

Technical Manual

pTARGET™ Mammalian Expression Vector System

INSTRUCTIONS FOR USE OF PRODUCT A1410.

PRINTED IN USA. Revised 2/06

Part# TM044

vww.promega.com

O Promega

pTARGET[™] Mammalian Expression Vector 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	1
II.	Vector Maps	2
	A. pTARGET TM Vector Multiple Cloning Region	2
	B. pTARGET TM Vector Map and Sequence Reference Points	3
III.	Product Components and Storage Conditions	4
IV.	Ligation and Transformation Protocols	4
	A. Ligations Using pTARGET TM Vector	4
	B. Transformations Using pTARGET [™] Vector	5
	C. Isolating Recombinant Plasmid DNA	7
	D. Transfecting Mammalian Cells with the pTARGET TM Vector	7
V.	General Considerations for Cloning PCR Products	8
	A. Generating PCR Product	8
	B. PCR Product Purity	8
	C. Optimizing Insert: Vector Molar Ratios	9
	D. Screening Transformants for Inserts	9
	E. Experimental Controls	10
VI.	Troubleshooting	11
VII.	Appendix	14
	A. Vector Components	14
	B. pTARGET TM Vector Restriction Enzyme Tables	17
	C. Composition of Buffers and Solutions	19
	D. References	20
	E. Related Products	23

I. Description

The pTARGETTM Mammalian Expression Vector System^(a,b,c) is a convenient system for cloning PCR^(d) products and for expression of cloned PCR products in mammalian cells. The vector is prepared by cutting the pTARGETTM Vector with *Eco*R V and adding a 3' terminal thymidine to both ends. These single 3'-T overhangs at the insertion site greatly improve the efficiency of ligation of a PCR product into the plasmid in two ways. First the overhangs prevent recircularization of the vector; second they provide a compatible overhang for PCR products generated by certain thermostable polymerases (1,2). As

summarized in Table 1, these polymerases often add a single deoxyadenosine in a template-independent fashion to the 3´-ends of amplified fragments (3,4). The pTARGETTM Vector also contains a modified version of the coding sequence of the α -peptide of β -galactosidase, allowing blue/white recombinant screening.

The pTARGETTM Vector carries the human cytomegalovirus (CMV) immediateearly enhancer/promoter^(c) region to promote constitutive expression of cloned DNA inserts in mammalian cells. This vector also contains the neomycin phosphotransferase gene, a selectable marker for mammalian cells. The pTARGETTM Vector can be used for transient expression or for stable expression by selecting transfected cells with the antibiotic G-418.

Table 1. Comparison of PCR Product Properties for Some Thermostable DNA Polymerases.

	Thermostable DNA Polymerase								
Characteristic	<i>Taq/</i> AmpliTaq	® Tfl	Tth	Vent _R ®/ (<i>Tli</i>)	Deep Vent _R ®	Pfu	Pwo		
Resulting DNA ends	3′ A	3´ A	3′ A	>95% Blunt	>95% Blunt	Blunt	Blunt		
5´→3´ exonuclease activit	y Yes	Yes	Yes	No	No	No	No		
3´→5´ exonuclease activit	y No	No	No	Yes	Yes	Yes	Yes		

II. Vector Maps

II.A. pTARGET[™] Vector Multiple Cloning Region



Figure 1. The promoter and multiple cloning region of the pTARGETTM Vector. The top strand of the sequence shown corresponds to the RNA synthesized by T7 RNA polymerase. The top strand shown is the same as the ssDNA produced by this vector. The codons of the *lacZ* α region within the multiple cloning region are indicated by the spacing of the nucleotides.

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison, WI
 53711-5399
 USA

 Toll Free in USA 800-356-9526
 Phone 608-274-4330
 Fax
 608-277-2516
 www.promega.com

 Part# TM044
 Printed in USA.
 Revised 2/06
 Revised 2/06



II.B. pTARGET™ Vector Map and Sequence Reference Points

Figure 2. Vector circle map and sequence reference points.

pTARGET[™] Vector sequence reference points:

Cytomegalovirus immediate early enhancer	1-659
Cytomegalovirus immediate early promoter	669–750
Chimeric intron	890-1022
pTargeT TM Sequencing Primer	1367-1344
<i>lac</i> Zα start codon	1377
<i>lac</i> Zα stop codon	1053
lac operon sequences	1066-1226, 1363-1499
lac operator	1397-1413
T7 promoter (-17 to +3)	1229-1248
Multiple cloning region	1250-1323
SV40 late polyadenylation signal	1535-1755
Phage f1 region	1798-2252
Neomycin selectable marker	
SV40 enhancer and early promoter	2260-2630
SV40 minimum origin of replication	2528-2593
coding region of neomycin phosphotransferase	2675-3469
synthetic polyadenylation signal	3533-3581
β-lactamase (Amp ^r) coding region	3978-4838



Product	Size	Cat.#
pTARGET™ Mammalian Expression Vector System	20 reactions	A1410
Includes:		

III. Product Components and Storage Conditions

- $1.2\mu g$ pTargeTTM Vector (60ng/µl)
- 12μl Control Insert DNA (4ng/μl)
- 100u T4 DNA Ligase
- 200µl T4 DNA Ligase 10X Buffer
- 1.2ml JM109 Competent Cells, High Efficiency (6 × 200µl)
- 1 Protocol

Storage Conditions: Store Competent Cells at -70°C. See the expiration date on the tube label. All other components can be stored at -20°C or -70°C. T4 DNA Ligase 10X Buffer contains ATP, which degrades during temperature fluctuations. Avoid multiple freeze-thaws by making smaller aliquots of the buffer.

IV. Ligation and Transformation Protocols

IV.A. Ligations Using pTARGET[™] Vector

- 1. Briefly centrifuge the pTARGETTM Vector and Control Insert DNA tubes to collect the contents at the bottom of the tube.
- 2. Set up ligation reactions as described below.

Note: Use 0.5ml tubes known to have low DNA-binding properties (e.g., Sarstedt Cat.# 72687005).

	Standard Reaction	Positive Control	Background Control
T4 DNA Ligase 10X Buffer	1µl	1µl	1µl
pTargeT TM Vector (60ng)	1µl	1µl	1µl
PCR product	Xµl*	-	-
Control Insert DNA (8ng)	-	2µ1	-
T4 DNA Ligase (3 Weiss units/μl)	1µl	1µl	1µl
deionized water to a final volume of	10µl	10µl	10µl

*Molar ratio of PCR product to pTARGETTM Vector may require optimization (see Section V.C).

3. Incubate overnight at 4°C. Shorter incubation times (3-hour minimum) may be used but may result in fewer colonies.

Notes:

- 1. Use only the T4 DNA Ligase supplied with this system in performing pTARGETTM Vector ligations. Other commercial preparations of T4 DNA ligase may contain exonuclease activity, which may remove the terminal thymidines from the vector.
- 2. T4 DNA Ligase 10X Buffer contains ATP, which can degrade with temperature fluctuations. Avoid multiple freeze-thaw cycles by making smaller aliquots of the buffer.
- 3. Vortex the thawed T4 DNA Ligase 10X Buffer before use.
- 4. When using T4 DNA Ligase 10X Buffer, low temperature ligations are necessary for annealing single-base overhangs. Ligation temperatures higher than 15°C may significantly reduce the number of recombinants.

IV.B. Transformations Using pTARGET[™] Vector

Use only high-efficiency competent cells ($\geq 1 \times 10^{8}$ cfu/µg DNA) for transformations. The ligation efficiency of fragments with a single-base overhang can be inefficient; use competent cells that have transformation efficiencies of 1×10^{8} cfu/µg DNA (or higher) in order to obtain a reasonable number of colonies (see Section V.E).

