



Promega

Technical Manual

Universal RiboClone[®] cDNA Synthesis System

INSTRUCTIONS FOR USE OF PRODUCT C4360.



www.promega.com

PRINTED IN USA.
Revised 10/05



AF9TM038 1005TM038

Part# TM038

Universal RiboClone[®] cDNA Synthesis System

All technical literature is available on the Internet at www.promega.com/tbs/
Please visit the web site to verify that you are using the most current version of this
Technical Manual. Please contact Promega Technical Services if you have questions on use
of this system. E-mail techserv@promega.com.

I. Description	2
II. Product Components	3
III. General Considerations	4
A. Methods of cDNA Synthesis.....	4
B. Choice of Primers	4
C. cDNA Cloning	5
D. Choice of Vector	7
E. RNA Preparation and Handling	7
IV. First-Strand Synthesis	8
A. Protocol	9
V. Second-Strand Synthesis	10
A. Protocol	10
VI. Incorporation Assays	12
A. TCA Precipitation.....	12
B. Calculation of First-Strand Yield	13
C. Calculation of Second-Strand Yield.....	13
VII. Gel Analysis	14
A. Labeling Size Markers	14
B. Electrophoresis.....	15
VIII. <i>EcoR</i> I Adaptor Ligation	16
A. Size Fractionation	16
B. Ligation of <i>EcoR</i> I Adaptors.....	16
C. Phosphorylation of DNA	17
IX. Removal of Excess Adaptors	17
A. Preparation of Spin Columns	17
B. Spin Column Protocol.....	18
C. Concentration of cDNA by Ethanol Precipitation.....	18
X. Ligation of cDNA Insert to <i>EcoR</i> I-Digested Plasmid Vector	19
A. Protocol	19
XI. Troubleshooting	20

XII. Composition of Buffers and Solutions	22
XIII. Related Products	22
XIV. References	23

I. Description

The Universal RiboClone® cDNA Synthesis System^(a,b,c) contains the reagents required for the synthesis of double-stranded cDNA from mRNA and subsequent ligation into a suitable vector. The system is based on the method described by Okayama and Berg (1) with modifications by Gubler and Hoffman (2). First-strand synthesis is driven by AMV (Avian Myeloblastosis Virus) Reverse Transcriptase and either Random Hexameric Primers or an Oligo(dT) Primer, followed directly by second-strand replacement synthesis using RNase H and DNA Polymerase I. After treatment with T4 DNA Polymerase to flush the ends, the double-stranded cDNA molecules are prepared for cloning by size fractionation and the addition of *EcoR* I Adaptors. The resulting cDNA preparation can then be cloned into a suitable vector.

Selected Citation using the Universal RiboClone® cDNA Synthesis System

- Krutovsky, K.V. *et al.* (2004) Comparative mapping in the Pinaceae. *Genetics* 168, 447–61.

Notes: To compare STS and ETS markers of a known conifer, loblolly pine, with Douglas fir, a cDNA library was made from 1-month-old Douglas fir seedlings. Total RNA was extracted from 5g tissue and double-stranded cDNA prepared using the Universal RiboClone® cDNA Synthesis System. The cDNA ligated to *EcoR* I adaptors were cloned into pGEM®-3Z, propagated and sequenced to give rise to 5031 EST sequences, several EST contigs and four Douglas fir STSs for comparison to the GenBank loblolly pine ESTs. Those regions that demonstrated homology between the two species were selected for further study.

For additional peer-reviewed articles that cite the use of the Universal RiboClone® cDNA Synthesis System, visit: www.promega.com/citations/

II. Product Components

Product	Cat.#
Universal RiboClone® cDNA Synthesis System	C4360

Each system contains sufficient reagents to process up to 40µg mRNA and to perform 10 adaptor ligation reactions. **C4360 consists of three systems:**

C4361 cDNA Synthesis Reagents

- 200µl First Strand 5X Buffer
- 20µg Oligo(dT)₁₅ Primer
- 20µg Random Hexameric Primers
- 50µl Sodium Pyrophosphate, 40mM
- 1,000u Recombinant RNasin® Ribonuclease Inhibitor^(a,b)
- 600u AMV Reverse Transcriptase (HC)
- 10µg 1.2kb Kanamycin Positive Control RNA
- 800µl Second Strand 2.5X Buffer
- 20u RNase H
- 500u DNA Polymerase I
- 100u T4 DNA Polymerase^(c)
- 1.25ml Nuclease-Free Water

C4362 EcoR I Adaptor Ligation Reagents

- 0.5ml T4 DNA Ligase 10X Buffer
- 100u T4 DNA Ligase
- 400µl Acetylated BSA, 1mg/ml
- 100u T4 Polynucleotide Kinase
- 0.5ml Kinase 10X Buffer
- 50µl ATP, 10mM
- 150pmol EcoR I Adaptors

C4363 Resin and Spin Columns

- 10 Spin Columns
- 10ml Sephacryl® S-400
- 1 Protocol

Note: C4362 contains the BSA needed for second-strand synthesis.

Storage and Stability: Store the Positive Control RNA at -70°C. Store the Sephacryl® S-400 Resin at 4°C and Spin Columns at room temperature. Store all other components at -20°C. Avoid multiple freeze-thaw cycles of the T4 DNA Ligase 10X Buffer.

III. General Considerations

III.A. Methods of cDNA Synthesis

One of the most challenging tasks in molecular biology is the synthesis and cloning of cDNA. A complex series of enzymatic steps is involved in copying mRNA into double-stranded cDNA and subsequently preparing the termini for vector ligation. Many approaches have been used to generate cDNA libraries, and these have tried to preserve as much of the original sequence as possible, to improve cloning efficiency and to facilitate screening and subsequent analysis (3).

Most cDNA molecules produced will lack a few nucleotides corresponding to the 5' end of the mRNA because second-strand replacement only proceeds from 3'-OH RNA primers. These primers are randomly generated by the addition of RNase H and the endogenous RNase H activity of AMV Reverse Transcriptase. The most 5' nick in the mRNA will generally occur several nucleotides from the end, and the remaining RNA oligonucleotide may be too short to remain hybridized. Thus, the 3'→5' exonuclease activity of DNA Polymerase I will remove the last few nucleotides of the cDNA first strand. However, since all eukaryotic mRNA molecules appear to have 5' noncoding leader sequences, which commonly range from 40–80 nucleotides (4), it is likely that the vast majority of double-stranded cDNA will contain all of the coding sequences present in the initial cellular mRNA molecules.

