



pcDNA[™] 6/His A, B, and C

Catalog no. V222-20

Rev. date: 28 October 2010

Manual part no. 25-0237

MAN0000079

User Manual

Table of Contents

Kit Contents and Storage	iv
Introduction	1
Product Overview	1
Methods	2
Cloning into pcDNA™6/His A, B, and C.....	2
Transfection and Analysis	7
Creating Stable Cell Lines.....	10
Appendix.....	14
pcDNA™6/His Vector	14
pcDNA™6/His/ <i>lacZ</i>	16
Accessory Products.....	17
Technical Support	19
Purchaser Notification	20
References	22

Kit Contents and Storage

Shipping and Storage

pcDNA™6/His vectors are shipped on wet ice. Upon receipt, store vectors at -20°C .

Contents

20 μg each of pcDNA™6/His A, B, and C are supplied at 0.5 $\mu\text{g}/\mu\text{L}$ in 10 mM Tris-HCl, 1 mM EDTA, pH 8.0 in a total volume of 40 μL .

20 μg of pcDNA™6/His/*lacZ* is supplied at 0.5 $\mu\text{g}/\mu\text{L}$ in 10 mM Tris-HCl, 1 mM EDTA, pH 8.0 in a total volume of 40 μL .

Introduction

Product Overview

Description of the System

pcDNA™6/His A, B, and C are 5.2 kb vectors designed for the overproduction of recombinant proteins in mammalian cell lines. Features of the vectors allow for the purification and detection of expressed proteins (see pages 14–15 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 an N-terminal peptide encoding the Xpress™ epitope and a polyhistidine (6xHis) metal-binding tag
- Blasticidin resistance gene (*bsd*) for the selection of stable cell lines (Kimura *et al.*, 1994)
- 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™6/His/*lacZ*, 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™6/His.

1. Consult the multiple cloning sites depicted on pages 3–5 to determine which vector (A, B, or C) should be used to clone your gene in-frame with the N-terminal Xpress™ epitope and the polyhistidine tag.
 2. Ligate your insert into the appropriate vector and transform into *E. coli*. Select transformants on 50 to 100 µg/mL ampicillin or 50 µg/mL blasticidin.
 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 N-terminal peptide.
 5. Transfect your construct into a 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 an antibody to the Xpress™ epitope, see page 17.
 7. To purify your recombinant protein, you may use a metal-chelating resin such as ProBond™. ProBond™ resin is available separately (see page 18 for ordering information).
-

Methods

Cloning into pcDNA™ 6/His A, B, and C

Before Starting

Diagrams are provided on pages 3–5 to help you ligate your gene of interest in-frame with the N-terminal peptide. General considerations for cloning and transformation are listed below.

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 growth of this vector including TOP10F', DH5αF', and INVαF' (see page 17 for ordering information). We recommend that you propagate vectors containing inserts in *E. coli* strains that are recombination deficient (*recA*) and endonuclease A deficient (*endA*).

Transformation Method

You may use any method of your choice for transformation. Chemical transformation is the most convenient for many researchers. Electroporation is the most efficient and the method of choice for large plasmids.

Maintaining pcDNA™ 6/His

To propagate and maintain the pcDNA™ 6/His vectors, use the supplied 0.5 µg/µL stock solution in TE, pH 8.0 to transform a *recA*, *endA* *E. coli* strain like TOP10F' or INVαF'. Select transformants on LB plates containing 50 to 100 µg/mL ampicillin or 50 µg/mL blasticidin. Prepare a glycerol stock of each plasmid for long-term storage (see page 6).



Note

The pcDNA™ 6/His vectors are fusion vectors. To ensure proper expression of your recombinant fusion protein, you must clone your gene in-frame with the ATG at base pairs 920-922. This will create a fusion with the N-terminal polyhistidine tag, Xpress™ epitope, and the enterokinase cleavage site. The vector is supplied in three reading frames to facilitate cloning. See pages 3–5 to develop a cloning strategy.

If you wish to clone as close as possible to the enterokinase cleavage site, follow the guidelines below:

1. Digest pcDNA™ 6/His A, B, or C with *Kpn* I.
2. Create blunt ends with T4 DNA polymerase and dNTPs. See (Ausubel *et al.*, 1994) for a detailed protocol.
3. Clone your blunt-ended insert in-frame with the lysine codon (AAG) of the enterokinase recognition site.

Following enterokinase cleavage, no vector-encoded amino acid residues will be present in your protein.

