

# **TOPO<sup>®</sup> Reporter Kits**

#### Five-minute Cloning of PCR Products for Analysis of Promoter Function in Mammalian Cells

Catalog nos. K4830-01, K4831-01

**Version K** 23 February 2006 25-0235

A Limited Label License covers this product (see Purchaser Notification). By use of this product, you accept the terms and conditions of the Limited Label License.

**User Manual** 

# **Table of Contents**

	Table of Contents	iii
	Kit Contents and Storage	iv
	Accessory Products	vi
Method	ls	1
	Overview	1
	General Cloning Considerations	
	Cloning into pBlue-TOPO <sup>®</sup>	4
	Cloning into pGlow-TOPO <sup>®</sup>	5
	Producing PCR Products	6
	TOPO <sup>®</sup> Cloning Reaction and Transformation	7
	Optimizing the TOPO <sup>®</sup> Cloning Reaction	11
	Transfection	
	Detection of β-Galactosidase Activity	
	Detection of GFP	14
	Creation of Stable Cell Lines	16
Append	dix	
••	TOPO <sup>®</sup> Reporter Control Reactions	
	Purifying PCR Products	
	Addition of 3' A-Overhangs Post-Amplification	
	pBlue-TOPO <sup>®</sup> Map	23
	pGlow-TOPO <sup>®</sup> Map	
	Features of the TOPO <sup>®</sup> Reporter Vectors	25
	Recipes	
	Technical Service	27
	Purchaser Notification	
	Product Qualification	
	References	

#### **Kit Contents and Storage**

	Each TOPO <sup>®</sup> Reporter Kit is shipped on dry ice. Each kit contains a box with TOPO <sup>®</sup> Reporter reagents (Box 1) and a box with One Shot <sup>®</sup> TOP10 chemically competent cells (Box 2). Store Box 1 at -20°C and <b>Box 2 at -80</b> °C.			
TOPO <sup>®</sup> Reporter Kits	Ordering information for the TOPO <sup>®</sup> Reporter Kits is provided below.			
	Kit	Vector	Reactions	Catalog no.
	pBlue TOPO <sup>®</sup> TA Expression Kit	pBlue-TOPO <sup>®</sup>	20	K4831-01
	pGlow TOPO <sup>®</sup> TA Expression Kit	pGlow-TOPO®	20	K4830-01

#### TOPO<sup>®</sup> Reporter Reagents

TOPO<sup>®</sup> Reporter reagents (Box 1) are listed below. **Please note that the user must supply** *Taq* **polymerase.** Store Box 1 at -20°C.

Item	Concentration	Amount
pBlue-TOPO <sup>®</sup>	10 ng/µl plasmid DNA in:	20 µl
or	50% glycerol	
pGlow-TOPO <sup>®</sup>	50 mM Tris-HCl, pH 7.4 (at 25°C)	
	1 mM EDTA	
	2 mM DTT	
	0.1% Triton X-100	
	100 µg/ml BSA	
	phenol red	
10X PCR Buffer	100 mM Tris-HCl, pH 8.3 (at 42°C)	100 µl
	500 mM KCl	
	25 mM MgCl <sub>2</sub>	
	0.01% gelatin	
dNTP Mix	12.5 mM dATP	10 µl
	12.5 mM dCTP	
	12.5 mM dGTP	
	12.5 mM dTTP	
	neutralized at pH 8.0 in water	
Salt Solution	1.2 M NaCl; 0.06 M MgCl <sub>2</sub>	50 µl
T7 Sequencing Primer	0.1 μg/μl in TE Buffer	20 µl
LacZ Reverse or GFP Reverse	0.1 μg/μl in TE Buffer	20 µl
Sequencing Primer		
Control PCR Template	0.05 µg/µl in TE Buffer	10 µl
Control PCR Primers	0.1 μg/μl <b>each</b> in TE Buffer	10 µl
Sterile Water		1 ml

#### Kit Contents and Storage, continued

#### One Shot<sup>®</sup> Reagents

The table below describes the items included in the One Shot® TOP10 chemically competent cell kit. Store at -80°C.

Item	Composition	Amount
SOC Medium	2% Tryptone	6 ml
(may be stored at +4°C	0.5% Yeast Extract	
or room temperature)	10 mM NaCl	
	2.5 mM KCl	
	10 mM MgCl <sub>2</sub>	
	10 mM MgSO <sub>4</sub>	
	20 mM glucose	
TOP10 cells		21 x 50 µl
pUC19 Control DNA	10 pg/µl in 5 mM Tris-HCl, 0.5 mM EDTA, pH 8	50 µl

#### Sequencing Primers

The table below provides the sequence and pmoles of the T7, *LacZ* Reverse, and GFP Reverse sequencing primers.

Primer	Sequence	Amount
Τ7	5'-TAATACGACTCACTATAGGG-3'	328 pmoles
LacZ Reverse	5'-CAGTCATGCTAGCCATACC-3'	350 pmoles
GFP Reverse	5'-GGGTAAGCTTTCCGTATGTAGC-3'	296 pmoles

# Genotype of TOP10 Cells

**TOP10**: Use this strain for general cloning. Please note that this strain cannot be used for single-strand rescue of DNA.

F- mcrA Δ(mrr-hsdRMS-mcrBC) Φ80lacZΔM15 ΔlacX74 recA1 araD139 Δ(ara-leu)7697 galU galK rpsL (Str<sup>R</sup>) endA1 nupG

## **Accessory Products**

#### Additional Products

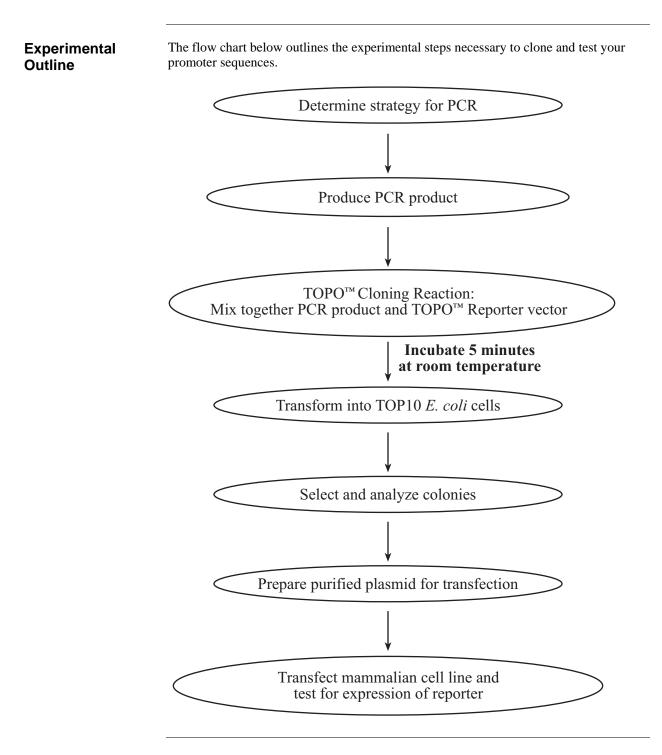
The table below lists additional products available from Invitrogen which you may use in conjunction with the  $TOPO^{\otimes}$  Reporter Kit.

Item	Amount	Catalog no.
One Shot <sup>®</sup> Kit	10 reactions	C4040-50
(TOP10 Electrocompetent Cells)	20 reactions	C4040-52
One Shot <sup>®</sup> Kit	10 reactions	C4040-10
(TOP10 Chemically Competent Cells)	20 reactions	C4040-03
	40 reactions	C4040-06
T7 Promoter Primer	2 µg (328 pmoles)	N560-02
S.N.A.P. <sup>™</sup> MiniPrep Kit	20 reactions	K1900-01
S.N.A.P. <sup>™</sup> MidiPrep Kit	20 reactions	K1910-01
Anti-Xpress <sup>™</sup> Antibody	25 Westerns	R910-25
β-galactosidase Antiserum	25 Westerns	R901-25
GFP Antiserum	25 Westerns	R970-01
β-Gal Assay Kit	100-400 reactions	K1455-01
β-Gal Staining Kit	1 kit	K1465-01

# Methods

Overview	
Introduction	<ul> <li>TOPO<sup>®</sup> Reporter Kits provide a highly efficient, 5 minute, one-step cloning strategy ("TOPO<sup>®</sup> Cloning") for the direct insertion of promoter sequences amplified by <i>Taq</i> polymerase into a reporter vector. Recombinant vectors can then be transfected into mammalian cells or transformed into <i>E. coli</i> (pGlow-TOPO only) and tested for promoter function <i>in vivo</i> or <i>in vitro</i>. No ligase, post-PCR procedures, or PCR primers containing specific sequences are required.</li> <li>Two types of TOPO<sup>®</sup> TA Expression Kit containing pBlue-TOPO<sup>®</sup> for insertion of promoter sequences upstream of the <i>E. coli</i> β-galactosidase gene (<i>lacZ</i>) for <i>in vitro</i> or <i>in vivo</i> assay.</li> <li>pGlow TOPO<sup>®</sup> TA Expression Kit containing pGlow-TOPO<sup>®</sup> for insertion of promoter sequences upstream of GFP for <i>in vitro</i> or <i>in vivo</i> assay of promoter function.</li> </ul>
How It Works	<text><list-item><text><text><text><text></text></text></text></text></list-item></text>
	Tyr-274

