

pcDNA[™]3.1/*myc*-His A, B, and C

Catalog no. V800-20

Rev. Date: 28 October 2010

Manual part no. 28-0137

MAN0000641

Table of Contents

Kit Contents and Storage	iv
Introduction	1
Overview	1
Methods	2
Cloning into pcDNA [™] 3.1/ <i>myc</i> -His A, B, and C	2
Transformation and Transfection	
Appendix	7
pcDNA [™] 3.1/ <i>myc</i> -His A, B, and C	7
pcDNA™3.1/ <i>myc</i> -His/lacZ	9
Accessory Products	10
Technical Support	11
Purchaser Notification	12
References	13

Kit Contents and Storage

Shipping and Storage

pcDNA[™]3.1/myc-His vectors are shipped on wet ice. Upon receipt, store vectors at -20°C.

Kit Contents

All vectors are supplied in aliquot detailed below. Store the vectors at -20°C.

Vector	Composition	Amount
pcDNA™3.1/ <i>myc</i> -His A, B, and C	40 μl of 0.5 μg/μl vector in 10 mM Tris-HCl, 1 mM EDTA, pH 8.0	20 μg
pcDNA™3.1/myc-His/lacZ	40 μl of 0.5 μg/μl vector in 10 mM Tris-HCl, 1 mM EDTA, pH 8.0	20 μg

Introduction

Overview

Description of the System

pcDNA $^{\text{M}}$ 3.1/myc-His A, B, and C are 5.5-kb vectors derived from pcDNA $^{\text{M}}$ 3.1(+) and designed for high-level expression, purification, and detection of recombinant proteins in mammalian hosts. High-level stable and non-replicative transient expression can be carried out in most mammalian cells. The vectors contain the following elements:

- Three reading frames to facilitate in frame cloning with a C-terminal tag encoding the *myc* (*c-myc*) epitope and a polyhistidine metal-binding peptide
- Human cytomegalovirus immediate-early (CMV) promoter for high-level expression in a wide range of mammalian cells
- Episomal replication in cell lines that are latently infected with SV40 or that express the SV40 large T antigen (*e.g.*, COS7)

The control plasmid, pcDNA $^{\text{\tiny{M}}}3.1/myc$ -His/lacZ, is the pcDNA $^{\text{\tiny{M}}}3.1/myc$ -His C vector with a 3.2-kb fragment containing the β -galactosidase gene cloned in frame with the C-terminal peptide (see page 9). It is included for use as a positive control for transfection, expression, purification, and detection in the cell line of choice.

Experimental Outline

Use the following outline to clone and express your gene of interest in pcDNA $^{\text{\tiny M}}$ 3.1/myc-His:

- Consult the multiple cloning sites described on pages 3-4 to determine which vector (A, B, or C) should be used to clone your gene in frame with the C-terminal *myc* epitope and the polyhistidine tag.
- Ligate your insert into the appropriate vector and transform into *E. coli*. Select transformants on $50-100 \mu g/ml$ ampicillin.
- Analyze your transformants for the presence of insert by restriction digestion.
- Select a transformant with the correct restriction pattern and confirm that your gene is in frame with the C-terminal peptide by sequencing.
- Transfect your construct into the cell line of choice using your own method of transfection.
- Test for expression of your recombinant gene by western blot analysis or functional assay. If you do not have an antibody to your protein, Invitrogen offers the Anti-*myc* antibody or the Anti-His(C-term) antibody to detect your recombinant protein (see page 10).
- To purify your recombinant protein, you may use a metal-chelating resin such as ProBond[™]. ProBond[™] resin is available separately (see page 10).

Methods

Cloning into pcDNA[™]3.1/*myc*-His A, B, and C

General Molecular Biology Techniques

For help with DNA ligations, *E. coli* transformations, restriction enzyme analysis, purification of single-stranded DNA, DNA sequencing, and DNA biochemistry, refer to *Molecular Cloning: A Laboratory Manual* (Sambrook *et al.*, 1989) or *Current Protocols in Molecular Biology* (Ausubel *et al.*, 1994).

E. coli Strain

Many *E. coli* strains are suitable for the propagation of pcDNA[™]3.1/myc-His vectors, including TOP10, TOP10F′, DH5[™]-T1^R. We recommend that you propagate vectors containing inserts in *E. coli* strains that are recombinant deficient (recA) and endonuclease A-deficient (endA).

For your convenience, TOP10F' is available as chemically competent or electrocompetent cells from Invitrogen (see page 10).