Continued on next page

Cloning into pcDNA™ 6/His A, B, and C, Continued

Kozak Sequence for Mammalian Expression

If you will be recombining your entry clone with a destination vector for mammalian expression, your insert should contain a Kozak consensus sequence with an ATG initiation codon for proper initiation of translation (Kozak, 1987; Kozak, 1990; Kozak, 1991). 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

Multiple Cloning Site of Version A

Below is the multiple cloning site for pcDNA™ 6/His A. Restriction sites are labeled to indicate the cleavage site. The boxed nucleotide indicates the variable region. **Note that there is a stop codon between the Xba I site and the Apa I site.** The multiple cloning site has been confirmed by sequencing and functional testing. **The sequence of pcDNA™ 6/His A is available at www.invitrogen.com or by request from Technical Support (see page 19).**

```

                                     T7 promoter priming site
839  CACTGCTTAC TGGCTTATCG AAATTAATAC GACTCACTAT AGGGAGACCC AAGCTGGCTA

                                     Polyhistidine Region
899  GCGTTTAAAC TTAAGCTTAC C ATG GGG GGT TCT CAT CAT CAT CAT CAT CAT
                                     Met Gly Gly Ser His His His His His His
                                     Xpress™ Epitope
950  GGT ATG GCT AGC ATG ACT GGT GGA CAG CAA ATG GGT CGG GAT CTG TAC
                                     Gly Met Ala Ser Met Thr Gly Gly Gln Gln Met Gly Arg Asp Leu Tyr

998  GAC GAT GAC GAT AAG GTA CCT AGG ATC CAG TGT GGT GGA ATT CTG CAG
     Asp Asp Asp Asp Lys Val Pro Arg Ile Gln Cys Gly Gly Ile Leu Gln
     Enterokinase recognition site ▲ EK cleavage site

     EcoR V      BstX I*      Not I      Xho I      Xba I      Apa I
1046 ATA TCC AGC ACA GTG GCG GCC GCT CGA GTC TAG AGGGCCCGTT TAAACCCGCT
     Ile Ser Ser Thr Val Ala Ala Ala Arg Val ***

                                     BGH reverse priming site
1099 GATCAGCCTC GACTGTGCCT TCTAGTTGCC AGCCATCTGT TGTTTGCCCC TCCCCCGTGC

1159 CTTCCCTTGAC CCTGGAAGGT GCCACTCCCA CTGTCCTTTC CTAATAAAAT GAGGAAATTG

```

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

Continued on next page

Cloning into pcDNA™ 6/His A, B, and C, Continued

Multiple Cloning Site of Version B

Below is the multiple cloning site for pcDNA™6/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 sequence of pcDNA™6/His B is available at www.invitrogen.com or by request from Technical Support (see page 19).**

```

                                     T7 promoter priming site
839  CACTGCTTAC TGGCTTATCG AAATTAATAC GACTCACTAT AGGGAGACCC AAGCTGGCTA

                                     Polyhistidine Region
899  GCGTTTAAAC TTAAGCTTAC C ATG GGG GGT TCT CAT CAT CAT CAT CAT CAT
                                     Met Gly Gly Ser His His His His His His

                                     Xpress™ Epitope
950  GGT ATG GCT AGC ATG ACT GGT GGA CAG CAA ATG GGT CGG GAT CTG TAC
                                     Gly Met Ala Ser Met Thr Gly Gly Gln Gln Met Gly Arg Asp Leu Tyr

998  GAC GAT GAC GAT AAG GTA CCT AAG GAT CCA GTG TGG TGG AAT TCT GCA
     Asp Asp Asp Asp Lys Val Pro Lys Asp Pro Val Trp Trp Asn Ser Ala
     Enterokinase recognition site ▲ EK cleavage site

     EcoR V      BstX I*   Not I     Xho I     Xba I           Apa I
1046  GAT ATC CAG CAC AGT GGC GGC CGC TCG AGT CTA GAG GGC CCG TTT AAA
     Asp Ile Gln His Ser Gly Gly Arg Ser Ser Leu Glu Gly Pro Phe Lys

                                     BGH reverse priming site
1094  CCC GCT GAT CAG CCT CGA CTG TGC CTT CTA GTT GCC AGC CAT CTG TTG
     Pro Ala Asp Gln Pro Arg Leu Cys Leu Leu Val Ala Ser His Leu Leu

1142  TTT GCC CCT CCC CCG TGC CTT CCT TGA CCCTGGAAGG TGCCACTCCC
     Phe Ala Pro Pro Pro Cys Leu Pro ***