If retaining these nucleotides is necessary, alternative procedures involving tailing after first-strand synthesis can be used (3,5). Homopolymeric tailing is best accomplished by adding dG residues (6), and the resulting clones will contain a long stretch (about 20bp) of dG/dC residues. These regions may interfere with screening using GC-rich oligomers, with dideoxy sequencing of clones, and possibly with their expression as part of a fusion protein (6). In addition, these regions can interfere with differential hybridization procedures in the construction of subtraction libraries (7).

III.B. Choice of Primers

The classical method of cDNA synthesis uses the Oligo(dT) Primer to prime first-strand synthesis (Figure 1). This method is suitable in most cases where poly(A)⁺ RNA of high quality can be prepared from the cell line or tissue of interest. Random Hexameric Primers (hexadeoxyribonucleotides) provide an alternative procedure (Figure 2) by which first-strand cDNA synthesis is initiated from internal sites within the mRNA molecule (8). Random Primers can be used to prime cDNA synthesis from mRNA molecules that do not possess a poly(A)⁺ tail or for RNA isolated from prokaryotic sources. Random Primers also provide a scheme by which cDNA can be synthesized representing mRNA with strong 5' secondary structure.

III.C. cDNA Cloning

Adaptors eliminate the *EcoR* I methylation and restriction enzyme steps necessary with conventional linker addition. The adaptor, a duplex DNA molecule, consists of one blunt end (for ligation to the cDNA) and an *EcoR* I sticky end. This feature of the Universal RiboClone® System allows cDNA to be cloned into unique *EcoR* I sites of vectors (Figure 3).

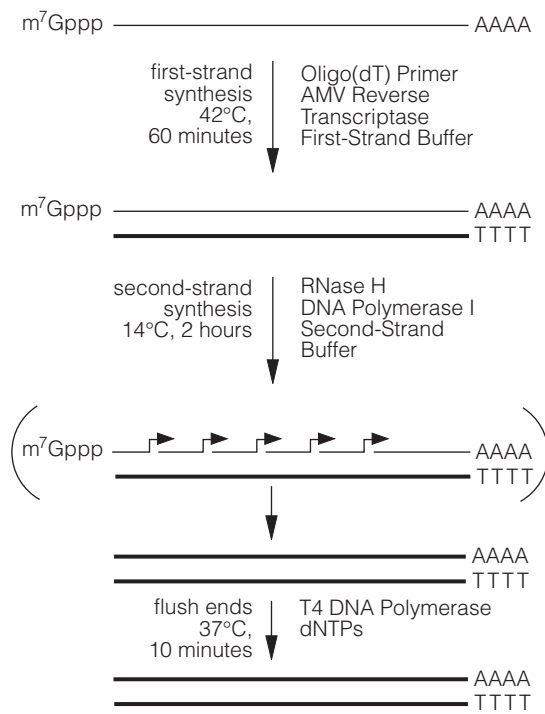
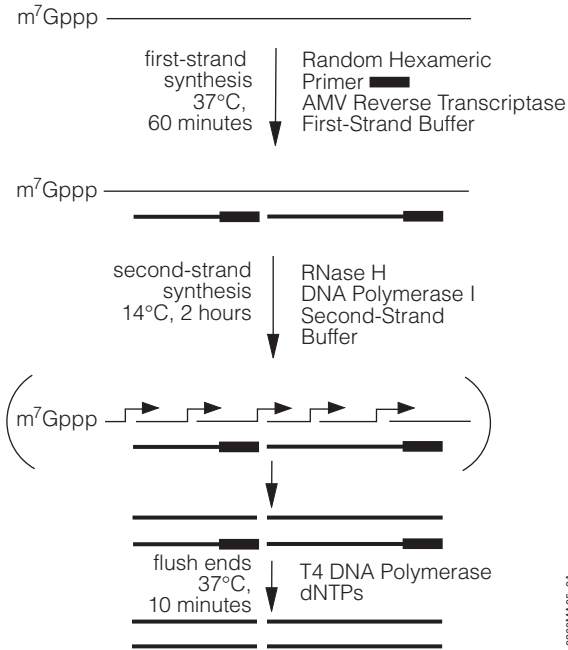


Figure 1. cDNA synthesis using the Universal RiboClone® cDNA Synthesis System and Oligo(dT) Primer.



0360MA05_3A

Figure 2. cDNA synthesis using the Universal RiboClone® cDNA Synthesis System and Random Hexameric Primers.

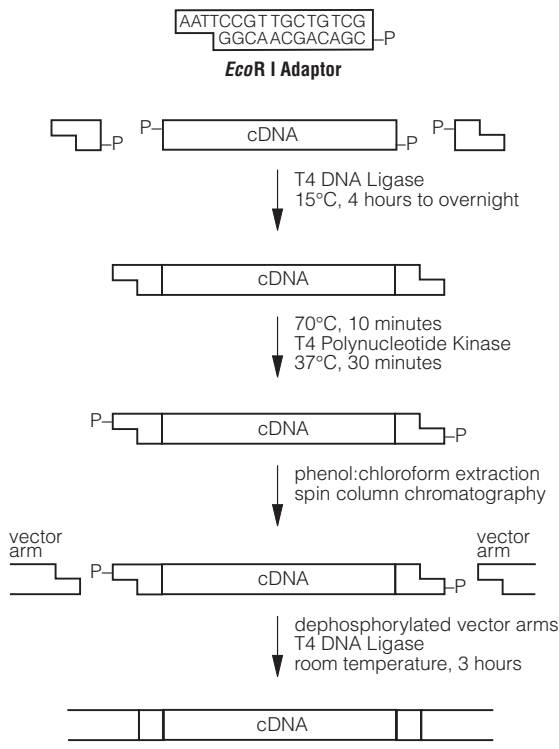


Figure 3. Schematic representation of EcoR I Adaptor ligation.

III.D. Choice of Vector

Many different plasmid and lambda vectors are available for cDNA cloning. The type of vector selected will depend on the possible applications for the cDNA library being constructed and other factors, such as blue-white cloning selection.

