

pcDNA[™]4/*myc*-His A, B, and C

Catalog no. V863-20

Rev. Date: 27 October 2010 Manual part no. 25-0236 MAN0000078

User Manual

Table of Contents

Kit Contents and Storage	iv
Introduction	1
Product Overview	1
Methods	2
Cloning into pcDNA [™] 4/ <i>myc</i> -His A, B, and C	
Transfection and Analysis	
Creating Stable Cell Lines	8
Appendix	
pcDNA [™] 4/ <i>myc</i> -His Vector pcDNA [™] 4/ <i>myc</i> -His/ <i>lac</i> Z	11
pcDNA [™] 4/ <i>myc-</i> His Vector	11
pcDNA [™] 4/ <i>myc</i> -His Vector pcDNA [™] 4/ <i>myc</i> -His/ <i>lac</i> Z Zeocin [™] Recipes	11
pcDNA [™] 4/ <i>myc</i> -His Vector pcDNA [™] 4/ <i>myc</i> -His/ <i>lac</i> Z Zeocin [™]	11
pcDNA [™] 4/ <i>myc</i> -His Vector pcDNA [™] 4/ <i>myc</i> -His/ <i>lac</i> Z Zeocin [™] Recipes	
pcDNA [™] 4/ <i>myc</i> -His Vector pcDNA [™] 4/ <i>myc</i> -His/ <i>lac</i> Z Zeocin [™] Recipes Accessory Products	

Kit Contents and Storage

Shipping and
Storage $pcDNA^{TM}4/myc$ -His vectors are shipped on wet ice. Upon receipt, store vectors
at $-20^{\circ}C$.

Kit Contents All vectors are supplied as detailed below. **Store the vectors at –20°C.**

Item	Composition	Amount
pcDNA ^{M} 4/ <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 [™] 4/ <i>myc</i> -His/ <i>lacZ</i>	40 μL of 0.5 μg/μL vector in 10 mM Tris-HCl, 1 mM EDTA, pH 8.0	20 µg

Introduction

Product Overview

Description of the System	pcDNA [™] 4/ <i>myc</i> -His A, B, and C are 5.1 kb vectors designed for overproduction of recombinant proteins in mammalian cell lines. Features of the vectors allow purification and detection of expressed proteins (see pages 11-12 for more information). High-level stable and transient expression can be carried out in most mammalian cells. The vectors contain the following elements:
	• Human cytomegalovirus immediate-early (CMV) promoter for high-level expression in a wide range of mammalian cells
	• Three reading frames to facilitate in-frame cloning with a C-terminal peptide encoding the <i>myc</i> (<i>c-myc</i>) epitope and a polyhistidine (6xHis) metal-binding tag
	• Zeocin [™] resistance gene for selection of stable cell lines (Mulsant <i>et al.</i> , 1988) (see page 14 for more information).
	• 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, $pcDNA^{M}4/myc$ -His/ <i>lacZ</i> is included for use as a positive control for transfection, expression, and detection in the cell line of choice.
Experimental Outline	Use the following outline to clone and express your gene of interest in pcDNA [™] 4/ <i>myc</i> -His:
	1. Consult the multiple cloning sites described on pages 3-4 to determine which vector (A, B, or C) to use for cloning your gene in frame with the C-terminal <i>myc</i> epitope and the polyhistidine tag.
	 Ligate your insert into the appropriate vector and transform into <i>E. coli</i>. Select transformants on 50 to 100 µg/mL ampicillin or 25 to 50 µg/mL Zeocin[™] in Low Salt LB. For more information, see page 16.
	3. Analyze your transformants for the presence of insert by restriction digestion.
	4. Select a transformant with the correct restriction pattern and use sequencing to confirm that your gene is cloned in-frame with the C-terminal peptide.
	5. Transfect your construct into the cell line of choice using your own method of transfection. Generate a stable cell line, if desired.
	6. Test for expression of your recombinant gene by western blot analysis or functional assay. For antibodies to the <i>myc</i> epitope or the C-terminal polyhistidine tag, see page 18.
	7. To purify your recombinant protein, you may use metal-chelating resin such as ProBond [™] . ProBond [™] resin is available separately (see page 17).