```

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

Continued on next page

Cloning into pcDNA™ 6/His A, B, and C, Continued

Multiple Cloning Site of Version C

Below is the multiple cloning site for pcDNA™6/His C. Restriction sites are labeled to indicate the cleavage site. The multiple cloning site has been confirmed by sequencing and functional testing. **The sequence of pcDNA™6/His C is available at www.invitrogen.com or by request from Technical Support (see page 19).**

```

      T7 promoter priming site
839  CACTGCTTAC TGGCTTATCG AAATTAATAC GACTCACTAT AGGGAGACCC AAGCTGGCTA

      Polyhistidine Region
899  GCGTTTAAAC TTAAGCTTAC C ATG GGG GGT TCT CAT CAT CAT CAT CAT CAT
      Met Gly Gly Ser His His His His His His His

      Xpress™ Epitope
950  GGT ATG GCT AGC ATG ACT GGT GGA CAG CAA ATG GGT CGG GAT CTG TAC
      Gly Met Ala Ser Met Thr Gly Gly Gln Gln Met Gly Arg Asp Leu Tyr

      Asp718 I Kpn I BamHI BstX I* EcoR I Pst I
998  GAC GAT GAC GAT AAG GTA CCA GGA TCC AGT GTG GTG GAA TTC TGC AGA
      Asp Asp Asp Asp Lys Val Pro Gly Ser Ser Val Val Glu Phe Cys Arg
      Enterokinase recognition site ▲ EK cleavage site

      EcoR V BstX I* Not I Xho I Xba I Apa I
1046 TAT CCA GCA CAG TGG CGG CCG CTC GAG TCT AGA GGG CCC GTT TAA
      Tyr Pro Ala Gln Trp Arg Pro Leu Glu Ser Arg Gly Pro Val ***

      BGH reverse priming site
1091 ACCCGCTGAT CAGCCTCGAC TGTGCCTTCT AGTTGCCAGC CATCTGTTGT TTGCCCTCC

1151 CCCGTGCCTT CCTTGACCCT GGAAGGTGCC ACTCCCCTG TCCTTTCCTA ATAAAATGAG
  
```

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

Continued on next page

Cloning into pcDNA™ 6/His A, B, and C, Continued

E. coli Transformation

Transform the ligation mixtures into a competent *recA*, *endA* *E. coli* strain (e.g. TOP10F', INVαF') and select on LB plates containing 50–100 µg/mL ampicillin or 50 µg/mL blasticidin. Select 10–20 clones and analyze for the presence and orientation of your insert.



Sequence your construct with the T7 Forward and BGH Reverse primers (see page 17 for ordering information) to confirm that your gene is fused in-frame with the Xpress™ epitope and the N-terminal polyhistidine tag. Refer to the diagrams on pages 3–5 for the sequences and locations of the priming sites. For your convenience, Invitrogen offers a custom primer service. For more information, visit www.invitrogen.com or call Technical Support (see page 19).

Primer	Sequence
T7 Forward	5'-TAATACGACTCACTATAGGG-3'
BGH Reverse	5'-TAGAAGGCACAGTCGAGG-3'

Preparing a Glycerol Stock

After identifying the correct clone, 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 on an LB plate containing 50 µg/mL ampicillin or 50 µg/mL blasticidin. Incubate the plate at 37°C overnight.
2. Isolate a single colony and inoculate into 1–2 mL of LB containing 50 µg/mL ampicillin (or 50 µg/mL blasticidin).
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.

Applying Selective Pressure

Take some (if not all) of the following precautions to prevent your clone from being “overrun” by background contaminants:

- **Use carbenicillin instead of ampicillin.** Carbenicillin is more stable than ampicillin, and allows for a longer period of selective pressure.
- **Increase the antibiotic concentration.** More antibiotic means that your clones will not be overwhelmed by β-lactamase buildup.
- **Periodically refresh plate media.** If you suspect that tubes/plates may be beginning to fail, spin them down, remove the old media, and replenish the wells with fresh LB media plus glycerol and antibiotic.

Streak clones on selective (preferably carbenicillin) LB agar plates. After about 12 hours, isolate colonies for downstream usage. This will isolate your desired clones from potential background contaminants.

Transfection and Analysis

Introduction

Once you have confirmed that your construct is in the correct orientation and fused in-frame with the N-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 lipid complexing, decreasing transfection efficiency. We recommend isolating DNA using the PureLink™ HiPure Miniprep Kit, the PureLink™ HiPure Midiprep Kit (see page 17 for ordering information), or CsCl gradient centrifugation.