Note: The JM109 competent cells supplied with the pTARGET[™] System are ready for blue/white selection. No further selection on M9 plates is necessary.

We highly recommend using JM109 High Efficiency Competent Cells (Cat.# L2001). In our experience, other strains such as DH5 α^{TM} and XL1-Blue have a much higher frequency of background white colonies than JM109. If you make your own competent cells, grow and maintain JM109 cells on M9 minimal medium plates supplemented with thiamine hydrochloride prior to the preparation of competent cells. This will ensure that the cells retain the F' episome. The episome contains the genetic information for proline biosynthesis, a nutritional requirement for growth on minimal medium. The F' episome also includes *lac*IqZ Δ M15, which is required in the blue/white color screening process. Selection for transformants should be on LB/ampicillin/IPTG/X-Gal plates (see Section VII.C). For best results, do not use plates more than 30 days old.

The genotype of JM109 is recA1, endA1, gyrA96, thi, hsdR17 ($r_{K}-,m_{K}+$), relA1, supE44, Δ (*lac-proAB*), [F', *traD36*, proAB, *lacI* $qZ\Delta$ M15] (5).

IV.B. Transformations Using pTARGET[™] Vector (continued)

Materials to Be Supplied by the User

(Solution compositions are provided in Section VII.C.)

- LB plates with ampicillin/IPTG/X-Gal
- SOC medium
- 0.1-1.0ng uncut plasmid for determination of transformation efficiency
- 1. Prepare two LB/ampicillin/IPTG/X-Gal plates for each ligation reaction, plus two plates for determining transformation efficiency (see Section V.E). Equilibrate the plates to room temperature prior to plating (Step 10).
- Remove frozen JM109 High Efficiency Competent Cells from -70°C storage and place on ice until just thawed (about 5 minutes). Mix the cells by gently flicking the tube. Prechill tubes on ice for each transformation reaction.
- 3. For each ligation reaction and control, carefully transfer 50µl of cells into each tube prepared in Step 2 (100µl of cells for control to determine transformation efficiency). Avoid excessive pipetting because the cells are very fragile.
- 4. Centrifuge the tubes containing the ligation reactions to collect the contents at the bottom of the tube. Add 2µl of each ligation reaction to a sterile 1.5ml microcentrifuge tube on ice. Add 0.1–1.0ng uncut plasmid to the tube containing the 100µl cells for determining the transformation efficiency of the competent cells (see Section V.E).

Note: In our experience, using larger 17 × 100mm polypropylene tubes (e.g., Falcon Cat.# 2059) can increase transformation efficiency. Tubes from some manufacturers bind DNA. These tubes should be avoided.

- 5. Gently flick the tubes to mix and place them on ice for 20 minutes.
- 6. Heat shock the cells for 45–50 seconds at exactly 42°C in a water bath. Do Not Shake.
- 7. Immediately return the tubes to ice for 2 minutes.
- Add 950µl of room temperature SOC medium to each sample tube and 900µl SOC medium to the transformation control tubes (LB broth may be substituted, but colony number may be lower).
- 9. Incubate for 1.5 hours at 37°C with shaking (~150rpm).
- 10. Plate 100 μ l of each transformation reaction onto duplicate antibiotic plates. If a higher number of colonies is desired, the cells may be pelleted by centrifugation at 1,000 × *g* for 10 minutes, resuspended in 100 μ l of SOC medium and plated on one plate. For the transformation control, we recommend plating 100 μ l a 1:10 dilution with SOC medium.
- 11. Incubate the plates overnight (16–24 hours) at 37°C. We routinely see approximately 100 colonies per plate when using competent cells that are 1 × 10%cfu/µg DNA if a 100µl aliquot is plated. Longer incubations or

 Promega Corporation
 2800
 Woods
 Hollow
 Road
 Madison,
 WI
 53711-5399
 USA

 Toll Free in USA 800-356-9526
 Phone 608-274-4330
 Fax
 608-277-2516
 www.promega.com

 Part# TM044
 Printed in USA.
 Revised 2/06

storing plates at 4°C (after 37°C overnight incubation) may facilitate blue/white screening. White colonies generally contain inserts, but inserts may also be present in blue colonies. Please see Section V.D for more information.

IV.C. Isolating Recombinant Plasmid DNA

A convenient and reliable method for isolating plasmid DNA is the Wizard® *Plus* SV Minipreps DNA Purification System (Cat.# A1330). Alternatively, a standard plasmid miniprep procedure, such as the one described in the *Protocols and Applications Guide* (6), may be used.

The Wizard[®] *Plus* SV Minipreps DNA System, a silica membrane-based spin or vacuum-driven system, isolates plasmid DNA that is ready for direct use in molecular biology applications, including restriction enzyme digestion and automated fluorescent sequencing.

IV.D. Transfecting Mammalian Cells with the pTARGET[™] Vector

The pTARGETTM Vector can be used for both transient and stable expression of genes. Transfection of DNA into mammalian cells may be mediated by cationic lipids (7,8), calcium phosphate (9,10), DEAE-dextran (11-13), polybrene-DMSO (14,15) or electroporation (16,17). Transfection systems based on the cationic lipids (TransFast™ Transfection Reagent, Transfectam® Reagent and Tfx[™] Reagents), calcium phosphate and DEAE-dextran (ProFection® Mammalian Transfection Systems) are available from Promega. For more information on using TransFast™ Transfection Reagent, please request the TransFast[™] Transfection Reagent Technical Bulletin (#TB260). For more information and a protocol for the Transfectam® Reagent, please request the Transfectam[®] Reagent for the Transfection of Eukaryotic Cells Technical Bulletin (#TB116). For more information on transfection using the Tfx[™] Reagents, please request the *Tfx*TM-20 and *Tfx*TM-50 Reagents for the Transfection of Eukaryotic Cells Technical Bulletin (#TB216). For transfection procedures using calcium phosphate or DEAE-dextran, please request the ProFection® Mammalian Transfection Systems Technical Manual (#TM012).

Note: Visit the Transfection Assistant for help in selecting a transfection reagent: www.promega.com/transfectionasst/

For stable expression, the transfected cells must be selected with the antibiotic G-418 (Cat.# V7981). Following transfection, seed the cells at a low density and apply the G-418 antibiotic to the medium at a concentration between $100\mu g/ml$ and 1mg/ml. The concentration of G-418 required to select and maintain drug resistance depends on the cell type and growth rate. In general, mammalian cells require a concentration of $400-600\mu g/ml$ of active G-418 for selection and $200-400\mu g/ml$ of active G-418 for maintenance of stable transformants (18). A stock solution of 100mg/ml of G-418 can be made up in 40mM HEPES (pH 7.3), PBS or water. (Store the stock solution of G-418 at -20° C.) For effective selection, the cells should be subconfluent, since confluent, nongrowing cells are very resistant to the effects of G-418. Change

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison, WI
 53711-5399
 USA

 Toll Free in USA 800-356-9526
 Phone 608-274-4330
 Fax
 608-277-2516
 www.promega.com

 Printed in USA.
 Part# TM044

 Revised 2/06
 Page 7

the medium every 3 days until drug-resistant clones appear (2–5 weeks, depending on the cell type). Death of cells not expressing neomycin phosphotransferase may require 3–9 days following addition of G-418. Cell strains vary in the level of resistance to antibiotics, so the level of resistance of a particular cell strain must be tested before attempting stable selection. A "kill curve" will determine the minimum concentration of antibiotic needed to kill nontransfected cells.

Note: Prepare a control plate for all selection experiments by treating untransfected cells with G-418 in medium under the experimental conditions. This control plate will indicate whether the conditions for antibiotic selection were sufficiently stringent to eliminate cells not expressing neomycin phosphotransferase.

