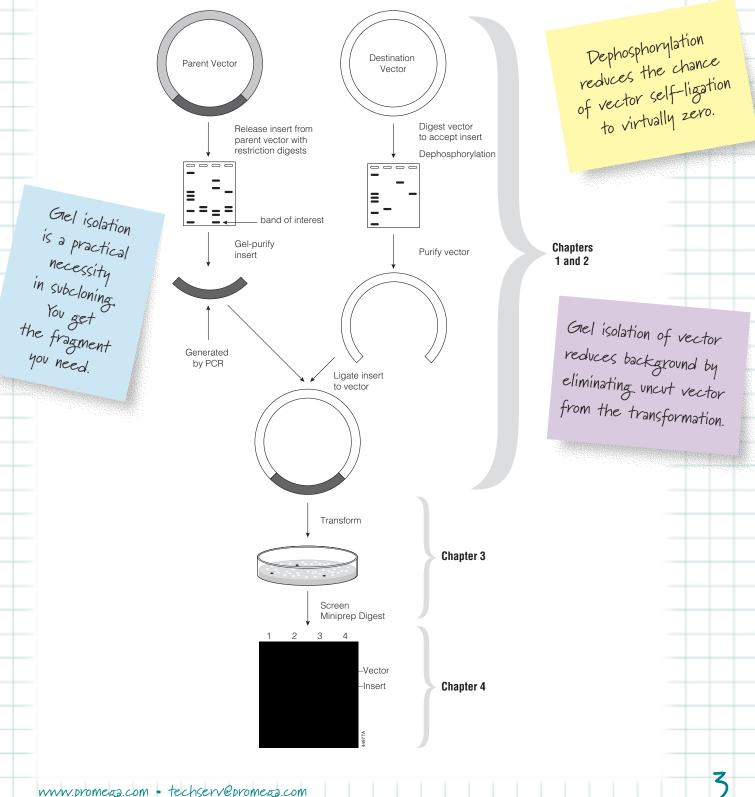
Basic Steps for Subcloning

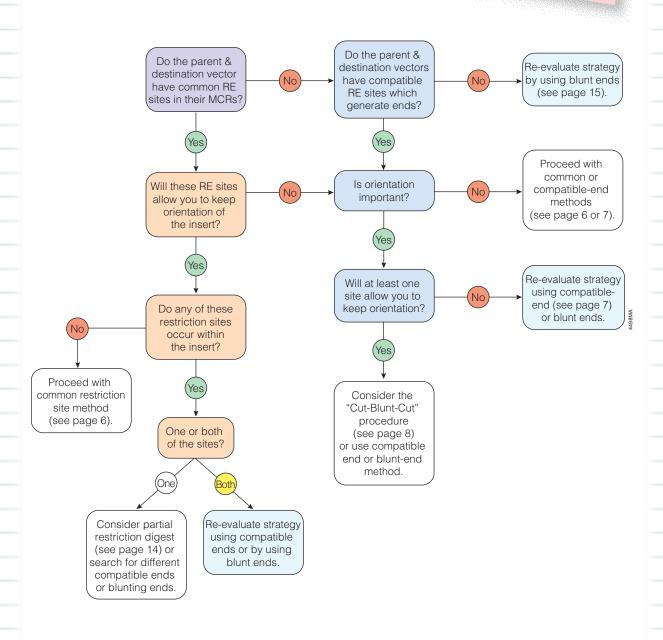
Subcloning is a basic procedure in molecular biology required to move inserts from one vector to another to gain the desired functionality to study your insert. Essentially all subcloning reactions proceed the same way as illustrated in the figure below. You release and purify your insert from the parent 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.



Subcloning Strategy

4

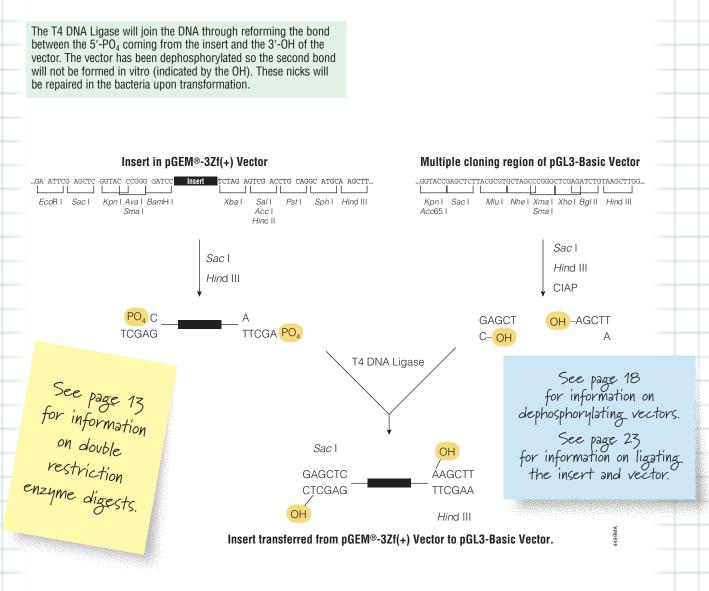
Before you begin your subcloning, you need to know: The restriction enzyme (RE) sites available for subcloning in your parent vector multiple cloning region (or in the insert if you need to digest the insert); the RE sites available in the destination vector multiple cloning region (MCR); and if these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