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. Precisely follow 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* (Reference section, page 22).

For high efficiency transfection in a broad range of mammalian cells, use Lipofectamine™ 2000 Reagent available from Invitrogen (see page 17). For more information on Lipofectamine™ 2000 and other transfection reagents, visit our web site at www.invitrogen.com or contact Technical Support (see page 19).

Positive Control

pcDNA™6/His/*lacZ* is provided as a positive control vector for mammalian cell transfection and expression (see page 16) 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.

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 for ordering information).

Continued on next page

Transfection and Analysis, Continued

Detecting Fusion Proteins

The Anti-Xpress™ Antibody is available from Invitrogen (see page 17) and can be used to detect expression of your fusion protein from pcDNA™6/His. To detect the fusion protein by western blot, prepare a cell lysate from transfected cells. We recommend that you perform a time course to optimize expression of the fusion protein (*e.g.* 24, 48, 72 hours, etc. after transfection). To lyse cells:

1. Wash cell monolayers (~10⁶ cells) once with phosphate-buffered saline (PBS).
2. Scrape cells into 1 mL PBS and pellet the cells at 1,500 × *g* for 5 minutes.
3. Resuspend in 50 µL Cell Lysis Buffer (see below) or other suitable lysis buffer.
4. Incubate cell suspension at 37°C for 10 minutes to lyse the cells.
5. Centrifuge at 10,000 × *g* for 10 minutes to pellet nuclei and transfer the post-nuclear lysate to a new 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 µg of lysate onto an SDS-PAGE gel and electrophorese. Use the appropriate percentage of acrylamide to resolve your fusion protein.

Cell Lysis Buffer

50 mM Tris, pH 7.8
150 mM NaCl
1% Nonidet P-40

1. Prepare the solution from the following common stock solutions.

For 100 mL, combine:

Stock Solution	Volume
1 M Tris base	5 mL
5 M NaCl	3 mL
Nonidet P-40	1 mL

2. Bring the volume to 90 mL with deionized water and adjust the pH to 7.8 with HCl.
3. Bring the volume to 100 mL. Store at room temperature.

Note: Protease inhibitors may be added at the following concentrations:

1 mM PMSF
1 µg/mL Pepstatin
1 µg/mL Leupeptin

Continued on next page

Transfection and Analysis, Continued



Note

The N-terminal peptide containing the Xpress™ epitope and the polyhistidine tag will add approximately 3.4 kDa to your protein.

Purification

You will need 5×10^6 to 1×10^7 of **transfected** cells for purification of your protein on a 2 mL ProBond™ column (or other metal-chelating column). Refer to the manufacturer's instructions before attempting to purify your fusion protein. To prepare cells for lysis, see the protocol on page 13.

Creating Stable Cell Lines

Introduction

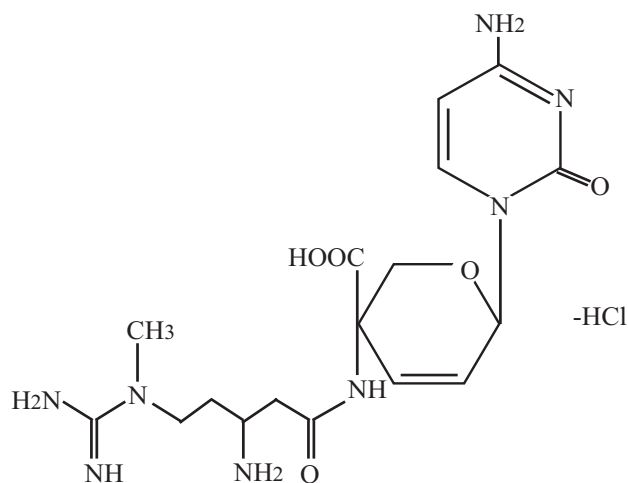
The pcDNA™6/His vectors contain the blasticidin resistance gene for selection of stable cell lines using blasticidin. Test the sensitivity of your mammalian host cell to blasticidin as natural resistance varies among cell lines. General information and guidelines are provided below for your convenience.

Blasticidin

Blasticidin S HCl is a nucleoside antibiotic isolated from *Streptomyces griseochromogenes* which inhibits protein synthesis in both prokaryotic and eukaryotic cells (Takeuchi *et al.*, 1958; Yamaguchi *et al.*, 1965). Resistance is conferred by expression of either one of two blasticidin S deaminase genes: *bsd* from *Aspergillus terreus* (Kimura *et al.*, 1994) or *bsr* from *Bacillus cereus* (Izumi *et al.*, 1991). These deaminases convert blasticidin S to a non-toxic deaminohydroxy derivative (Izumi *et al.*, 1991).

