bacterium (remember most common laboratory strains are <i>recA</i> minus, but there are other recombinases present). You can transform the plasmid into an <i>E. coli</i> strain deficient in more recombinases than just the <i>recA</i> . Some strains like SURE® cells from Stratagene are deficient in recombinases <i>recB</i> and <i>recJ</i> and may allow you to propagate the unstable insert. Also try growing at a lower temperature.
	Insert is toxic to the host strain	Certain inserts may produce toxic gene products. The bacteria responds by deleting a portion of either the plasmid, the insert or both. High copy-number plasmids will tend to produce more toxin and thus be more prone to deletion or rearrangement. Most cloning plasmids carry the modified ColE1 origin of replication derived from pUC vectors, which maintain the copy number of the plasmid as high as 100–400 copies per bacterium. Low-level, leaky transcription in this high copy number plasmid can yield significant quantities of the toxic product. One solution is to transfer the insert to a different vector with a lower copy number. Promega has some vectors with lower copy numbers available like the pALTER®-1 (based on the pBR322 ColE1 origin; as few as 25 copies per cell) and the pALTER®-Ex2 (based on the pACYC origin of replication; ~10 copies per cell). This solution requires moving to a new vector. If you need to stay with the same vector (e.g., a mammalian expression vector) there are <i>E. coli</i> strains that have mutations that limit the copy number of pUC-based ColE1 origin of replication. The ABLE® strains from Stratagene will reduce copy number four- to ten-fold compared to more common laboratory strains.

Screening for Recombinants: Ordering Information

Competent Bacteria

Product	Size	Cat. #
Select96™ Competent Cells (>10 ⁸ cfu/μg)	1 × 96 reactions	L3300
JM109 Competent Cells, >10 ⁸ cfu/μg*	5 × 200μl	L2001
JM109 Competent Cells, >10 ⁷ cfu/μg	5 × 200μl	L1001

*For Laboratory Use.

Chemicals for Blue/White Screening

Product	Size	Conc.	Cat. #
IPTG, Dioxane-Free	1g	—	V3955
	5g	—	V3951
	50g	—	V3953
X-Gal	100mg	50mg/ml	V3941

For Laboratory Use.

Reagents for Colony PCR

Product	Conc.	Size	Cat. #
GoTaq® DNA Polymerase	5u/μl	100u	M3001
	5u/μl	500u	M3005
	5u/μl	2,500u	M3008
PCR Nucleotide Mix	10mM	200μl	C1141
	10mM	1,000μl	C1145
Set of dATP, dCTP, dGTP and dTTP	100mM	10μmol	U1330
	100mM	25μmol	U1420
	100mM	40μmol	U1240
	100mM	200μmol	U1410

For Laboratory Use.

Plasmid DNA Purification

Product	Size	Cat. #
Wizard® <i>Plus</i> SV Minipreps DNA Purification System ^{(L,M)*}	50 preps	A1330
	250 preps	A1460

Ready for use as a spin prep requiring only a microcentrifuge.

Product	Size	Cat. #
Wizard® <i>Plus</i> SV Minipreps DNA Purification System plus Vacuum Adapters ^{(L,M)*}	50 preps	A1340
	250 preps	A1470

Ready for use as a spin prep or a vacuum prep. Spin protocol requires only a microcentrifuge. Vacuum protocol requires Vac-Man® Laboratory Vacuum Manifold and a microcentrifuge.

Product	Size	Cat. #
Vac-Man® Laboratory Vacuum Manifold, 20-sample capacity	1 each	A7231

*For Laboratory Use.

Technical Appendix

Restriction Enzyme Activity in Promega 10X Buffers, Reaction Temperature and Heat Inactivation.

The 10X Reaction Buffer supplied with each restriction enzyme is optimized to give 100% activity. In many cases good activity is also obtained using one of Promega's 4-CORE® 10X Buffers. Many commonly used cloning enzymes (i.e. restriction sites are found in vector multiple cloning sites) have buffers E and H as their optimal buffer and so we have determined the activity of many of our other restriction enzymes in those buffers as well. This table may be used to select the best buffer for digestion with multiple restriction enzymes. Enzyme activity is expressed as a percent of the activity obtained with the optimal buffer for each enzyme in a one-hour digest. Enzymes with 100% activity (green) perform as well as the optimal buffer. Enzymes with 50–75% or 75–100% (yellow) will give acceptable activity in that buffer. Enzymes with <10%, 10–25%, or 25–50% (pink) activity generally do not have acceptable activity in that buffer. Also, buffers leading to star activity of the enzyme (*; pink) should be avoided. If compatible buffers cannot be identified with acceptable activity for both enzymes, each digest should be performed separately in the optimal buffer for each enzyme.

Promega Enzyme	Cat. #	Buffer Supplied with Enzyme	Activity in							MULTI-CORE™	Heat Inactivation	Enzyme Assay Temperature
			A	B	C	D	E	H				
Aat II	R6541	J	50–75%	10–25%	<10%	<10%	10–25%	<10%	<10%	+	37°C	
Acc I	R6411	G	50–75%	25–50%	25–50%	10–25%	<10%	<10%	25–50%	–	37°C	
Acc III	R6581	F	<10%	10–25%	25–50%	25–50%	n.d.	n.d.	<10%	–	65°C	
Acc65 I	R6921	D	10–25%	50–75%	75–100%	100%	75–100%	100–125%**	100%	+	37°C	
AccB7 I	R7081	E	10–25%	50–75%	100%*	<10%	100%	n.d.	100%	+	37°C	
Age I	R7251	K	25–50%	25–50%	25–50%	50–75%	n.d.	n.d.	100%	+	37°C	
Alu I	R6281	B	75–100%	100%	75–100%	10–25%	n.d.	n.d.	10–25%	+	37°C	
Alw26 I	R6761	C	10–25%	25–50%	100%	10–25%	n.d.	n.d.	75–100%	+	37°C	
Alw44 I	R6771	C	<10%	25–50%	100%	25–50%	n.d.	n.d.	100%	+	37°C	
Apa I	R6361	A	100%	50–75%	50–75%	<10%	10–25%	<10%	75–100%	+	37°C	
Ava I	R6091	B	10–25%	100%	50–75%	25–50%	100%	10–25%	<10%	+/–	37°C	
Ava II	R6131	C	50–75%	50–75%	100%	25–50%	n.d.	n.d.	25–50%	+	37°C	
Bal I	R6691	G	10–25%	<10%	<10%	<10%	n.d.	n.d.	<10%	+	37°C	
BamH I	R6021	E	75–100%*	75–100%	75–100%	50–75%	100%	50–75%	75–100%	+	37°C	
Ban I	R6891	G	25–50%	25–50%	10–25%	<10%	n.d.	n.d.	100%	–	50°C	
Ban II	R6561	E	75–100%	75–100%	75–100%	25–50%	n.d.	n.d.	100%	+	37°C	
Bbu I	R6621	A	100%	75–100%	75–100%	<10%	10–25%	10–25%	100%	+	37°C	
Bcl I	R6651	C	10–25%	75–100%	100%	50–75%	50–75%	50–75%	10–25%	–	50°C	
Bgl I	R6071	D	10–25%	25–50%	75–100%	100%	25–50%	75–100%	100%	+	37°C	
Bgl II	R6081	D	25–50%	75–100%	75–100%	100%	n.d.	n.d.	<10%	–	37°C	
BsaM I	R6991	D	10–25%	25–50%	50–75%	100%	n.d.	n.d.	25–50%	–	65°C	
Bsp1286 I	R6741	A	100%	50–75%	25–50%	10–25%	n.d.	n.d.	75–100%	+	37°C	
BsrS I	R7241	D	10–25%	25–50%	10–25%	100%	n.d.	n.d.	100%	–	65°C	
BssH II	R6831	H	75–100%	50–75%	75–100%	50–75%	n.d.	100%	75–100%	–	50°C	
Bst98 I	R7141	D	<10%	10–25%	10–25%	100%	n.d.	n.d.	25–50%	–	37°C	
BstE II	R6641	D	25–50%	50–75%	50–75%	100%	n.d.	n.d.	100%	–	60°C	
BstO I	R6931	C	10–25%	25–50%	100%	25–50%	n.d.	n.d.	<10%	–	60°C	
BstX I	R6471	D	<10%	10–25%	25–50%	100%	100%	75–100%	10–25%	+/–	50°C	
BstZ I	R6881	D	<10%	<10%	10–25%	100%	10–25%	75–100%	10–25%	–	50°C	
Bsu36 I	R6821	E	<10%	25–50%	50–75%	25–50%	100%	n.d.	50–75%	–	37°C	
Cfo I	R6241	B	75–100%	100%	75–100%	25–50%	n.d.	n.d.	100%	+/–	37°C	
Cla I	R6551	C	75–100%	75–100%	100%	75–100%	100%	50–75%	100%	+	37°C	
Csp I	R6671	K	<10%	10–25%	25–50%	50–75%	100%	100–125%**	10–25%	+	30°C	
Csp45 I	R6571	B	25–50%	100%	50–75%	25–50%	100%	25–50%	50–75%	+	37°C	
Dde I	R6291	D	25–50%	25–50%	50–75%	100%	n.d.	n.d.	25–50%	+/–	37°C	
Dpn I	R6231	B	50–75%	100%	75–100%	50–75%	n.d.	n.d.	100%	+	37°C	
Dra I	R6271	B	75–100%	100%	75–100%	50–75%	n.d.	n.d.	25–50%	+	37°C	
EclHK I	R7111	E	<10%	<10%	75–100%	10–25%	100%	n.d.	50–75%	+	37°C	
Eco47 III	R6731	D	<10%	25–50%	50–75%	100%	n.d.	n.d.	25–50%	+	37°C	
Eco52 I	R6751	L	<10%	<10%	10–25%	25–50%	25–50%	50–75%	<10%	+	37°C	
EcoICR I	R6951	B	10–25%	100%	75–100%	<10%	25–50%	n.d.	100%	+	37°C	
EcoR I	R6011	H	25–50%	50–75%	50–75%	50–75%	75–100%	100%	100%*	+	37°C	
EcoR V	R6351	D	10–25%	25–50%	50–75%	100%	25–50%	50–75%	100%	+	37°C	
Fok I	R6781	B	75–100%	100%	75–100%	25–50%	n.d.	n.d.	50–75%	+	37°C	
Hae II	R6661	B	50–75%	100%	50–75%	10–25%	n.d.	n.d.	100%	–	37°C	
Hae III	R6171	C	75–100%	75–100%	100%	50–75%	n.d.	n.d.	100%	–	37°C	
Hha I	R6441	C	50–75%	75–100%	100%	50–75%	n.d.	n.d.	75–100%	+	37°C	
Hinc II	R6031	B	25–50%	100%	25–50%	50–75%	75–100%	50–75%	100%	+	37°C	
Hind III	R6041	E	25–50%	100%	75–100%	10–25%	100%	25–50%	50–75%	+	37°C	
Hint I	R6201	B	50–75%	100%	75–100%	75–100%	n.d.	n.d.	50–75%	–	37°C	

Technical Appendix

Restriction Enzyme Activity in Promega 10X Buffers, Reaction Temperature and Heat Inactivation (continued).