III.E. RNA Preparation and Handling

The quality and quantity of cDNA synthesized by any method is critically dependent on the integrity of the mRNA used as the template. Extremely stable RNases are ubiquitous in the laboratory environment (including the surface of fingers of the researcher) and in all cells. Therefore, certain precautions should be taken to eliminate the risk of RNase contamination whenever possible.

Procedures for creating an RNase-free laboratory environment have been described in detail (9). Sterile disposable plasticware should be used whenever possible, and disposable gloves should be worn at all times and changed often.

All glassware, nondisposable plasticware and electrophoresis apparatus used should be kept separately from other labware. Glassware can be soaked and rinsed in a solution of 0.1% diethyl pyrocarbonate (DEPC) and autoclaved. Proper safety precautions should be taken when using DEPC, which is a powerful acylating agent. This compound should always be used in a fume hood and never added to aqueous solutions containing ammonia, which would result in the formation of ethyl carbamate, a potent carcinogen (10).

Many procedures exist for isolating intact RNA from various cells and tissues. A widely used method that has been successful with RNase-rich sources was first described by Chirgwin *et al.* (11). The Promega RNAgents® Total RNA Isolation System provides a rapid and convenient means to isolate high-quality total RNA based on this method.

The standard method of preparing poly(A)⁺ RNA has been by chromatography on oligo(dT)-cellulose (12). Even after two cycles of oligo(dT) selection, most preparations still contain varying amounts of ribosomal RNA sequences. In general, these sequences will not interfere with cDNA synthesis using oligo(dT)-containing primers. However, reduced yields of cDNA may be observed with Random Hexameric Primers, and quantitation of the mRNA will be affected.

The PolyATtract® mRNA Isolation Systems use Promega MagneSphere® technology to eliminate the need for oligo(dT) cellulose and its associated problems. With total RNA as the starting material, the poly(A)⁺ mRNA fraction can be isolated free of all other nucleic acid contamination in approximately 45 minutes. The system uses a biotinylated oligo(dT) primer to hybridize at high efficiency in solution to the 3' poly(A) region present in most mature eukaryotic mRNA species. The hybrids are captured and washed at high stringency using streptavidin coupled to paramagnetic particles and a magnetic separation stand. The mRNA is eluted from the solid phase by the simple addition of ribonuclease-free deionized water. This procedure yields mature polyadenylated mRNA after only a single round of magnetic separation. Alternatively, the PolyATtract® System 1000 may be used to isolate mRNA directly from crude cell or tissue lysates and eliminates the need for total RNA isolations.

IV. First-Strand Synthesis

The standard reaction described below is performed in a total volume of 25 μ l, which is recommended for **up to 2 μ g** of mRNA. For each additional microgram of mRNA, increase the reaction volume by 10 μ l. For example, the reaction should be scaled up to 55 μ l when using 5 μ g of mRNA. Other components should be scaled up proportionally. After all of the components are added, a 5 μ l sample is removed to a separate tube containing a small amount of [α -³²P]dCTP as a tracer to follow first-strand synthesis. The remainder of the reaction is carried through second-strand synthesis.

IV.A. Protocol

Materials to Be Supplied by the User

(Solution compositions are provided in Section XII.)

- EDTA (50mM and 200mM)
- [α - 32 P]dCTP (>400Ci/mmol, 10mCi/ml)

1. Prepare the mRNA sample and control reactions as described below. The standard reactions described are optimized for **2 μ g** of mRNA template.

In a sterile, RNase-free microcentrifuge tube, add the appropriate primer to the mRNA or Control RNA sample. For mRNA samples, use 0.5 μ g primer/ μ g mRNA in a total volume of up to 15 μ l in water with either the Oligo(dT) Primer or Random Hexameric Primers.

! Do not alter the ratio of primer to mRNA.

mRNA sample

mRNA	2 μ g
Primer, 0.5mg/ml	2 μ l
Nuclease-Free Water to final volume	15 μ l

! The Control RNA uses a 1:1 ratio of primer to RNA. This ratio has been optimized for the 1.2kb Kanamycin Positive Control RNA and is not the same ratio used for mRNA.

control reaction

1.2kb Kanamycin Positive Control RNA,	
0.5mg/ml	2 μ l
Primer, 0.5mg/ml	2 μ l
Nuclease-Free Water	11 μ l
total volume	15 μ l

2. Heat to 70°C for 5–10 minutes. Chill the tube on ice for 5 minutes and spin briefly to collect the solution at the bottom of the tube.
3. Add the following components in the order shown.

sample or control reaction	15 μ l
First Strand 5X Buffer	5 μ l
RNasin® Ribonuclease Inhibitor	40u

Heat at 42°C (37°C for Random Primers) for 3–5 minutes, then add:

Sodium Pyrophosphate, 40mM	2.5 μ l
AMV Reverse Transcriptase	30u
Nuclease-Free Water to final volume	25 μ l

Note: First Strand 5X Buffer and Sodium Pyrophosphate may form a precipitate when frozen. If this occurs, warm the tube to 37°C to dissolve the precipitate before addition.

- Mix gently by flicking the tube, and transfer 5 μ l of each reaction mixture to another tube containing 2–5 μ Ci of [α -³²P]dCTP (>400Ci/mmol, 10mCi/ml). These are the sample and control tracer reactions, which will be used to measure first-strand synthesis by incorporation and alkaline agarose gel electrophoresis (Sections VI and VII).

Note: We recommend using [α -³²P]dCTP (aqueous form) with a specific activity of >400Ci/mmol, 10mCi/ml (such as Amersham Biosciences Cat.# PB10165) that is less than 1 week old.

- Incubate both labeled and unlabeled reactions for 60 minutes at 42°C for the Oligo(dT) Primer or 37°C for Random Hexameric Primers.

Note: The reaction mixture may appear cloudy at this point, but subsequent steps will not be adversely affected.

- Place reactions on ice.
- Add 95 μ l of 50mM EDTA to the tracer reactions (to a total volume of 100 μ l) and store on ice. The tracer reactions will be used later for incorporation assays (Section VI.A) and gel analysis (Section VII).
- Proceed to second-strand synthesis (Section V) using the larger, unlabeled first-strand reactions. No phenol extractions or ethanol precipitations are necessary.