#### **Overview**, continued



# **General Cloning Considerations**

Introduction	<ul><li>In general, promoter reporter vectors can be used to analyze</li><li>Tissue and cell-specific promoter function</li></ul>
	Transcriptional enhancers in a known promoter
	• Deletions within a promoter
	Please note that each TOPO <sup>®</sup> Reporter vector contains not only a TOPO <sup>®</sup> Cloning site but also additional unique restriction sites upstream of the TOPO <sup>®</sup> Cloning site. These may be exploited to analyze promoter function.
Note	When analyzing promoters in a reporter vector, it is important to realize that sequences within the native gene can influence regulation of its own promoter. In addition, sequences within the reporter gene can also affect transcription from the promoter under study. We recommend that you verify any observations of transcriptional control of the fusion gene with expression of the native gene. S1 mapping ( <i>Current Protocols in Molecular Biology</i> , pages 4.6.1 to 4.6.13) can be used to confirm that the subcloned promoter initiates transcription at the correct site.
<b>Q</b> Important	Since initiation of translation in eukaryotes occurs at the first available AUG codon, it is important that there are no AUG codons between the start of transcription and the AUG of the reporter gene.
PCR Primer Design	Use the diagrams on pages 4 and 5 and the sequence of your promoter to design PCR primers. Unique restriction sites may be included in the 5' and 3' primers to excise the fragment or facilitate analysis once it is TOPO <sup>®</sup> Cloned.
	For analysis of promoter function in <i>E. coli</i> , please use pGlow-TOPO and read page 5 before designing your primers.
Note	Once you have decided on the sequence of your PCR primers, do not add 5 <sup>'</sup> phosphates to your primers. Phosphates will inhibit topoisomerase I and the synthesized PCR product will not ligate into the TOPO <sup>®</sup> Reporter vectors. Please note that cloning efficiencies may vary depending on the 5 <sup>'</sup> nucleotide in the primers (see page 19).

# Cloning into pBlue-TOPO<sup>®</sup>

Important		There is a cryptic prokaryotic promoter upstream of the <i>lacZ</i> reporter gene. <i>E. coli</i> transformants may appear to be light blue when screened on plates containing X-Gal. We do not recommend using pBlue-TOPO <sup>®</sup> to evaluate promoter function in <i>E. coli</i> . Please use pGlow-TOPO <sup>®</sup> for these studies (next page). Please note that background expression of $\beta$ -galactosidase does not occur in mammalian cells.	
TOPO <sup>®</sup> Cloning Site of pBlue- TOPO <sup>®</sup>		Restriction sites are labeled to indicate the actual cleavage site. The vector is supplied linearized between base pair 116 and 117. This is the TOPO <sup>®</sup> Cloning site. <b>Please note that the full sequence of pBlue-TOPO<sup>®</sup> may be downloaded from our Web site (www.invitrogen.com) or requested from Technical Service (see page 27).</b> A map of pBlue-TOPO <sup>®</sup> is provided on page 23.	_
	Bgl II*	T7 promoter/priming site Afl II Hind III* BamH I	
1	GACGGATCGG GAGAT	T7 promoter/priming site Afl II Hind III* BamH I	
81	Spe I TCGGATCCAC TAGTO	Pst I* Hind III* Nco I* CAGTG TGGTGGAATT GCCCTT PCR AAGG GCAATTCTGC AGAAAGCTTA CC ATG GGG CGGGAA Product TTCC Met Gly	n
	Po	yhistidine region LacZ Reverse priming site	
149		CAT CAT CAT GGT ATG GCT AGC ATG ACT GGT GGA CAG CAA ATG GGT CGG GAT His His His Gly Met Ala Ser Met Thr Gly Gly Gln Gln Met Gly Arg Asp	
	Xpress™ ep	itope	
215		GAC GAT AAG GTA CCT AAG GAT CAG CTT GGA GTT GAT CCC GTT TTA CAA CGT Asp Asp Lys Val Pro Lys Asp Gln Leu Gly Val Asp Pro Val Val Leu Gln Arg	
		ecognition site	
281	CGT GAC TGG GAA		
*TL	Arg Asp Trp Glu	ue, but they may be used to excise the promoter sequence after TOPO <sup>®</sup> Cloning.	
. 111	ese snes are not unit		_

# Cloning into pGlow-TOPO<sup>®</sup>

Using pGlow- TOPO in <i>E. coli</i>		To use pGlow-TOPO as a reporter in <i>E. coli</i> , you must engineer your PCR product to ensure expression of GFP in the event that the sequences you are testing contain a promoter. Please note that there is no prokaryotic ribosomal binding site upstream of the GFP initiation codon. Your reverse PCR primer must include a ribosomal binding site (-AGGA-) and an initiation codon (ATG) in frame with the GFP initiation codon. Allow 8-12 nucleotides between the ribosomal binding site and the initiation codon to ensure proper spacing. There should not be any palindromic sequences within this region. Successful expression of GFP will result in additional amino acids at the N-terminus. This has been shown not to affect fluorescence. We recommend that you use a known promoter as a positive control and DNA sequences that do not contain a promoter as a negative control.	
Site	PO <sup>®</sup> Cloning of pGlow- PO <sup>®</sup>	linearized between base pair 116 an that the full sequence of pGlow-T (www.invitrogen.com) or requested	eate the actual cleavage site. The vector is supplied ad 117. This is the TOPO <sup>®</sup> Cloning site. <b>Please note</b> <b>OPO<sup>®</sup> may be downloaded from our Web site</b> <b>ed from Technical Service (see page</b> Error! pGlow-TOPO <sup>®</sup> is provided on page 24.
1	Bgl II*	T7 promoter/priming site	Pme I* Afl II Asp718 I Kpn I
1			
81	Spe I TCGGATCCAC TAGTC	BstX I* CAGTG TGGTGGAATT GCCCTT PCR CGGGAA Produc	Bgl II* Pst I* Xba I* GFP ORF AAGG GCAATTCTGC AGATCTAGA ATG GCT AGC TTCC Met Ala Ser
149			TT CTT GTT GAA TTA GAT GGT GAT GTT AAT GGG le Leu Val Glu Leu Asp Gly Asp Val Asn Gly GFP Reverse priming site
215			AT GCT ACA TAC GGA AAG CTT ACC CTT AAA TTT sp Ala Thr Tyr Gly Lys Leu Thr Leu Lys Phe
	ATT TGC ACT ACT Ile Cys Thr Thr	Gly Lys	
*The	ese sites are not uniq	ie, but they may be used to excise the	e promoter sequence after TOPO <sup>®</sup> Cloning.

## **Producing PCR Products**

Introduction	Once you have decided on a PCR strat ready to produce your PCR product.		ategy and have synthesized the primers you are	
Materials Supplied by the User	Yo • •	u will need the following reagents <i>Taq</i> polymerase Thermocycler DNA template and primers for P		
Polymerase Mixtures	If you wish to use a mixture containing $Taq$ polymerase and a proofreading polymerase, $Taq$ must be used in excess of a 10:1 ratio to ensure the presence of 3' A-overhangs on the PCR product.			
	-	1 1	do not have enough <i>Taq</i> polymerase or a proof- d 3' A-overhangs using the method on page 22.	
Producing PCR Products	1.	DNA as a template and more DN Use the cycling parameters suita	reaction. Use less DNA if you are using plasmid NA if you are using genomic DNA as a template. ble for your primers and template. Be sure to on at 72°C after the last cycle to ensure that all PCR denylated.	
		DNA Template	10-100 ng	
		10X PCR Buffer	5 µl	
		50 mM dNTPs	0.5 µl	
		Primers	100-200 ng each	
		Sterile water	add to a final volume of 49 $\mu$ l	
		Taq Polymerase (1 unit/µl)	<u>1 µl</u>	
		Total Volume	50 µl	
	2.		ose gel electrophoresis. You should see a single, a single band, please refer to the <b>Note</b> below.	



If you do not obtain a single, discrete band from your PCR, you may gel-purify your fragment before using the TOPO<sup>®</sup> Reporter Kits (see page 20). Take special care to avoid sources of nuclease contamination and long exposure to UV light. Alternatively, you may optimize your PCR to eliminate multiple bands and smearing (Innis *et al.*, 1990). The PCR Optimizer<sup>TM</sup> Kit (Catalog no. K1220-01) from Invitrogen can help you optimize your PCR. Please call Technical Service for more information (page **Error! Bookmark not defined.**).