Promega Enzyme	Cat.#	Buffer Supplied with Enzyme	Activity in							MULTI-CORE™	Heat Inactivation	Enzyme Assay Temperature
			A	B	C	D	E	H				
<i>Hpa</i> I	R6301	J	25–50%	50–75%	25–50%	10–25%	n.d.	n.d.	100%	—	37°C	
<i>Hpa</i> II	R6311	A	100%	50–75%	50–75%	10–25%	n.d.	n.d.	100%	—	37°C	
<i>Hsp92</i> I	R7151	F	10–25%	75–100%	50–75%	25–50%	n.d.	n.d.	10–25%	+	37°C	
<i>Hsp92</i> II	R7161	K	10–25%	25–50%	25–50%	<10%	n.d.	n.d.	<10%	+	37°C	
<i>I-Ppo</i> I	R7031	NA	10–25%	25–50%	25–50%	25–50%	n.d.	n.d.	—	+	37°C	
<i>Kpn</i> I	R6341	J	100%*	25–50%	25–50%	<10%	25–50%	<10%	75–100%	+/–	37°C	
<i>Mbo</i> I	R6711	C	10–25%	75–100%	100%	50–75%	n.d.	n.d.	<10%	+	37°C	
<i>Mbo</i> II	R6723	B	10–25%	100%	50–75%	75–100%	n.d.	n.d.	100%	+	37°C	
<i>Mlu</i> I	R6381	D	10–25%	25–50%	50–75%	100%	25–50%	100–125%**	10–25%	+/–	37°C	
<i>Msp</i> I	R6401	B	75–100%	100%	75–100%	25–50%	n.d.	n.d.	25–50%	+	37°C	
<i>MspA1</i> I	R7021	C	25–50%	100%*	100%	10–25%	n.d.	n.d.	100%	+	37°C	
<i>Nae</i> I	R7131	A	100%	50–75%	25–50%	<10%	n.d.	n.d.	50–75%	+	37°C	
<i>Nar</i> I	R6861	G	75–100%	50–75%	75–100%	25–50%	n.d.	n.d.	50–75%	+	37°C	
<i>Nci</i> I	R7061	B	100%*	100%	25–50%	25–50%	n.d.	n.d.	50–75%	+	37°C	
<i>Nco</i> I	R6513	D	50–75%	75–100%	75–100%	100%	100%	100–125%**	75–100%	+	37°C	
<i>Nde</i> I	R6801	D	<10%	<10%	25–50%	100%	n.d.	n.d.	25–50%	+	37°C	
<i>Nde</i> II	R7291	D	<10%	<10%	10–25%	100%	n.d.	n.d.	25–50%	+	37°C	
<i>NgoM</i> IV	R7171	MULTI-CORE™	100%*	100%*	100%*	<10%	n.d.	n.d.	100%	+	37°C	
<i>Nhe</i> I	R6501	B	75–100%	100%	75–100%	10–25%	75–100%	10–25%	100%	+	37°C	
<i>Not</i> I	R6431	D	<10%	10–25%	25–50%	100%	25–50%	100–125%**	25–50%	+	37°C	
<i>Nru</i> I	R7091	K	<10%	<10%	<10%	50–75%	n.d.	n.d.	10–25%	+	37°C	
<i>Nsi</i> I	R6531	D	10–25%	50–75%	50–75%	100%	25–50%	>125%**	10–25%	+/–	37°C	
<i>Pst</i> I	R6111	H	10–25%	50–75%	50–75%	50–75%	25–50%	100%	25–50%	+	37°C	
<i>Pvu</i> I	R6321	D	10–25%	25–50%	50–75%	100%	n.d.	n.d.	<10%	–	37°C	
<i>Pvu</i> II	R6331	B	25–50%	100%	50–75%	25–50%	n.d.	n.d.	50–75%	+	37°C	
<i>Rsa</i> I	R6371	C	75–100%	75–100%	100%	<10%	n.d.	n.d.	<10%	+	37°C	
<i>Sac</i> I	R6061	J	75–100%	25–50%	25–50%	<10%	100%	25–50%	100%	+	37°C	
<i>Sac</i> II	R6221	C	100%	50–75%	100%	50–75%	25–50%	>125%**	<10%	+	37°C	
<i>Sal</i> I	R6051	D	<10%	10–25%	25–50%	100%	25–50%	25–50%	<10%	+	37°C	
<i>Sau3A</i> I	R6191	B	25–50%	100%	75–100%	<10%	n.d.	n.d.	100%	+	37°C	
<i>Sca</i> I	R6211	K	<10%	100%*	50–75%	75–100%	n.d.	n.d.	10–25%	+	37°C	
<i>Sfi</i> I	R6391	B	75–100%	100%	75–100%	25–50%	75–100%	50–75%	75–100%	–	50°C	
<i>Sgf</i> I	R7103	C	25–50%	25–50%	100%	<10%	n.d.	n.d.	<10%	+/–	37°C	
<i>Sin</i> I	R6141	A	100%	75–100%	50–75%	10–25%	n.d.	n.d.	100%	+	37°C	
<i>Sma</i> I	R6121	J	<10%	<10%	<10%	<10%	<10%	<10%	100%	+	25°C	
<i>SnaB</i> I	R6791	B	50–75%	100%	50–75%	<10%	n.d.	n.d.	100%	–	37°C	
<i>Spe</i> I	R6591	B	75–100%	100%	75–100%	75–100%	100%	25–50%	100%	+	37°C	
<i>Sph</i> I	R6261	K	75–100%	75–100%	100%*	75–100%	100%	>125%**	10–25%	+	37°C	
<i>Ssp</i> I	R6601	E	10–25%	50–75%	50–75%	75–100%	100%	100–125%**	50–75%	+	37°C	
<i>Stu</i> I	R6421	B	75–100%	100%	75–100%	50–75%	n.d.	n.d.	50–75%	+	37°C	
<i>Sty</i> I	R6481	F	25–50%	75–100%	75–100%	75–100%	10–25%	50–75%	<10%	+	37°C	
<i>Taq</i> I	R6151	E	10–25%	25–50%	50–75%	50–75%	100%	n.d.	100%	–	65°C	
<i>Tru9</i> I	R7011	F	75–100%	50–75%	75–100%	25–50%	n.d.	n.d.	25–50%	–	65°C	
<i>Tth111</i> I	R6841	B	50–75%	100%	75–100%	25–50%	n.d.	n.d.	100%	–	65°C	
<i>Vsp</i> I	R6851	D	<10%	25–50%	75–100%	100%	n.d.	n.d.	<10%	+	37°C	
<i>Xba</i> I	R6181	D	50–75%	75–100%	75–100%	100%	100%	100–125%**	100%	–	37°C	
<i>Xho</i> I	R6161	D	25–50%	75–100%	75–100%	100%	25–50%	100–125%**	10–25%	+	37°C	
<i>Xho</i> II	R6811	C	25–50%	25–50%	100%	10–25%	n.d.	n.d.	<10%	+	37°C	
<i>Xma</i> I	R6491	B	50–75%	100%	25–50%	<10%	25–50%	<10%	50–75%	+	37°C	
<i>Xmn</i> I	R7271	B	75–100%	100%	75–100%	10–25%	n.d.	n.d.	75–100%	+	37°C	

* Not recommended due to potential star activity.

** Unit activity is based on recommended buffer. In Buffer H, some enzymes have enhanced activity.

n.d. = Not determined.

Heat Inactivation Key:

+	=greater than 95% inactivation (DNA is undigested)
—	=less than 95% inactivation (DNA digest is complete, i.e., ≥5% of the initial 20 activity units [≥1 unit] remains)
+/-	=partial inactivation (DNA is partially digested)

Heat Inactivation Conditions:

Twenty units of enzyme in 50µl of its optimal buffer were heated at 65°C for 15 minutes.

One microgram of DNA was added and incubated for 1 hour in accordance with the unit definition, then analyzed by agarose gel electrophoresis.

Technical Appendix

Isoschizomers. The enzymes in boldface type are available from Promega.

Enzyme	Isoschizomer(s)	Recognition Sequence
<i>mAat</i> I	Stu I, <i>Eco</i> 147 I, <i>Pme</i> 55 I, <i>Sse</i> B I	AGG▼CCT
Aaf II	—	GACGT▼C
Acc I	<i>Fbl</i> I, <i>Xmi</i> I	GT▼(A/C)(G/T)AC
Acc III	<i>Bsp</i> E II, <i>Mro</i> I	T▼CCGGA
Acc65 I	<i>Asp</i> 718 I Kpn I*	G▼GTACC GGTAC▼C
<i>Acc</i> B1 I	Ban I, <i>Bsh</i> N I, <i>Eco</i> 64 I	G▼G(C/T)(G/A)CC
AccB7 I	<i>Pfi</i> M I, <i>Van</i> 91 I	CCAN ₄ ▼NTGG
<i>Acc</i> N I	Spe I	A▼CTAGT
<i>Acc</i> W I	<i>Alw</i> I	GGATCNNNN▼
<i>Acy</i> I	<i>Bbi</i> II, <i>Hin</i> 1 I, Hsp92 I, <i>Bsa</i> H I, <i>Msp</i> 171 I	G(A/G)▼CG(T/C)C
<i>Acs</i> I	<i>Apo</i> I	(G/A)▼AATT(C/T)
<i>Afa</i> I	<i>Csp</i> 6 I*, Rsa I	GT▼AC
<i>Afe</i> I	Eco47 III	AGC▼GCT
<i>Afl</i> II	Bst 98 I	C▼TTAAGG
Age I	<i>Pin</i> A I	A▼CCGGT
<i>Aha</i> III	Dra I	TTT▼AAA
<i>Ahd</i> I	Ecl HK I	GACNNN▼NNGTC
Alu I	—	AG▼CT
<i>Alw</i> I	<i>Acc</i> W I	GGATCNNNN▼
Alw26 I ¹	<i>Bsm</i> A I	GTCTC(1/5)
Alw44 I	<i>Apa</i> L I	G▼TGCAC
<i>Aoc</i> I	Bsu 36 I, <i>Cvn</i> I	CC▼TNAGG
Apa I	<i>Bsp</i> 120 I	GGGCC▼C
<i>Apa</i> L I	Alw44 I, <i>Vne</i> I	G▼TGCAC
<i>Apo</i> I	<i>Acs</i> I	(G/A)▼AATT(C/T)
<i>Ase</i> I	Vsp I, <i>Asn</i> I	AT▼TAAT
<i>Asn</i> I	Vsp I, <i>Ase</i> I	AT▼TAAT
<i>Asp</i> I	Tih 111 I	GACN▼NNGTC
<i>Asp</i> E I	<i>Ahd</i> I, <i>Eam</i> 1105 I, Ecl HK I	GACNNN▼NNGTC
<i>Asp</i> 700 I	Xmn I	GAANN▼NNTTC
<i>Asp</i> 718 I	Acc65 I Kpn I*	G▼GTACC GGTAC▼C
<i>Asu</i> I	<i>Sau</i> 96 I, <i>Cfr</i> 13 I	G▼GNCC
<i>Asu</i> II	<i>Csp</i> 45 I, <i>Bst</i> B I	TT▼CGAA
<i>Asu</i> HP I	<i>Hph</i> I	GGTGAN ₈ ▼
Ava I	<i>Ama</i> 87 I, <i>Bco</i> I, <i>Bso</i> B I, <i>Eco</i> 88 I	C▼(C/T)CG(G/A)G
Ava II	Sin I, <i>Eco</i> 47 I, <i>Hgi</i> E I	G▼G(A/T)CC
<i>Axy</i> I	Bsu 36 I	CC▼TNAGG
Bal I	<i>Msc</i> I, <i>Mlu</i> N I	TGG▼CCA
Bam H I	—	G▼GATCC
Ban I	<i>Acc</i> B1, <i>Bsh</i> N I, <i>Eco</i> 64 I	G▼G(T/C)(A/G)CC
Ban II	<i>Eco</i> 24 I	G(A/G)GC(T/C)▼C
<i>Bbe</i> I	— Nar I*	GGCGC▼C GG▼CGCC
<i>Bbr</i> P I	<i>Eco</i> 72 I, <i>Pml</i> I	CAC▼GTG
<i>Bbs</i> I ¹	<i>Bsc</i> 91 I, <i>Bpi</i> I	GAAGAC(2/6)
Bbu I	<i>Pae</i> I, Sph I	GCATG▼C
Bcl I	<i>Bsi</i> Q I, <i>Fba</i> I	T▼GATCA
<i>Bcn</i> I	Nci I	CC▼(C/G)GG
<i>Bfr</i> I	Bst 98 I	C▼TTAAG
Bgl I	—	GCCNNNN▼NGGC
Bgl II	—	A▼GATCT
<i>Bmy</i> I	Bsp1286 I	G(G/A/T)GC(C/A/T)▼C
<i>Bpm</i> I	<i>Gsu</i> I	CTGGAG(16/14)
<i>Bsa</i> H I	Hsp92 I	G(A/G)▼CG(T/C)C
Bsa M I	<i>Bsm</i> I	GAATGC(1/–1)
<i>Bsa</i> O I	<i>Bsh</i> 1285 I, <i>Bsi</i> E I	CG(A/G)(T/C)▼C
<i>Bse</i> A I	Acc III	T▼CCGGA
<i>Bse</i> N I	Bsr S I, <i>Bsr</i> I	ACTGGN (1/–1)
<i>Bse</i> P I	Bss H II, <i>Pau</i> I	G▼CGCGC
<i>Bsh</i> 1285 I	<i>Bsa</i> O I	CG(A/G)(T/C)▼CG
<i>Bsh</i> N I	Ban I, <i>Acc</i> B1 I, <i>Eco</i> 64 I	G▼G(T/C)(A/G)CC
<i>Bsh</i> 1365 I	<i>Bsr</i> BR I	GATNN▼NATC
<i>Bsi</i> E I	<i>Bsa</i> O I	CG(A/G)(T/C)▼CG