V. General Considerations for Cloning PCR Products

V.A. Generating PCR Product

For efficient expression of the gene, the first ATG in the 5'-region of the mRNA transcript should be the translational start site, since upstream, out-of-frame ATGs can significantly reduce the rate of translation of genes in eukaryotes (19). Another consideration in designing PCR products is the inclusion of the consensus sequence ($-9GCCGCC(^A/_G)CCAUGG+4$) surrounding the initiation of translation. Kozak reported that efficient translation is obtained as long as the -3 position relative to the ATG contains a purine base (19). However, if the -3 position does not contain a purine, efficient translation can still be obtained if a guanine is at the +4 position.

The fidelity of the PCR depends upon the cycle number, the polymerase and the reaction conditions (20). Because successful expression of the inserted gene product depends upon integrity of the sequence, optimizing PCR fidelity is important (20). The DNA sequence cloned into the pTARGET[™] Vector should also be verified. To sequence upstream of the multiple cloning region toward the insert, use the T7 Promoter Primer (Cat.# Q5021). To sequence downstream of the multiple cloning region toward the insert, use the pTARGET[™] Sequencing Primer (Cat.# Q4461).

V.B. PCR Product Purity

Analyze an aliquot of the PCR amplification product on an agarose gel before using it in the ligation reaction. Several convenient methods exist for purifying the PCR products from an agarose gel. The Wizard® SV Gel and PCR Clean-Up System (Cat.# A9281) is designed to extract and purify DNA fragments of 100bp-10kb from standard or low-melt agarose gels. During gel purification, use longwave ultraviolet light (366nm) to visualize the PCR product so that the formation of pyrimidine dimers is minimized. If gel separation is not required, the Wizard® SV Gel and PCR Clean-Up System can be used to purify the bands of interest directly from the reaction mix.

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison, WI
 53711-5399
 USA

 Toll Free in USA 800-356-9526
 Phone 608-274-4330
 Fax
 608-277-2516
 www.promega.com

 Part# TM044
 Printed in USA.
 Revised 2/06

Even if distinct bands of the expected size are observed, primer-dimers should be removed. Using crude PCR product may also produce successful ligations; however, the number of white colonies containing the relevant insert may be reduced because of preferential incorporation of primer-dimers or other extraneous reaction products. Therefore, you may need to screen numerous colonies in order to identify clones that contain the insert of interest when you use crude product.

V.C. Optimizing Insert:Vector Molar Ratios

The pTARGETTM Vector System uses a 1.4:1 molar ratio of the Control Insert DNA to the pTARGETTM Vector. Ratios of 8:1 to 1:8 have been used successfully. If initial experiments with your PCR product are suboptimal, you may need to optimize the ratio of insert to vector. Ratios from 3:1 to 1:3 provide good initial parameters. Estimate the concentration of PCR product by comparing it to DNA mass standards on a gel or by using a fluorescent assay (21). The pTARGETTM Vector is 5.67kb in size and is supplied at a concentration of 60ng/µl. To calculate the appropriate amount of PCR product (insert) to include in the ligation reaction, use the following equation:

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

Sufficient $pTargeT^{\texttt{TM}}$ Vector is provided to vary insert:vector ratios as recommended and to perform control reactions.

Example of insert:vector ratio calculation:

How much 0.5kb PCR product should be added to a ligation in which 60ng of 5.67kb vector will be used if a 3:1 insert:vector molar ratio is desired?

 $\frac{60\text{ng vector} \times 0.5\text{kb insert}}{5.67\text{kb vector}} \times 3/1 = 15.9\text{ng insert}$

Note: Using the same parameters for a 1:1 insert:vector molar ratio, 5.3ng of a 0.5kb insert would be required.

V.D. Screening Transformants for Inserts

Successful cloning of an insert in the pTARGETTM Vector interrupts the coding sequence of the α -peptide of β -galactosidase. Recombinant clones can usually be identified by color screening on indicator plates. However, the characteristics of PCR products cloned into the pTARGETTM Vector can significantly affect the ratio of blue:white colonies obtained following transformation of competent cells. Clones that contain PCR products produce white colonies in most cases, but blue colonies can result from PCR fragments that are cloned in-frame with the *lacZ* gene. Such fragments are usually a multiple of 3 base pairs long (including the 3'-A overhangs) and do not contain in-frame stop codons.

V.D. Screening Transformants for Inserts (continued)

Even if your expected PCR product is not a multiple of 3 bases long, the amplification process can introduce mutations (e.g., deletions or point mutations) that may result in blue colonies when the fragment is inserted into the pTARGETTM Vector.

The Control Insert DNA supplied with the pTARGETTM Vector System is a 542bp sequence containing multiple stop codons in all six reading frames, which ensures a low background of blue colonies for the control reaction. Results obtained with the Control Insert DNA may not be representative of those achieved with your PCR product.

V.E. Experimental Controls

We strongly recommend performing the controls detailed below. These are necessary to accurately assess the performance of the pTARGETTM Vector System.

Positive Control

Set up a ligation reaction with the Control Insert DNA as described in the protocol (Section IV.A) and use it for transformations as described in Section IV.B. This control will allow you to determine the ligation efficiency. You will typically observe approximately 100 colonies, of which 10–30% are blue, if you use competent cells that have a transformation efficiency of 1 × 10⁸cfu/µg DNA. Greater than 60% of the colonies should be white and, therefore, recombinant, since the Control Insert DNA is designed to reduce the number of background blue colonies (discussed in Section V.D). Background blue colonies arise from non-T-tailed or undigested pTARGET[™] Vector. These blue colonies are a useful internal transformation control; if no colonies are obtained, the transformation has failed. If blue colonies are obtained, but no whites, the result suggests that the ligation reaction failed. If <50% white colonies are seen in this positive control reaction, then the ligation conditions were probably suboptimal. Also, fewer total colonies can indicate inefficient ligation.

Note: The Control Insert DNA should only be used as a ligation control and should not be used as a control for expression in mammalian cells.

Background Control

Set up a ligation reaction with 60ng of pTARGETTM Vector and no insert as described in the protocol (Section IV.A) and use it for transformations as described in Section IV.B. This ligation will allow you to determine the number of background blue colonies resulting from non-T-tailed or undigested pTARGETTM Vector.

If you follow the recommendations in Section IV.B closely, you should observe 20-60 colonies from competent cells with a transformation efficiency of 1×10^8 cfu/µg DNA. Of these, approximately 4–15 colonies will be white. The background white colonies seen with the pTARGETTM Vector result from

 Promega Corporation
 2800
 Woods
 Hollow
 Road
 Madison,
 WI
 53711-5399
 USA

 Toll
 Free in
 USA
 800-356-9526
 Phone
 608-274-4330
 Fax
 608-277-2516
 www.promega.com

 Part#
 TM044
 Printed in
 USA.

 Page 10
 Revised 2/06
 Revised 2/06

deletions generated in the $lacZ\alpha$ region during transformation. Deletions of plasmid DNA during the transformation of *E. coli* with linear plasmid molecules are a well-documented phenomenon (22–25).

Transformation Control

Check the transformation efficiency of the competent cells by transforming them with an uncut plasmid and calculating $cfu/\mu g$ DNA. If the transformation efficiency is lower than $1 \times 10^{8} cfu/\mu g$ DNA, prepare fresh cells.