Methods

Cloning into pcDNA[™]4/*myc*-His A, B, and C

General Molecular Biology Techniques	For help with DNA ligations, <i>E. coli</i> transformations, restriction enzyme analysis, purification of single-stranded DNA, DNA sequencing, and DNA biochemistry, refer to <i>Molecular Cloning: A Laboratory Manual</i> (Sambrook <i>et al.</i> , 1989) or <i>Current Protocols in Molecular Biology</i> (Ausubel <i>et al.</i> , 1994).
<i>E. coli</i> Strain	Many <i>E. coli</i> strains are suitable for the growth of this vector including TOP10F', DH5 α F', JM109, and INV α F'. We recommend that you propagate vectors containing inserts in <i>E. coli</i> strains that are recombination deficient (<i>rec</i> A) and endonuclease A deficient (<i>end</i> A).
	For your convenience, TOP10F' is available from Invitrogen as chemically competent or electrocompetent cells (see page 17).
Transformation Method	You may use any method of your choice for transformation. Chemical transformation is the most convenient for most researchers. Electroporation is the most efficient and the method of choice for large plasmids.
Maintaining pcDNA [™] 4/ <i>myc</i> -His	To propagate and maintain the pcDNA TM 4/ <i>myc</i> -His vectors, use a small amount of the supplied 0.5 μ g/ μ L stock solution in TE, pH 8.0 to transform a <i>recA</i> , <i>endA E. coli</i> strain like TOP10F', DH5 α , JM109, or equivalent. Select transformants on LB plates containing 50 to 100 μ g/mL ampicillin or 25 to 50 μ g/mL Zeocin TM in Low Salt LB. Be sure to prepare a glycerol stock of each plasmid for long-term storage (see page 5).
Cloning Considerations	Your insert should contain a Kozak translation initiation sequence for proper initiation of translation (Kozak, 1987; Kozak, 1991; Kozak, 1990). An example of a Kozak consensus sequence is provided below. Note that other sequences are possible, but the A at position -3 and the G at position +4 are the most critical for function (shown in bold). The ATG initiation codon is shown underlined. A NN <u>ATG</u> G
	To express your gene as a recombinant fusion protein, you must clone your gene in frame with the C-terminal peptide. The vector is supplied in three reading frames to facilitate cloning. See pages 3-4 to develop a cloning strategy. If you wish to express your protein WITHOUT the C-terminal peptide, be sure to include a stop codon.

Cloning into pcDNA[™]4/myc-His A, B, and C, Continued

Multiple Cloning Below is the multiple cloning site for pcDNA[™]4/myc-His A. Restriction sites are labeled to indicate the cleavage site. The boxed nucleotides indicate the variable region. Note that there is a stop codon between the *Bam*H I site and the *BstX* I site. The multiple cloning site has been confirmed by sequencing and functional testing. The vector sequence of pcDNA[™]4/myc-His A is available for downloading from our website (www.invitrogen.com) or from Technical Support (see page 19).

	_	Τ7	promo	oter/prir	ning sit	te	_					Hind I	I	Acc ₆₅	I Kp	nl		BamH I
861	ATT <i>I</i>	AATA	CGA (CTCAC	CTATA	AG GO	GAGA	CCCA	A GC	rggc	ΓAGT	TAA	GCT Ala	TGG Trp	TAC Tyr	0.011	GCT Ala	000
						Bst)	< I*	EcoR I			Pst I	EcoR V	<i>'</i>		BstX I	*	Not I	
922		CAC His		TCC Ser	AGT Ser	GTG Val	GTG Val	GAA Glu		TGC Cys		TAT Tyr	CCA Pro	0011	CÀG Gln	100	000	000
	Xho I		Xba I			Apa I	BstB	I				тус	epitop	е				
976	CTC	GAG	TCT	AGA Arg	GGG Gly	CCC] TTC	GAA				Myc ATC Ile	TCA	GAA	GAG Glu	-	~ - ~	AAT Asn
976	CTC	GAG Glu	TCT			CCC Pro	 TTC Phe	GAA	Gln			ATC Ile	TCA	GAA		-	~ - ~	

BGH Reverse priming site

1083 CTCGACTGTG CCTTCTAG

*Note that there are two *BstX* I sites in the polylinker.

Multiple Cloning
 Below is the multiple cloning site for pcDNA[™]4/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[™]4/myc-His B is available for downloading from our website (www.invitrogen.com) or from Technical Support (see page 19).

		Τ7	promo	oter/prir	ning si	te	_					<i>Hin</i> d II	I,	4 <i>cc</i> 65 I	Kp	n I		BamH I
861	ATTA	AATAG	CGA (CTCAC	CTATA	AG G	GAGAC	CCCAZ	A GCI	GGC	fagt	TAAC		r GGI 1 Gly		C GAC	- O - C	C GGA 1 Gly
						BstX	* Eco	RI		Pst	I Ecc	RV		E	ßstX I*	No	tl	
923		ACT Thr			GTG Val	TGG Trp	TGG Trp	AAT Asn	TCT Ser	GCA Ala	GAT Asp	ATC Ile		CAC His			GGC Gly	
Х	(ho I	Xb	al			Apa I	Sac II	<i>Bst</i> B	I				тус ер	itope				
977	TCG Ser			GAG Glu										TCA Ser				
					Age I			F	Polyhist	idine ta	ag			Pr	ne l			
1028	CTG Leu		ATG Met	CAT His	ACC Thr		CAT His				CAC His		TGA ***	GTTI	' AAA	ACCCC	GCTG	
			BGH F	Reverse	primir	ng site												
1081	ATCA	AGCCI	rcg A	ACTGI	[GCC]	FT C	ragt1	rgcc <i>i</i>	A									
	*Note	that	there	are tv	vo Bs	tX I si	tes in	the p	olylir	nker.								