Enzyme	Isoschizomer(s)	Recognition Sequence
<i>Bsm</i> I	Bsa M I	GAATGCN▼
<i>Bsm</i> A I ¹	Alw26 I	GTCTC(1/5)
<i>Bso</i> B I	Ava I, <i>Ama</i> 87 I, <i>Bco</i> I, <i>Eco</i> 88 I	C(C/T)CG(G/A)G
<i>Bsp</i> 19 I	Nco I	C▼CATGG
<i>Bsp</i> 68 I	Nru I	TCG▼CGA
<i>Bsp</i> 106 I	Cla I, <i>Bsa</i> D I	AT▼CGAT
<i>Bsp</i> 119 I	Csp45 I, <i>Nsp</i> V, <i>Bst</i> B I	TT▼CGAA
<i>Bsp</i> 120 I	Apa I	G▼GGCCC
<i>Bsp</i> 143 I	Mbo I, Sau3A I, Nde II	▼GATC
<i>Bsp</i> 143 II	Hae II	(A/G)GCGC▼(T/C)
Bsp1286 I	<i>Bmy</i> I, <i>Sdu</i> I	G(G/A/T)GC(C/A/T)▼C
<i>Bsp</i> C I	Pvu I	CGAT▼CG
<i>Bsp</i> D I	Cla I	AT▼CGAT
<i>Bsp</i> E I	Acc III	T▼CCGGA
<i>Bsr</i> I ¹	Bsr S I, <i>Bse</i> N I	ACTGGN(1/–1)
Bsr S I ¹	<i>Bse</i> N I, <i>Bsr</i> I	ACTGGN(1/–1)
Bss H II	<i>Bse</i> P I, <i>Pau</i> I	G▼CGCGC
Bst 98 I	<i>Afl</i> II, <i>Bfr</i> I	C▼TTAAG
<i>Bst</i> B I	Csp45 I, <i>Nsp</i> V, <i>Bsp</i> 119 I	TT▼CGAA
Bst E II	<i>Bst</i> P I, <i>Eco</i> 91 I, <i>Psp</i> E I	G▼GTNACC
<i>Bst</i> N I	Bst O I, <i>Mva</i> I, <i>Eco</i> R II	CC▼(A/T)GG
Bst O I	<i>Bst</i> N I, <i>Eco</i> R II, <i>Mva</i> I	CC▼(A/T)GG
Bst X I	—	CCANNNNN▼NTGG
<i>Bst</i> Y I	Xho II, <i>Mil</i> I	(A/G)▼GATC(T/C)
Bst Z I	Eco52 I, <i>Eag</i> I, <i>Xma</i> III, <i>Ecl</i> X I	C▼GGCCG
<i>Bsu</i> 15 I	Cla I	AT▼CGAT
Bsu 36 I	<i>Cvn</i> I, <i>Aoc</i> I, <i>Eco</i> 81 I	CC▼TNAGG
<i>Bsu</i> R I	Hae III, <i>Pal</i> I	GG▼CC
Cfo I	Hha I <i>Hin</i> 6 I <i>Hin</i> P1 I*	GCG▼C GCG▼C G▼CGC
<i>Cfr</i> 9 I	Xma I Sma I*	C▼CCGGG CCC▼GGG
<i>Cfr</i> 13 I	<i>Sau</i> 96 I	G▼GNCC
<i>Cfr</i> 42 I	Sac II	CCGC▼GG
Cla I	<i>Ban</i> III, <i>Bsp</i> 106 I, <i>Bsp</i> D I, <i>Bsu</i> 15 I	AT▼CGAT
<i>Cpo</i> I	Csp I, <i>Rsr</i> II	CG▼G(A/T)CCG
Csp I	<i>Cpo</i> I, <i>Rsr</i> II	CG▼G(A/T)CCG
<i>Csp</i> 6 I	Rsa I*, <i>Afa</i> I*	GT▼AC
Csp45 I	<i>Bst</i> B I, <i>Nsp</i> V, <i>Bsp</i> 119 I	TT▼CGAA
<i>Cvn</i> I	Bsu 36 I	CC▼TNAGG
Dde I	<i>Bst</i> DE I	C▼TNAG
Dpn I ²	<i>Dpn</i> II*	GmeA▼TC
<i>Dpn</i> II	Mbo I, Sau3A I, Nde II, Dpn I*	▼GATC
Dra I	—	TTT▼AAA
<i>Eag</i> I	Eco52 I, Bst Z I, <i>Ecl</i> X I, <i>Xma</i> III	C▼GGCCG
<i>Eam</i> 1105 I	Ecl HK I, <i>Ahd</i> I, <i>Asp</i> E I	GACNNN▼NNGTC
<i>Ecl</i> 136 II	Ecol CR I Sac I*	GAG▼CTC GAGCT▼C
Ecl HK I	<i>Ahd</i> I, <i>Eam</i> 1105 I, <i>Asp</i> E I	GACNNN▼NNGTC
<i>Ecl</i> X I	Bst Z I, <i>Eag</i> I, Eco52 I, <i>Xma</i> III	C▼GGCCG
<i>Eco</i> 24 I	Ban II, <i>Fri</i> O I	G(AG)GC(TC)▼C
<i>Eco</i> 32 I	Eco R V	GAT▼ATC
<i>Eco</i> 47 I	Ava II, Sin I	G▼G(A/T)CC
Eco47 III	<i>Afe</i> I	AGC▼GCT
Eco52 I	Bst Z I, <i>Xma</i> III, <i>Eag</i> I, <i>Ecl</i> X I	C▼GGCCG
<i>Eco</i> 64 I	Ban I, <i>Bsh</i> N I, <i>Eco</i> 64 I	G▼G(TC)(AG)CC
<i>Eco</i> 81 I	Bsu 36 I	CC▼TNAGG
<i>Eco</i> 88 I	Ava I	C▼(TC)CG(AG)G
<i>Eco</i> 91 I	Bst E II	G▼GTNACC
<i>Eco</i> 105 I	Sna B I	TAC▼CTA
<i>Eco</i> 130 I	Sty I	C▼C(A/T)(T/A)GG
<i>Eco</i> 147 I	Stu I	AGG▼CCT
Ecol CR I	<i>Ecl</i> 136 II	GAG▼CTC
	Sac I*	GAGCT▼C
	<i>Sst</i> I*	GAGCT▼C

Technical Appendix

Isoschizomers (continued). The enzymes in boldface type are available from Promega.