Example of transformation efficiency calculation:

After 100µl competent cells are transformed with 0.1ng uncut plasmid DNA, the transformation reaction is added to 900µl of SOC medium (0.1ng DNA/ml). From that volume, a 1:10 dilution with SOC medium (0.01ng DNA/ml) is made and 100µl plated on two plates (0.001ng DNA/100µl). If 200 colonies are obtained per plate (average of two plates), the transformation efficiency is calculated as shown below:

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

VI. 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
No colonies	Problem with the transformation reaction or cells not competent. Background undigested vector and religated non-T-tailed vector should yield 20–60 blue colonies independent of the presence of insert DNA. Check the background control (Section V.E). Test the efficiency of the cells by transforming them with an uncut plasmid that allows antibiotic selection, such as the pGEM®-5Zf(+) Vector. If the recommendations in Section IV.B are followed, cells that are 1 × 10 ⁸ cfu/µg DNA typically yield 100 colonies. Therefore, you would not see any colonies from cells that are 1 × 10 ⁶ cfu/µg DNA.
Less than 10% white colonies in the positive control ligation	Failed ligation. Ligase buffer may have low activity. T4 DNA Ligase 10X Buffer contains ATP, which degrades during temperature fluctuations. Avoid multiple freeze-thaw cycles by making smaller aliquots of the buffer. Use a fresh vial of buffer.

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison, WI
 53711-5399
 USA

 Toll Free in USA
 800-356-9526
 Phone 608-274-4330
 Fax
 608-277-2516
 www.promega.com

 Printed in USA.
 Part# TM044

 Revised 2/06
 Page 11



Symptoms	Causes and Comments
Less than 10% white colonies in the positive control ligation (continued)	Failed ligation. Ligase activity may be low. To test the activity of the ligase and buffer, set up a ligation with approximately 200ng of DNA markers (e.g., Lambda DNA/ <i>Hind</i> III Markers, Cat.# G1711). Compare ligated and nonligated DNA on a gel and check that the fragments have been religated into high molecular weight material.
	Loss of T-overhangs allowing blunt-ended ligation of vector. Avoid introducing nucleases that may degrade the T-overhangs. Use only the T4 DNA Ligase provided with the system that has been tested for minimal exonuclease activity.
Less than 60% white colonies in the positive control ligation	Ligation temperature too high. Low temperature ligations (i.e., 4°C) favor annealing of the A-overhang of the insert with the T-overhang of the vector. Higher temperatures (i.e., room temperature) favor blunt-ended ligations.
	Ligation temperature too high. There is a background of non-T-tailed vector that will favor blunt-ligated at higher ligation temperatures.
	Loss of T-overhangs allowing blunt-ended ligation of vector. Avoid introduction of nucleases that may degrade the T-overhangs. Use only the T4 DNA Ligase provided with the system that has been tested for minimal <u>exonuclease activity</u> .
	Ligation incubation too short. Overnight ligation yields optimal results.
High colony number but low percentage of white colonies in the positive control ligation	Failed ligation. If the ligation is suboptimal or fails, and the competent cells have a high transformation frequency $(1 \times 10^{\circ}$ cfu/µg), the total number of colonies will be high, but the amount of white colonies will be low.
Low number or no white colonies with PCR product	Failed ligation. See above comments.
-	Ligation incubation too short. Overnight ligation yields optimal results.

VI. Troubleshooting (continued)

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison,
 WI
 53711-5399
 USA

 Toll Free in USA 800-356-9526
 Phone 608-274-4330
 Fax
 608-277-2516
 www.promega.com

 Part# TM044
 Printed in USA.
 Revised 2/06
 Revised 2/06



Symptoms	Causes and Comments			
Low number or no white colonies with PCR product (continued)	Failed ligation due to inhibitory component in the PCR product. Mix some of the PCR product with the control ligation to see if it is inhibiting the reaction. If an inhibitor is suspected, repurify the PCR fragment.			
	No 3'-A overhangs to allow PCR product ligation. As summarized in Table 1, not all thermostable DNA polymerases create a 3'A overhang (3,4). Blunt-ended fragments may be subsequently A-tailed by treatment with an appropriate polymerase and dATP (26,27).			
	PCR fragment does not disrupt the <i>lacZ</i> gene. If there are a higher number of blue colonies resulting from the PCR fragment ligation than from the background control, some of these blue colonies may contain insert (especially with smaller insert fragments). Screen blue and pale blue colonies.			
	Nonoptimal insert:vector ratio. Check the integrity and quantity of your PCR fragment by gel analysis. Optimize the insert:vector ratio (see Section V.C).			
PCR product ligation reaction produces white colonies that do not contain insert (blue colonies are present)	Primer-dimers present in PCR fragment preparation. Primer-dimers will ligate into the pTARGET [™] Vector but may not be seen after gel analysis of recombinant plasmid DNA because of their small size. Gel purify the PCR fragment.			
	Multiple PCR products generated and cloned. Gel purify the PCR fragment of interest.			
	DNA rearranged. Check a number of clones to see whether rearrangement is random. If so, the clone of interest should be present and can be identified by screening several clones. Toxic, unstable or homologous elements may also be a factor. Grow cells at a lower temperature or use a recombination-deficient cell line.			
PCR product ligation reaction produces white colonies only (no blue colonies are present)	Inactive ampicillin. Check that ampicillin plates are made properly and used within two weeks. Test ampicillin by streaking plates with and without ampicillin with an ampicillin-sensitive strain of <i>E. coli</i> .			

VI. Troubleshooting (continued)

PromegaCorporation2800WoodsHollowRoadMadison,WI53711-5399USATollFreeinUSA800-356-9526Phone608-274-4330Fax608-277-2516•www.promega.comPrinted in USA.Part# TM044Page 13Page 13



Symptoms	Causes and Comments		
PCR product ligation reaction produces white colonies only (no blue colonies are present; continued)	No F' episome in bacterial strain. Check the background control. If these colonies are not blue, and fresh plates that contain ampicillin/ IPTG/X-Gal were used, the cells may have lost the F' episome. Be sure that the cells are prepared properly (see Section IV.B).		
	Plates incompatible with blue/white screening. Check the background control. If these colonies are not blue, check that the plates have ampicillin/IPTG/X-Gal and that they are fresh. If there is any question about the quality of the plates, repeat plating with fresh plates.		
Not enough clones contain the PCR product of interest	Insufficient A-tailing of of the PCR fragment. After purifying the PCR fragment, set up an A-tailing reaction (28,29). Clean up the sample and proceed with the protocol.		
Expression of recombinant protein not detected in the transfected cells	No ATG sequence present in insert. Insert DNA must contain an ATG for initiation of translation		
	Failed transfection. Check efficiency of transfection by cotransfecting with a reporter vector together with the pTARGET [™] Vector recombinant clone and measuring expression of reporter gene.		
	PCR-induced mutation in insert introducing a stop codon. Sequence insert to ensure no PCR-induced mutations are present.		

VI. Troubleshooting (continued)

VII. Appendix

VII.A. Vector Components

Enhancer/Promoter Regions

The CMV immediate-early enhancer/promoter region present in the pTARGETTM Vector allows strong, constitutive expression in a variety of cell types. The promiscuous nature of the CMV enhancer/promoter has been demonstrated in transgenic mice, where expression of the chloramphenicol acetyltransferase (CAT) gene regulated by the CMV enhancer/promoter was observed in 24 of the 28 tissues examined (28). The pTARGETTM Vector also contains the simian virus 40 (SV40) enhancer and early promoter region upstream of the neomycin phosphotransferase gene. The SV40 early promoter contains the SV40 origin of replication, which will induce transient, episomal replication of the pTARGETTM Vector in cells expressing the SV40 large T antigen such as COS-1 or COS-7 cells (29).