Maintenance of pcDNA[™]3.1/ *myc*-His

To propagate and maintain the pcDNA $^{\text{\tiny M}}3.1/myc$ -His vectors, use the supplied 0.5 $\mu g/\mu l$ stock solution in TE, pH 8.0 to transform a recA, endA E. coli strain like TOP10, TOP10F', DH5 α , JM109, or equivalent. Select transformants on LB plates containing 50–100 $\mu g/m l$ ampicillin. Be sure to prepare a glycerol stock of each plasmid for long-term storage.



Your insert should contain a Kozak consensus sequence with an ATG initiation codon for proper initiation of translation (Kozak, 1987; Kozak 1990). An example of a Kozak consensus sequence is provided below. Other sequences are possible, but the G or A at position –3 and the G at position +4 (shown in bold) illustrates the most commonly occurring sequence with strong consensus. Replacing one of the two bases at these positions provides moderate consensus, while having neither results in weak consensus. The ATG initiation codon is shown underlined.

(G/A)NNATGG

Continued on next page

Cloning into pcDNA[™]3.1/myc-His A, B, and C, Continued

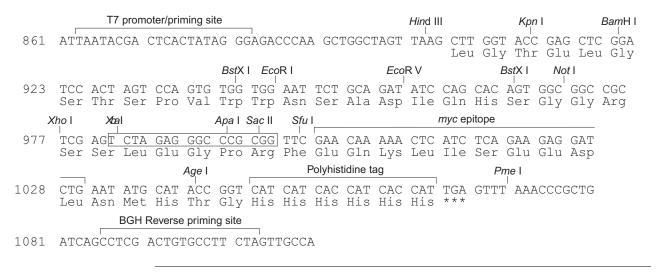
Multiple Cloning Site of Version A

Below is the multiple cloning site for pcDNA $^{\text{\tiny{M}}}$ 3.1/myc-His A. Restriction sites are labeled to indicate the cleavage site. **Note that there is a stop codon between the** BamH **I site and the** BstX **I site.** The boxed nucleotides indicate the variable region. The multiple cloning site has been confirmed by sequencing and functional testing. The vector sequence of pcDNA $^{\text{\tiny{M}}}$ 3.1/myc-His A is available for downloading from our website at www.invitrogen.com or from **Technical Support** (see page 11).

T7 promoter/priming site									Hind III Kpn I					BamH I				
861	ATTA	ATAC	CGA (CTCAC	CTATA	AG GO	; GAGA(CCCA	A GCT	rggc1	ragt	TAA	GCT Ala	TGG Trp	TAC Tyr	CGA Ara	GCT Ala	CGG Ara
	BstX I EcoR I						EcoR V				BstX I Not I			9				
922	ATC Ile	CAC His	TAG ***		AGT Ser	GTG Val	GTG Val	GAA Glu	TTC Phe	TGC Cys	AGA Arg	TAT Tyr	CCA Pro	GCA Ala	CAG Gln	TGG Trp	CGG Arg	CCG Pro
	Xho I Xba I Apa I Sfu <u>I</u>						<i>myc</i> epitope											
976	CTC Leu			AGA Arg	GGG Gly	CCC	TTC Phe					ATC Ile				GAT Asp	CTG Leu	AAT Asn
Age I Polyhistidine tag Pmel																		
1030	ATG Met	CAT His	ACC Thr	001	CAT His	CAT His	CAC His		CAC His	CAT His	TGA ***	GTTT	r'AAA(CCC (GCTGA	ATCA	GC .	
	BGH Reverse priming site																	
1083	1083 CTCGACTGTG CCTTCTAG																	

Multiple Cloning Site of Version B

Below is the multiple cloning site for pcDNA[™]3.1/*myc*-His B. Restriction sites are labeled to indicate the cleavage site. The boxed nucleotides indicate the variable region. The multiple cloning site has been confirmed by sequencing and functional testing. The vector sequence of pcDNA[™]3.1/*myc*-His B is available for downloading from our website at www.invitrogen.com or from **Technical Support** (see page 11).



Continued on next page

Cloning into pcDNA[™]3.1/myc-His A, B, and C, Continued

Multiple Cloning Site of Version C

Below is the multiple cloning site for pcDNA $^{\text{M}}3.1/myc$ -His C. Restriction sites are labeled to indicate the cleavage site. The boxed nucleotides indicate the variable region. The multiple cloning site has been confirmed by sequencing and functional testing. The vector sequence of pcDNA $^{\text{M}}3.1/myc$ -His C is available for downloading from our website at www.invitrogen.com or from **Technical Support** (see page 11).