Enzyme	Isoschizomer(s)	Recognition Sequence	Enzyme	Isoschizomer(s)	Recognition Sequence
EcoR I	—	G▼AATTC	Rsa I	<i>Afa I</i>	GT▼AC
<i>EcoR II</i>	BstO I , <i>BstN I</i> , <i>Mva I</i>	CC▼(A/T)GG	<i>Rsr II</i>	Csp I , <i>Cpo I</i>	CG▼G(A/T)CCG
EcoR V	<i>Eco32 I</i>	GAT▼ATC	Sac I	<i>Sst I</i>	GAGCT▼C
<i>EcoT14 I</i>	Sty I	C▼C(A/T)(A/T)GG		<i>Ec136 II*</i> , EcoICR I*	GAG▼CTC
<i>EcoT22 I</i>	Nsi I	ATGCA▼T	Sac II	<i>Sst II</i> , <i>Ksp I</i> , <i>Cfr42 I</i>	CCGC▼GG
<i>Ehe I</i>	Nar I*	GG▼CGCC	Sal I	—	G▼TCGAC
Fok I²	—	GGATG(9/13)	Sau3A I	Mbo I , Nde II , <i>Dpn II</i>	▼GATC
Hae II	<i>Bsp143 II</i>	(A/G)GCGC▼(T/C)	<i>Sau96 I</i>	<i>Cfr13 I</i>	G▼GNCC
Hae III	<i>BsuR I</i> , <i>Pal I</i>	GG▼CC	Sca I	—	AGT▼ACT
<i>Hap II</i>	Hpa II , Msp I	C▼CGG	<i>Sdu I</i>	Bsp1286 I	G(G/A/T)GC(C/A/T)▼C
<i>HgiE I</i>	<i>Eco47 I</i> , Sin I , Ava II	G▼G(A/T)CC	Sfi I	—	GGCCNNNN▼NGGCC
Hha I	Cfo I	GCG▼C	<i>Sfu I</i>	Csp45 I	TT▼CGAA
	<i>HinP1 I*</i> , <i>Hin6 I*</i>	G▼CGC	Sgf I	—	GCGAT▼CGC
<i>Hin1 I</i>	<i>Acy I</i> , Hsp92 I	G(A/G)▼CG(T/C)C	Sin I	Ava II , <i>Eco47 I</i>	G▼G(A/T)CC
Hinc II	<i>Hind II</i>	GT(T/C)▼(A/G)AC	Sma I	—	CCC▼GGG
<i>Hind II</i>	Hinc II	GT(T/C)▼(A/G)AC		Xma I* , <i>Cfr9 I*</i>	C▼CCGGG
Hind III	—	A▼AGCTT	SnaB I	<i>Eco105 I</i>	TAC▼GTA
Hinf I	—	G▼ANTC	Spe I	<i>AccI</i>	A▼CTAGT
<i>HinP1 I</i>	—	G▼CGC	Sph I	Bbu I , <i>Pae I</i>	GCATG▼C
	Hha I* , Cfo I*	GCG▼C	Ssp I	—	AAT▼ATT
Hpa I	<i>KspA I</i>	GTT▼AAC	<i>Sst I</i>	Sac I	GAGCT▼C
Hpa II³	Msp I , <i>Hap II</i>	C▼CGG		EcoICR I*	GAG▼CTC
Hsp92 I	<i>Acy I</i> , <i>BsaH I</i> , <i>Hin1 I</i>	G(A/G)▼CG(C/T)	<i>Sst II</i>	Sac II	CCGC▼GG
Hsp92 II	<i>Nla III</i>	CATG▼	Stu I	<i>Aat I</i> , <i>Eco147 I</i>	AGG▼CCT
I-Ppo I	—	CTCTCTAA▼GGTAGC	Sty I	<i>EcoT14 I</i>	C▼C(A/T)(A/T)GG
<i>Kas I</i>	Nar I*	GG▼CGCC	Taq I	<i>TthHB8 I</i>	T▼CGA
Kpn I	—	GGTAC▼C	Tru9 I	<i>Mse I</i>	T▼TAA
	Acc65 I* , <i>Asp718 I*</i>	G▼GTACC	Tth111 I	<i>Asp I</i>	GACN▼NNGTC
<i>Ksp I</i>	Sac II	CCGC▼GG	<i>TthHB8 I</i>	Taq I	T▼CGA
Mbo I	Sau3A I , Nde II , <i>Dpn II</i>	▼GATC	<i>Van91 I</i>	AccB7 I , <i>PfiM I</i>	CCAN ₄ ▼NTGG
Mbo II⁴	—	GAAGA(8/7)	<i>Vne I</i>	<i>ApaL I</i> , Alw44 I	G▼TGCAC
<i>Mil I</i>	Xho II	(A/G)▼GATC(T/C)	Vsp I	<i>Ase I</i> , <i>Asn I</i>	AT▼TAAT
Mlu I	—	A▼CGCGT	Xba I	—	T▼CTAGA
<i>MluN I</i>	Bal I , <i>Msc I</i>	TGG▼CCA	Xho I	<i>PaeR7 I</i>	C▼TCGAG
<i>Mro I</i>	Acc III	T▼CCGGA	Xho II	<i>BstY I</i> , <i>Mil I</i>	(A/G)▼GATC(T/C)
<i>Msc I</i>	Bal I , <i>MluN I</i>	TGG▼CCA	Xma I	<i>Cfr9 I</i> , <i>XmaC I</i> , Sma I*	C▼CCGGG
<i>Mse I</i>	Tru9 I	T▼TAA			CCC▼GGG
Msp I³	Hpa II , <i>Hap II</i>	C▼CGG	<i>Xma III</i>	Eco52 I , BstZ I , <i>Eag I</i> , <i>EcIX I</i>	C▼GGCCG
MspA1 I	<i>NspB II</i>	C(A/C)G▼C(G/T)G	<i>XmaC I</i>	Xma I	C▼CCGGG
<i>Mst II</i>	Bsu36 I	CC▼TNAGG		Sma I*	CCC▼GGG
<i>Mva I</i>	BstO I , <i>EcoR II</i> , <i>BstN I</i>	CC▼(A/T)GG	Xmn I	<i>Asp700 I</i>	GAANN▼NNTTC
Nae I	NgoM IV	G▼CCGGC	Key: N = A, C, G or T * = neoschizomer Notes: 1. The locations of cleavage sites falling outside the recognition site are indicated in parentheses. For example, GTCTC(1/5) indicates cleavage at: 5'...GTCTC▼...3' 3'...CAGAGNNNN▼...5'		
Nar I	—	GG▼CGCC			
	<i>Ehe I*</i>	GGC▼GCC			
	<i>Kas I*</i>	G▼GCGCC			
	<i>Bbe I*</i>	GGCGC▼C			
Nci I	<i>Bcn I</i>	CC▼(C/G)GG	Reference Roberts, R.J. (1991) <i>Nucl. Acids Res.</i> 19 (supp), 2077–109.		
Nco I	<i>Bsp19 I</i>	C▼CATGG			
Nde I	—	CA▼TATG			
Nde II	Mbo I , Sau3A I , <i>Dpn II</i>	▼GATC			
NgoM IV	Nae I	G▼CCGGC	2. <i>Dpn I</i> is unique among commercially available restriction enzymes in requiring methylation of a nucleotide (adenine) in its recognition sequence in order to cut. Therefore, <i>Dpn I</i> cannot be substituted for other enzymes recognizing the GATC sequence (e.g., <i>Mbo I</i> and <i>Sau3A I</i>). 3. Although <i>Hpa II</i> and <i>Msp I</i> recognize the same nucleotide sequence, <i>Hpa II</i> is sensitive to methylation of either cytosine in its recognition sequence, while <i>Msp I</i> is sensitive only to methylation of the external cytosine. These enzymes may not be interchanged for all applications.		
Nhe I	—	G▼CTAGC			
<i>Nla III</i>	Hsp92 II	CATG▼			
Not I	—	GC▼GGCCGC			
Nru I	<i>Bsp68 I</i>	TCG▼CGA	Reference Roberts, R.J. (1991) <i>Nucl. Acids Res.</i> 19 (supp), 2077–109.		
Nsi I	<i>EcoT22 I</i> , <i>Mph1103 I</i>	ATGCA▼T			
<i>Nsp V</i>	Csp45 I , <i>BstB I</i> , <i>Bsp119 I</i>	TT▼CGAA			
<i>NspB II</i>	MspA1 I	C(A/C)G▼C(G/T)G			
<i>Pae I</i>	Bbu I , Sph I	GCATG▼C	2. <i>Dpn I</i> is unique among commercially available restriction enzymes in requiring methylation of a nucleotide (adenine) in its recognition sequence in order to cut. Therefore, <i>Dpn I</i> cannot be substituted for other enzymes recognizing the GATC sequence (e.g., <i>Mbo I</i> and <i>Sau3A I</i>). 3. Although <i>Hpa II</i> and <i>Msp I</i> recognize the same nucleotide sequence, <i>Hpa II</i> is sensitive to methylation of either cytosine in its recognition sequence, while <i>Msp I</i> is sensitive only to methylation of the external cytosine. These enzymes may not be interchanged for all applications.		
<i>PaeR7 I</i>	Xho I	C▼TCGAG			
<i>Pal I</i>	Hae III , <i>BsuR I</i>	GG▼CC			
<i>PfiM I</i>	AccB7 I , <i>Vau91 I</i>	CCAN ₄ ▼NTGG			
<i>PinA I</i>	Age I	A▼CCGGT	Reference Roberts, R.J. (1991) <i>Nucl. Acids Res.</i> 19 (supp), 2077–109.		
Pst I	—	CTGCA▼G			
Pvu I	<i>BspC I</i>	CGAT▼CG			
Pvu II	—	CAG▼CTG			

Technical Appendix

Compatible Ends.

Promega Restriction Enzymes That Generate 5' Overhangs

Overhang	Definite Compatible Ends	Possible Compatible Ends
5'-N		<i>Tth</i> 111 I
5'-S		<i>Nci</i> I
5'-W		<i>Bst</i> O I
5'-AT	<i>Acc</i> I	
5'-CG	<i>Nar</i> I, <i>Msp</i> I, <i>Hsp</i> 92 I, <i>Taq</i> I, <i>Cla</i> I, <i>Csp</i> 45 I, <i>Hpa</i> II	
5'-GN		<i>Bsr</i> S I
5'-MK		<i>Acc</i> I
5'-TA	<i>Vsp</i> I, <i>Nde</i> I, <i>Tru</i> 9 I	
5'-ANT		<i>Hint</i> I
5'-GNC		<i>Sau</i> 96 I
5'-GWC		<i>Ava</i> II, <i>Csp</i> I, <i>Sin</i> I
5'-TNA		<i>Dde</i> I, <i>Bsu</i> 36 I
5'-AATT	<i>Eco</i> R I	
5'-AGCT	<i>Hind</i> III	
5'-CATG	<i>Nco</i> I	<i>Sty</i> I
5'-CCGG	<i>Age</i> I, <i>Xma</i> I, <i>Acc</i> III, <i>Ngo</i> M IV	<i>Ava</i> I
5'-CGCG	<i>Mlu</i> I, <i>Bss</i> H I	
5'-CTAG	<i>Spe</i> I, <i>Nhe</i> I, <i>Xba</i> I	<i>Sty</i> I
5'-CWWG		<i>Sty</i> I
5'-GATC	<i>Mbo</i> I, <i>Sau</i> 3A I, <i>Bam</i> H I, <i>Bgl</i> II, <i>Xho</i> II, <i>Bcl</i> I, <i>Nde</i> II	
5'-GCGC	<i>Ban</i> I	
5'-GGCC	<i>Not</i> I, <i>Bst</i> Z I, <i>Eco</i> 52 I	
5'-GTAC	<i>Acc</i> 65 I	<i>Ban</i> I
5'-GTNAC		<i>Bst</i> E II
5'-GYRC		<i>Ban</i> I
5'-TCGA	<i>Sal</i> I, <i>Xho</i> I	<i>Ava</i> I
5'-TGCA	<i>Alw</i> 44 I	
5'-TTAA	<i>Bst</i> 98 I	
5'-YCGR		<i>Ava</i> I

Promega Restriction Enzymes That Generate 3' Overhangs

Overhang	Definite Compatible Ends	Possible Compatible Ends
N-3'		<i>Ecl</i> HK I
AT-3'	<i>Sgf</i> I, <i>Pvu</i> I	
CG-3'	<i>Cfo</i> I, <i>Hha</i> I	
CN-3'		<i>Bsa</i> M I
GC-3'	<i>Sac</i> II	<i>Bsa</i> O I
NNN-3'		<i>Acc</i> B7 I, <i>Bgl</i> I, <i>Sfi</i> I
ACGT-3'	<i>Aat</i> II	
AGCT-3'	<i>Sac</i> I	<i>Ban</i> II, <i>Bsp</i> 1286 I
CATG-3'	<i>Hsp</i> 92 II, <i>Sph</i> I, <i>Bbu</i> I	
DGCH-3'		<i>Bsp</i> 1286 I
GCGC-3'	<i>Hae</i> II	
GGCC-3'	<i>Apa</i> I	<i>Ban</i> II, <i>Bsp</i> 1286 I
GTAC-3'	<i>Kpn</i> I	
NNNN-3'		<i>Bst</i> X I
RGCY-3'		<i>Ban</i> II, <i>Bsp</i> 1286 I
TGCA-3'	<i>Nsi</i> I, <i>Pst</i> I	<i>Bsp</i> 1286 I
TTAA-3'	<i>I-Ppo</i> I	

Key:

D = A or G or T
 H = A or C or T
 K = G or T
 M = A or C
 N = A or C or G or T
 R = A or G
 S = C or G
 W = A or T
 Y = C or T

Technical Appendix

Site-Specific Methylation Sensitivity of Promega Restriction Enzymes.