Molecular Weight, Formula, and Structure

The formula for blasticidin is C₁₇H₂₆N₈O₅-HCl, and the molecular weight is 458.9. The diagram below shows the structure of blasticidin.



Handling Blasticidin

Always wear gloves, mask, goggles, and protective clothing (e.g. a laboratory coat) when handling blasticidin. Weigh out blasticidin and prepare solutions in a hood.

Continued on next page

Creating Stable Cell Lines, Continued

Preparing and Storing Stock Solutions

Blasticidin may be obtained from Invitrogen in 50 mg aliquots (see page 17 for ordering information). Prepare stock solutions of 5 to 10 mg/mL of blasticidin in water as described below.

- Dissolve blasticidin in sterile water and filter-sterilize the solution.
 - Aliquot in small volumes suitable for one time use (see last point, below) and freeze at -20°C for long-term storage or store at 4°C for short term storage.
 - Aqueous stock solutions are stable for 1–2 weeks at 4°C and 6–8 weeks at -20°C .
 - The pH of the aqueous solution should not exceed 7 to prevent inactivation of blasticidin.
 - Do not subject stock solutions to freeze/thaw cycles (**do not store in a frost-free freezer**).
 - Upon thawing, use what you need and discard the unused portion.
-

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 likelihood that the vector will 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.** Be sure that your insert does not contain the restriction enzyme site you wish to use to linearize your vector.

Enzyme	Location	Supplier
<i>Bgl</i> II	Upstream of CMV promoter	Invitrogen*
<i>Mfe</i> I	Upstream of CMV promoter	New England Biolabs
<i>Bst</i> 1107 I	End of SV40 poly A	Roche
<i>Eam</i> 1105 I	Ampicillin gene	Roche
<i>Fsp</i> I	Ampicillin gene	New England Biolabs
<i>Sca</i> I	Ampicillin gene	Invitrogen*
<i>Ssp</i> I	Backbone	Invitrogen*

* see page 17 for ordering information

Continued on next page

Creating Stable Cell Lines, Continued

Selection in Mammalian Cell Lines

To generate a stable cell line expressing your protein, first determine the minimum concentration of blasticidin required to kill your untransfected host cell line. Typically, concentrations between 2 and 10 $\mu\text{g}/\text{mL}$ blasticidin are sufficient to kill the untransfected host cell line. Test a range of concentrations (see below) to determine the minimum concentration necessary for your cell line.

1. Seed cells (2×10^5 cells/60 mm plate) for each time point and allow cells to adhere overnight.
 2. The next day, substitute culture medium with medium containing varying concentrations of blasticidin (e.g., 0, 1, 3, 5, 7.5, and 10 $\mu\text{g}/\text{mL}$).
 3. Replenish the selective medium every 3–4 days. Cells sensitive to blasticidin will detach from the plate. Dead cells will accumulate in the medium.
 4. Count the number of viable cells at regular intervals to determine the appropriate concentration of blasticidin that prevents growth.
-

Selecting Stable Integrants

Once the appropriate blasticidin concentration is determined, 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 blasticidin 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 blasticidin-resistant colonies are detected.
 5. Pick and expand colonies.
-

Continued on next page

Creating Stable Cell Lines, Continued

Preparing Cells for Lysis

Use the procedure below to prepare cells for lysis prior to purifying your protein on ProBond™ (see page 18). You will need 5×10^6 to 1×10^7 cells for purification of your protein on a 2 mL ProBond™ column (see the ProBond™ Protein Purification 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 $250 \times g$ for 5 minutes. Resuspend the cell pellet in PBS.
 6. Centrifuge the cells at $250 \times g$ for 5 minutes. You may lyse the cells immediately or freeze in liquid nitrogen and store at -80°C until needed.
-

Lysing Cells

If you are using ProBond™ resin, refer to the ProBond™ Protein Purification manual for details about sample preparation for chromatography.