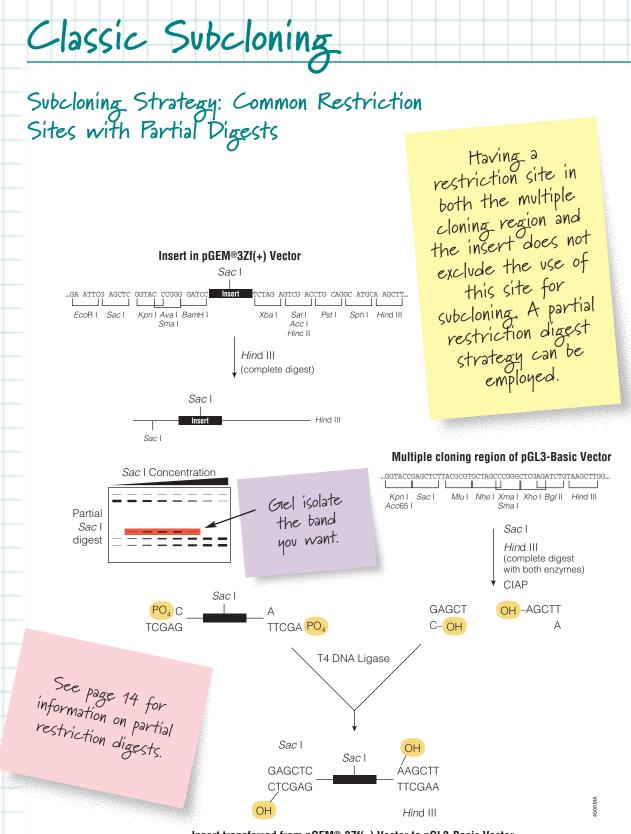
See the Compatible Ends Table on page 61 for a listing of overhangs compatible with Promega enzymes.

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.



5



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

6

Subcloning Strategy: Moving Inserts with Compatible Restriction Sites

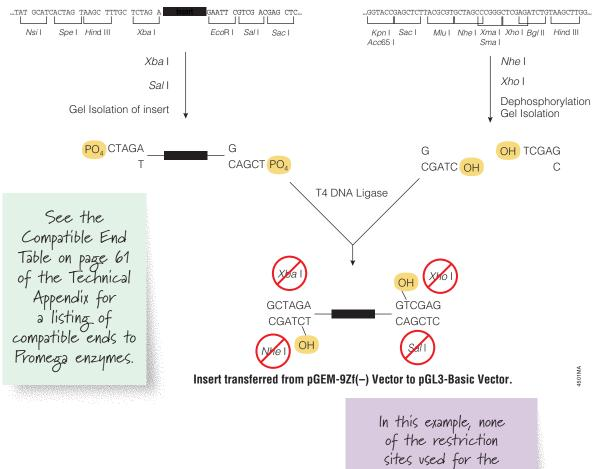
If you don't have common restriction sites in the parent and destination vector multiple cloning regions, you may have compatible restriction sites. Compatible restriction sites have the same overhang sequence and can be ligated together. In this example, *Xba* I and *Nhe* I both produce the same 5' 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 pGIL3-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 PGIL3-Basic Vector.

Multiple cloning region of pGL3-Basic Vector

Insert in pGEM®-9Zf(-) Vector

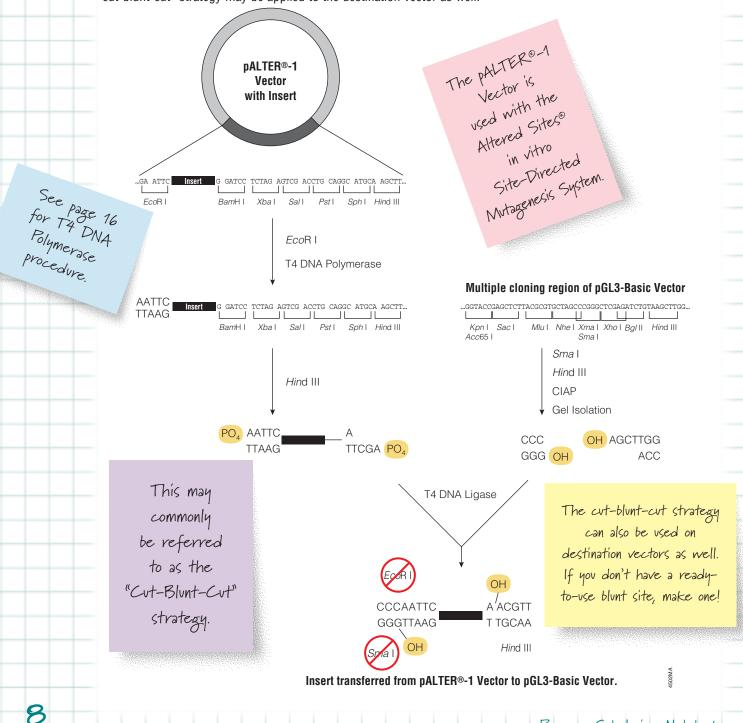


7

compatible-end subcloning are regenerated in the final ligation product.

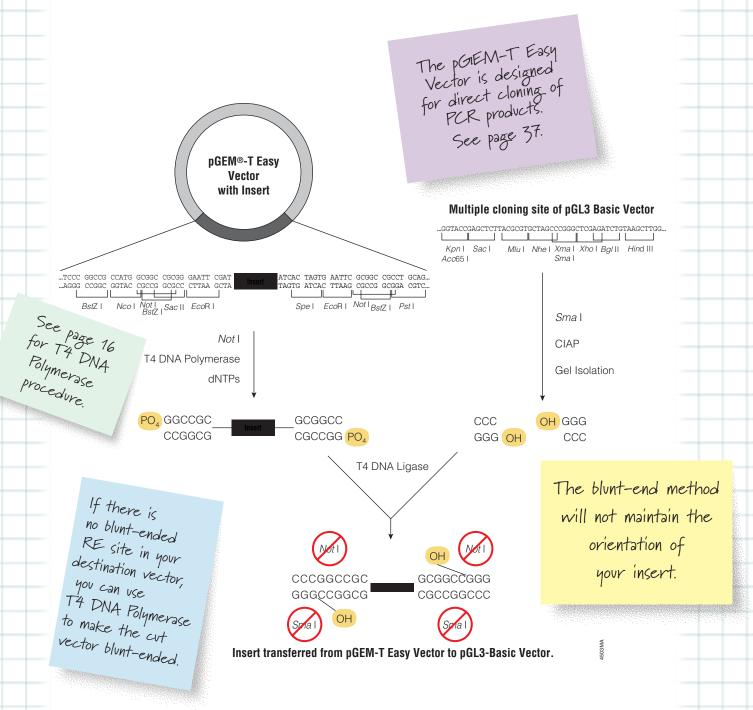
Subcloning Strategy: Moving Inserts with Only One Common Site

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.