V. Second-Strand Synthesis

The second-strand reaction volume is based on that of the first-strand reaction volume. Maintain a fivefold dilution of the first-strand reaction volume in the second-strand reaction so the component amounts are scaled proportionally.

V.A. Protocol

Materials to Be Supplied by the User

(Solution compositions are provided in Section XII.)

- [α -³²P]dCTP (>400Ci/mmol, 10mCi/ml)
- TE-saturated phenol:chloroform:isoamyl alcohol
- EDTA, 50mM and 200mM
- ammonium acetate, 7.5M (or sodium acetate, 2.5M [pH 5.2])
- ethanol (100% and 70%), –20°C
- TE buffer

1. Add the following components to the first-strand sample and Control RNA reactions in the order shown. Perform this step on ice.

second-strand sample

first-strand sample reaction	20 μ l
Second Strand 2.5X Buffer	40 μ l
Acetylated BSA, 1mg/ml	5 μ l
DNA Polymerase I	23u
RNase H	0.8u
Nuclease-Free Water to final volume	100μl

second-strand control

first-strand Control RNA reaction	20 μ l
Second Strand 2.5X Buffer	40 μ l
Acetylated BSA, 1mg/ml	5 μ l
DNA Polymerase I	23u
RNase H	0.8u
Nuclease-Free Water to final volume	100μl

2. Mix gently by flicking the tube. If a tracer reaction for the second-strand sample is desired, remove 10 μ l to another tube containing 2-5 μ Ci [α -³²P]dCTP (>400Ci/mmol, 10mCi/ml).
3. For the second-strand control tracer reaction, add 2-5 μ Ci [α -³²P]dCTP directly to the 100 μ l control reaction.
4. Incubate the reactions at 14°C for 2 hours.
Note: When cDNA longer than 3kb is desired, increase the second-strand incubation time to 3-4 hours.
5. Heat the unlabeled cDNA sample to 70°C for 10 minutes, collect the contents at the bottom of the tube by brief centrifugation and place on ice.
6. While the unlabeled samples are being heated, add 90 μ l of 50mM EDTA to the 10 μ l sample tracer reaction. Add 10 μ l of 200mM EDTA to the 100 μ l Control RNA tracer reaction. The tracer reactions will be used later for incorporation assays (Section VI) and gel analysis (Section VII). Store on ice.
7. Add 2 units of T4 DNA Polymerase/ μ g input RNA to the unlabeled reaction. Incubate at 37°C for 10 minutes.
8. Stop the reaction by adding 10 μ l of 200mM EDTA and place on ice.
9. To extract the cDNA, add an equal volume of TE-saturated phenol:chloroform:isoamyl alcohol, mix thoroughly, and centrifuge at top speed in a microcentrifuge for 2 minutes at room temperature.

10. Transfer the aqueous phase to a fresh tube, add 0.5 volume of 7.5M ammonium acetate (or 0.1 volume of 2.5M sodium acetate [pH 5.2]), mix and add 2 volumes of cold (-20°C) 100% ethanol. Place at -70°C for 30 minutes and centrifuge at top speed for 5 minutes.
11. Carefully remove the supernatant, add 0.5ml of cold (-20°C) 70% ethanol to the tube with the pellet. Centrifuge again at top speed for 2 minutes.
12. Carefully remove the supernatant and dry the pellet.
13. Dissolve the pellet in 10-50 μl of TE buffer. The cDNA is now ready for size fractionation prior to adaptor addition (Section VIII).

Note: The Promega Wizard[®] DNA Clean-Up System (Cat. # A7280) may be used instead of solvent extraction and ethanol precipitation in steps 9-13.

VI. Incorporation Assays

Trichloroacetic acid (TCA) assays of isotope incorporation are prone to variation (13), and thus some variability can be expected in assays of first- and second-strand conversion. In our experience, cDNA libraries can be successfully constructed from reactions yielding 12-50% first-strand conversion and at least 80% second-strand conversion values.

VI.A. TCA Precipitation

Materials to Be Supplied by the User

- 1mg/ml carrier DNA (e.g., herring sperm DNA, Cat.# D1811)
 - TCA 5% (w/v), ice-cold
 - glass fiber filters (e.g., Whatman[®] GF-A or GF-C)
 - acetone or ethanol, 100%
1. Spot 3 μl samples of the diluted first-strand tracer reaction (Section IV, Step 7) and the second-strand tracer reaction (Section V.A, Step 6) on glass fiber filters and let them air-dry. These samples represent the total cpm in the reactions.
 2. Add 3 μl samples of the same reactions to tubes containing 100 μl of a 1mg/ml solution of carrier DNA and mix. Add 0.5ml of 10% TCA (ice-cold) and vortex. Let sit on ice for 5-30 minutes.
 3. Filter the samples through glass fiber filters, washing 3 times each with 5ml of cold 5% TCA. Rinse the filters with 5ml of acetone or ethanol and let air-dry.
 4. Count both total and incorporated (TCA-precipitable) samples, either by Cerenkov radiation (without scintillant) or after adding an appropriate scintillation fluid.

VI.B. Calculation of First-Strand Yield

The first-strand yield is determined as follows:

$$\frac{\text{incorporated cpm}}{\text{total cpm}} \times 100\% = \% \text{ first-strand incorporation}$$

$$\frac{4 \text{ (nmol dNTP}/\mu\text{l}) \times \text{reaction vol. } (\mu\text{l}) \times \% \text{ first-strand incorp.}}{100} = \text{nmol dNTP incorp.}$$

$$\text{nmol dNTP incorporated} \times 330\text{ng/nmol} = \text{ng cDNA synthesized}$$

$$\frac{\text{ng cDNA synthesized}}{\text{ng mRNA in reaction}} \times 100\% = \% \text{ mRNA converted to cDNA}$$

Example: If there were 1,854cpm incorporated and 254,000 total cpm in the tracer reaction, and 2 μ g RNA was added to the original 25 μ l reaction, the calculations would be as follows:

$$\frac{1,854}{254,000} \times 100\% = 0.73\% \text{ first-strand incorporation}$$

$$\frac{4\text{nmol dNTP}/\mu\text{l} \times 25\mu\text{l} \times 0.73}{100} = 0.73\text{nmol dNTP incorporated}$$

$$0.73\text{nmol dNTP} \times 330\text{ng/nmol} = 240\text{ng cDNA synthesized}$$

$$\frac{240\text{ng cDNA synthesized}}{2,000\text{ng mRNA in reaction}} \times 100\% = 12\% \text{ mRNA converted to cDNA}$$

Of the original 2,000ng of mRNA, 20% (or 5 μ l/25 μ l) was removed for the tracer reaction, and 80% (or 20 μ l/25 μ l) remained in the main reaction. Therefore, 192ng (0.8 \times 240ng) first-strand cDNA was synthesized and 0.58nmol (0.8 \times 0.73nmol) dNTPs were incorporated in the reaction that was carried through to second-strand synthesis.