Cloning into pcDNA[™]4/myc-His A, B, and C, Continued

Multiple Cloning Site of Version C Below is the multiple cloning site for $pcDNA^{M}4/myc$ -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 $pcDNA^{M}4/myc$ -His C is available for downloading from our website (www.invitrogen.com) or from **Technical Support** (see page 19).

861	T7 pron ATTAATACGA	noter/priming site		CCCAA	A GCI	GGCI	TAGT		AGC 1		GTA (pnl CCG A Pro S	AGC Ser
	BamH I		E	ßstX I*	<i>Eco</i> R	I		Pst I	<i>Eco</i> R	V		BstX	I *
918	TCG GAT CCA Ser Asp Pro		CAG TGT Gln Cys	GGT Gly	GGA Gly	ATT Ile	CTG Leu	CAG Gln	ATA Ile	TCC Ser	AGC Ser	ACA Thr	GTG Val
	Not I Xho	ol <i>Bst</i> Ell		BstB I					<i>тус</i> ер	oitope			
969	GCG GCC GC Ala Ala Ala			TTC Phe			AAA Lys			TCA Ser			GAT Asp
		Age I		F	Polyhist	idine ta	ag			Pi	ne l		
1020	CTG AAT ATC Leu Asn Met		GGT CAT Gly His	CAT His	CAC His	CAT His	CAC His	CAT His	TGA ***	GTTI	' TAAA(CCC	
		BGH Reverse	priming site										
		BGH Reverse	priming site										

1069 GCTGATCAGC CTCGACTGTG CCTTCTAGTT GC

*Note that there are two *BstX* I sites in the polylinker.

(i.e., TOP10, DH5a, DH10, etc.).

E. coli
TransformationTransform your ligation mixtures into a competent *recA*, *endA E. coli* strain
(e.g., TOP10F', DH5 α) and select on LB plates containing 50–100 µg/mL
ampicillin or 25–50 µg/mL ZeocinTM in Low Salt LB (see page 16). Select 10–20
clones and analyze for the presence and orientation of your insert.**Q** ImportantAny *E. coli* strain that contains the complete Tn5 transposable element
(i.e., DH5 α F1Q, SURE, SURE2) encodes the *ble* (bleomycin resistance gene). These
strains will confer resistance to ZeocinTM. For the most efficient selection, we
recommend that you choose an *E. coli* strain that does not contain the Tn5 gene

Cloning into pcDNA[™]4/myc-His A, B, and C, Continued



Preparing a Glycerol Stock We recommend that you sequence your construct with the T7 Forward and BGH Reverse primers to confirm that your gene is fused in frame with the *myc* epitope and the C-terminal polyhistidine tag. For ordering primers, see page 17.

Once you have identified the correct clone, be sure to purify the colony and make a glycerol stock for long-term storage. It is also a good idea to keep a DNA stock of your plasmid at -20° C.

- 1. Streak the original colony out on an LB plate containing 50 μg/mL ampicillin or 25 μg/mL Zeocin[™] in Low Salt LB. Incubate the plate at 37°C overnight.
- 2. Isolate a single colony and inoculate into 1–2 mL of LB with 50 μ g/mL ampicillin.
- 3. Grow the culture to mid-log phase ($OD_{600} = 0.5-0.7$).
- 4. Mix 0.85 mL of culture with 0.15 mL of sterile glycerol and transfer to a cryovial.
- 5. Store at –80°C.

Transfection and Analysis

Introduction	Once you have confirmed that your construct is in the correct orientation and fused in frame with the C-terminal peptide, you are ready to transfect your cell line of choice. We recommend that you include the positive control vector and a mock transfection to evaluate your results.
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 lipids, decreasing transfection efficiency. We recommend isolating plasmid DNA using the PureLink [™] HiPure Miniprep Kit or the PureLink [™] HiPure Midiprep Kit (see page 17 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. We recommend 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> (see page 21). 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 Lipofectamine [™] 2000 Reagent for mammalian transfection For more details, call Technical Support (see page 19) or visit our website at www.invitrogen.com.
Positive Control	pcDNA TM 4/ <i>myc</i> -His/ <i>lacZ</i> is provided as a positive control vector for mammalian cell transfection and expression (see page 13) and may be used to optimize transfection conditions for your cell line. The gene encoding β -galactosidase is expressed in mammalian cells under the control of the CMV promoter. A successful transfection will result in β -galactosidase expression that can be easily assayed (see below).
Assay for β-galactosidase Activity	You may assay for β -galactosidase expression by activity assay using cell-free lysates (Miller, 1972) or by staining the cells for activity. Invitrogen offers the β -Gal Assay Kit and the β -Gal Staining Kit for fast and easy detection of β -galactosidase expression (see page 17).