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.



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:

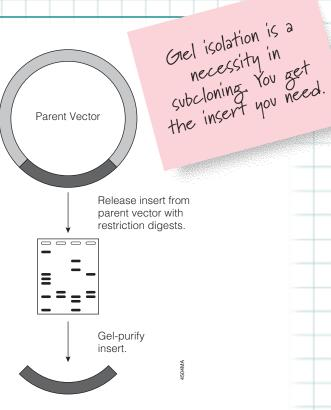
Nuclease-Free Water	14µl
10X Restriction Buffer	2µl
Acetylated BSA (1mg/ml)	2µl
DNA (~1µg)	1µl
Restriction Enzyme (10u)	1µl
Final Volume	20µl

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 4μ I 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).

10



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

Bard Pressage	Rendarita	Technical Reserves	CTTT:
school Reported	Restriction Enzy	mes Resource	-
lators		 that are meat useful for molecular blots requirement specificity is precision outto 	
Bolese,	Assource to an indepaction	e tool, designed to speed your research a	
adverse.		We offer four sourches, elon, as well as background it reference is	the state of the sector the
A()+	enzymes.		
	Searches		
605	SEARCH BY	DESCRIPTION	
Todoram	Enryme Name	Find information on recognition site, in resectionmer(s) and type of overhal	
Utilications	Reception Recentle	Liestly this that reception a specific	nucleolide sequence.
tours	Enzymes Generating Compatible Indo	Find compatible REs by the type of e blunt) and overhang sequence.	nd generated (overhand or
Fectores	Compatible Buffers	Identify a reaction buffer for annuita-	would, multiple restriction digests
aritani UAF Infranty Opeophica Look 4003-358-9526		(Promego erstytuns only).	
	Background & Ref	erence Information	



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.

Alexandreament Alexandr	ER ER	Usage Information
No. Image: Display in the second		 I. N. Operation of the second state of the second sta

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., *BstX* I and *BstZ* 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 DH5 α^{TM} 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?

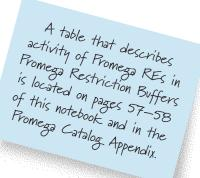
- Check to see if the enzyme has an isoschizomer or neoschizomer. The isoschizomer or neoschizomer may not be sensitive to the methylation.
- Transform the plasmid into a dam/dcm minus bacterial strain like JM110.

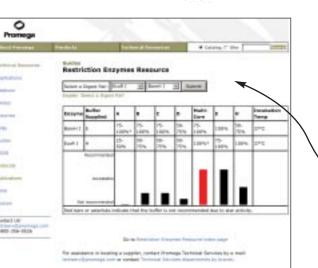
See the tables on pages 59-60 for listings of isoschizomers and neoschizomers.

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 *Eco*R I (optimal in Buffer H) and *Bam*H I (optimal in Buffer E) you would choose in Buffer E because the BamH 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.





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 15 μ l of water, and the buffer, enzyme and BSA can be added to bring the reaction to 20 μ l 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

Partial Restriction Digestion

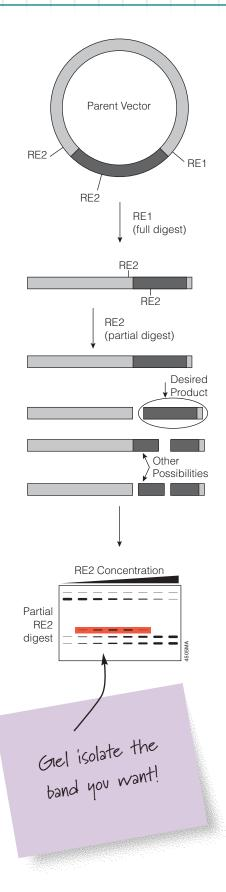
Controlling Cut Frequency in Restriction Digestion

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).
- Purify vector with the Wizard[®] SV Gel and PCR Clean-Up System directly from the reaction. Elute in 20µl nucleasefree water.
- 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 $2\mu I$ 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.

14



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.

Klenow Fragment or T4 DNA Polymerase

5'-A-3' → 5'-AGCTA-3' 3'-TCGAT-5' 3'-TCGAT-5'

Mg²⁺; dNTPs

5' Overhang Fill-In Reaction

T4 DNA Polymerase

5'-CTGCA-3'	5'-C-3'
3'-G-5'	 3'-G-5'

Mg²⁺; dNTPs

3' Overhang Blunting Reaction

1506MA

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-CORETM, 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.
- 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.
- 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 $0.1\mu g/\mu l$ 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-CORETM, 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.
- 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 Cət.# M4211 100v 5-10v/µl Cət.# M4215 500v 5-10v/µl

See the Product Information Sheet at: mmv.promega.com/tbs Note: With high concentrations of Note: With high concentrations of dNTPs (i.e., 100 µM), degradation of the DNA will stop at the duplex of the DNA will stop at the duplex active provided the highly active exonuclease DNA. However, if the difference than DNA. However, if the difference than DNA. However, if the difference than active the highly active exonuclease that of DNA polymerase i) of T4 DNA that of DNA polymerase i) of T4 DNA. Heat of DNA polymerase the dsDNA.