Note: The minimum quality control specification for first-strand conversion is 12%.

VI.C. Calculation of Second-Strand Yield

The second-strand yield is calculated in the same manner as the first-strand yield, except that the total dNTP in the reaction will be diminished by that which was incorporated during first-strand synthesis.

$$\frac{\text{incorporated cpm}}{\text{total cpm}} \times 100\% = \% \text{ second-strand incorporation}$$

$$[0.8 \text{ (nmol dNTP}/\mu\text{l}) \times \text{reaction vol. } (\mu\text{l}) - \text{nmol incorp. in first-strand reaction}] \times (\% \text{ second-strand incorporation}/100) = \text{nmol dNTP incorporated}$$

$$\text{nmol dNTP incorporated} \times 330\text{ng/nmol} = \text{ng second-strand cDNA synthesized}$$

$$\frac{\text{ng second-strand synthesized}}{\text{ng first-strand synthesized}} \times 100\% = \% \text{ converted to double-stranded cDNA}$$

Example: The first-strand reaction above was used for second-strand synthesis, which resulted in 1,390cpm incorporated out of a total of 235,000cpm.

$$\frac{1,390}{235,000} \times 100\% = 0.59\% \text{ second-strand incorporation}$$

$$[0.8\text{nmol}/\mu\text{l} \times 100\mu\text{l} - 0.58\text{nmol first-strand dNTP incorporated}] \times 0.59/100 \\ = 0.47\text{nmol dNTP incorporated}$$

$$0.47\text{nmol} \times 330\text{ng}/\text{nmol} = 155\text{ng second-strand cDNA synthesized}$$

$$\frac{155\text{ng second-strand cDNA}}{192\text{ng first-strand cDNA}} \times 100\% = 81\% \text{ conversion to ds cDNA}$$

Note: The minimum quality control specification for second-strand conversion is 80%.

VII. Gel Analysis

The size distribution of cDNA synthesized (usually 350–6,000 bases) in the first- and second-strand reactions should be checked by electrophoresis on a 1.4% alkaline agarose gel (14). The samples used for gel analysis are the first-strand and second-strand tracer reactions. Load approximately equivalent amounts of incorporated cpm for each strand so that comparisons can be made on the same autoradiograph. As a guideline, 5–10 μl of the control tracer reaction on a fixed and dried gel results in a dark, sharp signal in a 3–4 hour exposure with intensifying screens.

VII.A. Labeling Size Markers

DNA markers that have been ^{32}P -labeled may be used for estimating the sizes of the first- and second-strand products. A convenient way to prepare these markers is to use Lambda DNA/*Hind* III Markers in a fill-in reaction with Klenow DNA Polymerase.

Materials to Be Supplied by the User

(Solution compositions are provided in Section XII.)

- Klenow DNA Polymerase
- Klenow 10X Buffer
- dATP, dGTP (10mM)
- Lambda DNA/*Hind* III Markers (Cat.# G1711)
- [α - ^{32}P]dCTP (>400Ci/mmol)
- EDTA (200mM)
- sample 2X buffer

1. Assemble the following reaction:

Klenow 10X Buffer	2.5 μ l
dATP	0.2mM
dGTP	0.2mM
[α - ³² P]dCTP (>400Ci/mmol)	2 μ Ci
Lambda DNA/ <i>Hind</i> III Markers	1 μ g
Klenow DNA Polymerase	1u
Nuclease-Free Water to final volume	25μl

2. Incubate the reaction for 10 minutes at room temperature. Add 2.5 μ l of 200mM EDTA to stop the reaction. Add a sample directly to an equal volume of sample 2X buffer and store the remainder at -20°C. A 5 μ l sample routinely gives a sharp ladder of bands after a 1-hour exposure.

VII.B. Electrophoresis

Materials to Be Supplied by the User

(Solution compositions are provided in Section XII.)

- alkaline agarose gel
 - alkaline gel running buffer
 - sample 2X buffer
 - TCA, 7% (w/v)
1. Prepare a 1.4% alkaline agarose gel (see Section XII).
 2. Transfer the desired amount of each sample (usually 10,000–50,000 incorporated cpm) to separate tubes, adjust the volumes with TE buffer so that they are equal, and add an equal volume of sample 2X buffer to each.
 3. Load the samples and run the gel at up to 5V/cm for the desired time (usually until the tracking dye has migrated about 2/3 of the way through the gel).
 4. Soak the gel in several volumes of 7% TCA at room temperature for 30 minutes or until the dye changes from blue to yellow. Dry on a piece of Whatman® 3MM paper, either on a gel dryer or by placing the gel under a weighted stack of paper towels for several hours.
 5. Cover the dried gel with plastic wrap and expose to X-ray film at room temperature or at -70°C with an intensifying screen. Alternatively, scan the gel using phosphorimaging analysis.
 6. If the Positive Control mRNA was used with the Oligo(dT) Primer, the first- and second-strand reaction products should appear as sharp bands at 1.2kb in the gel. Random priming results in products that are 1.2kb and smaller.

VIII. *EcoR* I Adaptor Ligation

VIII.A. Size Fractionation

Materials to Be Supplied by the User

(Solution compositions are provided in Section XII.)