# **TOPO<sup>®</sup> Cloning Reaction and Transformation**

Introduction	TOPO <sup>®</sup> Cloning technology allows you to produce your PCR products, ligate them into either pBlue-TOPO <sup>®</sup> or pGlow-TOPO <sup>®</sup> , and transform the recombinant vector into <i>E. coli</i> all in one day. It is important to have everything you need set up and ready to use to ensure you obtain the best possible results. If this is the first time you have TOPO <sup>®</sup> Cloned, you may wish to perform the control reactions on pages 17-19 in parallel with your samples.
Note	Recent experiments at Invitrogen demonstrate that inclusion of salt (200 mM NaCl, 10 mM MgCl <sub>2</sub> ) in the TOPO <sup>®</sup> Cloning reaction increases the number of transformants 2-to 3-fold. We have also observed that in the presence of salt, incubation times of greater than 5 minutes can also increase the number of transformants. This is in contrast to earlier experiments <b>without salt</b> where the number of transformants decreases as the incubation time increases beyond 5 minutes.
	Inclusion of salt allows for longer incubation times because it prevents topoisomerase I from rebinding and potentially nicking the DNA after ligating the PCR product and dissociating from the DNA. The result is more intact molecules leading to higher transformation efficiencies.
Important	Because of the above results, we recommend adding salt to the TOPO <sup>®</sup> Cloning reaction. A stock salt solution is provided in the kit for this purpose. Please note that the amount of salt added to the TOPO <sup>®</sup> Cloning reaction varies depending on whether you plan to transform chemically competent cells (provided) or electrocompetent cells (see below). For this reason two different TOPO <sup>®</sup> Cloning reactions are provided to help you obtain the best possible results. Please read the following information carefully.
Chemically Competent <i>E. coli</i>	For TOPO <sup>®</sup> Cloning and transformation into chemically competent <i>E. coli</i> , adding sodium chloride and magnesium chloride to a final concentration of 200 mM NaCl, 10 mM MgCl <sub>2</sub> in the TOPO <sup>®</sup> Cloning reaction increases the number of colonies over time. A Salt Solution (1.2 M NaCl; 0.06 M MgCl <sub>2</sub> ) is provided to adjust the TOPO <sup>®</sup> Cloning reaction to the recommended concentration of NaCl and MgCl <sub>2</sub> .
Electrocompetent <i>E. coli</i>	For TOPO <sup>®</sup> Cloning and transformation of electrocompetent <i>E. coli</i> , salt must also be included in the TOPO <sup>®</sup> Cloning reaction, but the amount of salt <b>must be reduced</b> to 50 mM NaCl, 2.5 mM MgCl <sub>2</sub> to prevent arcing. The Salt Solution is diluted 4-fold to prepare a 300 mM NaCl, 15 mM MgCl <sub>2</sub> solution for convenient addition to the TOPO <sup>®</sup> Cloning reaction (see next page).
Materials Supplied by the User	<ul> <li>In addition to general microbiological supplies (i.e. plates, spreaders), you will need the following reagents and equipment.</li> <li>42°C water bath (or electroporator with cuvettes, optional)</li> <li>LB plates containing 50-100 μg/ml ampicillin (two for each transformation)</li> <li>Reagents and equipment for agarose gel electrophoresis</li> <li>37°C shaking and non-shaking incubator</li> </ul>

# **TOPO<sup>®</sup> Cloning Reaction and Transformation, continued**



There is no blue-white screening for the presence of inserts. Individual recombinant plasmids need to be analyzed by restriction analysis or sequencing for the presence and orientation of insert. Sequencing primers included in the kit can be used to sequence across an insert in the multiple cloning site to confirm orientation and reading frame.

#### Preparation for Transformation

For each transformation, you will need one vial of competent cells and two selective plates.

- Equilibrate a water bath to 42°C (for chemical transformation) or set up your electroporator if you are using electrocompetent *E. coli*.
- For electroporation, dilute a small portion of the Salt Solution 4-fold to prepare Dilute Salt Solution (e.g. add 5 µl of the Salt Solution to 15 µl sterile water)
- Warm the vial of SOC medium from Box 2 to room temperature.
- Warm selective plates at 37°C for 30 minutes.
- Thaw <u>on ice</u> 1 vial of One Shot<sup>®</sup> cells for each transformation.

# Setting Up the TOPO<sup>®</sup> Cloning Reaction

The table below describes how to set up your TOPO<sup>®</sup> Cloning reaction (6  $\mu$ l) for eventual transformation into either chemically competent TOP10 One Shot<sup>®</sup> *E. coli* (provided) or electrocompetent *E. coli*. Additional information on optimizing the TOPO<sup>®</sup> Cloning reaction for your needs can be found on page 11.

**Note**: The red or yellow color of the TOPO<sup>®</sup> vector solution is normal and is used to visualize the solution.

Reagent*	Chemically Competent E. coli	Electrocompetent E. coli
Fresh PCR product	0.5 to 4 µl	0.5 to 4 µl
Salt Solution	1 µl	
Dilute Salt Solution		1 μl
Sterile Water	add to a final volume of 5 $\mu$ l	Add to a final volume of 5 $\mu$ l
TOPO <sup>®</sup> vector	1 μl	1 µl

\*Store all reagents at -20°C when finished. Salt solutions and water can be stored at room temperature or +4°C.

# Performing the TOPO<sup>®</sup> Cloning Reaction

- 1. Mix reaction gently and incubate for 5 minutes at room temperature (22-23°C).
- **Note**: For most applications, 5 minutes will yield plenty of colonies for analysis. Depending on your needs, the length of the TOPO<sup>®</sup> Cloning reaction can be varied from 30 seconds to 30 minutes. For routine subcloning of PCR products, 30 seconds may be sufficient. For large PCR products (> 1 kb) or if you are TOPO<sup>®</sup> Cloning a pool of PCR products, increasing the reaction time will yield more colonies.
  - 2. Place the reaction on ice and proceed to **One Shot**<sup>®</sup> **Chemical Transformation** (next page) or **Transformation by Electroporation** (next page). **Note**: You may store the TOPO<sup>®</sup> Cloning reaction at -20°C overnight.

# **TOPO<sup>®</sup> Cloning Reaction and Transformation, continued**

One Shot <sup>®</sup> Chemical Transformation	1.	Add 2 $\mu$ l of the TOPO <sup>®</sup> Cloning reaction from Step 2 previous page into a vial of One Shot <sup>®</sup> TOP10 Chemically Competent <i>E. coli</i> and mix gently. <b>Do not mix by pipetting up and down.</b>
	2.	Incubate on ice for 5 to 30 minutes.
		<b>Note</b> : Longer incubations on ice seem to have a minimal effect on transformation efficiency. The length of the incubation is at the user's discretion (see page 11).
	3.	Heat-shock the cells for 30 seconds at 42°C without shaking.
	4.	Immediately transfer the tubes to ice.
	5.	Add 250 µl of room temperature SOC medium.
	6.	Cap the tube tightly and shake the tube horizontally (200 rpm) at 37°C for 1 hour.
	7.	Spread 25-200 $\mu$ l from each transformation on a pre-warmed selective plate and incubate overnight at 37°C. We recommend that you plate two different volumes to ensure that at least one plate will have well-spaced colonies.
	8.	An efficient TOPO <sup>®</sup> Cloning reaction will produce hundreds of colonies. Pick ~10 colonies for analysis (see <b>Analysis of Positive Clones</b> , next page).
Transformation by Electroporation	1.	Add 2 $\mu$ l of the TOPO <sup>®</sup> Cloning reaction into a 0.1 cm cuvette containing 50 $\mu$ l of electrocompetent <i>E. coli</i> and mix gently. <b>Do not mix by pipetting up and down. Avoid formation of bubbles.</b>
	2.	Electroporate your samples using your own protocol and your electroporator.
		Note: If you have problems with arcing, see below.
	3.	Immediately add 250 µl of room temperature SOC medium.
	4.	Transfer the solution to a 15 ml snap-cap tube (i.e. Falcon) and shake for at least 1 hour at 37°C to allow expression of the antibiotic resistance gene.
	5.	Spread 10-50 $\mu$ l from each transformation on a pre-warmed selective plate and incubate overnight at 37°C. To ensure even spreading of small volumes, add 20 $\mu$ l of SOC. We recommend that you plate two different volumes to ensure that at least one plate will have well-spaced colonies.
	6.	An efficient TOPO <sup>®</sup> Cloning reaction will produce hundreds of colonies. Pick ~10 colonies for analysis (see <b>Analysis of Positive Clones</b> , next page).
Note	cor res	Idition of the Dilute Salt Solution in the TOPO <sup>®</sup> Cloning Reaction brings the final ncentration of NaCl and MgCl <sub>2</sub> in the TOPO <sup>®</sup> Cloning reaction to 50 mM and 2.5 mM, spectively. To prevent arcing of your samples during electroporation, the volume of lls should be between 50 and 80 $\mu$ l (0.1 cm cuvettes) or 100 to 200 $\mu$ l (0.2 cm cuvettes)
	If y	you experience arcing during transformation, try one of the following suggestions:
	٠	Reduce the voltage normally used to charge your electroporator by 10%
	٠	Reduce the pulse length by reducing the load resistance to 100 ohms
		Ether all any similar the TODO <sup>®</sup> Classical production and assume and in suctain prior to

• Ethanol-precipitate the TOPO<sup>®</sup> Cloning reaction and resuspend in water prior to electroporation