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.
- Incubate at ambient room temperature for 10 minutes.
- 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



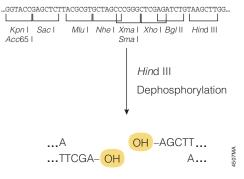
Dephosphorylating Vectors to Limit Self-Ligation

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 bluntends. 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.

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.

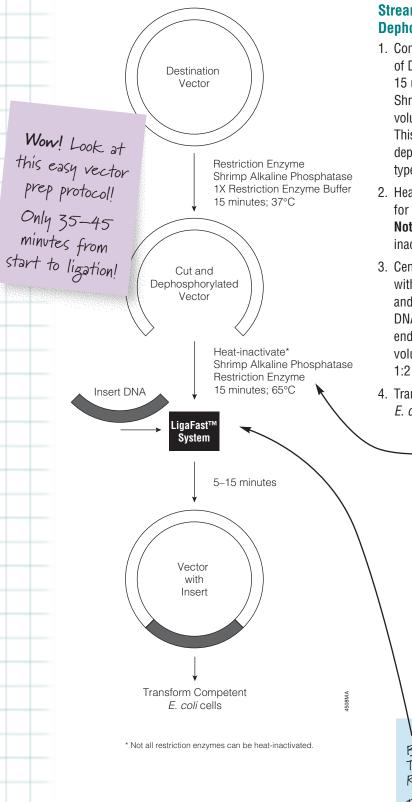
Multiple cloning region of pGL3-Basic Vector



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

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

Dephosphorylating Vectors: Shrimp Alkaline Phosphatase



Streamlined Restriction Digestion, Dephosphorylation and Ligation Procedure

- 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 30–50μl. 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).

- Centrifuge and remove 1–2µl 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 10–50µl. 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 Giel and PCR
 Clean-Up System for direct purification.
 Full purification in just 15 minutes, and you can elute the DNA in as little as 15µl 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.

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:

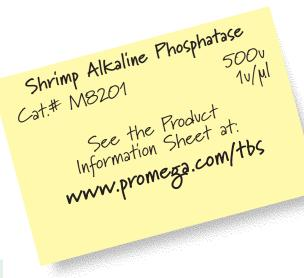
DNA (1–2µg)	ΧμΙ
10X SAP Buffer	3–5µl
SAP (1u/μl)	1µl/µg DNA
Nuclease-Free Water	to 30–50µl

- 3. Incubate at 37°C for 15 minutes (works for both 5' and 3' overhangs or blunt ends).
- 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

SAP

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



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: 0.015v SAP/pmol ends

Blunt Overhang: 0.031 SAP/pmol ends

3' Overhang: 0.4u SAP/pmol ends

Dephosphoylating 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:

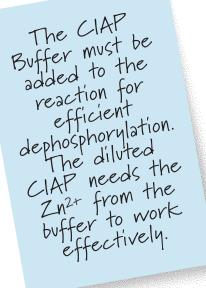
5' Overhangs: Incubate for 30 minutes at 37°C. Add another 0.01u CIAP/pmol ends and incubate an additional 30 minutes at 37°C.

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.

Calculating pmol of DNA from micrograms of DNA. $\mu g \text{ DNA} \times \frac{\text{pmol}}{660 \text{pg}} \times \frac{10^6 \text{pg}}{1 \mu g} \times \frac{1}{N} = \text{pmol DNA}$

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

Dephosphorylating Vectors: Calf Intestinal Alkaline Phosphatase

Dephosphorylation of Purified DNA

- 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 $(1u/\mu l)$ 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.

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

22

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.



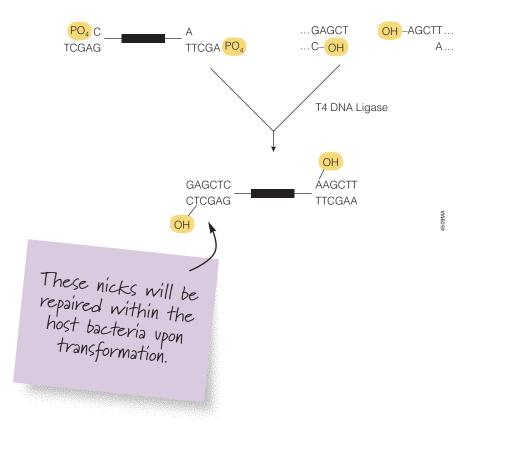
Ligation: Ligating Vector and Insert

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



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

ng of vector \times kb size of insert	× molar ratio of	insert	– na of insert
kb size of vector		vector	- ng or moore

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.

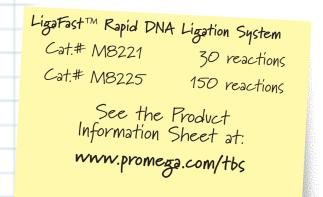
100ng vector \times 0.5kb insert	2	= 33.3ng insert
3kb vector	1	- 00.011g 11.0011

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.

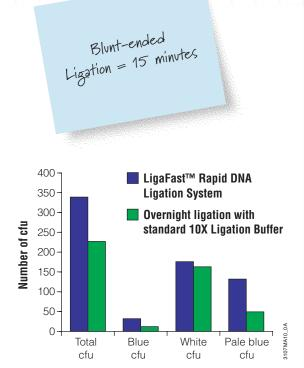
1. Assemble the following reaction in a sterile microcentrifuge tube:

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

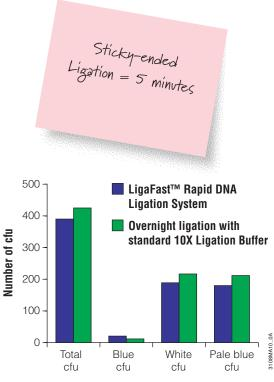
 Incubate the reaction at room temperature for 5 minutes for cohesive-ended ligations, or 15 minutes for blunt-ended ligations.