- spin columns

We recommend size fractionating cDNA samples before ligation with adaptors. Small-sized cDNA, less than 400bp, should be eliminated to ensure successful adaptor addition and cloning. This step can be accomplished by gel filtration using either column fractionation or spin columns. Alternatively, cDNA samples may be fractionated by agarose gel electrophoresis followed by selective elution of particular-sized cDNA molecules (14). We do not recommend precipitating DNA in the presence of ammonium ions after either of these fractionation procedures.

Note: The Sephacryl® S-400 Resin and Spin Columns in the Universal RiboClone® System are provided for the removal of adaptors. They may also be used for size fractionation of cDNA prior to ligation; however, the quantity of Resin and Spin Columns provided is sufficient for use in a total of 10 size fractionation and adaptor removal reactions. Additional Spin Columns and Resin may be purchased from Promega (see Section XIII).

! Do not freeze the Sephacryl® S-400 Resin.

Note: The size cutoff for the Sephacryl® S-400 Resin is 271bp.

VIII.B. Ligation of *EcoR* I Adaptors

1. Set up the following reaction (volumes can be scaled up proportionally, depending upon the amount of cDNA).

T4 DNA Ligase 10X Buffer	3µl
Acetylated BSA, 1mg/ml	3µl
cDNA (100ng/µl)	2.5µl
<i>EcoR</i> I Adaptors (20-fold molar excess:	
10pmol of adaptors)*	1µl
T4 DNA Ligase (Weiss units)	2.5u
Nuclease Free Water to final volume	30µl

***Example:** 250ng of a 1.5kb DNA (average size cDNA) =
 0.5pmol ends
 0.001 A₂₆₀ units of *EcoR* I Adaptors = 10pmol

2. Incubate at 15°C for 4 hours to overnight.
3. Inactivate the enzyme by heating the reaction to 70°C for 10 minutes.
4. Cool the reaction on ice and proceed to Section VIII.C.

VIII.C. Phosphorylation of DNA

For optimal cloning efficiency, the insert DNA should be phosphorylated and vector DNA should be dephosphorylated. The following protocol should be used for cDNA synthesized with the Oligo(dT) Primer or Random Hexameric Primers.

Materials to Be Supplied by the User

(Solution compositions are provided in Section XII.)

- TE buffer
- TE-saturated phenol:chloroform:isoamyl alcohol

1. Set up the following reaction.

ligation reaction (from Section VIII.B)	30 μ l
Kinase 10X Buffer	4 μ l
ATP, 0.1mM (1:100 dilution in water of 10mM stock)	2 μ l
T4 Polynucleotide kinase (5–10u/ μ l)	10u
Nuclease Free Water to final volume	40 μ l

2. Incubate at 37°C for 30 minutes.
3. Phosphorylated cDNA may be purified prior to ligation with vector. This can be accomplished using a standard DNA cleanup protocol (Steps 4–5) or using a resin-based purification system such as the Promega Wizard® DNA Clean-Up System (Cat.# A7280).
4. Add 1 volume of TE-saturated phenol:chloroform. Vortex for 30 seconds and centrifuge at top speed in a microcentrifuge for 3 minutes.
5. Transfer the upper, aqueous phase to a fresh tube. To increase recovery, the phenol:chloroform phase can be re-extracted with a small volume of TE buffer.
6. Proceed to removal of excess adaptors (Section IX).

IX. Removal of Excess Adaptors

IX.A. Preparation of Spin Columns

Materials to Be Supplied by the User

(Solution compositions are provided in Section XII.)

- TEN buffer

Sephacryl® S-400 Resin is provided in a solution of TEN buffer. It is a good idea to re-equilibrate with this buffer before use.

1. Thoroughly mix the Sephacryl® S-400 slurry and transfer 600 μ l to a Spin Column. Start the flow of buffer out of the bottom of the column by using a gentle push of air from a Pasteur pipette bulb. Allow the buffer to drain and add an additional 400 μ l for a total of 1ml.

2. Discard the collected fluid and add 600 μ l of TEN buffer, allow some to drain, then add an additional 400 μ l. Allow buffer to drain from the tube. This step should be repeated 2 times.
Note: Equilibration may also be achieved by repeated centrifugations using smaller volumes of buffer. If this method is used, be careful that the filtrate volume does not touch the frit of the column.
3. Place the column tip into the provided Wash Tube. Put the assembly inside a larger centrifuge tube (e.g., Falcon #25319), and centrifuge in a swinging bucket rotor at 800 \times g for 5 minutes.
4. The bed should have a dry appearance and may pull away from the sides of the tube. The column is now ready for immediate use.

IX.B. Spin Column Protocol

The optimal sample volume is 20–60 μ l. Do not exceed a 60 μ l sample volume. Spin Columns must be maintained in a vertical position during sample loading and subsequent handling procedures to ensure proper separation.

1. Pipet the reaction mixture from Section VIII.C. Hold the pipet tip within a few millimeters of the gel bed, and slowly add the sample dropwise to the center of the dried gel bed. Do not allow the sample to come in direct contact with the sides of the tube.
2. Place the Spin Column into the provided Collection Tube, carefully put this assembly inside a larger tube and centrifuge in a swinging bucket rotor at 800 \times g for 5 minutes.
3. Remove the assembly, discard the Spin Column and save the Collection Tube. The volume recovered should be approximately equal to the volume that was applied to the Spin Column.
4. The eluted cDNA can either be used directly or precipitated with ethanol to concentrate the sample.

IX.C. Concentration of cDNA by Ethanol Precipitation

Materials to Be Supplied by the User

(Solution compositions are provided in Section XII.)

- ammonium acetate, 7.5M (or sodium acetate, 2.5M [pH 5.2])
 - ethanol, 100% and 70% (-20°C)
 - TE buffer
1. To the sample from Section IX.B, add 0.5 volume of 7.5M ammonium acetate (or 0.1 volume of 2.5M sodium acetate [pH 5.2]) and 2.5 volumes of cold (-20°C) 100% ethanol. Mix and place at -70°C for 30 minutes.
 2. Centrifuge at top speed in a microcentrifuge for 15 minutes. Carefully remove and discard the supernatant.

3. Wash the pellet with 1ml of cold (-20°C) 70% ethanol and spin in microcentrifuge at top speed for 5 minutes. Carefully remove and discard the supernatant.
4. Dry the pellet briefly under vacuum. Resuspend the pellet in an appropriate volume of TE buffer for further reactions.