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison, WI
 53711-5399
 USA

 Toll Free in USA 800-356-9526
 Phone 608-274-4330
 Fax 608-277-2516
 www.promega.com

 Part# TM044
 Printed in USA.
 Revised 2/06



Chimeric Intron

Downstream of the CMV enhancer/promoter region is a chimeric intron composed of the 5'-donor site from the first intron of the human β -globin gene and the branch and 3'-acceptor site from the intron of an immunoglobulin gene heavy chain variable region (30). The sequences of the donor and acceptor sites, along with the branchpoint site, have been changed to match the consensus sequences for splicing (31). The intron is located upstream of the cDNA insert in order to prevent utilization of possible cryptic 5'-donor splice sites within the cDNA sequence (32).

Transfection studies have demonstrated that an intron flanking the cDNA insert frequently increases the level of gene expression (33–36). The increase in expression level due to the intron depends on the particular cDNA insert. For example, in transient transfections of 293 cells, the presence of this chimeric intron results in an approximate 20-fold increase in expression of the CAT gene (37). In contrast, the chimeric intron increases the gene expression level from luciferase cDNA by only threefold (37). In transgenic experiments, an intron is necessary to promote a high level of expression for virtually all cDNA inserts (38–40).

T7 RNA Polymerase Promoter

In the pTARGET[™] Vector, a T7 RNA polymerase promoter is located downstream of the intron. This promoter can be used to synthesize RNA transcripts in vitro using T7 RNA polymerase.

Multiple Cloning Region, Other Convenient Restriction Enzyme Sites and *lacZo* Region

The multiple cloning region in the pTARGETTM Vector is immediately downstream from the T7 promoter. Recognition sites for the restriction enzyme *Eco*R I flank the cloning site. This provides a convenient, single-enzyme digestion for release of the insert. Alternatively, a double-digestion may be used to release the insert.

Large hairpin structures in the 5´-end of untranslated mRNAs can reduce the level of in vitro and in vivo translation in higher eukaryotes (41–45). RNA transcribed from the multiple cloning region of the pTARGETTM Vector upstream of the cloning site does not contain hairpin structures that would interfere with translation (44).

Because there are no ATG sequences in either the multiple cloning region or between the transcriptional start site and the multiple cloning region, an ATG for the initiation of translation must be present in the inserted DNA.

Note: Additional requirements for efficient translation, including the Kozak consensus sequence (19), are described in Section V.A of this technical manual.

Unique restriction sites flank the CMV enhancer (*Bgl* II and *Sgf* I) and the CMV promoter (*Sgf* I and I-*Ppo* I), allowing easy replacement of these regulatory regions with other regulatory regions of interest.

VII.A. Vector Components (continued)

Multiple Cloning Region, Other Convenient Restriction Enzyme Sites and *lac*Zα Region (continued)

The multiple cloning region is within a $lacZ\alpha$ region, allowing identification of recombinant clones by color screening on indicator plates. The $lacZ\alpha$ region was modified to eliminate potential splice donor and acceptor sites and to eliminate an ATG sequence upstream of the cloning site. These modifications were made in an attempt to reduce any effect the $lacZ\alpha$ region may have on the expression level of the insert DNA. Recombinant clones may be identified by blue/white colony screening.

SV40 Late Polyadenylation Signal

Polyadenylation signals terminate transcription by RNA polymerase II and cause the addition of approximately 200 to 250 adenosine residues to the 3'-end of the RNA transcript (46). Polyadenylation enhances RNA stability and translation efficiency (47,48). To facilitate efficient processing of cloned DNA inserts not containing polyadenylation signals, the SV40 late polyadenylation signal has been positioned downstream from the multiple cloning region. The SV40 late polyadenylation signal is extremely efficient and can increase the steady-state level of RNA approximately fivefold more than the SV40 early polyadenylation signal (49).

Neomycin Phosphotransferase Selectable Marker

The neomycin phosphotransferase gene from Tn5 is present in the pTARGETTM Vector and is under the regulation SV40 enhancer and early promoter region. A synthetic polyadenylation signal based on the highly efficient polyadenylation signal of the rabbit β-globin gene (50) is located downstream of the neomycin phosphotransferase gene. To increase the translational efficiency of the neomycin phosphotransferase gene, the upstream, out-offrame ATG sequences present in the wildtype neomycin phosphotransferase gene have been eliminated. Additionally, the sequence upstream of the initiator ATG has been changed to match a sequence shown to improve the context for initiating translation (46).

Expression of the neomycin phosphotransferase gene in mammalian cells confers resistance to the antibiotic G-418 (18). G-418 is an aminoglycoside antibiotic produced by *Streptomyces*, which induces cytotoxicity by blocking translation. Neomycin phosphotransferase inactivates G-418 through phosphorylation, thereby blocking its toxic effects (51).

f1 Origin of Replication and Plasmid Replicon

The backbone for the pTARGETTM Vector is derived from the pGEM[®]-3Zf(+) Vector. As a result, this vector is a high-copy-number plasmid and contains the origin of replication of the filamentous phage f1. To generate single-stranded DNA (ssDNA) from the f1 origin, bacteria transformed with the pTARGETTM Vector carrying the cloned DNA of interest are infected with an appropriate

 Promega
 Corporation
 2800
 Woods
 Hollow
 Road
 Madison,
 WI
 53711-5399
 USA

 Toll
 Free in
 USA
 800-356-9526
 Phone
 608-274-4330
 Fax
 608-277-2516
 www.promega.com

 Part#
 TM044
 Printed in USA.
 Revised 2/06

helper phage. The plasmid then enters the f1 replication mode, and the resulting ssDNA is exported from the cell as an encapsidated virus particle. The ssDNA molecule exported corresponds to the top strand shown for the multiple cloning region (Figure 1). The ssDNA is purified from the supernatant by simple precipitation and extraction procedures, which are described in detail in the *Protocols and Applications Guide* (6).

Note: For further information about ssDNA purification, please contact your local Promega Branch Office or Distributor. In the U.S., contact Technical Services at 1-800-356-9526.

The "poison" sequence present in pBR322 that has been shown to inhibit replication of SV40 origin-containing vectors in COS cells has been deleted in the pTARGET[™] Vector (52). This results in more efficient expression of the cloned cDNAs in COS cells and other cells transformed with the SV40 large T antigen.

VII.B. pTARGET[™] Vector Restriction Enzyme Tables

The following restriction enzyme tables were constructed using DNASTAR[®] sequence analysis software with the pTARGETTM Vector prior to cutting with *Eco*R V and T-tailing. Please note that we have not verified this information by restriction digestion with each enzyme listed. The location given specifies the 3' end of the cut DNA (the base to the left of the cut site). For more information about the cut sites for these enzymes, or if you identify a discrepancy, please contact your local Promega Branch or Distributor. In the U.S., contact Promega Technical Services at 800-356-9526. The pTARGETTM Vector is supplied linearized with a "T" added to both 3' ends. The *Eco*R V site will not be recovered upon ligation of vector and insert. Vector sequences are available in the GenBank[®] database (GenBank[®]/EMBL Accession Number **AY540613**) and on the Internet at: www.promega.com/vectors/

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
Aat II	5	278, 331, 414,	Bal I	3	10, 64, 2885
		600, 3846	BamH I	1	1256
Acc I	1	1304	Ban II	3	729, 1956, 3168
Acc III	1	1775	Bbe I	1	2806
Acc65 I	1	1297	BbrP I	1	1849
Afl II	3	828, 847, 2657	Bbs I	2	961, 1540
Áfl III	1	1276	Bbu I	3	2357, 2429, 3208
Álw44 I	3	3596, 4093, 5339	Bgl II	1	5665
AlwN I	1	5244	Bsa I	2	915, 4699
Ava I	2	1270, 1291	BsaA I	4	493, 1849, 2027,
Ava II	3	3319, 4401, 4623			3107
Avr II	1	2609	BsaB I	1	1766

Table 2	Restriction	Enzymes	That (Cut the	pTargeT™	Vector	Between	1 and 5
Times.								