Transfection and Analysis, Continued

Detecting Fusion Proteins	fusi To from opt	eral antibodies are available from Invitrogen to detect expression of your ion protein from pcDNA ^{M} 4/ <i>myc</i> -His (see page 18). detect fusion protein by western blot, you will need to prepare a cell lysate m transfected cells. We recommend that you perform a time course to imize expression of the fusion protein (<i>e.g.</i> 24, 48, 72 hours, etc. after hsfection). To lyse cells:
		Wash cell monolayers (~10 ⁶ cells) once with phosphate-buffered saline (PBS).
	2.	Scrape cells into 1 mL PBS and pellet the cells at $1,500 \times g$ for 5 minutes.
	3.	Resuspend in 50 µL Cell Lysis Buffer (see page 16). Other lysis buffers may be suitable.
	4.	Incubate cell suspension at 37°C for 10 minutes to lyse the cells.
	5.	Centrifuge the cell lysate at $10,000 \times g$ for 10 minutes to pellet nuclei and transfer the supernatant to a fresh tube. Assay the lysate for protein concentration.
		Note: Do not use protein assays utilizing Coomassie [®] Blue or other dyes. NP-40 interferes with the binding of the dye with the protein.
	6.	Add SDS-PAGE sample buffer to a final concentration of 1X and boil the sample for 5 minutes.
	7.	Load 20 μ g of lysate onto an SDS-PAGE gel and electrophorese. Use the appropriate percentage of acrylamide to resolve your fusion protein.
Note		e C-terminal peptide containing the <i>myc</i> epitope and the polyhistidine tag will approximately 3 kDa to the size of your protein.
Purification	on a mai	a will need 5×10^6 to 1×10^7 transfected cells for purification of your protein a 2 mL ProBond [™] column (or other metal-chelating column). Refer to the nufacturer's instructions before attempting to purify your fusion protein. To pare cells for lysis, refer to the protocol on page 10.

Creating Stable Cell Lines

Introduction	The pcDNA [™] 4/ <i>myc</i> -His vectors contain the Zeocin [™] resistance gene for selection of stable cell lines using Zeocin [™] . We recommend that you test the sensitivity of your mammalian host cell to Zeocin [™] as natural resistance varies among cell lines. General information and guidelines are provided below for your convenience. For more information about Zeocin [™] , refer to page 14.
Effect of Zeocin [™] on Sensitive and Resistant Cells	 The method of killing with Zeocin[™] is quite different from neomycin and hygromycin. Cells do not round up and detach from the plate. Sensitive cells will exhibit the following morphological changes upon exposure to Zeocin[™]: Vast increase in size Abnormal cell shape
	• Presence of large empty vesicles in the cytoplasm (breakdown of the
	 endoplasmic reticulum and golgi apparatus or scaffolding proteins) Breakdown of plasma and nuclear membrane (appearance of many holes in these membranes). Eventually, these "cells" will completely break down and only "strings" of protein will remain.
	Zeocin [™] -resistant cells should continue to divide at regular intervals to form distinct colonies. There should not be any distinct morphological changes in Zeocin [™] -resistant cells when compared to cells not under selection with Zeocin [™] .
Selection in Mammalian Cell Lines	To generate a stable cell line expressing your protein, you need to determine the minimum concentration of Zeocin [™] required to kill your untransfected host cell line. Typically, concentrations between 50 and 1,000 µg/mL Zeocin [™] are sufficient to kill the untransfected host cell line. Test a range of concentrations (see below) to ensure that you determine the minimum concentration necessary for your cell line.
	1. Seed cells $(2 \times 10^5 \text{ cells}/60 \text{ mm plate})$ for each time point and allow cells to adhere overnight.
	 The next day, substitute culture medium with medium containing varying concentrations of Zeocin[™] (e.g., 0, 50, 125, 250, 500, 750, and 1,000 µg/mL).
	3. Replenish the selective medium every 3–4 days, and observe the percentage of surviving cells.
	 Count the number of viable cells at regular intervals to determine the appropriate concentration of Zeocin[™] that prevents growth.
	Continued on next page

Creating Stable Cell Lines, Continued

Possible Sites for Linearization

To obtain stable transfectants, you may choose to linearize your vector before transfection. While linearizing your vector may not improve the efficiency of transfection, it increases the chances that the vector does not integrate in a way that disrupts the gene of interest. The table below lists unique sites that may be used to linearize your construct prior to transformation. **Other restriction sites are possible. Note that the cleavage site is indicated for versions A, B, and C of pcDNA[™]4/***myc***-His. Be sure that your insert does not contain the restriction enzyme site you wish to use to linearize your vector.**