X. Ligation of cDNA Insert to *EcoR* I-Digested Plasmid Vector

Synthesized cDNA can be cloned into a plasmid vector, which offers versatility and ease of manipulation after cloning. The ratio of insert DNA to vector is critical to successful ligation and should be carefully optimized for each cDNA preparation following the protocol described below. As a general recommendation, an insert:vector molar ratio of 1:3 may be used for dephosphorylated vectors. We do not recommend using phosphorylated vector due to the low cloning efficiency obtained as a result of vector self-ligation.

In the example given below, the molar ratio of plasmid (3kb) to insert DNA is varied from 1:3 to 1:1. A tube containing no insert DNA is also used to determine background levels of religated vector.

X.A. Protocol

1. Set up sample reactions as described. The amounts of insert and vector specified are based on an average cDNA size of 1.5kb. Maintain a constant molar concentration of the vector while adjusting the amount of cDNA.

Sample Reaction: Plasmid Vector

	Tube			
	A	B	C	D
Vector DNA (10ng/μl; 0.005pmol/μl)	2μl	2μl	2μl	2μl
Insert DNA (10ng/μl; 0.01pmol/μl)	0μl	3μl	2μl	1μl
T4 DNA Ligase 10X Buffer	1μl	1μl	1μl	1μl
Nuclease-Free Water	6μl	3μl	4μl	5μl
T4 DNA Ligase	1μl	1μl	1μl	1μl

2. Incubate the reactions at room temperature for 3 hours. Alternatively, the incubation may be performed at 4°C for 16-18 hours.
3. Once optimal ligation conditions have been established, reactions may be scaled up accordingly for library construction.

Note: Synthesized cDNA can also be cloned into bacteriophage lambda vectors. These vectors may be preferred when high-efficiency cloning is desired. If substituting a lambda vector for plasmid in the sample reaction, the molar ratio of lambda arms (43kb) to DNA insert is varied from 3.6:1 to

1.2:1 (e.g., use the same volumes as for Plasmid Vector, but use 0.5µg/µl; 0.018pmol/µl lambda vector). Recombinant lambda vectors can be packaged using the Packagene® Systems (see Section XIII for a list of related products).

XI. Troubleshooting

For questions not addressed here, please contact your local Promega Branch Office or Distributor. Contact information available at: www.promega.com. E-mail: techserv@promega.com

Symptoms	Causes and Comments
Low yield of first- and second-strand cDNA, or no cDNA synthesized*	RNase contamination in synthesis reaction. Check control reaction for yield. Using the Oligo(dT) Primer, the Positive Control RNA should give at least 12% first-strand conversion, 80% second-strand conversion and clear bands of 1.2kb on alkaline agarose gels. Follow guidelines in Section III.E to minimize RNase contamination.
	Degraded mRNA template. Check mRNA preparation by electrophoresis using either denaturing or native conditions. If total RNA is run on a gel, bands corresponding to 28S and 18S ribosomal RNA should be apparent, and 28S RNA should stain with approximately twice the intensity of 18S RNA. Procedures outlined in Section III.E should be followed to ensure the isolation of undegraded mRNA template.
	mRNA template has high degree of secondary structure. Heat mRNA to 70°C for 5–10 minutes, then place immediately on ice to ensure denaturation of template.
	Incorrect primer:RNA ratio. Use 0.5µg primer/µg mRNA for the Oligo(dT) or Random Hexameric Primers and mRNA samples. The control RNA uses a 1:1 ratio of primer:RNA.
	Inhibitors in mRNA preparation. Inhibitors such as SDS, EDTA, polysaccharides or salts may affect first-strand synthesis. Add a small amount of the mRNA preparation to the Positive Control RNA first-strand reaction to check for inhibition of synthesis.

XI. Troubleshooting (continued)

Symptoms	Causes and Comments
Low cloning efficiency	<p>Poor ligation of adaptors to cDNA. Check ligation reactions with pure DNA preparations (e.g., purified plasmid DNA or Lambda DNA/<i>Hind</i> III Markers). Longer incubation time at a lower temperature may increase ligation efficiency (e.g., overnight at 4°C).</p>
	<p>Hairpin priming. First-strand synthesis should be performed in the presence of sodium pyrophosphate to suppress formation of hairpins.</p>
	<p>Insufficient phosphorylation of ligated <i>Eco</i>R I Adaptors. The efficiency of the kinase reaction may be monitored using [γ-³²P]ATP.</p>
	<p>Do not dissolve or precipitate DNA in the presence of ammonium ions after fractionation. Ammonium ions inhibit T4 polynucleotide kinase.</p>
	<p>Poor yield of cDNA from Sephacryl® S-400 columns. Resin may not be sufficiently equilibrated. Wash column 3 times with TEN buffer before use.</p>
	<p>Ratio of cDNA to vector not optimal for ligation. Determine optimal conditions for ligation by adjusting the amount of cDNA in small-scale reactions while maintaining a constant concentration of vector.</p>
<p>Poor ligation of cDNA to vector. Cloning vector should be dephosphorylated for optimal results.</p>	

*A low percent conversion for second-strand product may be due to a high percent conversion for first-strand synthesis. If a high yield is achieved for first-strand synthesis (as from extended incubation), the percent converted during second-strand synthesis may be lower, but the amount of DNA should be adequate for most applications.

XII. Composition of Buffers and Solutions

alkaline gel running buffer

30mM NaOH
1mM EDTA

Make up fresh for each use.

alkaline agarose gel

1.4% agarose
50mM NaOH
1mM EDTA

Equilibrate solidified gel for at least 30 minutes in alkaline gel running buffer prior to electrophoresis.

First Strand 5X Buffer

250mM Tris-HCl (pH 8.3 at 42°C)
250mM KCl
50mM MgCl₂
2.5mM spermidine
50mM DTT
5mM ea dATP, dCTP, dGTP, dTTP

Kinase 10X Buffer

700mM Tris-HCl (pH 7.6)
100mM MgCl₂
50mM DTT

Klenow 10X Buffer

500mM Tris-HCl (pH 7.2)
100mM MgSO₄
1mM DTT

sample 2X buffer

20mM NaOH
20% glycerol
0.025% bromophenol blue (added fresh for each use)

Second Strand 2.5X Buffer

100mM Tris-HCl (pH 7.2)
225mM KCl
7.5mM MgCl₂
7.5mM DTT

TE buffer

10mM Tris-HCl (pH 8.0)
1mM EDTA

TEN buffer

100mM NaCl in TE buffer

TE-saturated phenol: chloroform:isoamyl alcohol

Mix equal parts of TE buffer and phenol and allow the phases to separate. Then mix 1 part of the lower, phenol phase with 1 part of chloroform:isoamyl alcohol (24:1).