	T7 promoter/priming site								<i>Hin</i> d III					Kpn I		
861	ATTAATA	CGA (CTCA	CTATA	AG GC	GAGA(CCCA	A GCI	[GGC]	TAGT				GTA (Val 1		AGC Ser
	<i>Bam</i> H I						BstX I	<i>Eco</i> R	.1			<i>Eco</i> R	V		Bst	ΧI
918	TCG GAT Ser Asp	CCA Pro	0	0 - 0		TGT Cys	GGT Gly			-	-			AGC Ser		
	Not I Xho I BstE II					Sfu I				myc epitope						
969	GCG GCC Ala Ala						TTC Phe	GAA Glu	CAA Gln	AAA Lys	CTC Leu	ATC Ile	TCA Ser	GAA Glu	GAG Glu	GAT Asp
				Age I			F	Polyhist	idine ta	ag			P	me I		
1020	CTG AAT Leu Asn	ATG Met		ACC Thr			0	CAC His	OLIL	CAC His	OLIL	TGA ***	GTT	TAAA(CCC	
			BGH R	Reverse	primin	g site	_									
1069	GCTGATCA	AGC (CTCGA	ACTG	rg co	CTTCT	ragt:	r GC								

Transformation and Transfection

Introduction

If you need more details about the techniques discussed, refer to *Molecular Cloning: A Laboratory Manual* (Sambrook *et al.*, 1989) or *Current Protocols in Molecular Biology* (Ausubel *et al.*, 1994).

Method of Transformation

Transform your ligation mixtures into a competent recA, endA E. coli strain $(e.g. TOP10, TOP10F', DH5\alpha)$ and select on LB plates containing 50–100 $\mu g/ml$ ampicillin. Select 10–20 clones and analyze for the presence and orientation of your insert. For your convenience, TOP10F' is available as chemically competent or electrocompetent cells from Invitrogen (see page 10).



We recommend that you sequence your construct to confirm that your gene is fused in frame with the *myc* epitope and the C-terminal polyhistidine tag. We suggest using the T7 Promoter and BGH Reverse primer sequences. Refer to the diagrams on pages 3–4 for the sequence and location of the primer binding sites.

For your convenience, Invitrogen offers a custom primer synthesis service. For more information, visit www.invitrogen.com or contact **Technical Support** (page 11).

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 using the PureLink $^{\text{\tiny M}}$ HiPure Miniprep Kit or the PureLink $^{\text{\tiny M}}$ HiPure Midiprep Kit (see page 10 for ordering information).

Methods of Transfection

For established cell lines (*e.g.*, HeLa), 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 *Current Protocols in Molecular Biology*.

Methods of transfection include calcium phosphate (Chen & Okayama, 1987; Wigler *et al.*, 1977), lipid-mediated (Felgner *et al.*, 1989; Felgner & Ringold, 1989) and electroporation (Chu *et al.*, 1987; Shigekawa & Dower, 1988). For high efficiency transfection in a broad range of mammalian cells, we recommend using Lipofectamine[™] 2000 Reagent available from Invitrogen. For more information on Lipofectamine[™] 2000 and other transfection reagents available, visit our website at www.invitrogen.com or contact **Technical Support** (page 11).

Positive Control

pcDNATM3.1/myc-His/lacZ is provided as a positive control vector for mammalian transfection and expression (see page 9). It may be used to optimize transfection conditions for your cell line. The gene encoding β -galactosidase is expressed in mammalian cells under the CMV promoter. A successful transfection will result in β -galactosidase expression that can be easily assayed (see below).

Detection of Fusion Proteins

A number of antibodies are available from Invitrogen (see page 10) that can be used to detect expression of your fusion protein from pcDNA $^{\text{\tiny{TM}}}$ 3.1/myc-His.

Continued on next page

Transformation and Transfection, Continued

Assay for βgalactosidase Activity

Transform your ligation mixtures into a competent recA, endA E. coli strain $(e.g., TOP10, TOP10F', DH5<math>\alpha$) and select on LB plates containing 50–100 μ g/ml ampicillin. Select 10–20 clones and analyze for the presence and orientation of your insert.