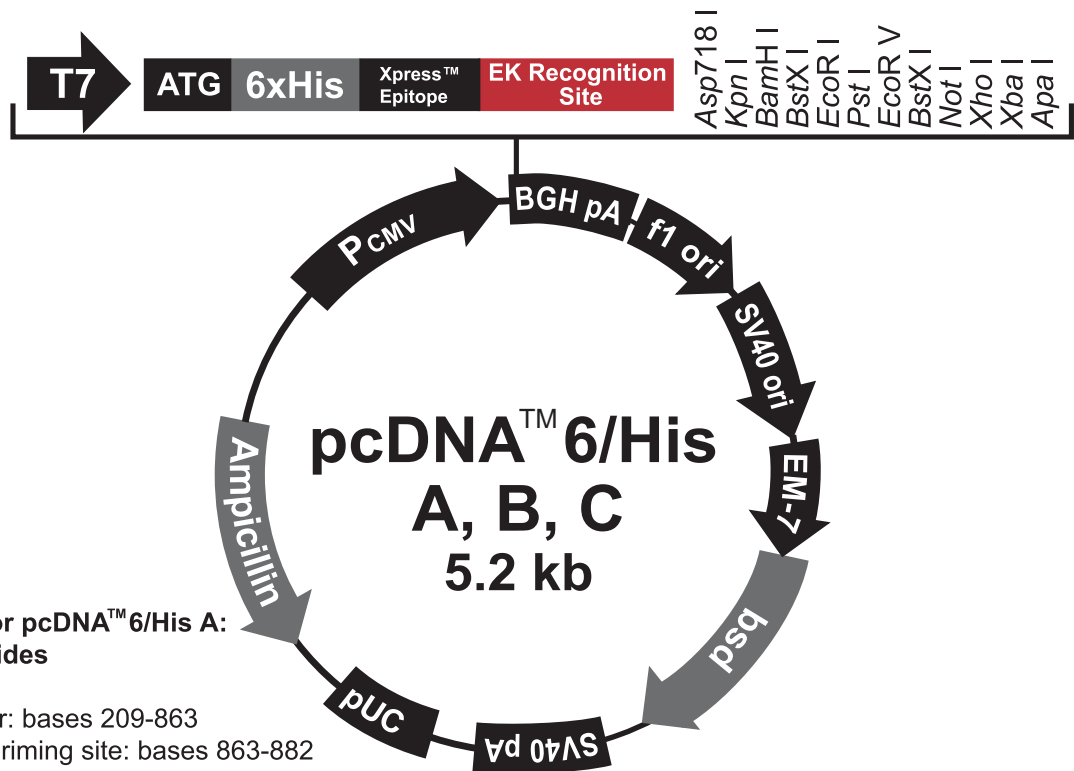
If you are using another metal-chelating resin, refer to the manufacturer's instructions for recommendations on sample preparation.

Appendix

pcDNA™ 6/His Vector

Map of pcDNA™ 6/His

The figure below summarizes the features of the pcDNA™ 6/His vectors. The sequences for pcDNA™ 6/His A, B, and C are available at www.invitrogen.com or by request from Technical Support (see page 19).



Comments for pcDNA™ 6/His A: 5149 nucleotides

- CMV promoter: bases 209-863
- T7 promoter/priming site: bases 863-882
- ATG initiation codon: bases 920-922
- Polyhistidine region: bases 932-949
- Xpress™ epitope: bases 989-1012
- Enterokinase recognition site: bases 998-1012
- Multiple cloning site: bases 1012-1085
- BGH reverse priming site: bases 1105-1122
- BGH polyadenylation signal: bases 1108-1335
- f1 origin: bases 1389-1809
- SV40 promoter and origin: bases 1837-2144
- EM-7 promoter: bases 2192-2247
- Blasticidin resistance gene: bases 2266-2664
- SV40 polyadenylation signal: bases 2822-2952
- pUC origin: bases 3335-4008
- bla* promoter bases: 5014-5112 (C)
- Ampicillin (*bla*) resistance gene: bases 4153-5013 (C)
- (C) = Complementary strand

Continued on next page

pcDNA™ 6/His Vector, Continued

Features of pcDNA™ 6/His pcDNA™ 6/His A (5,150 bp), pcDNA™ 6/His B (5,151 bp), and pcDNA™ 6/His C (5,149 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 <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.
N-terminal polyhistidine tag	Permits purification of recombinant protein on metal-chelating resin such as ProBond™.
Xpress™ epitope tag	Allows for the detection of an 8 amino acid epitope (Asp-Leu-Tyr-Asp-Asp-Asp-Lys) on the recombinant protein with the Anti-Xpress™ Antibody.
Enterokinase cleavage site	Allows for the removal of the N-terminal polyhistidine tag from the recombinant protein using an enterokinase such as EKMax™ Enterokinase (see page 17).
Multiple cloning site in three reading frames	Allows for the insertion of your gene and facilitates cloning in-frame with the Xpress™ epitope and N-terminal polyhistidine tag.
BGH reverse priming site	Permits sequencing through the insert.
Bovine growth hormone (BGH) polyadenylation signal	Efficient transcription termination and polyadenylation of mRNA (Goodwin & Rottman, 1992).
f1 origin	Allows rescue of single-stranded DNA.
SV40 early promoter and origin	Allows for efficient, high-level expression of the blasticidin 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 blasticidin resistance gene in <i>E. coli</i> .
Blasticidin resistance gene (<i>bsd</i>)	Selection of transformants in <i>E. coli</i> and stable transfectants in mammalian cells (Kimura <i>et al.</i> , 1994).
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 transformants in <i>E. coli</i> .