# **TOPO<sup>®</sup> Cloning Reaction and Transformation, continued**

Analysis of Positive Clones	<ol> <li>Pick 10 colonies and culture them overnight in LB medium containing 50-100 µg/ml ampicillin (3-5 ml). Note: In cells transformed with pGlow-TOPO, a hand-held UV light can be used to detect fluorescence.</li> <li>Isolate plasmid DNA using your method of choice. If you need ultra-pure plasmid DNA for automated or manual sequencing, we recommend the S.N.A.P.<sup>™</sup> MiniPrep Kit (Catalog no. K1900-01).</li> <li>Please note that PCR products will clone bidirectionally. Analyze the plasmids for insertion and orientation by restriction analysis or by sequencing. Use the sequencing primers included in the kit to help you sequence your insert. Please refer to the diagrams on page 4 and page 5 for restriction sites and sequence surrounding the TOPO Cloning<sup>®</sup> site. For the complete sequence of either vector, please see our Web site (www.invitrogen.com) or contact Technical Service (page 27). If you need help with setting up restriction enzyme digests or DNA sequencing, refer</li> </ol>
	to general molecular biology texts (Ausubel et al., 1994; Sambrook et al., 1989).
Alternative Method of Analysis	You may wish to use PCR to directly analyze positive transformants. Use a combination of either the T7 or <i>LacZ</i> Reverse (or GFP Reverse) sequencing primer and a primer that binds within your insert as PCR primers. You will have to determine the amplification conditions If this is the first time you have used this technique, we recommend that you perform restriction analysis in parallel to confirm that PCR gives you the correct result. Artifacts may be obtained because of mispriming or contaminating template.
	<ul> <li>The following protocol is provided for your convenience. Other protocols are suitable.</li> <li>Prepare a PCR cocktail consisting of PCR buffer, dNTPs, primers, and <i>Taq</i> polymerase. Use a 20 µl reaction volume. Multiply by the number of colonies to be analyzed (e.g. 10).</li> </ul>
	2. Pick 10 colonies and resuspend them individually in 20 µl of the PCR cocktail. (Don't
	<ol> <li>forget to make a patch plate to preserve the colonies for further analysis.)</li> <li>Incubate the reaction for 10 minutes at 94°C to lyse the cells and inactivate nucleases.</li> </ol>
	<ol> <li>Amplify for 20 to 30 cycles using parameters previously determined (see text, above)</li> <li>For the final extension, incubate at 72°C for 10 minutes. Store at +4°C.</li> </ol>
	6. Visualize by agarose gel electrophoresis.
<b>O</b> Important	If you have problems obtaining transformants or the correct insert, perform the control reactions described on pages 17-19. These reactions will help you troubleshoot your experiment.
Long-Term Storage	Once you have identified the correct clone, be sure to isolate a single colony and prepare a glycerol stock for long term storage. We recommend that you store a stock of plasmid DNA at -20°C.
	1. Streak the original colony on LB plates containing 50-100 $\mu$ g/ml ampicillin.
	<ol> <li>Isolate a single colony and inoculate into 1-2 ml of LB containing 50-100 μg/ml ampicillin. Grow until culture reaches stationary phase.</li> </ol>
	<ol> <li>Mix 0.85 ml of culture with 0.15 ml of sterile glycerol and transfer to a cryovial.</li> <li>Store at -80°C.</li> </ol>

# **Optimizing the TOPO<sup>®</sup> Cloning Reaction**

Introduction	The information below will help you optimize the TOPO <sup>®</sup> Cloning reaction for your particular needs.
Faster Subcloning	The high efficiency of TOPO <sup>®</sup> Cloning technology allows you to streamline the cloning process. If you routinely clone PCR products and wish to speed up the process, consider the following:
	• Incubate the TOPO <sup>®</sup> Cloning reaction for only 30 seconds instead of 5 minutes.
	<ul> <li>You may not obtain the highest number of colonies, but with the high efficiency of TOPO<sup>®</sup> Cloning, most of the transformants will contain your insert.</li> <li>After adding 2 µl of the TOPO<sup>®</sup> Cloning reaction to chemically competent cells, incubate on ice for only 5 minutes.</li> </ul>
	Increasing the incubation time to 30 minutes does not significantly improve transformation efficiency.
More Transformants	If you are TOPO <sup>®</sup> Cloning large PCR products, toxic genes, or cloning a pool of PCR products, you may need more transformants to obtain the clones you want. To increase the number of colonies:
	• Incubate the salt-supplemented TOPO <sup>®</sup> Cloning reaction for 20 to 30 minutes instead of 5 minutes.
	Increasing the incubation time of the salt-supplemented TOPO <sup>®</sup> Cloning reaction allows more molecules to ligate, increasing the transformation efficiency. Addition of salt appears to prevent topoisomerase from rebinding and nicking the DNA after it has ligated the PCR product and dissociated from the DNA.
Cloning Dilute	To clone dilute PCR products, you may:
PCR Products	Increase the amount of the PCR product
	• Incubate the TOPO <sup>®</sup> Cloning reaction for 20 to 30 minutes
	• Concentrate the PCR product

# Transfection

Introduction	Once you obtain the desired construct, you are ready to transfect the plasmid into the mammalian cells of choice. Please note the following guidelines for transfection.
	We recommend that you include a positive and a negative control to evaluate expression of the reporter genes. A negative control can be either a mock transfection, or TOPO <sup>®</sup> Clone a PCR product that does not contain a promoter (stuffer DNA) into the desired TOPO <sup>®</sup> Reporter vector. For a positive control, we recommend cloning a known promoter that is active in your cell line.
Plasmid Preparation	Plasmid DNA for transfection into eukaryotic cells must be very clean and free from phenol and sodium chloride. Contaminants will kill the cells and salt will interfere with lipid complexing decreasing transfection efficiency. We recommend isolating plasmid DNA (up to 200 µg) using the S.N.A.P. <sup>™</sup> MidiPrep Kit (Catalog no. K1910-01) or CsCl gradient centrifugation.
Methods of Transfection	For established cell lines (e.g. HeLa), please consult original references or the supplier of your cell line for the optimal method of transfection. It is recommended that you follow exactly the protocol for your cell line. Pay particular attention to medium requirements, when to pass the cells, and at what dilution to split the cells. Further information is provided in <i>Current Protocols in Molecular Biology</i> ( <b>Reference</b> section, page 31).
	Methods for transfection include calcium phosphate (Chen and Okayama, 1987; Wigler <i>et al.</i> , 1977), lipid-mediated (Felgner <i>et al.</i> , 1989; Felgner and Ringold, 1989) and electroporation (Chu <i>et al.</i> , 1987; Shigekawa and Dower, 1988). Invitrogen offers the Calcium Phosphate Transfection Kit (Catalog no. K2780-01) and a large selection of reagents for transfection. For more information on the reagents available, please visit our Web site (www.invitrogen.com) or call Technical Service (see page 27).
Detection of Reporter	Once you have transfected your cell line with one of the TOPO <sup>®</sup> Reporter vectors and the appropriate controls, you are ready to assay for reporter function. See the next page for information on how to assay for $\beta$ -galactosidase activity and page 14 for GFP assays.

# Detection of $\beta$ -Galactosidase Activity

Introduction	$\beta$ -galactosidase is one of the most versatile reporters available. It can be assayed both <i>in vitro</i> and <i>in vivo</i> and a wide variety of substrates are available for detection. A few assays and substrates are described below. Other assays and substrates may be used. In addition to its use as a reporter for uncharacterized promoters, constitutive promoters may be cloned upstream of the <i>lacZ</i> gene for use as an internal control to normalize variability with other promoter reporter assays (Alam and Cook, 1990).
Choosing an <i>In</i> <i>Vitro</i> Assay	$\beta$ -galactosidase activity can be detected using cell-free lysates and <i>o</i> -nitrophenyl- $\beta$ -D-galactopyranoside (ONPG). This colorimetric assay is easy to perform and useful for determining whether or not a promoter is active, but it lacks the sensitivity needed for promoter analysis (detects about 100 pg of $\beta$ -galactosidase). Invitrogen offers a $\beta$ -Gal Assay Kit (Catalog no. K1455-01) which contains ONPG and all the buffers necessary to assay for $\beta$ -galactosidase activity. Instructions are also included for a 96-well format (see page vi for ordering information).
	If you need greater sensitivity for promoter analysis, we recommend using chemi- luminescent 1,2-dioxetane substrates (i.e. Galacton, Tropix) (Beale <i>et al.</i> , 1992; Jain and Magrath, 1991). Use of these substrates increases the sensitivity of the assay and extends the range of detection (Bronstein <i>et al.</i> , 1994). If endogenous enzyme activity is minimized, sensitivity is enhanced (Young <i>et al.</i> , 1993). As little as 2 fg of $\beta$ -galacto- sidase can be detected using chemiluminescent substrates. For more information on this assay, please see the references cited above and <i>Current Protocols in Molecular Biology</i> , pages 9.7.15 to 9.7.21.
Choosing an <i>In</i> <i>Vivo</i> Assay	<i>In vivo</i> detection systems are defined as those in which the reporter gene is detected in live cells or tissues or in cells or tissues fixed for histochemical staining. This is a less quantitative approach but provides important information about cell-type specificity, temporal and tissue expression patterns, and distribution of transcription factors.
	The precipitating substrate X-Gal may be used to determine <i>in vivo</i> levels of $\beta$ -galactosidase in eukaryotic cells, tissue sections, and intact embryos (Alam and Cook, 1990). Please note that staining with X-Gal requires that the cells or tissue be fixed. Invitrogen offers the $\beta$ -Gal Staining Kit (Catalog no. K1465-01) to stain cells expressing $\beta$ -galactosidase (see page vi for ordering information).
	Alternatively, detection in live cultured cells may be achieved with the substrate fluorescein di- $\beta$ -D-galactopyranoside (FDG) (Jongkind <i>et al.</i> , 1986). Using hypotonic loading, FDG is introduced into the cell and cleaved by $\beta$ -galactosidase. The resulting fluorescein compound is trapped in the cell because of its hydrophobic nature and easily assayed using fluorescence.
Detection of β-galactosidase	If you do not detect activity of $\beta$ -galactosidase, check for expression by Western blot. You may use antibody to $\beta$ -galactosidase (see page vi for ordering information), or, since $\beta$ -galactosidase is expressed as a fusion to an N-terminal peptide containing the Xpress <sup>TM</sup> epitope, use the Anti-Xpress <sup>TM</sup> Antibody (see page vi for ordering information).