Geneticin[®] Selection Guidelines

Geneticin® is available separately from Invitrogen. Geneticin® blocks protein synthesis in mammalian cells by interfering with ribosomal function. It is an aminoglycoside, similar in structure to neomycin, gentamycin, and kanamycin. Expression of the bacterial aminoglycoside phosphotransferase gene (APH), derived from Tn5, in mammalian cells results in detoxification of Geneticin® (Southern and Berg, 1982). Use as follows:

- Prepare Geneticin[®] in a buffered solution (*e.g.*, 100 mM HEPES, pH 7.3).
- Use 100–1000 μg/ml of Geneticin® in complete medium.
- Calculate concentration based on the amount of active drug (check lot label).
- Test varying concentrations of Geneticin® on your cell line to determine the
 concentration that kills your cells (kill curve). Cells differ in their
 susceptibility to Geneticin®.

Cells will divide once or twice in the presence of lethal doses of Geneticin[®], so the effects of the drug take several days to become apparent. Complete selection can take from 3 to 6 weeks of growth in selective medium.

Preparing Cells for Lysis

Use the procedure below to prepare cells for lysis prior to purification of your protein on ProBond^{TM}. You will need 5×10^6 to 1×10^7 cells for purification of your protein on a 2-ml ProBond^{TM} column (see the ProBond^{TM} Purification manual).

- 1. Seed cells in five T-75 flasks or two to three T-175 flasks.
- 2. Grow the cells in selective medium until they are 80–90% confluent.
- 3. Harvest the cells by treating with trypsin-EDTA for 2–5 minutes or by scraping the cells in PBS.
- 4. Inactivate the trypsin by diluting with fresh medium (if necessary) and transfer the cells to a sterile microcentrifuge tube.
- 5. Centrifuge the cells at 1,500 rpm for 5 minutes. You may lyse the cells immediately or freeze in liquid nitrogen and store at -70°C until needed.

Lysis of Cells

If you are using ProBond™ resin, refer to the ProBond™ Purification manual for details about sample preparation for chromatography. If you are using another resin, refer to the manufacturer's instruction for recommendations on sample preparation.



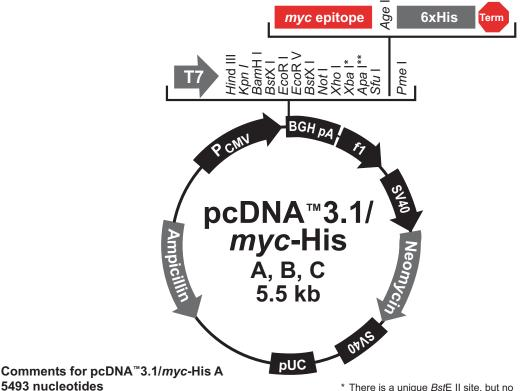
The C-terminal peptide containing the *myc* epitope and polyhistidine tag will add approximately 3 kDa to the size of your protein. The size of the lacZ/*myc*-His fusion protein is approximately 121 kDa.

Appendix

pcDNA[™]3.1/*myc*-His A, B, and C

Map of pcDNA[™]3.1/*myc*-His

The figure below summarizes the features of the pcDNA^m3.1/myc-His vectors. The nucleotide sequence for pcDNA m 3.1/myc-His A is available for downloading from www.invitrogen.com or from Technical Support (page 11). Details of the multiple cloning sites for pcDNA m 3.1/myc-His A, B, and C are shown on pages 3–4.



CMV promoter: bases 209-863

T7 promoter/priming site: bases 863-882 Multiple cloning site: bases 902-999

myc epitope: bases 997-1026 Polyhistidine tag: bases 1042-1059

BGH reverse priming site: bases 1082-1099 BGH polyadenylation signal: bases 1081-1295 f1 origin of replication: bases 1358-1771 SV40 promoter and origin: bases 1836-2160 Neomycin resistance gene: bases 2196-2990 SV40 polyadenylation signal: bases 3166-3296

pUC origin: bases 3679-4352

Ampicillin resistance gene: bases 4497-5357 (complementary strand)

- * There is a unique BstE II site, but no Xba I or Apa I sites in version C.
- ** There is a unique Sac II site between the Apa I site and the Sfu I site in version B only.

pcDNA[™]3.1/myc-His A, B, and C, Continued

Features of pcDNA[™]3.1/*myc*-His

pcDNA[™]3.1/*myc*-His A (5,493 bp), pcDNA[™]3.1/*myc*-His B (5,497 bp), and pcDNA[™]3.1/*myc*-His C (5,489 bp) contain the following elements. All features have been functionally tested.