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison, WI
 53711-5399
 USA

 Toll Free in USA
 800-356-9526
 Phone 608-274-4330
 Fax
 608-277-2516
 www.promega.com

 Printed in USA.
 Part# TM044

 Revised 2/06
 Page 17

VII.B. pTARGETTM Vector Restriction Enzyme Tables (continued)

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
BsaM I	2	1585, 1678	Kpn I	1	1301
Bsm I	2	1585, 1678	Mlu I	1	1276
BspH I	3	3820, 3925, 4933	Nar I	1	2803
BspM I	4	877, 2690, 3071,	Nco I	3	513, 2516, 3235
		3521	Nde I	1	387
BsrG I	1	96	NgoM IV	2	1922, 3303
BssH II	1	3200	Nhe I	1	1264
BssS I	4	3395, 3789, 4096,	Not I	1	1311
		5480	Nsi I	2	2359, 2431
Bst98 I	3	828, 847, 2657	Nsp I	4	2357, 2429, 3208,
BstX I	1	3524			3740
BstZ I	2	1311, 2709	PaeR7 I	1	1270
Cfr10 I	4	1922, 3122, 3303,	Pml I	1	1849
		4680	<i>Рри</i> 10 I	2	2355, 2427
Cla I	1	1770	PspA I	1	1291
Csp I	1	3319	Pst I	2	838, 2856
Csp45 I	1	3485	Pvu I	2	664, 4396
Dra I	4	1725, 4187, 4879,	Pvu II	3	1135, 2285, 2909
		4898	Rsr II	1	3319
Dra II	1	3785	Sac I	1	729
Dra III	1	2030	Sal I	1	1303
Dsa I	3	513, 2516, 3235	Sca I	1	4284
Eag I	2	1311, 2709	Sfi I	1	2562
Ear I	4	1123, 3147, 3357,	Sgf I	1	664
		3966	Sin I	3	3319, 4401, 4623
EclHK I	1	4765	Sma I	1	1293
<i>Eco</i> 52 I	2	1311, 2709	SnaB I	1	493
<i>Eco</i> 72 I	1	1849	Spe I	1	152
EcoICR I	1	727	Sph I	3	2357, 2429, 3208
EcoR I	2	1250, 1318	Ssp I	4	5, 52, 2235, 3960
Ehe I	1	2804	Stu I	1	2608
Fsp I	3	1085, 2905, 4542	Sty I	4	513, 2516, 2609,
Hae II	4	1872, 1880, 2806,			3235
		5413	Tfi I	4	1772, 2631, 3288,
Hinc II	3	677, 1305, 1664			3422
Hind II	3	677, 1305, 1664	<i>Tth</i> 111 I	1	2921
Hind III	3	756, 1325, 2625	Vsp I	3	160, 1485, 4590
Hpa I	1	1664	Xho I	1	1270
I-Ppo I	1	851	Xma I	1	1291
Kas I	1	2802	Xmn I	1	4165

Table 2. Restriction Enzymes That Cut the $pT_{\text{ARGET}^{\text{TM}}}$ Vector Between 1 and 5 Times (continued).

Note: The enzymes listed in boldface type are available from Promega.

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison, WI
 53711-5399
 USA

 Toll Free in USA 800-356-9526
 Phone 608-274-4330
 Fax 608-277-2516
 www.promega.com

 Part# TM044
 Printed in USA.
 Revised 2/06

VII.B. pTARGETTM	Vector Restriction	Enzyme	Tables	(continued)
------------------	--------------------	--------	--------	-------------

BsiW I	EcoN I	Pme I	Srf I
Bsp120 I	Fse I	РриМ I	Sse8387 I
Bst1107 I	Ksp I	PshA I	Swa I
BstE II	Nru I	Psp5 II	Xba I
Bsu36 I	Pac I	Sac II	Xcm I
Eco47 III	PflM I	SgrA I	
<i>Eco</i> 81 I	PinA I	Spl I	
	BsiW I Bsp120 I Bst1107 I BstE II Bsu36 I Eco47 III Eco81 I	BsiW I EcoN I Bsp120 I Fse I Bst1107 I Ksp I BstE II Nru I Bsu36 I Pac I Eco47 III PflM I Eco81 I PinA I	BsiW I EcoN I Pme I Bsp120 I Fse I PpuM I Bst1107 I Ksp I PshA I BstE II Nru I Psp5 II Bsu36 I Pac I Sac II Eco47 III PflM I SgrA I Eco81 I PinA I Spl I

Table 3. Restriction Enzymes That Do Not Cut the pTARGET[™] Vector.

Table 4. Restriction Enzymes That Cut the pTARGET[™] Vector 6 or More Times.

Aci I	Bsp1286 I	Eae I	Hsp92 II	Nla III
Acy I	Bsr I	Fnu4H I	Mae I	Nla IV
Alu I	BsrS I	Fok I	Mae II	Ple I
<i>Alw</i> 26 I	Bst71 I	Hae III	Mae III	Rsa I
AspH I	BstO I	Hga I	Mbo I	Sau3A I
Ban I	BstU I	Hha I	Mbo II	Sau96 I
Bbv I	Cfo I	HinP I	Mnl I	ScrF I
Bgl I	Dde I	Hinf I	Mse I	SfaN I
BsaH I	Dpn I	Hpa II	Msp I	Ťaq I
BsaJ I	Dpn II	Hph I	MspA1 I	Tru9 I
BsaO I	Drd I	Hsp92 I	Nci I	Xho II

VII.C. Composition of Buffers and Solutions

IPTG stock solution (0.1M)

1.2g IPTG (Cat.# V3951)

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

X-Gal (2ml)

100mg 5-bromo-4-chloro-3 indolyl-β-D-galactoside

Dissolve in 2ml N,N'-dimethylformamide. Cover with aluminum foil and store at -20° C.

LB medium (per liter)

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

Adjust pH to 7.0 with NaOH.

LB plates with ampicillin

Add 15g agar to 1L of LB medium. Allow the medium to cool to 50° C before adding ampicillin to a final concentration of 100μ g/ml. Pour 30–35ml of medium into 85mm petri dishes. Let the agar harden. Store at 4°C for up to one month or at room temperature for up to one week.

LB plates with ampicillin/IPTG/ X-Gal

Make the LB plates with ampicillin as above; then supplement with 0.5mM IPTG and 80μ g/ml X-Gal and pour the plates. Alternatively, 100µl of 100mM IPTG and 20µl of 50mg/ml X-Gal may be spread over the surface of an LB-ampicillin plate and allowed to absorb for 30 minutes at 37°C prior to use.

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison, WI
 53711-5399
 USA

 Toll Free in USA 800-356-9526
 Phone 608-274-4330
 Fax 608-277-2516
 www.promega.com

 Printed in USA.
 Part# TM044

 Revised 2/06
 Page 19

VII.C. Composition of Buffers and Solutions (continued)

SOC medium (100ml)

- 2.0g Bacto®-tryptone
- 0.5g Bacto®-yeast extract
- 1ml 1M NaCl
- 0.25ml 1M KCl
 - 1ml 2M Mg²⁺ stock, filter-sterilized (as prepared below)
 - 1ml 2M glucose, filtersterilized

Add Bacto[®]-tryptone, Bacto[®]-yeast extract, NaCl and KCl to 97ml deionized water. Stir to dissolve. Autoclave and cool to room temperature. Add 2M Mg²⁺ stock and 2M glucose, each to a final concentration of 20mM. Bring to 100ml with sterile deionized water. Filter the complete medium through a 0.2µm filter unit. The final pH should be 7.0.