## **Detection of GFP**

Introduction	Green fluorescent protein (GFP) is very usef function. <i>In vivo</i> assays, while less quantitati regarding cell-type specificity of promoters/ transcription factors. Use of pGlow-TOPO <sup>®</sup> in real time. Please note that detection of GF promoter. For low-level expression it may be a fluorimeter. For detection in <i>E. coli</i> , assay	ive than enhanc allows P <i>in vi</i> e neces	n <i>in vi</i> ers an you t yo wil sary t	<i>tro</i> ass d the t o mon l depe	says, provide information tissue distribution of spec itor transcriptional chang end on the strength of the	cific ges
GFP Gene Used in pGlow-TOPO <sup>®</sup>	The GFP gene used in pGlow-TOPO <sup>®</sup> is des usage was optimized for expression in <i>E. co.</i> to generate a collection of mutants. The GFF fluorescence in mammalian cells is utilized i has the following characteristics:	<i>li</i> and th P mutar	ree c t that	ycles o exhibi	of DNA shuffling were us ited the greatest	
	• Excitation and emission maxima that are 478 nm for primary and secondary excitemission)					
	• High solubility in <i>E. coli</i> for visual dete a promoter recognized by <i>E. coli</i> )	ction of	f trans	sforme	ed cells (if expressed from	1
	• >40-fold increase in fluorescent yield or	ver wild	l-type	GFP		
	This GFP protein will be subsequently referred wild-type GFP.		• •		FP to differentiate it from	1
<i>In Vivo</i> Detection of Cycle 3 GFP Fluorescence	To detect fluorescent cells, it is important detection. The primary excitation peak of C secondary excitation peak at 478 nm. Excita fluorescent emission peak with a maximum Please note that the quantum yield can va the wavelength of light that is used to excit	ycle 3 tion at at 507 i i <b>ry as n</b>	GFP i either nm, as nuch	s at 39 of the s show <b>as 5- t</b>	<ul><li>25 nm. There is a see wavelengths yields a on in the figure below.</li><li>36 10-fold depending on</li></ul>	
	Use of the best filter set will ensure that the	700 -		Excitatio	on and Emission Spectra for GFP	
	optimal regions of the Cycle 3 GFP spectra	700				
	are excited and passed (emitted). For best	600 -			507	
	results, use a filter set designed to detect fluorescence from wild-type GFP (e.g.	500 -			٨	
	XF76 filter from Omega Optical, www.	9 9 9 9 400 -			(	
	omegafilters.com). FITC filter sets can also	uə 400 -				
	be used to detect Cycle 3 GFP fluores- cence. For example, the FITC filter set that	aion]. 300 -				
	we use excites Cycle 3 GFP with light from	Relative Fluoresc				
	460 to 490 nm, which covers the secondary excitation peak. The filter set passes light from 515 to 550, allowing detection of most of the Cycle 3 GFP fluorescence.	200 - 22 100 -		395	5 478	
	For general information about GFP fluor- escence and detection, refer to <i>Current</i> <i>Protocols in Molecular Biology</i> , pages 9.7.22 to 9.7.28 (Ausubel <i>et al.</i> , 1994).	ŝ	330	370	04 Wavelength (nm)	630

# Detection of GFP, continued

Detection of Transformed <i>E. coli</i>	After transformation of <i>E. coli</i> , screen colonies using a hand-held UV light and select glowing cells. To quantitatively assay fluorescence, prepare cell lysates $(10^8 \text{ to } 10^9 \text{ cells/ml})$ from mid-log phase cells using your method of choice. Pellet cell debris and assay supernatant for fluorescence. Be sure to include positive and negative controls.
Detection of Transfected Cells	After transfection, allow the cells to recover and monitor the cells by fluorescence for expression of Cycle 3 GFP. Please note that the CMV promoter is a strong promoter and usually allows detection of Cycle 3 GFP by 24 hours posttransfection. If your promoter is not as strong as CMV, it will take longer to observe fluorescence.
Note	Most media fluoresce because of the presence of riboflavin (Zylka and Schnapp, 1996) and may interfere with detection of Cycle 3 GFP fluorescence. Medium can be removed and replaced with PBS during the assay to alleviate this problem. If cells will be cultured further after assaying, do not keep cells in PBS for a prolonged time. Remove PBS and replace with fresh medium prior to re-incubation.
<i>In Vitro</i> Detection of Cycle 3 GFP	If promoter activity is too low to be detected <i>in vivo</i> , you may prepare mammalian cell lysates and assay fluorescence in a fluorimeter if available. A sample protocol is provided below to prepare lysates.
	1. Wash cell monolayers ( $\sim 10^6$ cells) two times with PBS.
	2. Scrape cells into 1 ml PBS and pellet the cells at 1500 x g for 5 minutes.
	3. Resuspend in 100 µl Cell Lysis Buffer (see recipe on page 26).
	4. Incubate cell suspension on ice or at room temperature for 5 to 10 minutes to lyse the cells.
	5. Centrifuge the cell lysate at 10,000 x g for 10 minutes to pellet nuclei and transfer the supernatant to a fresh tube. Assay the lysate for protein concentration. <b>Note:</b> Do not use protein assays utilizing Coomassie Blue or other dyes. NP-40 interferes with the binding of the dye with the protein.
	6. Assay 20 to 100 µg protein in 0.5-1 ml of PBS. Excite at 395 nm and detect at 510 nm.
Detection of Cycle 3 GFP by Western	If you do not detect fluorescence activity, check for expression of Cycle 3 GFP by Western blot. Antiserum to Cycle 3 GFP is available from Invitrogen as a rabbit polyclonal antibody (see page vi for ordering information).

#### **Creation of Stable Cell Lines**

Introduction	If you wish to create stable cell lines, select for foci using Geneticin <sup>®</sup> Selective Antibiotic. General information and guidelines are provided below for your convenience.				
Geneticin <sup>®</sup> Selective Antibiotic	Geneticin <sup>®</sup> Selective Antibiotic blocks protein synthesis in mammalian cells by interfering with ribosomal function. It is an aminoglycoside, similar in structure to neomycin, gentamycin, and kanamycin. Expression in mammalian cells of the bacterial aminoglycoside phosphotransferase gene (APH), derived from Tn <i>5</i> , results in detoxification of Geneticin <sup>®</sup> Selective Antibiotic (Southern and Berg, 1982).				
Geneticin <sup>®</sup> Selection	Geneticin <sup>®</sup> Selective Antibiotic is available from Invitrogen (Catalog no. 11811-031). Use as follows:				
Guidelines	<ol> <li>Prepare Geneticin<sup>®</sup> Selective Antibiotic in a buffered solution (e.g. 100 mM HEPES, pH 7.3).</li> </ol>				
	2. Use 100 to 1000 $\mu$ g/ml of Geneticin <sup>®</sup> Selective Antibiotic in complete medium.				
	3. Calculate concentration based on the amount of active drug.				
	<ol> <li>Test varying concentrations of Geneticin<sup>®</sup> Selective Antibiotic on your cell line to determine the concentration that kills your cells (kill curve). Cells differ in their susceptibility to Geneticin<sup>®</sup> Selective Antibiotic.</li> </ol>				
	Note: Cells will divide once or twice in the presence of lethal doses of Geneticin <sup>®</sup>				
	Selective Antibiotic, so the effects of the drug take several days to become apparent.				
	Complete selection can take from 2 to 4 weeks of growth in selective medium.				

#### Possible Linearization Sites

To obtain stable transfectants, you may choose to linearize your vector before transfection. While linearizing your vector may not improve your chances of obtaining stable transfectants, it ensures that the vector does not integrate in a way that disrupts the gene of interest. The table below lists some unique sites that can be used to linearize your construct prior to transformation. Other sites are possible. Be sure that your insert does not contain the restriction enzyme site you wish to use to linearize your vector.

Enzyme	Site in pBlue-TOPO <sup>®</sup>	Site in pGlow-TOPO <sup>®</sup>	Location	Supplier
Aat II		5333	Backbone	Many
Afl II	63	63	Upstream of TOPO <sup>®</sup> Cloning site	Many
AlwN I		3934	pUC origin	Many
BamH I	84		Upstream of TOPO <sup>®</sup> Cloning site	Invitrogen, Cat. no. 15201-023
Bgl II	13		Upstream of TOPO <sup>®</sup> Cloning site	Invitrogen, Cat. no. 15213-010
Eam1105 I	6871	4411	Ampicillin gene	AGS*, Fermentas, Takara
Hind III	66		Upstream of TOPO <sup>®</sup> Cloning site	Invitrogen, Cat. no. 15207-012
Kpn I		76	Upstream of TOPO <sup>®</sup> Cloning site	Invitrogen, Cat. no. 15232-010
Pvu I		4781	Ampicillin gene	Invitrogen, Cat. no. 25420-019
Sap I	5862	3402	Backbone	New England Biolabs
Sca I	7351	4891	Ampicillin gene	Invitrogen, Cat. no. 15436-017
Spe I	90	90	Upstream of TOPO <sup>®</sup> Cloning site	Invitrogen, Cat. no. 15443-013

## Appendix

# **TOPO<sup>®</sup> Reporter Control Reactions**

Introduction	you the <i>lac</i>	We recommend performing the following control TOPO <sup>®</sup> Cloning reactions the first time you use the kit to help you evaluate your results. Performing the control reactions using the reagents included in the kit involves producing a control PCR product containing the <i>lac</i> promoter and the LacZ $\alpha$ protein. Successful TOPO <sup>®</sup> Cloning of the control PCR product will yield blue colonies on LB agar plates containing ampicillin and X-gal.					
Before Starting	Be	sure to prepare the following rea	• •	•	action:		
	•	<ul> <li>40 mg/ml X-gal in dimethylformamide (see page 26 for recipe)</li> <li>LB plates containing 100 µg/ml ampicillin and X-gal (two per transformation)</li> </ul>					
		To add X-gal to previously mac of the 40 mg/ml stock solution Protect plates from light.	le agar plates, w	varm the plate to 37°C.	Pipette 40 µl		
Producing the Control PCR	1.	To produce the 500 bp control PCR product containing the <i>lac</i> promoter and LacZ $\alpha$ , set up the following 50 µl PCR:					
Product		Control DNA Template (50 ng	()	1 µl			
		10X PCR Buffer		5 µl			
		50 mM dNTPs		0.5 µl			
		Control PCR Primers (0.1 µg/µ	ıl each)	2 µl			
		Sterile Water		40.5 µl			
		Taq Polymerase (1 unit/µl)		1 µl			
		Total Volume		50 µl			
	2.	Overlay with 70 $\mu$ l (1 drop) of	mineral oil.				
	3.	Amplify using the following c	ycling paramete	rs:			
		Step	Time	Temperature	Cycles		
		Initial Denaturation	2 minutes	94°C	1X		
		Denaturation	1 minute	94°C			
		Annealing	1 minute	60°C	25X		
		Extension	1 minute	72°C			
		Final Extension	7 minutes	72°C	1X		
	4.	Remove 10 µl from the reaction	n and analyze b	y agarose gel electroph	oresis. A		

Remove 10 μl from the reaction and analyze by agarose gel electrophoresis. A discrete 500 bp band should be visible. Proceed to the Control TOPO<sup>®</sup> Cloning Reactions, next page.