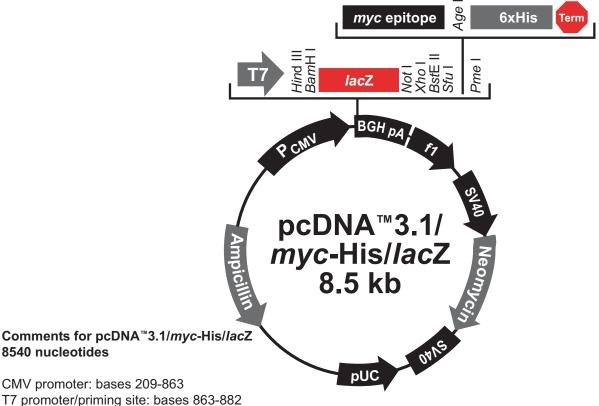
Feature	Benefit
Human cytomegalovirus (CMV) immediate-early promoter/enhancer	Allows efficient, high-level expression of your recombinant protein (Andersson <i>et al.</i> , 1989; Boshart <i>et al.</i> , 1985; Nelson <i>et al.</i> , 1987).
T7 promoter/priming site	Allows for <i>in vitro</i> transcription in the sense orientation and sequencing through the insert.
Multiple cloning site in three reading frames	Allows insertion of your gene and facilitates cloning in frame with the <i>myc</i> epitope and polyhistidine C-terminal tag.
<i>myc</i> epitope (c <i>-myc</i>) (Glu-Gln-Lys-Leu-Ile-Ser-Glu-Glu-Asp-Leu)	Allows detection of your recombinant protein with the Anti- <i>myc</i> Antibody or Anti- <i>myc</i> -HRP Antibody (Evan <i>et al.</i> , 1985).
C-terminal polyhistidine tag	Allows purification of your recombinant protein on metal-chelating resin such as $ProBond^{TM}$.
	In addition, the C-terminal polyhistidine tag is the epitope for the Anti-His (C-term) Antibody and the Anti-His (C-term)-HRP Antibody.
BGH reverse priming site	Allows 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 (Geneticin®) 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> .

pcDNA[™]3.1/*myc*-His/lacZ

Map of Control Vector

pcDNA[™]3.1/*myc*-His/*lacZ* is a 8,540-bp control vector containing the gene for β-galactosidase. pcDNA[™]3.1/myc-His C was digested with EcoR V and Not I. A 3.2-kb blunt-Not I fragment containing the β-galactosidase gene was then ligated into pcDNA[™]3.1/*myc*-His C in frame with the C-terminal peptide.

The figure below summarizes the features of the pcDNA[™]3.1/*myc*-His/*lac*Z vector. The nucleotide sequence for pcDNA™3.1/myc-His/lacZ is available by downloading it from our website (www.invitrogen.com) or by contacting Technical Support (page 11).



lacZ with C-terminal tag: 963-4106 lacZ ORF: bases 963-4019 myc epitope: bases 4044-4073 Polyhistidine tag: bases 4089-4106 BGH reverse priming site: bases 4129-4146 BGH polyadenylation signal: bases 4128-4342 f1 origin of replication: bases 4405-4818 SV40 promoter and origin: bases 4883-5207 Neomycin resistance gene: bases 5243-6037

SV40 polyadenylation signal: bases 6213-63473

pUC origin: bases 67267-7399

Ampicillin resistance gene: bases 7544-8404 (complementary strand)

Accessory Products

Introduction

The following additional products may be used with the pcDNA $^{\text{\tiny{TM}}}$ 3.1/myc-His vectors. For more information, $\underline{www.invitrogen.com}$ or contact **Technical Support** (see page 11).

Product	Amount	Catalog no.		
ProBond™ Purification System	6 purifications	K850-01		
ProBond™ Resin	50 ml	R801-01		
Flodolid Resili	150 ml	R801-15		
PureLink™ HiPure Plasmid Miniprep Kit	100 preps	K2100-03		
PureLink™ HiPure Plasmid Midiprep Kit	25 preps	K2100-04		
Electrocomp™ TOP10F′	$2 \times 20 \text{ rxns}$ $6 \times 20 \text{ rxns}$	C665-11 C665-24		
One Shot™ TOP10F′ (chemically competent cells)	$21 \times 50 \mu l$	C3030-03		

Detection of Fusion Proteins

A number of antibodies are available from Invitrogen that can be used to detect expression of your fusion protein from pcDNA $^{\text{\tiny{M}}}3.1/myc$ -His. The table below describes the antibodies available and ordering information. The amount supplied is sufficient for 25 Westerns.