This table lists the sensitivities of several Promega restriction enzymes to site-specific methylation at *dam*, *dcm*, CpG and CpNpG sites (p = phosphoryl group). These four modifications are frequently found in DNA of bacteria, eukaryotes or their viruses. Many strains of *E. coli* contain the site-specific *dam* and *dcm* DNA methylases. Higher eukaryotes contain the site-specific CpG and CpNpG DNA methylases. In mammalian genomes, methylation occurs mainly at the CG dinucleotide. In plant genomes, methylation may occur at both the CG and CNG sequences.

Prokaryotic Methylation

dcm Cytosine methylase mutation—methylates the C5 position of the internal cytosine residue in the sequence 5'...CCTGG...3'.

dam Adenine methylase mutation—methylates the N6 position of the adenine residue in the sequence 5'...GATC...3'.

Eukaryotic Methylation

CpG Methylates the C5 position of the cytosine residue in the dinucleotide recognition sequence 5'...CG...3'.

CpNpGp Methylates the C5 position of the cytosine residue in the trinucleotide recognition sequence 5'...CNG...3' (N = any base).

For further information regarding site-specific methylation, refer to McClelland, M., Nelson, M. and Raschke, E. (1994) *Nucl. Acids Res.* **22**, 3640–59.

Key:

- s = sensitive to this methylation
- i = insensitive to this methylation
- s(ol) = overlapping (sensitive when restriction site overlaps methylation sequence)
- n/a = information not available

Enzyme	Recognition Sequence	<i>dam</i>	<i>dcm</i>	CpG	CpNpG	Enzyme	Recognition Sequence	<i>dam</i>	<i>dcm</i>	CpG	CpNpG
<i>Aat</i> II	GACGTC	i	i	s	i	<i>Kpn</i> I	GGTACC	i	i	i	i
<i>Acc</i> B7 I	CCANNNNTGG	i	s(ol)	i	i	<i>Mbo</i> II	GAAGA(8/7)	s(ol)	i	i	i
<i>Acc</i> III	TCCGGA	s(ol)	i	i	i	<i>Mlu</i> I	ACGCGT	i	i	s	i
<i>Acc</i> 65 I	GGTACC	i	s(ol)	i	i	<i>Msp</i> I	CCGG	i	i	i	s
<i>Apa</i> I	GGGCCC	i	s(ol)	s(ol)	i	<i>Nae</i> I	GCCGGC	i	i	s	s
<i>Ava</i> I	CYCGRG	i	i	s	i	<i>Nar</i> I	GGCGCC	i	i	s	i
<i>Ava</i> II	GGWCC	i	s(ol)	s(ol)	s(ol)	<i>Nde</i> II	GATC	s	i	i	i
<i>Bal</i> I	TGGCCA	i	s(ol)	i	s(ol)	<i>Ngo</i> M IV	GCCGGC	i	i	s	s
<i>Bam</i> H I	GGATCC	i	i	i	s(ol)	<i>Nhe</i> I	GCTAGC	i	i	s(ol)	s(ol)
<i>Ban</i> II	GRGCTC	i	i	i	i	<i>Not</i> I	GCGGCCGC	i	i	s	s
<i>Bbu</i> I	GCATGC	i	i	i	i	<i>Nru</i> I	TCGCGA	s(ol)	i	s	i
<i>Bcl</i> I	TGATCA	s	i	i	i	<i>Pst</i> I	CTGCAG	i	i	i	s
<i>Bgl</i> I	GCCNNNNNGGC	i	i	s(ol)	s(ol)	<i>Pvu</i> I	CGATCG	i	i	s	s(ol)
<i>Bgl</i> II	AGATCT	i	i	i	s(ol)	<i>Pvu</i> II	CAGCTG	i	i	i	s
<i>Bsp</i> 1286 I	GDGCHC	i	i	i	i	<i>Sac</i> I	GAGCTC	i	i	i	i
<i>Bss</i> H II	GCGCGC	i	i	s	i	<i>Sac</i> II	CCGCGG	i	i	s	s
<i>Bst</i> E II	GGTNACC	i	i	i	i	<i>Sal</i> I	GTCGAC	i	i	s	n/a
<i>Bst</i> O I	CCWGG	i	i	i	n/a	<i>Sau</i> 3A I	GATC	i	i	s(ol)	s(ol)
<i>Bst</i> X I	CCANNNNNNTGG	i	i	i	i	<i>Sau</i> 96 I	GGNCC	i	s(ol)	s(ol)	s(ol)
<i>Bst</i> Z I	CGGCCG	i	i	s(ol)	s(ol)	<i>Sca</i> I	AGTACT	i	i	i	i
<i>Cfo</i> I	GCGC	i	i	s	n/a	<i>Sfi</i> I	GGCCNNNNNGGCC	i	s(ol)	s(ol)	s(ol)
<i>Cla</i> I	ATCGAT	s(ol)	i	s	i	<i>Sgf</i> I	GCGATCGC	i	i	s	n/a
<i>Csp</i> I	CGGWCCG	i	i	i	s	<i>Sin</i> I	GGWCC	i	i	i	s(ol)
<i>Csp</i> 45 I	TTCGAA	i	i	s	i	<i>Sma</i> I	CCCGGG	i	i	s	s
<i>Dde</i> I	CTNAG	i	i	i	s(ol)	<i>Sna</i> B I	TACGTA	i	i	s	i
<i>Eco</i> 47 III	AGCGCT	i	i	s	i	<i>Sph</i> I	GCATGC	i	i	i	i
<i>Eco</i> 52 I	CGGCCG	i	i	s	i	<i>Stu</i> I	AGGCCT	i	s(ol)	i	s(ol)
<i>Eco</i> R I	GAATTC	i	i	s(ol)	i	<i>Taq</i> I	TCGA	s(ol)	i	i	i
<i>Fok</i> I	GGATC	i	i	i	i	<i>Xba</i> I	TCTAGA	s(ol)	i	i	i
<i>Hae</i> III	GGCC	i	i	i	s(ol)	<i>Xho</i> I	CTCGAG	i	i	s	i
<i>Hha</i> I	GCGC	i	i	s	s(ol)	<i>Xho</i> II	RGATCY	i	i	i	s(ol)
<i>Hinc</i> II	GTYRAC	i	i	i	i	<i>Xma</i> I	CCCGGG	i	i	i	n/a
<i>Hind</i> III	AAGCTT	i	i	i	i	<i>Xmn</i> I	GAANNNN	i	i	n/a	n/a
<i>Hpa</i> II	CCGG	i	i	s	s						

Technical Appendix

Restriction Enzyme Buffer Composition.

Buffer	pH (at 37°C)	Tris-HCl (mM)	MgCl ₂ (mM)	NaCl (mM)	KCl (mM)	DTT (mM)
A	7.5	6	6	6	—	1
B	7.5	6	6	50	—	1
C	7.9	10	10	50	—	1
D	7.9	6	6	150	—	1
E	7.5	6	6	100	—	1
F	8.5	10	10	100	—	1
G	8.2	50	5	—	—	—
H	7.5	90	10	50	—	—
J	7.5	10	7	—	50	1
K	7.4	10	10	—	150	—
L	9.0	10	3	100	—	—

MULTI-CORE™ Buffer (1X) = 25mM Tris-acetate (pH 7.5 at 37°C), 100mM potassium acetate, 10mM magnesium acetate, 1mM DTT.

Notes:

- For each 10°C rise in temperature between 0°C and 25°C, the pH of Tris buffers decreases 0.31 pH units.
- For each 10°C rise in temperature between 25°C and 37°C, the pH of Tris buffers decreases 0.25 pH units.
- All of Promega's Restriction enzymes are supplied with 10mg/ml Acetylated BSA. Although BSA is not absolutely required for activity, it has been shown to enhance activity of many restriction enzymes. We recommend adding BSA to all restriction digests at a final concentration of 0.1mg/ml.

Copy Number of Commonly Used Plasmids.

Plasmid	Plasmid Size (approx.)	Origin of Replication*	Copy Number	**Yield per ml of Culture	Reference
pGEM®	2,700bp	mutated pMB1	300–700	1.8–4.1µg	1
pUC	2,700bp	mutated pMB1	500–700	2.9–4.1µg	1
pBR322	4,400bp	pMB1	>25	>0.23µg	2
ColE1	4,500bp	ColE1	>15	>0.15µg	3
pACYC	4,000bp	p15A	~10	~0.09µg	4
pSC101	9,000bp	pSC101	~6	~0.12µg	5
pGL Series	5,000bp	mutated pMB1	300–700	3.3–7.6µg	1
pRL Series	4,000bp	mutated pMB1	300–700	2.7–6.0µg	1
phRL Series	4,000bp	mutated pMB1	300–700	2.7–6.0µg	1
phRG Series	4,000bp	mutated pMB1	300–700	2.7–6.0µg	1
pGEM®-T/ T easy	3,000bp	mutated pMB1	300–700	2.0–4.6µg	1
psiLentGene™ Series	4,000bp	mutated pMB1	300–700	2.7–6.0µg	1
psiCHECK™ 1/2	3,500bp	mutated pMB1	300–700	2.7–5.3µg	1
psiSTRIKE™ Series	4,000bp	mutated pMB1	300–700	2.7–6.0µg	1
pALTER®-1/Ex1	5,800bp	pMB1	>25	>0.3µg	3
pALTER®-Ex2	5,800bp	p15A	~10	~0.13µg	4
pSP	2,500bp	mutated pMB1	300–700	1.6–3.8µg	1
pCI, pSI	3,600bp	mutated pMB1	300–700	2.4–5.5µg	1

* Plasmids carrying the pMB1, mutated pMB1 and ColE1 belong to the same incompatibility group, so they are not compatible with one another, but they are fully compatible with those carrying p15A and pSC101 replicons.

** Theoretical plasmid yields were calculated from the reported copy number and size of each plasmid assuming 2.0×10^9 cells per milliliter of culture grown for 16 hours at 37°C.

References

- Summerton, J., Atkins, T. and Bestwick, R. (1983) *Anal. Biochem.* **133**, 79–84.
- Holmes, D.S. and Quigley, M. (1981) *Anal. Biochem.* **114**, 193–7.
- Jansz, H.S., Pouwels, P.H. and Schiphorst, J. (1966) *Biochem. Biophys. Acta* **123**, 626.
- Birnboim, H.C. and Doly, J. (1979) *Nucl. Acids Res.* **7**, 1513–23.
- Birnboim, H.C. (1983) *Meth. Enzymol.* **100**, 243–55.

Star Activity.

Restriction enzymes, under nonstandard conditions, can demonstrate the ability to cleave DNA at sequences different from their defined recognition sites. The term "star activity" has been given to this nonsequence-specific cleavage of DNA under nonoptimal reaction conditions. The most common types of altered activity are single-base substitutions, truncation of the outer bases in the recognition sequence and single-strand nicking (1). In general, star activity is not a concern if restriction endonucleases are used in the recommended buffers at the appropriate temperatures. Star activity is evident with a number of restriction enzymes when the following parameters are altered in the reaction environment (2):

- High enzyme concentration (generally >100 units/µg).
- High glycerol content (>5% v/v).
- Substitution of Mn²⁺ for Mg²⁺ (or substitution of other divalent cations).
- Low salt concentration (generally <25mM).
- Extremes of pH, especially pH>8.0.
- Presence of DMSO, ethanol or other organic solvents.