# **TOPO<sup>®</sup> Reporter Control Reactions, continued**

#### Control TOPO<sup>®</sup> Cloning Reactions

Using the control PCR product produced on the previous page and either the pBlue-TOPO<sup>®</sup> or the pGlow-TOPO<sup>®</sup> vectors set up two 6 µl TOPO<sup>®</sup> Cloning reactions as described below.

1. Set up control TOPO<sup>®</sup> Cloning reactions:

	Reagent	"Vector Only"	"Vector + PCR Insert"		
	Sterile Water	4 μl	3 µl		
	Salt Solution or Dilute Salt Solution	1 μl	1 μl		
	Control PCR Product		1 µl		
	TOPO <sup>®</sup> vector	1 µl	1 µl		
2.	Incubate at room temperature for <b>5 minutes</b> and place on ice.				
3.	Transform 2 $\mu$ l of each reaction into separate vials of TOP10 One Shot <sup>®</sup> cells (page 9).				
4.	Spread 10-50 µl of each transformation mix onto LB plates containing 50-100 µg/ml				

- 4. Spread 10-50 μl of each transformation mix onto LB plates containing 50-100 μg/ml ampicillin and X-Gal (see page 26). Be sure to plate two different volumes to ensure that at least one plate has well-spaced colonies. For plating small volumes, add 20 μl of SOC to allow even spreading.
- 5. Incubate overnight at 37°C.

Analysis of Hundreds of colonies from the vector + PCR insert reaction should be produced. Greater than 85% of these will be blue and contain the 500 bp insert. Very few colonies (<10% Results of the vector + PCR insert) will be present on the vector only plate. These colonies should be white. Note: pBlue-TOPO<sup>®</sup> will yield dark blue colonies on the "vector + PCR insert" plate and light blue colonies on the "vector only" plate. This is apparently because of a cryptic prokaryotic promoter upstream of the reporter. Please note that no expression of β-galactosidase has been detected in mammalian cells without a promoter. pUC19 plasmid is included to check the transformation efficiency of the One Shot® Transformation competent E. coli. Transform one vial of One Shot® TOP10 cells with 10 pg of pUC19 Control using the protocol on page 9 Plate 10 µl of the transformation mixture plus 20 µl SOC on LB plates containing 100 µg/ml ampicillin. Transformation efficiency should be ~1 x  $10^9$  cfu/µg DNA.

# **TOPO<sup>®</sup> Reporter Control Reactions, continued**

#### Factors Affecting Cloning Efficiency

Please note that lower transformation and/or cloning efficiencies will result from the following variables. Most of these are easily corrected, but if you are cloning large inserts, you may not obtain the expected 85% (or more) cloning efficiency.

Variable	Solution	
pH>9	Check the pH of the PCR amplification reaction and adjust with 1 M Tris-HCl, pH 8.	
Incomplete extension during PCR	Be sure to include a final extension step of 7 to 30 minutes during PCR. Longer PCR products will need a longer extension time.	
Cloning large inserts (>3 kb)	Gel-purify as described on page 20.	
Excess (or overly dilute) PCR product	Reduce (or concentrate) the amount of PCR product. Please note that you can add up to $4 \mu l$ of the PCR to the TOPO <sup>®</sup> Cloning reaction.	
Cloning blunt-ended fragments	Add 3' A-overhangs by incubating with <i>Taq</i> polymerase (page 22).	
PCR cloning artifacts ("false positives")	TOPO <sup>®</sup> Cloning is very efficient for small fragments (< 100 bp) present in certain PCR reactions. Gel-purify your PCR product (page 20).	
PCR product does not contain sufficient 3' A-overhangs even	Increase the time of the final extension to ensure that the 3 <sup>-</sup> ends are adenylated.	
though you used <i>Taq</i> polymerase	Please note that $Taq$ polymerase is less efficient at adding a nontemplate 3' A next to another A. $Taq$ is most efficient at adding a nontemplate 3' A next to a C. You may have to redesign your primers so that they contain a 5' G instead of a 5' T (Brownstein <i>et al.</i> , 1996).	
Size of promoter sequences cloned	For large plasmids, you may have to use electroporation to transform into <i>E. coli</i> . <b>Do</b> <b>not use the chemically competent TOP10</b> <b>cells included in the kit for electroporation.</b> Use electrocompetent TOP10 cells (see page vi for ordering information).	

# **Purifying PCR Products**

Introduction	nece care isol Mol	earing, multiple banding, primer-dimer artifacts, or large PCR products (>1 kb) may essitate gel purification. If you intend to purify your PCR product, be extremely eful to remove all sources of nuclease contamination. There are many protocols to ate DNA fragments or remove oligonucleotides. Please refer to Current Protocols in lecular Biology, Unit 2.6 (Ausubel <i>et al.</i> , 1994) for the most common protocols. ee simple protocols are provided below for your convenience.		
Note	Please note that cloning efficiency may decrease with purification of the PCR product. You may wish to optimize your PCR to produce a single band (see <b>Producing PCR</b> <b>Products</b> , page 6).			
Using the S.N.A.P. <sup>™</sup> Gel		S.N.A.P. <sup>™</sup> Gel Purification Kit (Catalog no. K1999-25) allows you to rapidly purify R products from regular agarose gels.		
Purification Kit	1.	Electrophorese amplification reaction on a 1 to 5% regular TAE agarose gel.		
		Note: Do not use TBE. Borate will interfere with the NaI step (Step 2.)		
	2.	Cut out the gel slice containing the PCR product and melt it at 65°C in 2 volumes of 6 M NaI.		
	3.	Add 1.5 volumes of Binding Buffer.		
	4.	Load solution (no more than 1 ml at a time) from Step 3 onto a S.N.A.P. <sup>TM</sup> column. Centrifuge 1 minute at 3000 x g in a microcentrifuge and discard the supernatant.		
	5.	If you have solution remaining from Step 3, repeat Step 4.		
	6.	Add 900 µl of the Final Wash Buffer.		
	7.	Centrifuge 1 minute at full speed in a microcentrifuge and discard the flow-through.		
	8.	Repeat Step 7.		
	9.	Elute the purified PCR product in 40 $\mu$ l of TE or sterile water. Use 4 $\mu$ l for the TOPO <sup>®</sup> Cloning reaction and proceed as described on page 8.		
Quick S.N.A.P. <sup>™</sup> Method	it oı 1-2	even easier method is to simply cut out the gel slice containing your PCR product, place n top of the S.N.A.P. <sup><math>M</math></sup> column bed, and centrifuge at full speed for 10 seconds. Use $\mu$ l of the flow-through in the TOPO <sup>®</sup> Cloning reaction (page 8). Be sure to make the gel e as small as possible for best results.		

# Purifying PCR Products, continued

Low-Melt Agarose Method	Please note that gel purification will result in a dilution of your PCR product. Use chemically competent cells for transformation.		
	1.	Electrophorese as much as possible of your PCR reaction on a low-melt agarose gel (0.8 to 1.2%) in TAE buffer.	
	2.	Visualize the band of interest and excise the band.	
gel slice melts.		Place the gel slice in a microcentrifuge tube and incubate the tube at 65°C until the gel slice melts.	
		Place the tube at 37°C to keep the agarose melted.	
	5.	Add 4 µl of the melted agarose containing your PCR product to the TOPO <sup>®</sup> Cloning reaction as described on page 8.	
	6.	Incubate the TOPO <sup>®</sup> Cloning reaction at 37°C for 5 to 10 minutes. This is to keep the agarose melted.	
	7.	Transform 2 to 4 $\mu$ l directly into chemically competent TOP10 One Shot <sup>®</sup> <i>E. coli</i> using the method on page 9.	