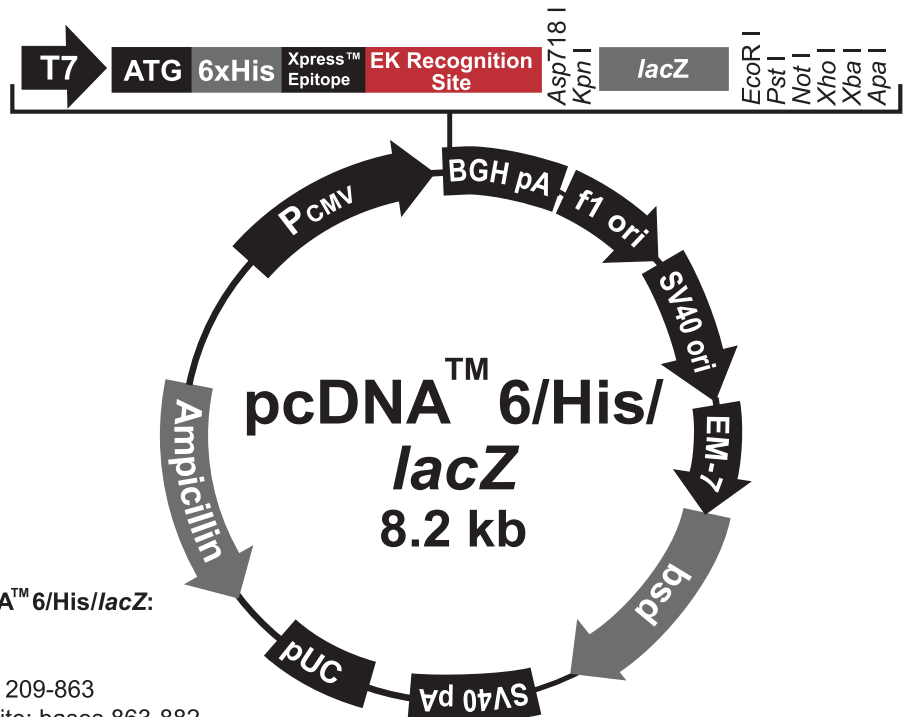
pcDNA™ 6/His/lacZ

Description

pcDNA™ 6/His/lacZ is a 8,213 bp control vector containing the gene for β-galactosidase. This vector was constructed by ligating a 3,082 bp *Kpn* I-*Eco*R I fragment containing the *lacZ* gene from pcDNA4™/His/lacZ to a 5,131 bp *Kpn* I-*Eco*R I fragment containing the CMV promoter, Xpress™ epitope, and polyhistidine tag from pcDNA™ 6/His B.

Map of Control Vector

The figure below summarizes the features of the pcDNA™ 6/His/lacZ vector. The nucleotide sequence for pcDNA™ 6/His/lacZ is available at www.invitrogen.com or by request from Technical Support (see page 19).



Comments for pcDNA™ 6/His/lacZ: 8212 nucleotides

- CMV promoter: bases 209-863
- T7 promoter/priming site: bases 863-882
- ATG initiation codon: bases 920-922
- Polyhistidine region: bases 932-949
- Xpress™ epitope: bases 989-1012
- Enterokinase recognition site: bases 998-1012
- LacZ* ORF: bases 1037-4087
- BGH reverse priming site: bases 4168-4185
- BGH polyadenylation signal: bases 4171-4398
- f1 origin: bases 4444-4872
- SV40 promoter and origin: bases 4900-5207
- EM-7 promoter: bases 5255-5310
- Blasticidin resistance gene: bases 5329-5727
- SV40 polyadenylation signal: bases 5885-6015
- pUC origin: bases 6398-7071
- bla* promoter bases: 8077-8175 (C)
- Ampicillin (*bla*) resistance gene: bases 7216-8076 (C)
- (C)= Complementary strand

Accessory Products

Additional Products

The following additional products may be used with the pcDNATM6/His vectors. For more information, visit www.invitrogen.com or contact Technical Support (see page 19).