T4 DNA Ligase 10X Buffer

300mM Tris-HCl (pH 7.8)
100mM MgCl₂
100mM DTT
5mM ATP

XIII. Related Products

cDNA Synthesis and Cloning Reagents

Product	Size	Cat.#
Oligo(dT) ₁₅ Primer	20µg	C1101
Random Primers*	20µg	C1181
EcoR I Adaptors	150pmol	C1291
Spin Columns	10 each	C1281
Sephacryl® S-400*	10ml	V3181
1.2kb Kanamycin Positive Control RNA*	5µg	C1381

*For Laboratory Use

XIII. Related Products (continued)

Product	Size	Cat.#
Packagene® Lambda DNA Packaging System	6 extracts	K3154
	11 extracts	K3152

RNA Isolation Systems

Product	Size	Cat.#
RNAgents® Total RNA Isolation System	Scalable	Z5110

For Laboratory Use. Each system contains sufficient reagents for six RNA isolations, each from 1g of tissue or 10⁸ cultured cells.

Product	Cat.#
PolyAtract® mRNA Isolation System I	Z5210
PolyAtract® mRNA Isolation System II	Z5200

For Laboratory Use. These systems contain sufficient reagents to perform three separate mRNA isolations, each from 1–5mg total RNA. System II includes the Magnetic Separation Stand.

Product	Cat.#
PolyAtract® mRNA Isolation System III	Z5300
PolyAtract® mRNA Isolation System IV	Z5310

For Laboratory Use. These systems contain sufficient reagents to perform fifteen separate mRNA isolations, each from approximately 100–1,000µg of total RNA. System III includes the Magnetic Separation Stand.

Product	Size	Cat.#
PolyAtract® System 1000	Scalable	Z5400
PolyAtract® System 1000 with Magnetic Separation Stand	Scalable	Z5420

For Laboratory Use. These systems contain sufficient reagents to isolate poly(A) RNA from up to 2g of tissue.

Product	Size	Cat.#
PolyAtract® System 1000 Magnetic Separation Stand	1 each	Z5410

XIV. References

1. Okayama, H. and Berg, P. (1982) High-efficiency cloning of full-length cDNA. *Mol. Cell. Biol.* **2**, 161–70.
2. Gubler, U. and Hoffman, B.J. (1983) A simple and very efficient method for generating cDNA libraries. *Gene* **25**, 263–69.
3. Kimmel, A.R. and Berger, S.L. (1987) Preparation of cDNA and the generation of cDNA libraries: Overview. *Meth. Enzymol.* **152**, 307–16.
4. Kozak, M. (1983) Comparison of initiation of protein synthesis in procaryotes, eucaryotes, and organelles. *Microbiol. Rev.* **47**, 1–45.

5. Han, J.H., Stratowa, C. and Rutter, W.J. (1987) Isolation of full-length putative rat lysophospholipase cDNA using improved methods for mRNA isolation and cDNA cloning. *Biochemistry* **26**, 1617-25.
6. Eschenfeldt, W.H., Puskas, R.S. and Berger, S.L. (1987) Homopolymeric tailing. *Meth. Enzymol.* **152**, 337-42.
7. Palazzolo, M.J. and Meyerowitz, E.M. (1987) A family of lambda phage cDNA cloning vectors, lambda SWAJ, allowing the amplification of RNA sequences. *Gene* **52**, 197-206.
8. Koike, S., Sakai, M. and Muramatsu, M. (1987) Molecular cloning and characterization of rat estrogen receptor cDNA. *Nucl. Acids Res.* **15**, 2499-513.
9. Blumberg, D.D. (1987) Creating a ribonuclease-free environment. *Meth. Enzymol.* **152**, 20-4.
10. Ehrenberg, L., Fedorcsak, I. and Solymosy, F. (1976) Diethyl pyrocarbonate in nucleic acid research. *Prog. Nucleic Acid Res. and Mol. Biol.* **16**, 189-262.
11. Chirgwin, J.M. *et al.* (1979) Isolation of biologically active ribonucleic acid from sources enriched in ribonuclease. *Biochemistry* **18**, 5294-99.
12. Aviv, H. and Leder, P. (1972) Purification of biologically active globin messenger RNA by chromatography on oligothymidylic acid-cellulose. *Proc. Natl. Acad. Sci. USA* **69**, 1408-12.
13. Krug, M. and Berger, S. (1987) First-strand cDNA synthesis primed with oligo(dT). *Meth. Enzymol.* **152**, 316-25.
14. Sambrook, J., Fritsch, E.F. and Maniatis, T. (1989) *Molecular Cloning: A Laboratory Manual*, 2nd ed., Cold Spring Harbor Laboratory, Cold Spring Harbor, New York.

©U.S. Pat. No. 5,552,302, European Pat. No. 0 422 217, Australian Pat. No. 646803 and Japanese Pat. Nos. 3009458 and 3366596 have been issued to Promega Corporation for the methods and compositions for production of human recombinant placental ribonuclease inhibitor.

©U.S. Pat. Nos. 4,966,964, 5,019,556 and 5,266,687, Australian Pat. Nos. 616881 and 641261 and other pending and issued patents, which claim vectors encoding a portion of human placental ribonuclease inhibitor, are exclusively licensed to Promega Corporation.

©Licensed under U.S. Pat. No. 4,935,361.

© 1996-2005 Promega Corporation. All Rights Reserved.

MagneSphere, Packagene, pGEM, PolyATtract, RiboClone, RNasin, RNAagents and Wizard are registered trademarks of Promega Corporation.

Sephacryl is a registered trademark of Amersham Biosciences Ltd. Whatman is a registered trademark of Whatman Paper Company, Ltd.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All prices and specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.