# Addition of 3<sup>´</sup> A-Overhangs Post-Amplification

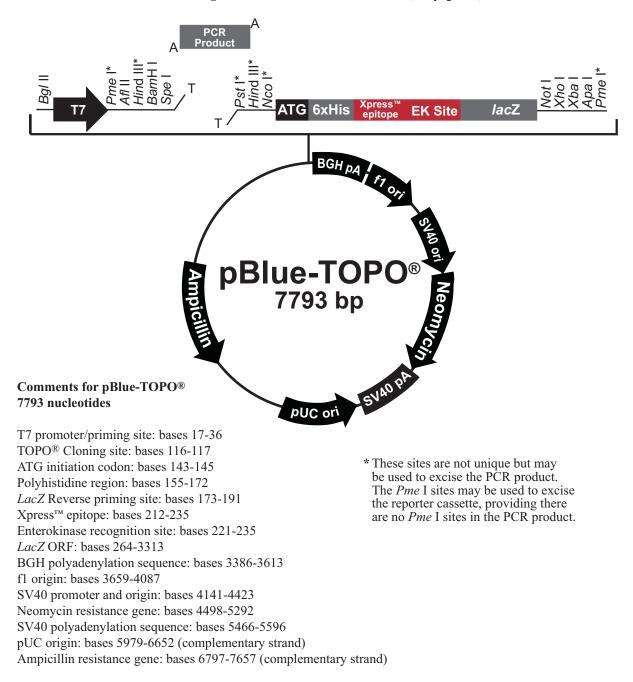
Introduction	Direct cloning of DNA amplified by <i>Vent</i> <sup>®</sup> or <i>Pfu</i> polymerases into TOPO TA Cloning <sup>®</sup> vectors is often difficult because of very low cloning efficiencies. These low efficiencies are caused by the 3' to 5' exonuclease activity associated with proofreading polymerases which removes the 3' A-overhangs necessary for TA Cloning <sup>®</sup> . A simple method is provided below to clone these blunt-ended fragments.			
Before Starting	You will need the following items:			
•	• <i>Taq</i> polymerase			
	• A heat block equilibrated to 72°C			
	• Phenol-chloroform (optional)			
	• 3 M sodium acetate (optional)			
	• 100% ethanol (optional)			
	• 80% ethanol (optional)			
	• TE buffer (optional)			
Procedure	This is just one method for adding 3 <sup>-</sup> adenines. Other protocols may be suitable.			
	1. After amplification with <i>Vent</i> <sup>®</sup> or <i>Pfu</i> polymerase, place vials on ice and add 0.7-1 unit of <i>Taq</i> polymerase per tube. Mix well. It is not necessary to change the buffer.			
	2. Incubate at 72°C for 8-10 minutes (do not cycle).			
	3. Place the vials on ice. The DNA amplification product is now ready for ligation into pBlue-TOPO <sup>®</sup> or pGlow-TOPO <sup>®</sup>			
	<b>Note:</b> If you plan to store your sample(s) overnight before proceeding with TOPO <sup>®</sup> Cloning, you may want to extract your sample(s) with phenol-chloroform to remove the polymerases. After phenol-chloroform extraction, precipitate the DNA with ethanol and resuspend the DNA in TE buffer to the starting volume of the amplification reaction.			
Note	You may also gel-purify your PCR product after amplification with $Vent^{\text{(B)}}$ or $Pfu$ (see previous page). After purification, add <i>Taq</i> polymerase buffer, dATP, and 0.5 unit of <i>Taq</i> polymerase and incubate 10-15 minutes at 72°C. Use 4 µl in the TOPO <sup>(B)</sup> Cloning reaction.			

Vent<sup>®</sup> is a registered trademark of New England Biolabs.

# pBlue-TOPO<sup>®</sup> Map

Мар

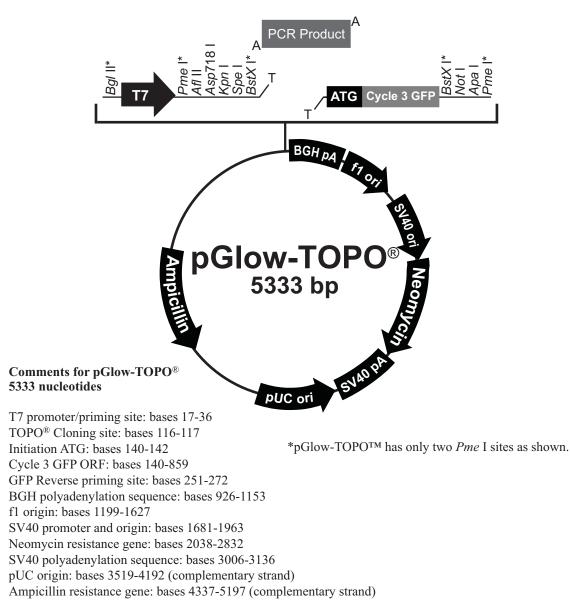
The figure below summarizes the features of the pBlue-TOPO<sup>®</sup> vector. The vector is supplied linearized between base pairs 116 and 117. This is the TOPO<sup>®</sup> Cloning site. **The complete nucleotide sequence is available for downloading from our Web site** (www.invitrogen.com) or from Technical Service (see page 27.).



# pGlow-TOPO<sup>®</sup> Map

Мар

The figure below summarizes the features of the pGlow-TOPO<sup>®</sup> vector. The vector is supplied linearized between base pairs 116 and 117. This is the TOPO<sup>®</sup> Cloning site. **The complete nucleotide sequence is available for downloading from our Web site** (www.invitrogen.com) or from Technical Service (see page 27.).



# Features of the TOPO<sup>®</sup> Reporter Vectors

#### Features

The TOPO<sup>®</sup> Reporter vectors pBlue-TOPO<sup>®</sup> (7793 bp) and pGlow-TOPO<sup>®</sup> (5333 bp) contain the following elements. All features have been functionally tested.

Feature	Benefit
T7 promoter/priming site	Allows for <i>in vitro</i> transcription in the sense orientation and sequencing through the insert
TOPO <sup>®</sup> Cloning site	Allows insertion of your PCR product containing the promoter of interest upstream of the reporter gene.
β-galactosidase (pBlue-TOPO <sup>®</sup> ) or Cycle 3 GFP (pGlow-TOPO <sup>®</sup> )	Allows assay of promoter function either <i>in vitro</i> or <i>in vivo</i> .
<i>LacZ</i> Reverse priming site (pBlue-TOPO <sup>®</sup> ) or GFP Reverse priming site (pGlow-TOPO <sup>®</sup> )	Permits sequencing through the insert
Bovine growth hormone (BGH) polyadenylation signal	Efficient transcription termination and polyadenylation of mRNA (Goodwin and Rottman, 1992)
f1 origin	Allows rescue of single-stranded DNA
SV40 early promoter and origin	Allows efficient, high-level expression of the neomycin resistance gene and episomal replication in cells expressing the SV40 large T antigen
Neomycin resistance gene	Selection of stable transfectants in mammalian cells (Southern and Berg, 1982)
SV40 polyadenylation signal	Efficient transcription termination and polyadenylation of mRNA
pUC origin	High-copy number replication and growth in <i>E. coli</i>
Ampicillin resistance gene (β-lactamase)	Selection of vector in <i>E. coli</i>

# Recipes

LB (Luria-Bertani) Medium and Plates	1.0% 0.5% 1.0%	Composition: 1.0% Tryptone 0.5% Yeast Extract 1.0% NaCl pH 7.0		
	1.	<ol> <li>For 1 liter, dissolve 10 g tryptone, 5 g yeast extract, and 10 g NaCl in 950 ml deionized water.</li> </ol>		
	2.	Adjust the pH of the solution to 7.0 with NaOH and bring the volume up to 1 lit		
	3.	Autoclave on liquid cycle for 20 minutes at 15 psi. Allow solution to cool to $55^{\circ}$ C and add antibiotic (100 µg/ml ampicillin) if needed.		
	4.	Store at room temperature or	at $+4^{\circ}$ C.	
	LB	agar plates		
	1.	Prepare LB medium as above	e, but add 15 g/L agar before autoclaving.	
	2.			
	3.			
	4.	Let harden, then invert and st	ore at $+4^{\circ}$ C, in the dark.	
X-Gal Stock Solution	<ol> <li>To make a 40 mg/ml stock solution, dissolve 400 mg X-Gal in 10 ml dimethylformamide.</li> </ol>			
	2.	Protect from light by storing in a brown bottle at -20°C.		
	3.	To add to previously made agar plates, warm the plate to $37^{\circ}$ C. Pipette 40 µl of the 40 mg/ml stock solution onto the plate, spread evenly, and let dry 15 minutes. Protect plates from light.		
Cell Lysis Buffer	150	50 mM Tris-HCl, pH 7.8 150 mM NaCl 1% Nonidet P-40		
	1.	This solution can be prepared 100 ml, combine:	from the following common stock solutions. For	
		1 M Tris base	5 ml	
		5 M NaCl	3 ml	
		Nonidet P-40	1 ml	
	2.	Bring the volume up to 90 m	with deionized water and adjust the pH to 7.8 with HCl.	
	3.		nl. Store at room temperature.	
	<b>Note:</b> Protease inhibitors may be added at the following concentrations: 1 mM PMSF; 1 $\mu$ g/ml pepstatin; 1 $\mu$ g/ml leupeptin.			