24



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 *Eco*R 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.



Comparison of overnight ligations and the LigaFast[™] Rapid DNA Ligation System using a DNA insert with 5' overhangs. Experiment performed with bluntend insert ligated into an *Sal* 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.

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	1µI
T4 DNA Ligase (3u/µl)	1µI
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 aliguot.

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

T4 DNA Ligase Blue/White Cloning Qualified Cat.# M1801 1-3u/µl Cat.# M1804 1-3u/µl Cat.# M1794 500u Cat.# M1794 10-20u/µl See the Product Information Sheet at: Information Sheet at:



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 μ g of a DNA digest marker (e.g., Lambda DNA *Hin*d 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.

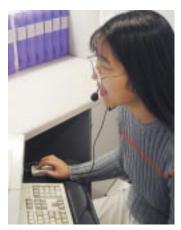
Another quick test is to cut a plasmid with a single restriction enzyme. Add this vector to a ligation reaction and transform.

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.

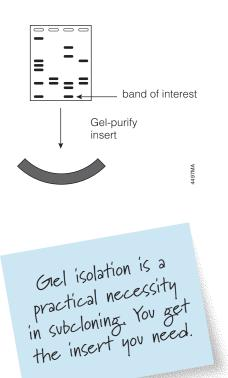
Contact Promega Technical Services directly or through your Branch Office or by email at: techserv@promega.com

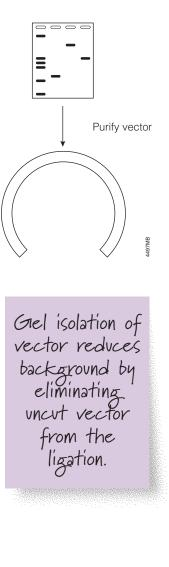
26



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.





Purifying Vector and Insert

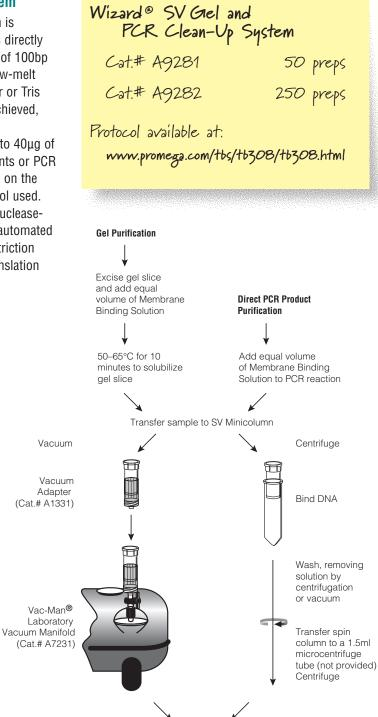
Wizard® SV Gel and PCR Clean-Up System

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 nucleasefree 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 40 µg of DNA on a single column!

From start to purified DNA in 15 minutes!

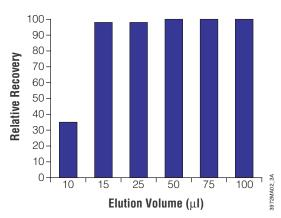


Flow chart of DNA fragment gel purification or direct PCR product purification using the Wizard SV Gel and PCR Clean-Up System. 3792MA07_2A

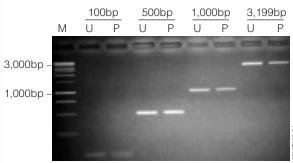
Elute DNA fragment

or PCR product

Purifying Vector and Insert



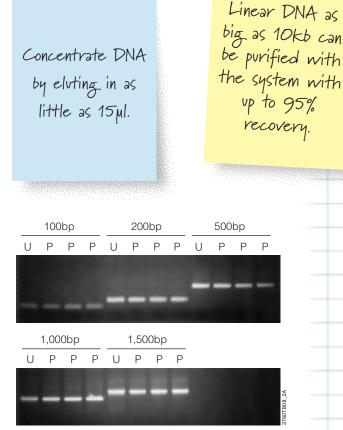
Elution volume versus recovery for a 700bp PCR product. One hundred percent is based on recovery with 50µl elution. Adapted from Table 4 in Betz, N. and Strader, T. (2002) Clean Up with Wizard® SV for Gel and PCR. Promega Notes 82, 2-5.



U: Unpurified P: Purified

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



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

> Wizard SV Giel 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.

Giel 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.
- 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.
- 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.
- 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.

 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])

Э

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 Ingof dSDNA in a band. Brown, T.A. (1998) In: Molecular Biology LABFAX II: Giene Analysis. 2nd ed. Academic Press, 101.

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 Conventional Cat.# G1741

pGEM[®] DNA Markers BenchTop Cat.# G7521

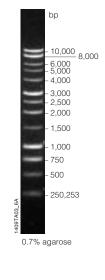
bp

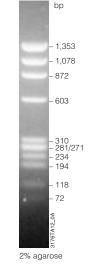
2.645

ΦX174 DNA/ Hae III Markers BenchTop Cat.# G7511 Conventional Cat.# G1761

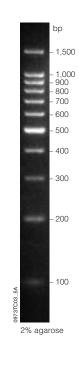
PCR Markers BenchTop Cat.# G7531 Conventional Cat.# G3161