Product	Purpose	Catalog no.		
Anti-myc	Detects 10 amino acid epitope derived from c-myc	R950-25		
Anti-myc-HRP	See above. Provided as an HRP conjugate for time-saving detection.	R951-25		
Anti-His(C-term)	Detects the C-terminal polyhistidine tag (requires the free carboxyl group for detection)	R930-25		
Anti-His(C-term)-HRP	See above. Provided as an HRP conjugate for time-saving detection.	R931-25		

Technical Support

Web Resources



Visit the Invitrogen website at <u>www.invitrogen.com</u> for:

- Technical resources, including manuals, vector maps and sequences, application notes, MSDSs, FAQs, formulations, citations, handbooks, etc.
- Complete technical support contact information
- Access to the Invitrogen Online Catalog
- Additional product information and special offers

Contact Us

For more information or technical assistance, call, write, fax, or email. Additional international offices are listed on our website (www.invitrogen.com).

Corporate Headquarters:

5791 Van Allen Way Carlsbad, CA 92008 USA Tel: 1 760 603 7200

Tel (Toll Free): 1 800 955 6288

Fax: 1 760 602 6500

E-mail: tech_support@invitrogen.com

Japanese Headquarters:

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:

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

MSDS

Material Safety Data Sheets (MSDSs) are available on our website at www.invitrogen.com/msds.

Certificate of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to www.invitrogen.com/support and search for the Certificate of Analysis by product lot number, which is printed on the box.

Limited Warranty

Invitrogen (a part of Life Technologies Corporation) 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 Support Representatives. All Invitrogen products are warranted to 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. This warranty limits the Company's liability to only the price of the product. 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. The Company reserves the right to select the method(s) used to analyze a product unless the Company 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 the Company 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 Support Representatives.

Life Technologies Corporation shall have 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.

Purchaser Notification

Limited Use Label License No. 22: Vectors and Clones Encoding Histidine Hexamer This product is licensed under U.S. Patent Nos. 5,284,933 and 5,310,663 and foreign equivalents from Hoffmann-LaRoche, Inc., Nutley, NJ and/or Hoffmann-LaRoche Ltd., Basel, Switzerland and is provided only for use in research. Information about licenses for commercial use is available from QIAGEN GmbH, Max-Volmer-Str. 4, D-40724 Hilden, Germany.

References

- Andersson, S., Davis, D. L., Dahlbäck, H., Jörnvall, H., and Russell, D. W. (1989). Cloning, Structure, and Expression of the Mitochondrial Cytochrome P-450 Sterol 26-Hydroxylase, a Bile Acid Biosynthetic Enzyme. J. Biol. Chem. 264, 8222-8229.
- 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).
- Boshart, M., Weber, F., Jahn, G., Dorsch-Häsler, K., Fleckenstein, B., and Schaffner, W. (1985). A Very Strong Enhancer is Located Upstream of an Immediate Early Gene of Human Cytomegalovirus. Cell 41, 521-530.
- 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.
- Evan, G. I., Lewis, G. K., Ramsay, G., and Bishop, V. M. (1985). Isolation of Monoclonal Antibodies Specific for *c-myc* Proto-oncogene Product. Mol. Cell. Biol. *5*, 3610-3616.
- Felgner, P. L., Holm, M., and Chan, H. (1989). Cationic Liposome Mediated Transfection. Proc. West. Pharmacol. Soc. 32, 115-121.
- Felgner, P. L., 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.
- Kozak, M. (1987). An Analysis of 5'-Noncoding Sequences from 699 Vertebrate Messenger RNAs. Nucleic Acids Res. 15, 8125-8148.
- Kozak, M. (1990). Downstream Secondary Structure Facilitates Recognition of Initiator Codons by Eukaryotic Ribosomes. Proc. Natl. Acad. Sci. USA 87, 8301-8305.
- Miller, J. H. (1972). Experiments in Molecular Genetics (Cold Spring Harbor, New York: Cold Spring Harbor Laboratory).
- Nelson, J. A., Reynolds-Kohler, C., and Smith, B. A. (1987). Negative and Positive Regulation by a Short Segment in the 5′-Flanking Region of the Human Cytomegalovirus Major Immediate-Early Gene. Molec. Cell. Biol. 7, 4125-4129.
- 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.
- 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.

©2009, 2010 Life Technologies Corporation. All rights reserved.

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

Notes



Corporate Headquarters

Invitrogen Corporation 5791 Van Allen Way Carlsbad, CA 92008

T: 1 760 603 7200

F: 1 760 602 6500

E: tech_support@invitrogen.com

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