Enzyme	Restriction Site (bp)	Location	Supplier
	(A,B,C)		
Bgl II	13	Upstream of CMV promoter	Many
Mfe I	161	Upstream of CMV promoter	New England Biolabs
Nru I	209	Upstream of CMV promoter	Many
Mlu I	229	5' end of CMV promoter	Many
Bst1107 I	2881 (A), 2885 (B), 2877 (C)	End of SV40 poly A	AGS*, Fermentas, Takara, Boehringer-Mannhiem
Eam1105 I	4153 (A), 4157 (B), 4149 (C)	Ampicillin gene	AGS*, Fermentas, Takara
Fsp I	4375 (A), 4379 (B), 4371 (C)	Ampicillin gene	Many
Pvu I	4523 (A), 4527 (B), 4519 (C)	Ampicillin gene	Many
Sca I	4633 (A), 4637 (B), 4629 (C)	Ampicillin gene	Many
Ssp I	4957 (A), 4961 (B), 4953 (C)	Ampicillin gene	Many

*Angewandte Gentechnologie Systeme

Selecting Stable Integrants

Once the appropriate Zeocin[™] concentration is determined, you can generate a stable cell line with your construct.

- 1. Transfect your cells using the appropriate protocol for your cell line. Include a sample of untransfected cells as a negative control.
- 2. After transfection, wash the cells once with 1X PBS and add fresh medium to the cells.
- 3. 48 hours after transfection, split the cells into fresh medium containing Zeocin[™] at the appropriate concentration for your cell line. Split the cells such that they are no more than 25% confluent.
- 4. Replenish selective medium every 3–4 days until Zeocin[™]-resistant colonies are detected.
- 5. Pick and expand colonies.

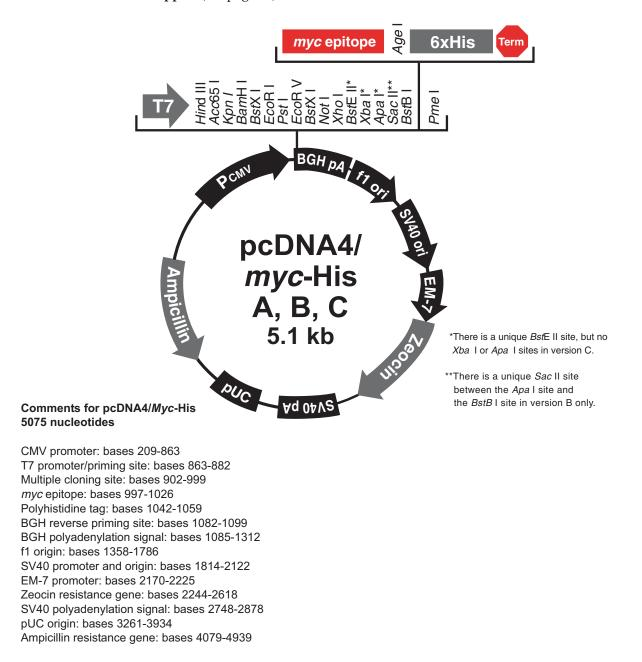
Creating Stable Cell Lines, Continued

Preparing Cells for Lysis	Use the procedure below to prepare cells for lysis prior to purification of your protein on ProBond [™] . You will need 5 × 10 ⁶ to 1 × 10 ⁷ cells for purification of your protein on a 2 mL ProBond [™] column (see ProBond [™] Purification System manual).							
	1.	Seed cells in five T-75 flasks or 2 to 3 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 to 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 240 × g for 5 minutes. Resuspend the cell pellet in PBS.						
	6.	Centrifuge the cells at 240 × g for 5 minutes. You may lyse the cells immediately or freeze in liquid nitrogen and store at –80°C until needed.						
Lysis of Cells	ma	you are using ProBond [™] resin, refer to the Probond [™] Purification System nual for details about sample preparation for chromatography. you are using other metal-chelating resin, refer to the manufacturer's						
		truction for recommendations on sample preparation.						

Appendix

pcDNA[™]4/*myc*-His Vector

Map of pcDNA[™]4/*myc*-His The figure below summarizes the features of the pcDNATM4/*myc*-His vectors. The vector sequences for pcDNATM4/*myc*-His A, B, and C are available for downloading from our website (www.invitrogen.com) or from **Technical Support** (see page 19).