100bp DNA Ladder BenchTop Cat.# G8291 Conventional Cat.# G2101





bp 1,000 750 500 300 150 50 2% agarose



www.promega.com - techserv@promega.com

2% agarose

31

Classic Subcloning: Ordering Information

Enzyme lr	Heat nactivated	Ruffor	Recognition Site	Size (u)	Conc. (u/µl)	Cat.#		Enzyme	Heat Inactivated	Buffor	Recognition Site	Size (u)	Conc. (u/µl)	Cat.#	
Aat II	+	J	GACGT▼C	<u>(u)</u> 50	<u>(α/μι)</u> 3–5	R6541	\$	Csp45 I	+	B	TTVCGAA	2,500	10	R6571	
nat II	Ŧ	J		250	3-5	R6545		Dde I	+/-	D	C▼TNAG	200	10	R6291	
Accl	_	G	GT▼(A/C)(T/G)AC	100	3–10	R6411		Duoi	17	D	0 • 110/10	1,000	10	R6295	
71001		ŭ		500	3–10	R6415		Dpn I	+	В	G™A▼TC	200	10	R6231	
Acc III	_	F	T▼CCGGA	200	10	R6581		(<i>Sau</i> 3A I)							
Acc65 I	+	D	G▼GTACC	1,500	10	R6921		Dra I	+	В	TTT▼AAA	2,000	10	R6271	
(<i>Kpn</i> I)							.0	Ec/HKI	+		GACNNN▼NNGTC	100	10	R7111	
AccB7 I	+		CCANNNNVNTGG	200	10	R7081	୍ବର ଜୁନ	Eco47 III	+	D	AGCVGCT	50	2-5	R6731	
Agel	+	K	A▼CCGGT	100	3-10	R7251	J 6-	<i>Eco</i> 52 (<i>Bst</i> Z)	+	L	C▼GGCCG	50	1–5	R6751	
Alul	+	B	AGVCT	500	10	R6281		EcolCR I	+	В	GAG▼CTC	1,000	10	R6951	
<i>Alw</i> 26 I	+	С	GTCTC(N) ₁ ▼	100	8–12 8–12	R6761		(Sac I)				5,000	40-80	R4954	
<i>Alw</i> 44 I		С	GTCTC(N)₅ G▼TGCAC	500 1,000	8 <u>-12</u> 10	R6765 R6771		<i>Eco</i> R I	+	Н	G▼AATTC	5,000	12	R6011	
Apa I	+ +	A		5,000	10	R6361						15,000	12	R6017	
πρατ	Ŧ	~	uuuco • c	25,000	40-80	R4364						25,000	40-80	R4014	
Aval	+/	В	C▼(T/C)CG(A/G)G	20,000	8–12	R6091						50,000	40-80	R4017	
/1/21	17	D	0 • (1/0)00(/(0)0	1,000	8–12	R6095		<i>Eco</i> R V	+	D	GAT▼ATC	2,000	10	R6351	
Ava II	+	С	G▼G(A/T)CC	100	1-10	R6131						10,000	10	R6355	
(Sin I)				1,000	1–10	R6135						10,000	40-80		
Bal I	+	G	TGG▼CCA	50	2–10	R6691		Fokl	+	В	GGATG(N)9 GGATG(N)(13)▼	100	2–10	R6781	
				250	2–10	R6695		Hae II	_	В	$(A/G)GCGC \mathbf{\nabla}(T/C)$	1,000	10	R6661	
<i>Bam</i> H I	+	Е	G▼GATCC	2,500	10	R6021		Hae III	_	C	GG▼CC	2,500	10	R6171	
				12,500	10	R6025						10,000	10	R6175	
				12,500	40-80	R4024						12,500	40-80	R4174	
				50,000	40-80	R4027		Hha I	+	С	GCG▼C	1,000	10	R6441	
Ban I	-	G	G▼G(T/C)(A/G)CC	200	8–12	R6891		(Cfol)							
Ban II	+	Е	G(A/G)GC(T/C)▼C	1,000	8–12	R6561		<i>Hin</i> c II	+	В	GT(T/C)▼(A/G)AC	200	10	R6031	
Bbu I	+	А	GCATG▼C	200	10	R6621						1,000	10	R6035	
(Sph I)				1,000	40-80	R4624						5,000	10 40-80	R6037	
Bcll	-	С	T▼GATCA	1,000	10	R6651		<i>Hin</i> d III	+	E	A▼AGCTT	1,000 5,000	40-60	R4034 R6041	
				5,000	40-80	R4654	•	<i>Timu</i> III	÷	L	A V AGUTT	15,000	10	R6045	
Bg/ I	+	D	GCCNNNN▼NGGC		10	R6071						25,000		R4044	
				5,000	10 40-80	R6077						50,000	40-80	R4047	
Bg/ II		D	A▼GATCT	5,000 500	40-60	R4074 R6081		<i>Hin</i> f I	_	В	G▼ANTC	1,000	10	R6201	
Dyr II	-	D	A V GATCT	2,500	10	R6085						5,000	10	R6205	
				10,000	10	R6087						5,000	40-80	R4204	
				2,500	40-80	R4084		Hpa I	-	J	GTT▼AAC	100	3–10	R6301	
<i>Bsa</i> M I	_	D	GAATGCN▼	500	10	R6991						500	3–10	R6305	
<i>Bsp</i> 1286 I	+	AG(G/A/T)GC(C/A/T)▼		10	R6741		Hpa II	-	A	C▼CGG	1,000	10	R6311	
BsrS I	-	D	ACTGGN	500	10	R7241		(Mspl)				5,000	10	R6315	
<i>Bss</i> H II	-	Н	G▼CGCGC	100	10	R6831		Hsp92	+		G(A/G)▼CG(T/C)C	500	10	R7151	
				500	10	R6835		Hsp92 II	+	K	CATG	1,000	10	R7161	
<i>Bst</i> 98 I	-	D	C▼TTAAG	500	8–12	R7141		I- <i>Ppo</i> I (Intron-En	+ coded Endon		CTCTCTTAA▼GGTAGC I- <i>Ppo</i> I	; 10,000	100-200	18/031	
<i>Bst</i> E II	-	D	G▼GTNACC	2,000	10	R6641		Kpn I ^(b)	+/	J	GGTAC▼C	2,500	8–12	R6341	
Bst0 I	-	С	CC▼(A/T)GG	2,000	10	R6931		(<i>Acc</i> 65 I)				10,000	8–12	R6345	
BstX I	+/	DC	CANNNNN▼NTGG		8–12	R6471						12,500	40-80	R4344	
				1,000	8–12	R6475		Mbo II	+	В	GAAGA(N) ₈	100	2-10	R6723	
BstZ I	-	D	C▼GGCCG	500	10	R6881		MAL: 1	1	r.	GAAGA(N)7	1.000	10	DCOOT	
<i>Bsu</i> 36 I	_	E	CC▼TNAGG	500	10	R6821	\$	Man	+/	D	A▼CGCGT	1,000	10	R6381	
Cfo I (Hha I)	+/	В	GCG▼C	3,000	10	R6241	9 0	Msp I (Hpa II)	+	В	C▼CGG	2,000	10 10	R6401 R6405	
Clal	+	С	AT▼CGAT	500	10	R6551		(·				10,000		R6405	
	·	Ŭ		2,500	10	R6555		MspA1 I	+	С	C(A/C)G▼C(G/T)G	1,000	40-60	R7021	
Csp I	+	K	CG▼G(A/T)CCG	100	10	R6671		Turbo™ N		 Turbo™		250	4	R7231	
				500	10	R6675	ତ		ome Qualified						
										o					