References

- Barany, F. (1988) *Gene* **65**, 149–65.
- Brown, T.A., Hames, B.D. and Rickwood, D. (1991) In: *Molecular Biology Lab Fax*, BIOS Scientific Publishers Limited, Oxford, United Kingdom.

Technical Appendix

Genotypes of Frequently Used Bacterial Strains.

All genes in the bacterium are presumed to be in the wildtype state, except for those listed, which are mutant alleles carried by that bacterium. Genes listed on the F' episome, however, represent wildtype alleles unless specified otherwise. Strains are λ^- unless specified otherwise. *Strains available from Promega as competent cells are indicated by an asterisk. Strains shown in **bold** are available from Promega as glycerol freezer stocks.

Strain	Genotype
BL21(DE3)	F ⁻ , <i>ompT</i> , <i>hsdS_B</i> (<i>r_B⁻</i> , <i>m_B⁻</i>), <i>dcm</i> , <i>gal</i> , λ (DE3)
*BL21(DE3)pLysS	F ⁻ , <i>ompT</i> , <i>hsdS_B</i> (<i>r_B⁻</i> , <i>m_B⁻</i>), <i>dcm</i> , <i>gal</i> , λ (DE3), pLysS (Cm ^r)
*BMH 71-18 mutS	<i>thi</i> , <i>supE</i> , Δ (<i>lac-proAB</i>), [<i>mutS</i> ::Tn10(<i>tet^r</i>)] [F', <i>tra</i> D36, <i>proAB</i> , <i>laqI^a</i> Δ M15]
C600 (1)	<i>thi-1</i> , <i>thr-1</i> , <i>leuB6</i> , <i>lacY1</i> , <i>tonA21</i> , <i>supE44</i>
C600 <i>hfl</i> (1)	<i>thi-1</i> , <i>thr-1</i> , <i>leuB6</i> , <i>lacY1</i> , <i>tonA21</i> , <i>supE44</i> , <i>hflA150</i> ::Tn10(<i>tet^r</i>)
DH1 (2)	<i>recA1</i> , <i>endA1</i> , <i>gyrA96</i> , <i>thi-1</i> , <i>hsdR17</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>supE44</i> , <i>relA1</i>
DH10B	F ⁻ , <i>mcrA</i> Δ (<i>mrr-hsdRMS-mcrBC</i>) ϕ 80 <i>lacZ</i> Δ M15, Δ <i>lacX74</i> , <i>deoR</i> , <i>recA1</i> , <i>endA1</i> , <i>araD139</i> , Δ (<i>ara, leu</i>)7697, <i>galU</i> , <i>galK</i> , λ^- , <i>rpsL</i> (<i>str^r</i>), <i>nupG</i>
DH5 α TM	ϕ 80d <i>lacZ</i> Δ M15, <i>recA1</i> , <i>endA1</i> , <i>gyrA96</i> , <i>thi-1</i> , <i>hsdR17</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>supE44</i> , <i>relA1</i> , <i>deoR</i> , Δ (<i>lacZYA-argF</i>) U169, <i>phoA</i>
DM1 (3)	F', <i>dam-13</i> ::Tn9(Cm ^r) <i>dcm</i> , <i>mcrB</i> , <i>hsdR-M^r</i> , <i>gal1</i> , <i>gal2</i> , <i>ara-</i> , <i>lac-</i> , <i>thr-</i> , <i>leu-</i> , <i>ton^R</i> , <i>tsx^R</i> , <i>Su^o</i>
ES1301 mutS	<i>lacZ53</i> , <i>thyA36</i> , <i>rha-5</i> , <i>metB1</i> , <i>deoC</i> , IN(<i>rrnD-rrnE</i>), [<i>mutS201</i> ::Tn5]
*HB101 (4)	<i>thi-1</i> , <i>hsdS20</i> (<i>r_B⁻</i> , <i>m_B⁻</i>), <i>supE44</i> , <i>recA13</i> , <i>ara-14</i> , <i>leuB6</i> , <i>proA2</i> , <i>lacY1</i> , <i>galK2</i> , <i>rpsL20</i> (<i>str^r</i>), <i>xyl-5</i> , <i>mtl-1</i>
JM101 (5)	<i>supE</i> , <i>thi</i> , Δ (<i>lac-proAB</i>), F' (<i>traD36</i> , <i>proAB</i> , <i>lacI^a</i> Δ M15)
*JM109 (5)	<i>endA1</i> , <i>recA1</i> , <i>gyrA96</i> , <i>thi-1</i> , <i>hsdR17</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>relA1</i> , <i>supE44</i> , Δ (<i>lac-proAB</i>), [F', <i>traD36</i> , <i>proAB</i> , <i>lacI^a</i> Δ M15]
JM109(DE3) (5)	<i>endA1</i> , <i>recA1</i> , <i>gyrA96</i> , <i>thi-1</i> , <i>hsdR17</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>relA1</i> , <i>supE44</i> , Δ (<i>lac-proAB</i>), [F', <i>traD36</i> , <i>proAB</i> , <i>lacI^a</i> Δ M15], λ (DE3)
JM110 (5)	<i>rpsL</i> (<i>str^r</i>), <i>thr</i> , <i>leu</i> , <i>thi</i> , <i>hsdR17</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>lacY</i> , <i>galK</i> , <i>galT</i> , <i>ara</i> , <i>tonA</i> , <i>tsx</i> , <i>dam</i> , <i>dcm</i> , <i>supE44</i> , Δ (<i>lac-proAB</i>), [F', <i>traD36</i> , <i>proAB</i> , <i>lacI^a</i> Δ M15]
KW251	<i>supE44</i> , <i>galK2</i> , <i>galT22</i> , <i>metB1</i> , <i>hsdR2</i> , <i>mcrB1</i> , <i>mcrA</i> , [<i>argA81</i> ::Tn10(<i>tet^r</i>)], <i>recD1014</i>
LE392 (6)	<i>hsdR514</i> , (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>supE44</i> , <i>supF58</i> , <i>lacY1</i> or Δ (<i>lacIZY</i>)6, <i>galK2</i> , <i>galT22</i> , <i>metB1</i> , <i>trpR55</i>
NM522 (7)	<i>supE</i> , <i>thi</i> , Δ (<i>lac-proAB</i>), Δ <i>hsd5</i> (<i>r_K⁻</i> , <i>m_K⁻</i>), [F', <i>proAB</i> , <i>lacI^a</i> Δ M15]
NM538 (8)	<i>supF</i> , <i>hsdR</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>trpR</i> , <i>lacY</i>
NM539 (8)	<i>supF</i> , <i>hsdR</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>lacY</i> , (P2)
*Select96 TM	<i>mcrA</i> , Δ (<i>mrr-hsdRMS-mcrBC</i>), ϕ 80 <i>lacZ</i> Δ M15, Δ <i>lacX74</i> , <i>recA1</i> , <i>araD139</i> (<i>ara-leu</i>)7697, <i>galU</i> , <i>galK</i> , <i>rpsL</i> , <i>endA1</i> , <i>nupG</i>
Stb12 TM	F ⁻ , <i>mcrA</i> , Δ (<i>mcrBC-hsdRMS-mrr</i>), <i>recA1</i> , <i>endA1</i> , <i>gyrA96</i> , <i>thi-1</i> , <i>supE44</i> , <i>relA1</i> , λ^- , Δ (<i>lac-proAB</i>)
Stb14 TM	<i>mcrA</i> , Δ (<i>mcrBC-hsdRMS-mrr</i>), <i>recA1</i> , <i>endA1</i> , <i>gyrA96</i> , <i>thi-1</i> , <i>supE44</i> , <i>relA1</i> , λ^- , Δ (<i>lac-proAB</i>), <i>gal</i> , F' { <i>proAB⁺</i> , <i>lacI^a</i> , Δ M15, <i>Tn10</i> (<i>tet^R</i>)}
SURE [®]	<i>e14-</i> , (<i>mcrA-</i>) Δ (<i>mcrCB-hsdSMR-mrr</i>)171, <i>endA1</i> , <i>supE44</i> , <i>thi-1</i> , <i>gyrA96</i> , <i>relA1</i> , <i>lac</i> , <i>recB</i> , <i>recJ</i> , <i>sbcC</i> , <i>umuC</i> ::Tn5 (<i>kan^r</i>), <i>uvrC</i> , [F' <i>proAB</i> , <i>lacI^a</i> Δ M15::Tn10 (<i>tet^r</i>)]
TOP10	F ⁻ , <i>mcrA</i> , Δ (<i>mrr-hsdRMS-mcrBC</i>), ϕ 80 <i>lacZ</i> Δ M15, Δ <i>lacX74</i> , <i>deoR</i> , <i>recA1</i> , <i>araD139</i> , Δ (<i>ara, leu</i>)7697, <i>galU</i> , <i>galK</i> , <i>rpsL</i> (<i>str^R</i>), <i>endA1</i> , <i>nupG</i>
TOP10F'	F' { <i>lacI^a</i> Tn10 (<i>tet^R</i>)}, <i>mcrA</i> , Δ (<i>mrr-hsdRMS-mcrBC</i>), ϕ 80 <i>lacZ</i> Δ M15, Δ <i>lacX74</i> , <i>deoR</i> , <i>recA1</i> , <i>araD139</i> , Δ (<i>ara-leu</i>)7697, <i>galU</i> , <i>galK</i> , <i>rpsL</i> (<i>str^r</i>), <i>endA1</i> , <i>nupG</i>
XL1-Blue	<i>recA1</i> , <i>endA1</i> , <i>gyrA96</i> , <i>thi-1</i> , <i>hsdR17</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>supE44</i> , <i>relA1</i> , <i>lac</i> , [F', <i>proAB</i> , <i>lacI^a</i> Δ M15::Tn10(<i>tet^r</i>)]
Y1089 (9)	Δ (<i>lacU169</i>), <i>proA⁺</i> , Δ (<i>lon</i>), <i>araD139</i> , <i>strA</i> , <i>hflA150</i> , [<i>chr</i> ::Tn10(<i>tet^r</i>)], (pMC9)
Y1090 (9)	Δ (<i>lacU169</i>), <i>proA⁺</i> , Δ (<i>lon</i>), <i>araD139</i> , <i>strA</i> , <i>supF</i> , <i>rpsL</i> (<i>str^r</i>), [<i>trpC22</i> ::Tn10 (<i>tet^r</i>)], (pMC9), <i>hsdR</i> (<i>r_K⁻</i> , <i>m_K⁺</i>)

Miscellaneous

F' Host contains an F' episome with the stated features

λ (DE3) Bacteriophage λ carrying the gene for T7 RNA polymerase is integrated into the host genome.

pMC9 is pBR322 with *lacI^a* inserted and confers amp and tet resistance.