2M Mg²⁺ stock

20.33g MgCl₂ • 6H₂O 24.65g MgSO₄ • 7H₂O

Add distilled water to 100ml. Filter sterilize.

T4 DNA Ligase 10X Buffer (provided)

 300mM
 Tris-HCl (pH 7.8)

 100mM
 MgCl₂

 100mM
 DTT

 10mM
 ATP

Store in single-use aliquots at -20°C. Avoid multiple freeze-thaw cycles.

TYP broth (per liter)

16g	Bacto [®] -tryptone
()	

- 16g Bacto®-yeast extract
- 5g NaCl
- 2.5g K₂HPO₄

VII.D. References

- Mezei, L.M. and Storts, D.R. (1994) In: *PCR Technology: Current Innovations*, Griffin, H.G. and Griffin, A.M., eds., CRC Press, Boca Raton, FL, 21–8.
- Robles, J. and Doers, M. (1994) pGEM®-T Vector Systems Troubleshooting Guide. Promega Notes 45, 19–20.
- Clark, J.M. (1988) Novel non-templated nucleotide addition reactions catalyzed by procaryotic and eucaryotic DNA polymerases. *Nucl. Acids Res.* 16, 9677–86.
- Newton, C.R. and Graham, A. (1994) In: PCR, BIOS Scientific Publishers, Ltd., Oxford, UK, 13.
- Messing, J. et al. (1981) A system for shotgun DNA sequencing. Nucl. Acids Res. 9, 309–21.
- 6. Protocols and Applications Guide, Third Edition (1996) Promega Corporation.
- 7. Behr, J.P. *et al.* (1989) Efficient gene transfer into mammalian primary endocrine cells with lipopolyamine-coated DNA. *Proc. Natl. Acad. Sci. USA* **86**, 6982–6.
- Loeffler, J.P. et al. (1990) Lipopolyamine-mediated transfection allows gene expression studies in primary neuronal cells. J. Neurochem. 54, 1812–5.
- Graham, F.L. and van der Eb, A.J. (1973) A new technique for the assay of infectivity of human adenovirus 5 DNA. *Virology* 52, 456–67.
- 10. Wigler, M. et al. (1977) Transfer of purified herpes virus thymidine kinase gene to cultured mouse cells. Cell 11, 223–32.

 Promega Corporation
 2800
 Woods
 Hollow
 Road
 Madison,
 WI
 53711-5399
 USA

 Toll
 Free in
 USA
 800-356-9526
 Phone
 608-274-4330
 Fax
 608-277-2516
 www.promega.com

 Part#
 TM044
 Printed in
 USA.

 Page 20
 Revised 2/06
 2/06

- McCutchan, J.H. and Pagano, J.S. (1968) Enchancement of the infectivity of simian virus 40 deoxyribonucleic acid with diethylaminoethyl-dextran. J. Natl. Cancer Inst. 41, 351-7.
- Al-Moslih, M.I. and Dubes, G.R. (1973) The kinetics of DEAE-dextran-induced cell sensitization to transfection. J. Gen. Virol. 18, 189–93.
- Luthman, H. and Magnusson, G. (1983) High efficiency polyoma DNA transfection of chloroquine treated cells. *Nucl. Acids Res.* 11, 1295–8.
- Kawai, S. and Nishizawa, M. (1984) New procedure for DNA transfection with polycation and dimethyl sulfoxide. *Mol. Cell. Biol.* 4, 1172-4.
- Aubin, R.J., Weinfeld, M. and Paterson, M.C. (1988) Factors influencing efficiency and reproducibility of polybrene-assisted gene transfer. Son. Cell Mol. Genet. 14, 155–67.
- Andreason, G.L. and Evans, G.A. (1988) Introduction and expression of DNA molecules in eukaryotic cells by electroporation. *BioTechniques* 6, 650–60.
- Neumann, E. et al. (1982) Gene transfer into mouse lyoma cells by electroporation in high electric fields. EMBO J. 1, 841–5.
- Southern, P.J. and Berg, P. (1982) Transformation of mammalian cells to antibiotic resistance with a bacterial gene under control of the SV40 early region promoter. *J. Molec. Appl. Genet.* 1, 327–41.
- Kozak, M. (1989) The scanning model for translation: an update. J. Cell Biol. 108, 229-41.
- Eckert, K.A. and Kunkel, T.A. (1992) In: PCR: A Practical Approach, McPherson, M.J., Quirke, P. and Taylor, G.R., eds., IRL Press, NY.
- Sambrook, J. and Russell, W.D., eds. (2001) Molecular Cloning: A Laboratory Manual Cold Spring Harbor Laboratory Press, Cold Spring Harbor, NY, A8.22–8.23.
- Conley, E.C. and Saunders, J.R. (1984) Recombination-dependent recircularization of linearized pBR322 plasmid DNA following transformation of *Escherichia coli*. *Mol. Gen. Genet.* 194, 211–8.
- Conley, E.C. et al. (1986) Mechanism of intramolecular recyclization and deletion formation following transformation of *Escherichia coli* with linearized plasmid DNA. *Nucl. Acids Res.* 14, 8919–32.
- Conley, E.C. et al. (1986) Deletion and rearrangement of plasmid DNA during transformation of *Escherichia coli* with linear plasmid molecules. *Nucl. Acids Res.* 14, 8905–17.
- Chua, K.L. and Oliver, P. (1992) Intramolecular homologous recombination of linearized plasmids in *Escherichia coli* K12. *Mol. Gen. Genet.* 232, 199–205.
- Kobs, G. (1995) pGEM®-T Vector: Cloning of modified blunt-ended DNA fragments. Promega Notes 55, 28–9.
- Zhou, M.Y., Clark, S.E. and Gomez-Sanchez, C.E. (1995) Universal cloning method by TA strategy. *BioTechniques* 19, 34–5.
- Schmidt, E.V. et al. (1990) The cytomegalovirus enhancer: a pan-active control element in transgenic mice. Mol. Cell. Biol. 10, 4406–11.

VII.D. References (continued)

- Gluzman, Y. (1981) SV40-transformed simian cells support the replication of early SV40 mutants. *Cell* 23, 175–83.
- Bothwell, A.L.M. *et al.* (1981) Heavy chain variable region contribution to the NPb family of antibodies: somatic mutation evident in a gamma 2a variable region *Cell* 24, 625–37.
- Senapathy, P., Shapiro, M.B. and Harris, N.L. (1990) Splice junctions, branch point sites, and exons: sequence statistics, identification, and applications to genome project. *Methods Enzymol.* 183, 252–78.
- Huang, M.T.F. and Gorman, C.M. (1990) The simian virus 40 small-t intron, present in many common expression vectors, leads to aberrant splicing. *Mol. Cell. Biol.* 10, 1805–10.
- Gross, M.K., Kainz, M.S. and Merrill, G.F. (1987) Introns are inconsequential to efficient formation of cellular thymidine kinase mRNA in mouse L cells. *Mol. Cell. Biol.* 7, 4576–81.
- Buchman, A.R. and Berg, P. (1988) Comparison of intron-dependent and intronindependent gene expression. *Mol. Cell. Biol.* 8, 4395–405.
- Evans, M.J. and Scarpulla, R.C. (1989) Introns in the 3'-untranslated region can inhibit chimeric CAT and beta-galactosidase gene expression. *Gene* 84, 135–42.
- Huang, M.T.F. and Gorman, C.M. (1990) Intervening sequences increase efficiency of RNA 3' processing and accumulation of cytoplasmic RNA. Nucl. Acids Res. 18, 937–9.
- Brondyk, B. (1994) pCI and pSI Mammalian Expression Vectors. Promega Notes 49, 7–11.
- Brinster, R.L. et al. (1988) Introns increase transcriptional efficiency in transgenic mice. Proc. Natl. Acad. Sci. USA 85, 836–40.
- Choi, T. *et al.* (1991) A generic intron increases gene expression in transgenic mice. *Mol. Cell. Biol.* 11, 3070–4.
- Palmiter, R.D. et al. (1991) Heterologous introns can enhance expression of transgenes in mice. Proc. Natl. Acad. Sci. USA 88, 478–82.
- Fu, L. et al. (1991) Translational potentiation of messenger RNA with secondary structure in Xenopus. Science 251, 807–10.
- Kim, S.J. et al. (1992) Post-transcriptional regulation of the human transforming growth factor-beta 1 gene. J. Biol. Chem. 267, 13702–7.
- Kozak, M. (1986) Influences of mRNA secondary structure on initiation by eukaryotic ribosomes. Proc. Natl. Acad. Sci. USA 83, 2850–4.
- Kozak, M. (1989) Circumstances and mechanisms of inhibition of translation by secondary structure in eucaryotic mRNAs. *Mol. Cell. Biol.* 9, 5134–42.
- Rao, C.D. et al. (1988) The 5' untranslated sequence of the c-sis/platelet-derived growth factor 2 transcript is a potent translational inhibitor. Mol. Cell. Biol. 8, 284–92.
- 46. Proudfoot, N.J. (1991) Poly(A) signals. Cell 64, 671-4.