Indicates Blue/White Cloning Qualified.

e olonning Quanneu.		
Picana -	Subcloning	Matchart
IVONERA	SUCCIONING	NOTEDOOR
0	0	

	ACC III	-	F	I ▼CCGGA	200	10	K65
	<i>Асс</i> 65 I (<i>Крп</i> I)	+	D	G ▼ GTACC	1,500	10	R69
	AccB7 I	+	E	CCANNNN ▼ NTGG	200	10	R70
	Age I	+	Κ	A▼CCGGT	100	3–10	R72
	Alu I	+	В	AG▼CT	500	10	R62
	<i>Alw</i> 26 I	+	С	GTCTC(N)₁▼	100	8–12	R67
				GTCTC(N) ₅	500	8–12	R67
_	<i>Alw</i> 44 I	+	С	G▼TGCAC	1,000	10	R67
	Apa I	+	А	GGGCC▼C	5,000	10	R63
					25,000	40-80	R43
	Ava I	+/	В	$C \mathbf{\nabla}(T/C)CG(A/G)G$	200	8–12	R60
_					1,000	8–12	R60
	Ava II	+	С	G▼G(A/T)CC	100	1–10	R61
_	(<i>Sin</i> I)				1,000	1–10	R61
	Ball	+	G	TGG▼CCA	50	2–10	R66
					250	2–10	R66
	<i>Bam</i> H I	+	E	G▼GATCC	2,500	10	R60
					12,500	10	R60
					12,500	40-80	R40
_					50,000	40-80	R40
	Ban I	_	G	G▼G(T/C)(A/G)CC	200	8–12	R68
_	Ban II	+	Ε	G(A/G)GC(T/C)▼C	1,000	8–12	R65
	Bbu I	+	А	GCATG▼C	200	10	R66
	(Sph I)				1,000	40-80	R46
SO	Bc/ I	-	С	T▼GATCA	1,000	10	R66
					5,000	40-80	R46
SO.	Bgl I	+	D	GCCNNNN▼NGGC	1,000	10	R60
					5,000	10	R60
					5,000	40-80	R40
	Bg/ II	-	D	A▼GATCT	500	10	R60
					2,500	10	R60
					10,000	10	R60
					2,500	40-80	R40
	<i>Bsa</i> M I	-	D	GAATGCN▼	500	10	R69
	<i>Bsp</i> 1286 I	+	AC	G(G/A/T)GC(C/A/T)▼	C 500	10	R67
	<i>Bsr</i> S I	-	D	ACTGGN	500	10	R72
SO	<i>Bss</i> H II	-	Н	G▼CGCGC	100	10	R68
					500	10	R68
	<i>Bst</i> 98 I	-	D	C▼TTAAG	500	8–12	R714
	<i>Bst</i> E II	-	D	G▼GTNACC	2,000	10	R66
	Bst0 I	-	С	CC▼(A/T)GG	2,000	10	R69
	BstX I	+/	D	CCANNNNN▼NTGG	250	8–12	R64
					1,000	8–12	R64
_	BstZ I	-	D	C▼GGCCG	500	10	R68
୍ଦ୍ର ତ	Bsu36 I	-	E	CC▼TNAGG	500	10	R68
L	Cfo I (Hha I)	+/-	В	GCG▼C	3,000	10	R624
	Clal	+	С	AT▼CGAT	500	10	R65
					2,500	10	R65
ço	Csp I	+	Κ	CG▼G(A/T)CCG	100	10	R66
					F00	40	DCC