pcDNA[™]4/myc-His Vector, Continued

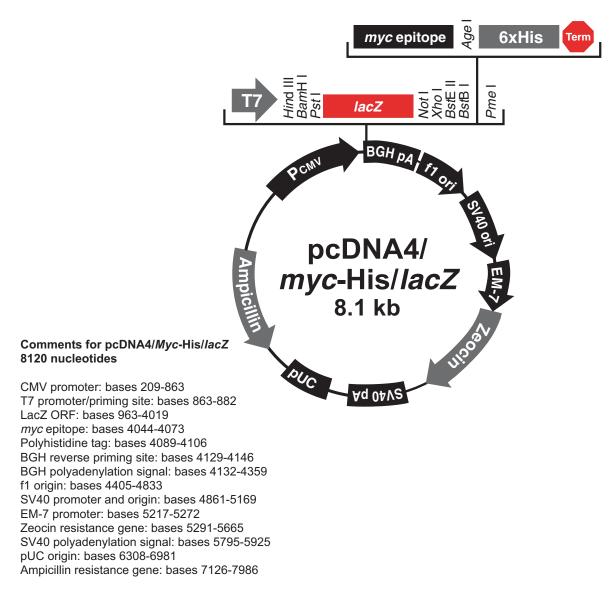
Features of pcDNA[™]4/*myc*-His A (5075 bp), pcDNA[™]4/*myc*-His B (5079 bp), and pcDNA[™]4/*myc*-His C (5071 bp) contain the following elements. All features have been functionally tested.

Feature	Benefit
Human cytomegalovirus (CMV) immediate-early promoter/enhancer	Permits efficient, high-level expression of your recombinant protein (Andersson et al., 1989; Boshart et al., 1985; Nelson et al., 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 (Glu-Gln-Lys-Leu-Ile-Ser-Glu-Glu-Asp- Leu)	Allows detection of your recombinant protein with the Anti- <i>myc</i> Antibody the Anti- <i>myc</i> -HRP Antibody, or the Anti- <i>myc</i> -AP Antibody (Evans et al., 1985) (see page 18 for ordering).
C-terminal polyhistidine (6xHis) tag	Permits purification of your recombinant protein on metal- chelating resin such as ProBond [™] . In addition, the C-terminal polyhistidine tag is the epitope for the Anti-His(C-term) Antibody, the Anti-His (C-term)-HRP Antibody and the Anti-His(C-term)-AP (Lindner et al., 1997) (see page 18).
BGH reverse priming site	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 Zeocin [™] resistance gene and episomal replication in cells expressing the SV40 large T antigen.
EM-7 promoter	Synthetic promoter based on the bacteriophage T7 promoter for expression of the Zeocin [™] resistance gene in <i>E. coli</i> .
Zeocin [™] resistance gene	Selection of transformants in <i>E. coli</i> and stable transfectants in mammalian cells (Drocourt et al., 1990; Mulsant et al., 1988).
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	Selection of transformants in <i>E. coli</i> .
(β-lactamase)	

pcDNA[™]4/*myc*-His/*lac*Z

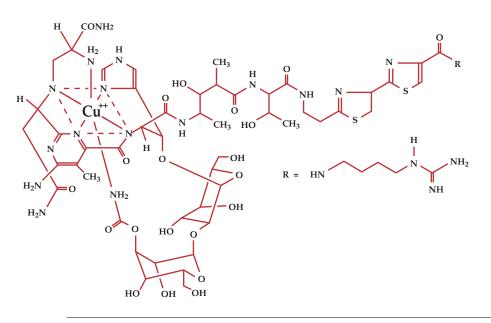
Map of Control Vector

pcDNATM4/*myc*-His/*lacZ* is a 8120 bp control vector containing the gene for β -galactosidase. This vector was constructed by ligating a 3,880 bp *Bam*H I-*Stu* I fragment containing the CMV promoter and the ZeocinTM resistance gene from pcDNATM4/*myc*-His B to a 4,240 bp *Bam*H I-*Stu* I fragment containing the *lacZ* gene, *myc* epitope, and polyhistidine tag from pcDNATM3.1/*myc*-His/*lacZ*. The figure below summarizes the features of the pcDNATM4/*myc*-His/*lacZ* vector. The vector sequence for pcDNATM4/*myc*-His/lacZ is available for downloading from our website (www.invitrogen.com) or by contacting **Technical Support** (see page 19).



Zeocin[™]

Introduction	The pcDNA [™] 4/ <i>myc</i> -His vectors contain the Zeocin [™] resistance gene for selection of stable cell lines using Zeocin [™] . We recommend that you test the sensitivity of your mammalian host cell to Zeocin [™] as natural resistance varies among cell lines. General information and guidelines are provided in this section for your convenience.
Zeocin [™]	Zeocin [™] is a member of the bleomycin/phleomycin family of antibiotics isolated from <i>Streptomyces</i> . Antibiotics in this family are broad spectrum antibiotics that act as strong anti-bacterial and anti-tumor drugs. They show strong toxicity against bacteria, fungi (including yeast), plants, and mammalian cells. The Zeocin [™] resistance protein has been isolated and characterized (Calmels <i>et al.</i> ,
	1991; Drocourt <i>et al.</i> , 1990). This protein, the product of the <i>Sh ble</i> gene (<i>Streptoalloteichus hindustanus</i> bleomycin gene), is a 13.7 kDa protein that binds Zeocin [™] in a stoichiometric manner to inhibit its DNA strand cleavage activity. Expression of this protein in eukaryotic and prokaryotic hosts confers resistance to Zeocin [™] .
Molecular Weight, Formula, and Structure	The formula for Zeocin TM is $C_{55}H_{86}O_{21}N_{20}S_2Cu$ -HCl and the molecular weight is 1,527.5 daltons. Zeocin TM is an HCl salt. The diagram below shows the structure of Zeocin TM .