# **Technical Service**

Web Resources	Visit the Invitrogen Web site at <u>www.invitrogen.com</u> for:			
	• Technical resources, including manuals, vector maps and sequences, application notes, MSDSs, FAQs, formulations, citations, handbooks, etc.			
	• Comple	te technical service contact inform	nation	
	Access t	to the Invitrogen Online Catalog		
	Addition	nal product information and speci	al offers	
Contact Us	For more information or technical assistance, call, write, fax, or email. Additional international offices are listed on our Web page ( <u>www.invitrogen.com</u> ).			
Corporate Headquarters Invitrogen Corporation 1600 Faraday Avenue Carlsbad, CA 92008 USA Tel: 1 760 603 7200 Tel (Toll Free): 1 800 955 Fax: 1 760 602 6500 E-mail: tech_service@invitrogen.co	6288	Japanese Headquarters: Invitrogen Japan LOOP-X Bldg. 6F 3-9-15, Kaigan Minato-ku, Tokyo 108-0022 Tel: 81 3 5730 6509 Fax: 81 3 5730 6519 E-mail: jpinfo@invitrogen.com	European Headquarters: Invitrogen Ltd Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF, UK Tel: +44 (0) 141 814 6100 Tech Fax: +44 (0) 141 814 6117 E-mail: eurotech@invitrogen.com	
Material Data Safety Sheets (MSDSs)	MSDSs are available on our Web site at <u>www.invitrogen.com</u> . On the home page, click on <b>Technical Resources</b> and follow instructions on the page to download the MSDS for your product.			
Limited Warranty	Invitrogen is committed to providing our customers with high-quality goods and services. Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about an Invitrogen product or service, contact our Technical Service Representatives. Invitrogen warrants that all of its products will perform according to specifications stated on the certificate of analysis. The company will replace, free of charge, any product that does not meet those specifications. <u>This warranty limits Invitrogen Corporation's liability only to the cost of the product</u> . No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. Invitrogen reserves the right to select the method(s) used to analyze a product unless Invitrogen agrees to a specified method in writing prior to acceptance of the order. Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore Invitrogen makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, please report it to our Technical Service Representatives. <b>Invitrogen assumes no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive. No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.</b>			

#### **Purchaser Notification**

#### Limited Use Label License No: 5 Invitrogen Technology

The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of patents owned or controlled by Invitrogen Corporation which cover this product based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500. Email: outlicensing@invitrogen.com

# Limited Use Label<br/>LicenseThis product is licensed under U.S. Patent Nos. 5,284,933 and 5,310,663 and<br/>foreign equivalents from Hoffmann-LaRoche, Inc., Nutley, NJ and/or<br/>Hoffmann-LaRoche Ltd., Basel, Switzerland and is provided only for use in<br/>research. Information about licenses for commercial use is available from<br/>QIAGEN GmbH, Max-Volmer-Str. 4, D-40724 Hilden, Germany.

# **Purchaser Notification, continued**

Limited Use Label License No: 55 Cycle 3 GFP	The 'cycle 3' mutant GFP was produced by Maxygen, Inc. using DNA shuffling technology. Commercial licensing inquiries should be directed to: Affymax Research Institute, 4001 Miranda Avenue, Palo Alto, CA 94304, U.S.A.	
Limited Use Label License No: 127 GFP with Heterologous Promoter	This product and its use is the subject of one or more of U.S. Patent Nos. 5,491,084 and 6,146,826, and foreign equivalents. This product is sold under license from Columbia University. Rights to use this product are limited to research use only, and expressly exclude the right to manufacture, use, sell or lease this product for use for measuring the level of toxicity for chemical agents and environmental samples in cells and transgenic animals. No other rights are conveyed. Not for human use or use in diagnostic or therapeutic procedures. Inquiry into the availability of a license to broader rights or the use of this product for commercial purposes should be directed to Columbia Innovation Enterprise, Columbia University, Engineering Terrace-Suite 363, New York, New York 10027.	

# **Product Qualification**

Introduction	Invitrogen qualifies the TOP	O <sup>®</sup> Reporter Kits as descri	bed below.	
Restriction Digest	The parental supercoiled pBlue and pGlow vectors are qualified by restriction digest prior to adaptation with topoisomerase. The table below lists the restriction enzymes and the expected fragments.			
	Restriction Enzyme	pBlue	pGlow	
	EcoR I	7777 bp (linearizes)	5317 bp (linearizes)	
	BamH I	7777 bp (linearizes)		
	Xho I	7777 bp (linearizes)	323, 4994 bp	
	Kpn I		5317 bp (linearizes)	
TOPO <sup>®</sup> Cloning Efficiency	Once the vectors have been adapted with topoisomerase I, they are lot-qualified using the control reagents included in the kit. Under conditions described on pages 17-19, a 500 bp control PCR product was TOPO <sup>®</sup> Cloned into each vector and subsequently transformed into the One Shot <sup>®</sup> competent <i>E. coli</i> included with the kit. Each lot of vector should yield greater than 85% cloning efficiency.			
Primers	Both primers have been lot-qualified by DNA sequencing experiments using the dideoxy chain termination technique.			
One Shot <sup>®</sup> Chemically Competent <i>E. coli</i>	All competent cells are tested for transformation efficiency using the control plasmid included in the One Shot <sup>®</sup> kit. Transformed cultures are plated on LB plates containing 100 $\mu$ g/ml ampicillin and the transformation efficiency is calculated. Test transformations are performed in duplicate. Transformation efficiency should be greater than 1 x 10 <sup>9</sup> cfu/ $\mu$ g plasmid DNA. In addition, untransformed cells are tested for the appropriate antibiotic sensitivity and lack of phage contamination.			

#### References

Alam, J., and Cook, J. L. (1990). Reporter Genes: Application to the Study of Mammalian Gene Transcription. Anal. Biochem. *188*, 245-254.

Ausubel, F. M., Brent, R., Kingston, R. E., Moore, D. D., Seidman, J. G., Smith, J. A., and Struhl, K. (1994). Current Protocols in Molecular Biology (New York: Greene Publishing Associates and Wiley-Interscience).

Beale, E. G., Deeb, E. A., Handley, R. S., Akhaven-Tafti, H., and Schaap, A. P. (1992). A Rapid and Simple Chemiluminescent Assay for *Escherichia coli* β-galactosidase. BioTechniques *12*, 320-322.

Bronstein, I., Fortin, J., Stanley, P. E., Stewart, G. S., and Kricka, L. J. (1994). Chemiluminescent and Bioluminescent Reporter Gene Assays. Anal. Biochem. 219, 169-181.

Brownstein, M. J., Carpten, J. D., and Smith, J. R. (1996). Modulation of Non-Templated Nucleotide Addition by *Taq* DNA Polymerase: Primer Modifications that Facilitate Genotyping. BioTechniques 20, 1004-1010.

Chen, C., and Okayama, H. (1987). High-Efficiency Transformation of Mammalian Cells by Plasmid DNA. Molec. Cell. Biol. 7, 2745-2752.

Chu, G., Hayakawa, H., and Berg, P. (1987). Electroporation for the Efficient Transfection of Mammalian Cells with DNA. Nucleic Acids Res. *15*, 1311-1326.

Crameri, A., Whitehorn, E. A., Tate, E., and Stemmer, W. P. C. (1996). Improved Green Fluorescent Protein by Molecular Evolution Using DNA Shuffling. Nature Biotechnology *14*, 315-319.

Felgner, P. L., Holm, M., and Chan, H. (1989). Cationic Liposome Mediated Transfection. Proc. West. Pharmacol. Soc. 32, 115-121.

Felgner, P. L. a., and Ringold, G. M. (1989). Cationic Liposome-Mediated Transfection. Nature 337, 387-388.

Goodwin, E. C., and Rottman, F. M. (1992). The 3'-Flanking Sequence of the Bovine Growth Hormone Gene Contains Novel Elements Required for Efficient and Accurate Polyadenylation. J. Biol. Chem. 267, 16330-16334.

Innis, M. A., Gelfand, D. H., Sninsky, J. J., and White, T. S. (1990) PCR Protocols: A Guide to Methods and Applications. Academic Press, San Diego, CA.

Jain, V. K., and Magrath, I. T. (1991). A Chemiluminescent Assay for Quantitation of  $\beta$ -galactosidase in the Femtogram Range: Application to Quantitation of  $\beta$ -galactosidase in *lacZ*-transfected Cells. Anal. Biochem. *199*, 119-124.

Jongkind, J. F., Verkerk, A., and Sernetz, M. (1986). Detection of Acid-β-galactosidase Activity in Viable Human Fibroblasts by Flow Cytometry. Cytometry 7, 463-466.

Sambrook, J., Fritsch, E. F., and Maniatis, T. (1989). Molecular Cloning: A Laboratory Manual, Second Edition (Plainview, New York: Cold Spring Harbor Laboratory Press).

Shigekawa, K., and Dower, W. J. (1988). Electroporation of Eukaryotes and Prokaryotes: A General Approach to the Introduction of Macromolecules into Cells. BioTechniques *6*, 742-751.

Shuman, S. (1994). Novel Approach to Molecular Cloning and Polynucleotide Synthesis Using Vaccinia DNA Topoisomerase. J. Biol. Chem. 269, 32678-32684.

#### **References**, continued

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. Gen. 1, 327-339.

Wigler, M., Silverstein, S., Lee, L.-S., Pellicer, A., Cheng, Y.-C., and Axel, R. (1977). Transfer of Purified Herpes Virus Thymidine Kinase Gene to Cultured Mouse Cells. Cell *11*, 223-232.

Young, D. C., Kingsley, S. D., Ryan, K. A., and Dutko, F. J. (1993). Selective Inactivation of Eukaryotic  $\beta$ -galactosidase in Assays for Inhibitors of HIV-1 TAT using Bacterial  $\beta$ -galactosidase as a Reporter Enzyme. Anal. Biochem. 215, 24-30.

Zylka, M. J., and Schnapp, B. J. (1996). Optimized Filter Set and Viewing Conditions for the S65T Mutant of GFP in Living Cells. BioTechniques 21, 220-226.

©1999-2006 Invitrogen Corporation. All rights reserved.

For research use only. Not intended for any animal or human therapeutic or diagnostic use.

# invitrogen®

Corporate Headquarters Invitrogen Corporation 1600 Faraday Avenue Carlsbad, CA 92008 T: 1 760 603 7200 F: 1 760 602 6500 E: tech.service@invitrogen.com

For country-specific contact information visit our web site at **www.invitrogen.com**