32

Classic Subcloning: Ordering Information

Enzym	e	Heat Inactivated	Buffer	Recognition Site	Size (u)	Conc. (u/µl)	Cat.#
	Nae I	+	А	GCC▼GGC	250	4	R7131
	(<i>Ngo</i> M I	V)			1,000	4	R7135
6 ⁰	Turbo™ /	Var I ^(c) +	Turbo™	GG▼CGCC	200	10	R7261
^{co}	Narl	+	G	GG▼CGCC	200	10	R6861
	Nci I	+	В	CC▼(C/G)GG	1,000	10	R7061
0	Nco I	+	D	C▼CATGG	200	10	R6513
					1,000	10	R6515
	Nde I	+	D	CA▼TATG	500	10	R6801
	Nde II	+	D	▼GATC	200	10	R7291
	(Dpn I, Sa	au3A I)			1,000	10	R7295
	Ngo M IV (Nae I)		JLTI-COF	RE™ G▼CCGGC	500	10	R7171
^{co}	Nhel	+	В	G▼CTAGC	250	10	R6501
	11101		D		1,250	10	R6505
🖉 ୍ଚତ	Not I	+	D	GC▼GGCCGC	200	10	R6431
•••	10011	т	D	400000	1,000	10	R6435
					1,000	40-80	R4434
e G	Nru I	.1	K	TCG▼CGA	200	40-60	R7091
	Nsil	+ +/-	D	ATGCAVT	200	10	R6531
-	-						
-	Pst I	+/	Н	CIGCAVG	3,000	10	R6111
					15,000	10	R6115
					15,000	40-80	R4114
					50,000	40-80	R4117
	Pvu I	-	D	CGAT▼CG	100	2-10	R6321
			-		500	2-10	R6325
	Pvu II	+	В	CAG▼CTG	1,000	8–12	R6331
					5,000	8–12	R6335
	Rsa I	+	С	GT▼AC	1,000	10	R6371
					5,000	40-80	R4374
	Sacl	+	J	GAGCT▼C	1,000	10	R6061
	(<i>Eco</i> ICR I)			5,000	10	R6065
					5,000	40-80	R4064
	Sac II	+	С	CCGC▼GG	500	10	R6221
🖉 େ	Sal I	+	D	G▼TCGAC	2,000	10	R6051
					10,000	10	R6055
					10,000	40-80	R4054
	Sau3A I	+	В	▼GATC	100	3–10	R6191
	(Dpn I, N	de II)			500	3–10	R6195
	Scal	+	Κ	AGT▼ACT	1,000	8–12	R6211
					5,000	40-80	R4214
🖉 🕹	Sfi I(d)	+	B G(GCCNNNN▼NGGC		10	R6391
					1,250	40-80	R4394
^{co}	Sgfl	+/	С	GCGAT▼CGC	250	8–12	R7103
	0		-		1,250	40-80	R5104
	Sin I	+	A	G▼G(A/T)CC	200	8–12	R6141
	(Ava II)			2. 2. 4. 100	1,000	40-80	R4144
🖉 📀	Smal	+	J	CCC▼GGG	1,000	8–12	R6121
	(Xma I)		U	000 ¥ uuu	5,000	8–12	R6125
	/				5,000	40-80	R4124
	SnaB I		В	TAC▼GTA	100	2–10	
	JIIAD I	_	D		-		R6791
	Snal		В	A▼CTAGT	500 200	<u>2–10</u> 10	R6795
-	Spe I	+	D	A▼UIAGI			R6591
					1,000	10	R6595

Enzym	e	Heat Inactivated	Buffer	Recognition Site	Size (u)	Conc. (u/µl)	Cat.#
6 ⁰	Sph I	+	Κ	GCATG▼C	200	10	R6261
	(Bbu I)				1,000	10	R6265
	Ssp I	+	Е	AATVATT	500	10	R6601
					2,500	40-80	R4604
	Stu I	+	В	AGG▼CCT	400	10	R6421
	Sty I	+	F	C▼C(A/T)(T/A)GG	2,000	10	R6481
	Taq I	—S	Е	T▼CGA	1,000	10	R6151
					10,000	10	R6155
					5,000	40-80	R4154
	Tru9 I	-	F	T▼TAA	200	8–12	R7011
	<i>Tth</i> 111	-	В	GACN▼NNGTC	500	8–12	R6841
60 C	Vsp I	+	D	AT▼TAAT	500	8–12	R6851
🖉 ୍ଦ୍ରେ	Xba I	-	D	T▼CTAGA	2,000	8–12	R6181
					10,000	8–12	R6185
					10,000	40-80	R4184
	Xho I	+	D	C▼TCGAG	3,000	10	R6161
					10,000	10	R6165
					15,000	40-80	R4164
	Xho II	+	С	$(A/G) \mathbf{\nabla} GATC(T/C)$	100	5–10	R6811
					500	5–10	R6815
	Xma I	+	В	C▼CCGGG	50	1–5	R6491
	(<i>Sma</i> I)				250	1–5	R6495
	Xmn I	+	В	GAANN▼NNTTC	500	10	R7271
					2,500	10	R7273
					Cor		
Produc	t			Size	(mg/	/ml)	Cat.#
BSA, (B	ovine Serı	um Albumin) A	cetylated	d <u>400µ</u> l	1		R9461
				1ml	1()	R3961
MULTI-	CORE [™] Bu	Iffer Pack		3 × 1ml	_	-	R9991
4-CORE	Buffer P	ack (1 each A-	D)	4×1 ml		-	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-HCI (mM)	MgCl₂ (mM)	NaCI (mM)	KCI (mM)	DTT (mM)
А	7.5	6	6	6	_	1
В	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	_	_	_
Н	7.5	90	10	50	_	_
J	7.5	10	7	_	50	1
K	7.4	10	10	_	150	_
L	9.0	10	3	100	_	_
				a 0500) 100-	Marshare turns	

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

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.

Indicates Genome Qualified.

Indicates Blue/White Cloning Qualified.

Classic Subcloning: Ordering Information

Enzymes	Size	Conc.	Cat.#
T4 DNA Polymerase ^(d)	100u	5—10u/µl	M4211
		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*	50 preps	A9281
(ready for spin protocol)	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/ <i>Hae</i> 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/ <i>Hae</i> 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*	400µI	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 Loboratory Llos			

*For Laboratory Use.