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison, WI
 53711-5399
 USA

 Toll Free in USA 800-356-9526
 Phone 608-274-4330
 Fax 608-277-2516
 •
 www.promega.com

 Part# TM044
 Printed in USA.
 Revised 2/06

Size

Cat #

- Bernstein, P. and Ross, J. (1989) Poly(A), poly(A) binding protein and the regulation of mRNA stability. *Trends Biochem. Sci.* 14, 373–7.
- Jackson, R.J. and Standart, N. (1990) Do the poly(A) tail and 3´ untranslated region control mRNA translation? *Cell* 62, 15–24.
- Carswell, S. and Alwine, J.C. (1989) Efficiency of utilization of the simian virus 40 late polyadenylation site: effects of upstream sequences. *Mol. Cell. Biol.* 9, 4248–58.
- Levitt, N. et al. (1989) Definition of an efficient synthetic poly(A) site. Genes and Dev. 3, 1019–25.
- Davies, J. and Jimenez, A. (1980) A new selective agent for eukaryotic cloning vectors. Am. J. Trop. Med. Hyg. 29, 1089–92.
- Lusky, M. and Botchan, M. (1981) Inhibition of SV40 replication in simian cells by specific pBR322 DNA sequences. *Nature* 293, 79–81.

VII.E. Related Products

Basic PCR Cloning Systems

Product

Todate	OILC	Cuttin
pGEM®-T Vector System I	20 reactions	A3600
pGEM®-T Vector System II	20 reactions	A3610
pGEM®-T Easy Vector System I	20 reactions	A1360
pGEM®-T Easy Vector System II	20 reactions	A1380

For Laboratory Use.

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

*For Laboratory Use.

Gel and PCR Purification Products

Product	Size	Cat.#
Wizard [®] SV Gel and PCR Clean-Up System	50 preps	A9281
	250 preps	A9282

For Laboratory Use.

Amplification Products

Product	Size	Cat.#
PCR Master Mix	10 reactions*	M7501
GoTaq® DNA Polymerase	100 units*	M3001

For Laboratory Use. *Available in additional sizes.



VII.E. Related Products (continued)

Reverse Transcription

Product	Size	Cat. #
Access RT-PCR System	100 reactions*	A1250
ImProm-II [™] Reverse Transcription System	100 reactions	A3800
AccessQuick™ RT PCR System	20 reactions*	A1701

For Laboratory Use. *Available in additional sizes.

Transfection Products

Product	Size	Cat.#
TransFast [™] Transfection Reagent	1.2mg	E2431
Tfx TM -20 Reagent	4.8mg	E2391
Tfx TM -50 Reagent	2.1mg	E1811
Tfx™ Reagents Transfection Trio	5.4mg	E2400
Transfectam [®] Reagent for the Transfection of		
Eukaryotic Cells	0.5mg	E1232
	1.0mg	E1231
ProFection [®] Mammalian Transfection System –		
Calcium Phosphate	1 system	E1200
ProFection [®] Mammalian Transfection System –		
DEAE-Dextran	1 system	E1210

Plasmid DNA Purification Systems

Product	Size	Cat.#
PureYield™ Plasmid Midiprep System	25 preps	A2492
	100 preps	A2495
Wizard [®] Plus SV Minipreps DNA Purification System	50 preps*	A1330
(ready for spin protocols using a microcentrifuge)		
Wizard [®] Plus SV Minipreps DNA Puification System	50 preps*	A1340
(with vacuum adapters: ready for spin or vacuum protocols**)		

For Laboratory Use. *Available in additional sizes.

**Vacuum protocol requires Vac-Man® Laboratory Vacuum Manifold.

Product	Capacity	Cat.#
Vac-Man [®] Laboratory Vacuum Manifold	20 samples	A7231

Sequencing Primers

Product	Size	Cat.#
T7 Promoter Primer	2μg	Q5021
pTargeT™ Sequencing Primer	2µg	Q4461

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison,
 WI
 53711-5399
 USA

 Toll Free in USA 800-356-9526
 Phone 608-274-4330
 Fax
 608-277-2516
 www.promega.com

 Part# TM044
 Printed in USA.
 Revised 2/06
 Revised 2/06

Reagents

Product	Size	Cat.#
X-Gal	100mg	V3941
IPTG, Dioxane Free	1g	V3955
		V3951
	50g	V3953
Antibiotic G-418 Sulfate	100mg	V7981
	1g	V7982
		V7983
Antibiotic G-418 Sulfate Solution	20ml	V8091

(a)Licensed under U.S. Pat. No. 5,075,430.

(b)Licensed under one or more of U.S. Pat. Nos. 5,487,993 and 5,827,657 and European Pat. No. 0 550 693.

^(c)The CMV promoter and its use are covered under U.S. Pat. Nos. 5,168,062 and 5,385,839 owned by the University of Iowa Research Foundation, Iowa City, Iowa, and licensed FOR RESEARCH USE ONLY. Research Use includes contract research for which monetary or other consideration may be received. Other commercial users must obtain a license to these patents directly from the University of Iowa Research Foundation.

(d)The PCR process is covered by patents issued and applicable in certain countries*. Promega does not encourage or support the unauthorized or unlicensed use of the PCR process.

*In Europe, effective March 28, 2006, European Pat. Nos. 201,184 and 200,362 will expire. In the U.S., the patents covering the foundational PCR process expired on March 29, 2005.

© 1996-2006 Promega Corporation. All Rights Reserved.

GoTaq, pGEM, ProFection, Transfectam, Vac-Man and Wizard are registered trademarks of Promega Corporation. AccessQuick, ImProm-II, pTARGET, PureYield, Select%, Tfx and TransFast are trademarks of Promega Corporation.

AmpliTaq is a registered trademark of Roche Molecular Systems, Inc. Bacto is a registered trademark of Difco Laboratories, Detroit, MI. Deep Vent_R and Vent_R are registered trademarks of New England Biolabs, Inc. DH5 α is a trademark of Life Technologies. DNASTAR is a registered trademark of DNASTAR, Inc. GenBank is a registered trademark of the U.S. Department of Health and Human Services.

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.

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison, WI
 53711-5399
 USA

 Toll Free in USA
 800-356-9526
 Phone 608-274-4330
 Fax
 608-277-2516
 www.promega.com

 Printed in USA.
 Part# TM044

 Revised 2/06
 Page 25