References

- Jendrisak, J., Young, R.A. and Engel, J. (1987) In: *Guide to Molecular Cloning Techniques*, Berger, S. and Kimmel, A., eds., Academic Press, San Diego, CA.
- Hanahan, D. (1983) *J. Mol. Biol.* **166**, 557–80.
- Lorow-Murray, D. and Bloom, F. (1991) *Focus* **13**, 20.
- Lacks, S. and Greenberg, J.B. (1977) *J. Mol. Biol.* **114**, 153–60.
- Yanisch-Perron, C., Vieira, J. and Messing, J. (1985) *Gene* **33**, 103–19.
- Murray, N. et al. (1977) *Mol. Gen. Genet.* **150**, 53–61.
- Gough, J. and Murray, N. (1983) *J. Mol. Bio.* **166**, 1–19.
- Frischauf, A. et al. (1983) *J. Mol. Biol.* **170**, 827–42.
- Huynh, T., Young, R.A. and Davis, R. (1985) In: *DNA Cloning*, Vol. 1, Glover, D., ed., IRL Press Ltd., Oxford, UK.

Technical Appendix

Genetic Markers in *E. coli*.

Symbol	Description	Effect of Mutation
<i>ara-14</i>	Mutation in arabinose metabolism	Blocks arabinose catabolism.
<i>araD</i>	L-ribulose phosphate 4-epimerase mutation; part of an inducible operon <i>araBAD</i> repressed by L-arabinose	Blocks arabinose catabolism.
<i>argA</i>	N-Acetylglutamate synthase mutation; inhibited by the presence of arginine	Arginine required from growth in minimal media.
<i>cycA</i>	Involved in D-alanine, glycine, D-serine and D-cycloserine transport, and an L-alanine carrier	Mutants cannot use D-alanine as a carbon source.
<i>dam</i>	DNA adenine methylase mutation	Blocks methylation of adenine residues in the sequence 5'...G ^m ATC...3'.
<i>dapD</i>	Succinyl-diaminopimelate aminotransferase mutation	Mutant reflects impaired synthesis of succinyl CoA and needs to be supplemented with succinate or lysine + methionine.
<i>dcm</i>	DNA cytosine methylase mutation	Blocks methylation of cytosine in the sequence 5'...C ^m CAGG...3' or 5'...C ^m CTGG...3'.
<i>deoC</i>	Deoxyribose-phosphate aldolase mutation	
<i>deoR</i>	Regulatory gene mutation allowing constitutive expression of genes for deoxyribose synthesis	Allows efficient propagation of large plasmids.
<i>dut1</i>	Mutation of deoxyuridine triphosphatase, which catalyzes dUTP the conversion to dUMP and PPI	Mutants are impaired in conversion of dUTP to dUMP, leading to higher dUTP pools that can lead to misincorporation of uracil instead of thymidine. Stable incorporation of dUTP needs mutation in <i>ung</i> gene.
<i>endA1</i>	DNA-specific endonuclease I mutation	Improves quality of plasmid DNA isolations.
<i>galE</i>	Part of the <i>galETK</i> operon that encodes UDP galactose-4-epimerase	Mutant is more resistant to bacteriophage P1 infection.
<i>galK</i>	Galactokinase mutation	Blocks catabolism of galactose.
<i>galT</i>	Galactose-1-phosphate uridylyltransferase mutation	Blocks catabolism of galactose.
<i>gyrA96</i>	DNA gyrase mutation	Confers resistance to nalidixic acid.
<i>hflA150</i>	Protease mutation that leads to stabilization of <i>cll</i> gene products	Leads to high frequency of lysogeny by λ phages (1).
<i>hflB</i>	Gene encodes a possible protease component	Mutations lead to high frequency of bacteriophage lambda lysogenization.
<i>hsdR</i> (r_K^- , m_K^+)	Host DNA restriction and methylation system mutation: Restriction minus, modification positive for the <i>E. coli</i> K strain methylation system	Allows cloning without cleavage of transformed DNA by endogenous restriction endonucleases. DNA prepared from this strain can be used to transform r_K^+ <i>E. coli</i> strains.
<i>hsdS20</i> (r_B^- , m_B^-)	Mutation of specificity determinant for host DNA restriction and methylation system. Restriction minus, modification minus for the <i>E. coli</i> B strain methylation system	Allows cloning without cleavage of transformed DNA by endogenous restriction endonucleases. DNA prepared from this strain is unmethylated by the <i>hsdS20</i> methylases.
<i>lacI^q</i>	Overproduction of the <i>lac</i> repressor protein	Leads to high levels of the <i>lac</i> repressor protein, inhibiting transcription from the <i>lac</i> promoter.
<i>lacY</i>	Galactoside permease mutation	Blocks lactose utilization.
<i>lacZΔM15</i>	Partial deletion of β -D-galactosidase gene	Allows complementation of β -galactosidase activity by α -complementation sequence in pGEM [®] -Z Vectors. Allows blue/white selection for recombinant colonies when plated on X-Gal.
<i>leuB</i>	β -isopropylmalate dehydrogenase mutation	Requires leucine for growth on minimal media.
Δ (<i>lon</i>)	Deletion of <i>lon</i> protease	Reduces proteolysis of expressed proteins.
<i>LysS</i>	pLysS plasmid is integrated into the host genome	Strains carrying this plasmid will be tet resistant and produce T7 lysozyme, a natural inhibitor of T7 RNA polymerase, thus lowering background transcription of sequences under the control of the T7 RNA polymerase promoter (2).
<i>mcrA</i>	Mutation in methylcytosine restriction system	Blocks restriction of DNA methylated at the sequence 5'...G ^m CGC...3'.
<i>mcrB</i>	Mutation in methylcytosine restriction system	Blocks restriction of DNA methylated at the sequence 5'...AG ^m CT...3'.
<i>metB</i>	Cystathionine γ -synthase mutation	Requires methionine for growth on minimal media.
<i>metC</i>	Cystathionine beta-lyase mutation; involved in methionine biosynthesis	Methionine required from growth in minimal media.
<i>mtl</i>	Mutation in mannitol metabolism	Blocks catabolism of mannitol.
<i>mutS</i>	Methyl-directed mismatch repair mutation	Prevents repair of the newly synthesized, unmethylated strand.
<i>ompT</i>	Mutation of protease VII, an outer membrane protein	Reduces proteolysis of expressed proteins.
P2	P2 bacteriophage lysogen present in host	λ phages containing the <i>red</i> and <i>gam</i> genes of λ are growth inhibited by P2 lysogens (3).
<i>proA</i>	γ -glutamyl phosphate reductase mutation	<i>proA/argD</i> mutant will not block proline synthesis, but will be repressed by arginine. Mutants excrete proline on minimal media and are resistant to proline analogs. <i>proA/argD/argR</i> triple mutant grows slowly on minimal media + arginine.
<i>proAB</i>	Mutations in proline metabolism	Requires proline for growth in minimal media.

Technical Appendix

Genetic Markers in *E. coli* (continued).

Symbol	Description	Effect of Mutation
<i>recA1</i> , <i>recA13</i>	Mutation in recombination	Minimizes recombination of introduced DNA with host DNA, increasing stability of inserts. Inserts are more stable in <i>recA1</i> than <i>recA13</i> hosts.
<i>recB</i> , <i>recC</i> <i>recD</i>	Exonuclease V mutations The Rec BCD trimer (exonuclease V) progressively degrades ssDNA and dsDNA in an ATP-dependent manner to form oligonucleotides; implicated in homologous recombination	Reduces general recombination and affects repair of radiation damage. Allows easier propagation of sequences with inverted repeats.
<i>recF</i>	Recombination and repair mutation	Mutant cannot repair daughter strand gaps (post-replicative repair).
<i>relA</i>	ppGpp synthetase I mutation, a novel nucleotide guanosine 5'-diphosphate-3'-diphosphate produced in response to starvation by <i>relA</i> ribosomal protein sensing uncharged tRNA	Allows RNA synthesis in the absence of protein synthesis.
<i>rha</i>	Utilization of L-rhamnose, a methylpentose	Blocks rhamnose catabolism.
<i>rpsL</i>	Mutation in subunit S12 of 30S ribosome	Confers resistance to streptomycin.
<i>sbcB</i>	Exonuclease I mutation	Allows general recombination in <i>recBC</i> mutant strains.
<i>strA</i>	Mutant alters ribosome protein S12	Confers resistance to streptomycin
<i>supB</i> , <i>supC</i> , <i>supG</i> , <i>supL</i> , <i>supM</i> , <i>supN</i> , <i>supO</i>	Suppressor mutations	Suppresses ochre (UAA) and amber (UAG) mutations.
<i>supD</i> , <i>supE</i> , <i>supF</i>	Suppressor mutations	Suppresses amber (UAG) mutations.
<i>thi-1</i>	Mutation in thiamine metabolism	Thiamine required for growth in minimal media.
<i>thr</i>	Threonine biosynthesis mutation	Mutants are obligate threonine auxotrophs.
<i>thyA</i>	Thymidylate synthase; dTTP biosynthesis	Mutants are obligate thymidine auxotrophs.
Tn5	Transposon	Encodes resistance to kanamycin.
Tn10	Transposon	Encodes resistance to tetracycline.
<i>tonA</i>	Mutation in outer membrane protein	Confers resistance to bacteriophage T1.
<i>traD36</i>	Transfer factor mutation	Prevents transfer of F' episome.
<i>trpC</i>	Phosphoribosyl anthranilate isomerase mutation; part of tryptophan biosynthesis pathway	
<i>trpR</i>	<i>trpR</i> aporepressor; regulates the biosynthesis of tryptophan and its transport	
<i>tsx</i>	T6 and colicin K phage receptor; outer membrane protein involved in specific diffusion of nucleosides; transports the antibiotic albicidin	Resistant to bacteriophage T6 and colicin K.
<i>ung1</i>	Uracil-DNA N-glycosylase	Allows uracil to exist in plasmid DNA.
<i>xyl-5</i>	Mutation in xylose metabolism	Blocks catabolism of xylose.

References

1. Hoyt, M.A. *et. al.* (1982) *Cell* **31**, 565–73.
2. Studier, F.W. (1991) *J. Mol. Biol.* **219**, 37–44.
3. Kaiser, K. and Murray, N. (1985) In: *DNA Cloning*, Vol. 1, Glover, D., ed., IRL Press Ltd., Oxford, UK.
4. Neidhardt, F. ed. (1996) *Escherichia coli and Salmonella Cellular and Molecular Biology* 2nd ed, ASM Press, Washington, D.C.

Technical Appendix

Nucleic Acids and Proteins: Calculations.

An online calculator for these values is available in the "tools" section of Promega's Web site at: www.promega.com/techserv/tools/

Metric Prefixes

Prefix	Symbol	Factor
kilo	k	10 ³
centi	c	10 ⁻²
milli	m	10 ⁻³
micro	μ	10 ⁻⁶
nano	n	10 ⁻⁹
pico	p	10 ⁻¹²
femto	f	10 ⁻¹⁵
atto	a	10 ⁻¹⁸
zepto	z	10 ⁻²¹

Spectrophotometric Conversions

- 1 A₂₆₀ unit of double-stranded DNA = 50 μg/ml
- 1 A₂₆₀ unit of single-stranded DNA = 33 μg/ml
- 1 A₂₆₀ unit of single-stranded RNA = 40 μg/ml

DNA Molar Conversions

- 1 μg of 1,000bp DNA = 1.52 pmol (3.03 pmol of ends)
- 1 μg of pBR322 DNA = 0.36 pmol DNA
- 1 pmol of 1,000bp DNA = 0.66 μg
- 1 pmol of pBR322 DNA = 2.8 μg

Formulas for DNA Molar Conversions

For dsDNA:

To convert pmol to μg:

$$\text{pmol} \times N \times \frac{660 \text{ pg}}{\text{pmol}} \times \frac{1 \mu\text{g}}{10^6 \text{ pg}} = \mu\text{g}$$

To convert μg to pmol:

$$\mu\text{g} \times \frac{10^6 \text{ pg}}{1 \mu\text{g}} \times \frac{\text{pmol}}{660 \text{ pg}} \times \frac{1}{N} = \text{pmol}$$

where N is the number of nucleotide pairs and 660 pg/pmol is the average MW of a nucleotide pair.