Zeocin[™], Continued

Applications of Zeocin[™]

ZeocinTM is used for selection in mammalian cells (Mulsant *et al.*, 1988); plants (Perez *et al.*, 1989); yeast (Baron *et al.*, 1992); and prokaryotes (Drocourt *et al.*, 1990). Suggested concentrations of ZeocinTM for selection in mammalian cell lines and *E. coli* are listed below:

Organism	Zeocin [™] Concentration and Selective Medium
E. coli	25–50 μg/mL in low salt LB medium* (see page 16 for recipe)
Mammalian Cells	50–1,000 μ g/mL (varies with cell line)

^{*}Efficient selection requires that the concentration of NaCl be no more than 5 g/liter (< 90 mM).

Handling Zeocin[™]

- High salt and acidity or basicity inactivates Zeocin[™]. Therefore, we recommend that you reduce the salt in bacterial medium and adjust the pH to 7.5 to keep the drug active (see page 16).
- Store Zeocin[™] at –20°C and thaw on ice before use.
- Zeocin[™] is light sensitive. Store drug, plates, and medium containing drug in the dark.
- Wear gloves, a laboratory coat, and safety glasses or goggles when handling solutions containing Zeocin[™].
- Zeocin[™] is toxic. Do not ingest or inhale solutions containing the drug.

Recipes

Low Salt LB Medium with Zeocin [™]	(<90 prep cont) mM) and the pH must be 7 pare LB broth and plates usi	t concentration of the medium must remain low 7.5. For selection in <i>E. coli</i> , it is imperative that you ng the following recipe. Note the lower salt to use low salt LB medium will result in n of the drug.	
	Low Salt LB Medium:			
	10 g Tryptone 5 g NaCl 5 g Yeast Extract			
		 Combine the dry reagents above and add deionized, distilled water to 950 n Adjust pH to 7.5 with 5 M NaOH. Bring the volume up to 1 liter. For plates, add 15 g/L agar before autoclaving. 		
	2.			
	3. Thaw Zeocin ^{TM} on ice and vortex before removing an aliquot.			
	5. Store plates at 4°C in the dark. Plates containing Zeocin [™] are stable for 1-2 weeks.			
Cell Lysis Buffer	50 mM Tris, pH 7.8 150 mM NaCl 1% Nonidet P-40			
	1.	1. This solution can be prepared from the following common stock solutions For 100 mL, combine:		
		1 M Tris base	5 mL	
		5 M NaCl	3 mL	
		Nonidet P-40	1 mL	
	2.	Bring the volume up to 90 mL with deionized water and adjust the pH to 7.8 with HCl.		
	3. Bring the volume up to 100 mL. Store at room temperature.		mL. Store at room temperature.	
		Note: Protease inhibitors may be added at the following concentrations: 1 mM PMSF		
		g/mL pepstatin g/mL leupeptin		

Accessory Products

Introduction

The following products may be used with the pcDNA[™]4/*myc*-His vectors. For details, visit www.invitrogen.com or contact **Technical Support** (page 19).

Item	Amount	Catalog no.
ProBond [™] Purification System	6 × 2 mL precharged, prepacked ProBond [™] resin columns and buffers for native and denaturing purification	K850-01
ProBond [™] Resin	50 mL	R801-01
rrodonu kesin	150 mL	R801-15
Electrocomp [™] TOP10F′	$5 \times 80 \ \mu L$	C665-55
One Shot [®] TOP10F´ (chemically competent cells)	21 × 50 μL	C3030-03
EKMax [™] Enterokinase	250 units	E180-01
PureLink [™] HiPure Plasmid Miniprep Kit	100 preps	K2100-03
PureLink [™] HiPure Plasmid Midiprep Kit	25 preps	K2100-04
β-Gal Assay Kit	80 mL	K1455-01
β-Gal Staining Kit	1 kit	K1465-01
Zeocin™	1 gram	R250-01
Zeocin	5 grams	R250-05
Lipofectamine [™] 2000 Reagent	0.75 mL	11668-027

Primers

For your convenience, Invitrogen offers a custom primer synthesis service. Visit www.invitrogen.com for more details.

Accessory Products, Continued

Antibodies

If you do not have an antibody specific to your protein, Invitrogen offers the Anti-*myc*, or Anti-His(C-term) antibodies to detect your recombinant fusion protein. Horseradish peroxidase (HRP)- and alkaline phosphatase (AP)– conjugated antibodies are available for convenient one-step detection.