Item	Quantity	Cat. no.
Electrocomp TM Kit (TOP10F')	2 × 20 reactions	C665-11
	6 × 20 reactions	C665-24
One Shot [®] TOP10F' Chemically Competent <i>E. coli</i>	20 × 50 µL	C3030-03
One Shot [®] INVαF' Chemically Competent <i>E. coli</i>	20 × 50 µL	C2020-03
	40 × 50 µL	C2020-06
Ampicillin	200 mg	11593-027
Blasticidin	50 mg	R210-01
Carbenicillin	5 g	10177-012
T7 promoter primer	2 µg	N560-02
BGH Reverse primer	2 µg	N575-02
PureLink TM HiPure Plasmid Miniprep Kit	100 preps	K2100-03
PureLink TM HiPure Plasmid Midiprep Kit	25 preps	K2100-04
Lipofectamine TM 2000 Reagent	1.5 mL	11668-019
β-Gal Assay Kit	1 kit	K1455-01
β-Gal Staining Kit	1 kit	K1465-01
Anti-Xpress TM Antibody	50 µL	R910-25
EKMax TM Enterokinase	250 units	E180-01

Restriction Enzymes

For your convenience, Invitrogen offers an extensive selection of restriction enzymes, including the following:

- *Bgl* II
- *Ssp* I
- *Sca* I

Visit www.invitrogen.com for more details.

Continued on next page

Accessory Products, Continued

ProBond™ Resin

Ordering information for ProBond™ resin is provided below.

Item	Quantity	Cat. no.
ProBond™ Purification System	12 mL precharged ProBond™ resin, 6 columns, and buffers for native and denaturing purification	K850-01
ProBond™ Resin	50 mL	R801-01
	150 mL	R801-15

Other Mammalian Expression Vectors

Invitrogen offers a wide variety of mammalian expression vectors utilizing the CMV or EF-1 α promoters. Vectors are available with the Xpress™ (N-terminal), *c-myc* (C-terminal), V5 (C-terminal), or polyhistidine epitopes for detection and either the neomycin, blasticidin, or Zeocin™ resistance genes. All vectors utilize the polyhistidine tag for purification. For more information on the mammalian expression vectors available, visit www.invitrogen.com or contact Technical Support (see page 19).

Technical Support

Web Resources



Visit the Invitrogen website at www.invitrogen.com 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. 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.

Continued on next page

Purchaser Notification, Continued

**Limited Use Label
License No. 22:
Vectors and
Clones Containing
Sequences
Coding for
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.

**Limited Use Label
License No. 51:
Blasticidin and the
Blasticidin
Selection Marker**

Blasticidin and the blasticidin resistance gene (bsd) are the subject of U.S. Patent No. 5,527,701 sold under patent license for research purposes only. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500.

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*, Greene Publishing Associates and Wiley-Interscience, New York
- 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
- 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
- Izumi, M., Miyazawa, H., Kamakura, T., Yamaguchi, I., Endo, T., and Hanaoka, F. (1991) Blastocidin S-Resistance Gene (*bsr*): A Novel Selectable Marker for Mammalian Cells. *Exp. Cell Res.* 197, 229-233
- Kimura, M., Takatsuki, A., and Yamaguchi, I. (1994) Blastocidin S Deaminase Gene from *Aspergillus terreus* (*BSD*): A New Drug Resistance Gene for Transfection of Mammalian Cells. *Biochim. Biophys. ACTA* 1219, 653-659
- 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
- Kozak, M. (1991) An Analysis of Vertebrate mRNA Sequences: Intimations of Translational Control. *J. Cell Biology* 115, 887-903
- Miller, J. H. (1972) *Experiments in Molecular Genetics*, Cold Spring Harbor Laboratory, Cold Spring Harbor, New York
- 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 Ed., Cold Spring Harbor Laboratory Press, Plainview, New York
- Takeuchi, S., Hirayama, K., Ueda, K., Sakai, H., and Yonehara, H. (1958) Blastocidin S, A New Antibiotic. *The Journal of Antibiotics, Series A* 11, 1-5
- Yamaguchi, H., Yamamoto, C., and Tanaka, N. (1965) Inhibition of Protein Synthesis by Blastocidin S. I. Studies with Cell-free Systems from Bacterial and Mammalian Cells. *J. Biochem (Tokyo)* 57, 667-677

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

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



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