For ssDNA:

To convert pmol to μg:

$$\text{pmol} \times N \times \frac{330 \text{ pg}}{\text{pmol}} \times \frac{1 \mu\text{g}}{10^6 \text{ pg}} = \mu\text{g}$$

To convert μg to pmol:

$$\mu\text{g} \times \frac{10^6 \text{ pg}}{1 \mu\text{g}} \times \frac{\text{pmol}}{330 \text{ pg}} \times \frac{1}{N} = \text{pmol}$$

where N is the number of nucleotides and 330 pg/pmol is the average MW of a nucleotide.

Dalton (Da) is an alternate name for the atomic mass unit, and kiloDalton (kDa) is 1,000 Daltons. Thus a protein with a mass of 64 kDa has a molecular weight of 64,000 grams per mole.

- ^(a) The PCR process is covered by patents issued and applicable in certain countries*. Promega does not encourage or support the unauthorized or unlicensed use of the PCR process. *In the U.S., effective March 29, 2005, U.S. Pat. Nos. 4,683,195, 4,965,188 and 4,683,202 will expire. In Europe, effective March 28, 2006, European Pat. Nos. 201,184 and 200,362 will expire.
- ^(b) Purchase of this product is accompanied by a limited license under U.S. Pat. Nos. 5,082,784 and 5,192,675 for the internal research use of the computer.
- ^(c) Turbo™ Nae I and Turbo™ Nar I are the subjects of U.S. Pat. Nos. 5,248,600 and 5,418,150 and are licensed exclusively to Promega Corporation, as well as a license under DD 264 231.
- ^(d) Licensed using U.S. Pat. No. 4,935,361.
- ^(e) Licensed under U.S. Pat. No. 5,7075.
- ^(f) Licensed using one or more of U.S. Pat. Nos. 5,487,993 and 5,827,657 and European Pat. No. 0 550 693.
- ^(g) U.S. Pat. No. 4,766,072.
- ^(h) The PCR process, which is the subject of European Pat. Nos. 201,184 and 200,362 and U.S. Pat. Nos. 4,683,195, 4,965,188 and 4,683,202 owned by Hoffmann-LaRoche*, is covered by patents issued and applicable in certain countries. Promega does not encourage or support the unauthorized or unlicensed use of the PCR process. Use of this product is recommended for persons that either have a license to perform PCR or are not required to obtain a license. *In the U.S., effective March 29, 2005, the above primary U.S. Pat. Nos. 4,683,195, 4,965,188 and 4,683,202 will expire. In Europe, effective March 28, 2006, the above primary European Pat. Nos. 201,184 and 200,362 will expire.
- ⁽ⁱ⁾ Certain applications of this product are covered by patents issued and applicable in certain countries. Because purchase of this product does not include a license to perform any patented application, users of this product may be required to obtain a patent license depending upon the particular application and country in which the product is used.
- ^(j) Australian Pat. No. 730718 and other patents and patents pending.
- ^(k) U.S. Pat. No. 5,981,235, Australian Pat. No. 729932 and other patents pending.

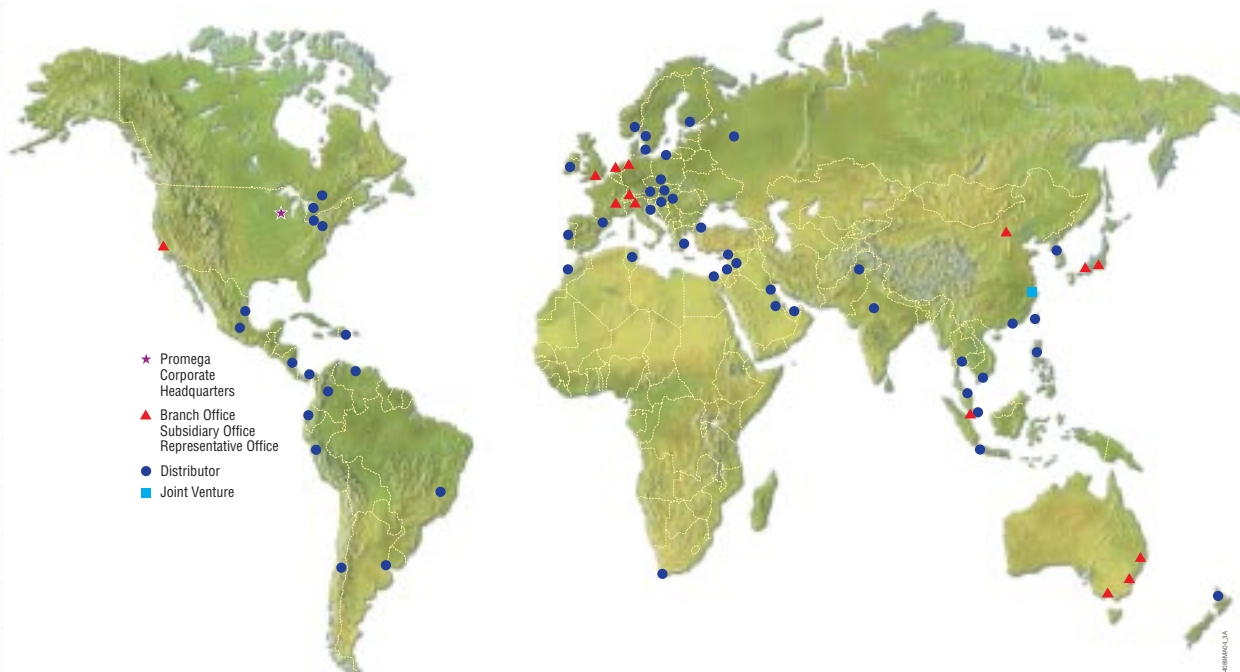
©2004 Promega Corporation. All Rights Reserved.

All prices and specifications are subject to change without prior notice. Product claims are subject to change. Please contact Promega Technical Services or access Promega online for the most up-to-date information on Promega products.

GeneEditor, LigaFast, MULTICORE, Select96, TagBead, and Turbo are trademarks of Promega Corporation. 4-CORE, Altered Sites, GoTaq, pALTER, pGEM, VacMan and Wizard are trademarks of Promega Corporation and are registered with the U.S. Patent and Trademark Office.

ABLE and SURE are registered trademarks of Stratagene. Bacto is a registered trademark of Difco Laboratories, Detroit, Michigan. DH5α is a trademark of Life Technologies, Inc. Ficoll is a registered trademark of Amersham Biosciences Ltd. Stb12 adn Stb14 are trademarks of Invitrogen Corporation.

Contact Us



Promega Corporation

2800 Woods Hollow Road
Madison, WI 53711-5399 USA
Tel: 608-274-4330
Fax: 608-277-2516
Toll-Free: 800-356-9526
Toll-Free Fax: 800-356-1970
Internet: www.promega.com

Promega Biosciences, Inc.

A Division of Promega Corporation
San Luis Obispo, California

Australia, Sydney

Tel: 02 9565 1100
Fax: 02 9550 4454
Freecall: 1800 225 123
Freefax: 1800 626 017
E-mail: aus_custserv@au.promega.com

China, Beijing

Tel: 10 6849 8287
Fax: 10 6849 8390
E-mail: promega@promega.com.cn

France, Lyon

Tel: 04 37 22 50 00
Fax: 04 37 22 50 10
Numero Vert: 0 800 48 79 99
E-mail: fr_custserv@fr.promega.com

Germany/Austria, Mannheim

Tel: (+49) (0) 621-8501-0
Fax: (+49) (0) 621-8501-222
Free Phone: 00800-77663422
Free Fax: 00800-77663423
E-mail: de_custserv@de.promega.com

Italy, Milan

Tel: 02 54 05 01 94
Fax: 02 55 18 56 64
Numero Verde: 800 69 18 18
E-mail: it_custserv@it.promega.com

Japan, Tokyo

Tel: 03-3669-7981
Fax: 03-3669-7982
E-mail: prometec@jp.promega.com

Belgium/Luxembourg/ The Netherlands, Leiden

Tel: (+31) (0) 71-5324244
Fax: (+31) (0) 71-5324907
Free Tel BE: 0800-18098
Free Fax BE: 0800-16971
Free Tel NL: 0800-0221910
Free Fax NL: 0800-0226545
E-mail: bnl_custserv@nl.promega.com

Pacific Asia Region, Singapore

Tel: 65 6254 5265
Fax: 65 6254 8645
E-mail: nicholas.ng@promega.com

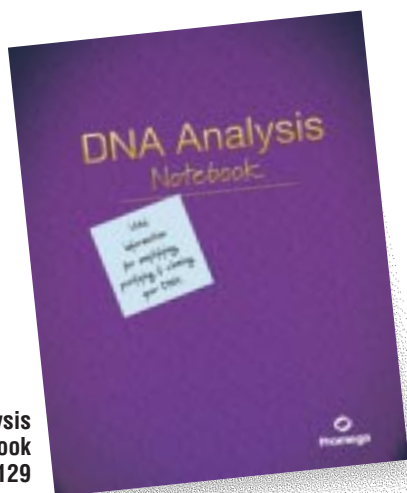
Switzerland, Wallisellen

Tel: 044 878 90 00
Fax: 044 878 90 10
Technical Service: 044 878 90 20
E-mail: catalys_custserv@promega.com

United Kingdom, Southampton

Tel: 023 8076 0225
Fax: 023 8076 7014
Free Phone: 0800 378994
Free Fax: 0800 181037
E-mail: ukcustserve@promega.com

Learn More. Ask Promega!



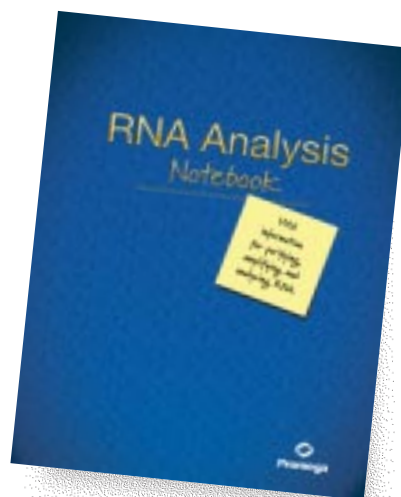
DNA Analysis
Notebook
#BR129

DNA Analysis Notebook

Get the latest information
to help you purify genomic DNA,
amplify and clone your targets.

RNA Analysis Notebook

Isolate RNA and analyze by
RT-PCR, microarrays
and even RNAi.



RNA Analysis
Notebook
#BR120

**Call your local Promega
Branch Office or Distributor
to get your copies.**

www.promega.com

Toll Free in USA: 800-356-5926



Promega Corporation • 2800 Woods Hollow Road • Madison, WI 53711-5399 USA • Telephone 608-274-4330 • Fax 608-277-2601

©2004 Promega Corporation. All Rights Reserved.
Prices and specifications subject to change
without prior notice.

Printed in USA 08/04
11469-BR-PU
Part #BR152



Promega
www.promega.com