Antibody	Epitope	Catalog no.
Anti-myc	Detects a 10 amino acid epitope	R950-25
Anti-myc-HRP	derived from <i>c-myc</i> (Evan <i>et al.</i> , 1985):	R951-25
Anti-myc-AP	EQKLISEEDL	R952-25
Anti-His(C-term)	Detects the C-terminal polyhistidine	R930-25
Anti-His(C-term)-HRP	tag (requires the free carboxyl group for detection) (Lindner et al., 1997):	R931-25
Anti-His(C-term)-AP	HHHHHH-COOH	R932-25

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 (<u>www.invitrogen.com</u>).

Corporate Headquarte 5791 Van Allen Way Carlsbad, CA 92008 US Tel: 1 760 603 7200 Tel (Toll Free): 1 800 99 Fax: 1 760 602 6500 E-mail: <u>tech_support@</u>	SA 55 6288	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: <u>eurotech@invitrogen.com</u>		
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	with high-quality satisfied with our about an Invitrog All Invitrogen pr- certificate of anal meet those specif <u>product.</u> No warn warranty is appli instructions. The product unless th the order. Invitrogen makes occasional typog warranty of any I discover an error Representatives. Life Technologie incidental, indire	to f Life Technologies Corporation) is committed to providing our customers goods and services. Our goal is to ensure that every customer is 100% products and our service. If you should have any questions or concerns en product or service, contact our Technical Support Representatives. oducts are warranted to perform according to specifications stated on the ysis. The Company will replace, free of charge, any product that does not ications. This warranty limits the Company's liability to only the price of the anty is granted for products beyond their listed expiration date. No cable unless all product components are stored in accordance with Company reserves the right to select the method(s) used to analyze a e Company agrees to a specified method in writing prior to acceptance of every effort to ensure the accuracy of its publications, but realizes that the aphical or other error is inevitable. Therefore the Company makes no tind regarding the contents of any publications or documentation. If you in any of our publications, please report it to our Technical Support s Corporation shall have no responsibility or liability for any special, ect or consequential loss or damage whatsoever. The above limited and exclusive. No other warranty is made, whether expressed or implied, arranty of merchantability or fitness for a particular purpose.			

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. For products that are subject to multiple limited use label licenses, the terms of the most restrictive limited use label license shall control. Life Technologies Corporation will not assert a claim against the buyer of infringement of patents owned or controlled by Life Technologies 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, Life Technologies is willing to accept return of the product with a full refund. For information about purchasing a license to use this product or the technology embedded in it for any use other than for research use please contact Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008; Phone (760) 603-7200 or e-mail: outlicensing@lifetech.com.

Limited Use LabelTheLicense No. 22:forVectors andHadClones EncodingresHistidine HexamerQI

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).
- Baron, M., Reynes, J. P., Stassi, D., and Tiraby, G. (1992). A Selectable Bifunctional b-Galactosidase::Phleomycin-resistance Fusion Protein as a Potential Marker for Eukaryotic Cells. Gene 114, 239-243.
- 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.
- Calmels, T., Parriche, M., Burand, H., and Tiraby, G. (1991). High Efficiency Transformation of Tolypocladium geodes Conidiospores to Phleomycin Resistance. Curr. Genet. 20, 309-314.
- 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.
- Drocourt, D., Calmels, T. P. G., Reynes, J. P., Baron, M., and Tiraby, G. (1990). Cassettes of the Streptoalloteichus hindustanus ble Gene for Transformation of Lower and Higher Eukaryotes to Phleomycin Resistance. Nucleic Acids Res. 18, 4009.
- Evans, 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. (1991). An Analysis of Vertebrate mRNA Sequences: Intimations of Translational Control. J. Cell Biology 115, 887-903.
- Kozak, M. (1990). Downstream Secondary Structure Facilitates Recognition of Initiator Codons by Eukaryotic Ribosomes. Proc. Natl. Acad. Sci. USA 87, 8301-8305.
- Lindner, P., Bauer, K., Krebber, A., Nieba, L., Kremmer, E., Krebber, C., Honegger, A., Klinger, B., Mocikat, R., and Pluckthun, A. (1997). Specific Detection of His-tagged Proteins With Recombinant Anti-His Tag scFv-Phosphatase or scFv-Phage Fusions. BioTechniques 22, 140-149.
- Miller, J. H. (1972). Experiments in Molecular Genetics (Cold Spring Harbor, New York: Cold Spring Harbor Laboratory).

- Mulsant, P., Tiraby, G., Kallerhoff, J., and Perret, J. (1988). Phleomycin Resistance as a Dominant Selectable Marker in CHO Cells. Somat. Cell Mol. Genet. 14, 243-252.
- 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.
- Perez, P., Tiraby, G., Kallerhoff, J., and Perret, J. (1989). Phleomycin Resistance as a Dominant Selectable Marker for Plant Cell Transformation. Plant Mol. Biol. 13, 365-373.
- 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.
- 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.

invitrogen

Corporate Headquarters 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