

USER GUIDE

invitrogen™
by *life* technologies™

pMT/V5-His A, B, and C

Catalog number V4120-20

Revision date 20 February 2012

Publication Part number 28-0177

MAN0000658

**For Research Use Only. Not intended for any animal or human
therapeutic or diagnostic use.**

life
technologies™

Contents

Kit Contents and Storage.....	iv
Methods	1
Product Overview	1
Cloning into pMT/V5-His A, B, and C	3
Transfection and Analysis.....	9
Appendix.....	10
pMT/V5-His Vector.....	10
pMT/V5-His/ <i>lacZ</i> Vector	12
Accessory Products	13
Technical Support.....	15
Purchaser Notification.....	16
References.....	17

Kit Contents and Storage

Kit Contents

20 µg each of pMT/V5-His A, B, and C are supplied at 0.5 µg/µL in 10 mM Tris-HCl, 1 mM EDTA, pH 8.0 in a total volume of 40 µL.

20 µg of pMT/V5-His/*lacZ* is supplied at 0.5 µg/µL in 10 mM Tris-HCl, 1 mM EDTA, pH 8.0 in a total volume of 40 µL.

Shipping/Storage

Vectors are shipped at room temperature. Upon receipt, store at -30°C to -10°C .

Product Use

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

Methods

Product Overview

Introduction

pMT/V5-His is a 3.5 kb expression vector designed for use with the *Drosophila* Inducible Expression System (DES[®]; Catalog nos. K4120-01, K5120-01). Upon transfection, the vector allows transient, inducible expression of your protein of interest in *Drosophila* cells. When cotransfected with the selection vectors, pCoHygro or pCoBlast, included with the appropriate DES[®] Inducible Kit, pMT/V5-His allows selection of stable cell lines exhibiting inducible expression of the protein of interest. The pMT/V5-His vector contains the following elements:

- The *Drosophila* metallothionein (MT) promoter for high-level, metal-inducible expression of the gene of interest in S2 cells (Angelichio *et al.*, 1991; Bunch *et al.*, 1988; Maroni *et al.*, 1986; Olsen, 1992)
- Multiple cloning site to facilitate cloning the gene of interest
- C-terminal peptide containing the V5 epitope and polyhistidine (6xHis) tag for detection and purification of your protein of interest (if desired)
- Three reading frames to facilitate in-frame cloning with the C-terminal peptide
- Ampicillin resistance gene for selection of transformants in *E. coli*

The control plasmid, pMT/V5-His/*lacZ*, is included for use as a positive control for transfection and expression.

For more information about the DES[®] Inducible Kits, pCoHygro, and pCoBlast, refer to the *Drosophila* Expression System manual. The *Drosophila* Expression System manual is supplied with each DES[®] Inducible Kit, but is also available from www.lifetechnologies.com or by contacting Technical Support (see page 15).

Description of MT Promoter

The *Drosophila* MT promoter allows high-level, inducible expression of the gene of interest in *Drosophila* S2 (or D.Mel-2) cells. When used to express heterologous proteins, the promoter is extremely efficient and tightly regulated, even at high copy number (Johansen *et al.*, 1989). The MT promoter is well characterized (Angelichio *et al.*, 1991; Bunch *et al.*, 1988; Maroni *et al.*, 1986; Olsen, 1992), with regulatory elements and the start of transcription well defined.

The MT promoter is inducible by addition of copper sulfate or cadmium chloride to the culture medium (Bunch *et al.*, 1988). Copper sulfate is generally the preferred inducer due to its reduced toxicity as compared to cadmium. While cadmium is an effective inducer, it also induces a heat-shock response in S2 cells.

Continued on next page

Product Overview, Continued

Experimental Outline

The table below describes the general steps needed to clone and express your gene of interest. For more details, refer to the manual and pages indicated.

Step	Action	Source
1	Develop a cloning strategy to ligate your gene of interest into pMT/V5-His A, B, or C in frame with the C-terminal peptide encoding the V5 epitope and the polyhistidine tag (if desired).	Pages 3–7, this manual
2	Transform your ligation reactions into a <i>recA</i> , <i>endA</i> <i>E. coli</i> strain (e.g. TOP10). Select on LB agar plates containing 50–100 µg/mL ampicillin.	Page 8, this manual
3	Analyze your transformants for the presence of insert.	Page 8, this manual
4	Select a transformant with the correct restriction pattern and sequence it to confirm that your gene is cloned in frame with the C-terminal peptide.	Page 8, this manual
5	Transfect your pMT/V5-His construct into S2 cells and induce expression of the gene of interest with copper sulfate.	Page 9, this manual and DES [®] manual
6	Assay for transient expression of your recombinant protein.	Page 9, this manual and DES [®] manual
7	To generate stable cell lines, cotransfect your pMT/V5-His construct and pCoHygro or pCoBlast into S2 cells and select for hygromycin or blasticidin resistant clones, as appropriate.	DES [®] manual
8	Scale up expression for purification.	DES [®] manual
9	Purify your recombinant protein by chromatography on metal-chelating resin (i.e. ProBond [™]).	DES [®] manual

Cloning into pMT/V5-His A, B, and C

Introduction

Diagrams are provided on pages 5–7 to help you clone your gene of interest into pMT/V5-His. General considerations for cloning and transformation are discussed below.

General Molecular Biology Techniques

For help with DNA ligations, *E. coli* transformations, restriction enzyme analysis, 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 and maintenance of the pMT/V5-His vectors including TOP10, DH5 α TM-T1^R, and JM109. We recommend that you propagate the vectors in *E. coli* strains that are recombination deficient (*recA*) and endonuclease A deficient (*endA*).

For your convenience, the TOP10 and DH5 α TM-T1^R strains are available as chemically competent cells (see page 13 for ordering information). TOP10 cells are also available as electrocompetent cells.

Transformation Method

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

Maintenance of Plasmids

To propagate and maintain the pMT/V5-His and pMT/V5-His/*lacZ* vectors, we recommend that you use the supplied 0.5 $\mu\text{g}/\mu\text{L}$ stock solution in TE buffer to transform a *recA*, *endA* *E. coli* strain like TOP10, DH5 α TM-T1^R, JM109, or equivalent. Select transformants on LB agar plates containing 50 to 100 $\mu\text{g}/\text{mL}$ ampicillin. Be sure to prepare a glycerol stock of each plasmid for long-term storage (see page 8).

Continued on next page

Cloning into pMT/V5-His A, B, and C, Continued

Cloning Considerations

Consider the following points when designing a strategy to clone your gene of interest into pMT/V5-His.

- Your insert should contain a Kozak translation initiation sequence with an ATG initiation codon 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 G or 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.

(G/A)NNATGG

- It is possible to clone and express a secreted protein in pMT/V5-His if your protein includes a native signal sequence. If your protein does not have a secretion signal, you may wish to consider using the pMT/BiP/V5-His vector for secreted expression. For more information, refer to the DES[®] manual.
- If you wish to use the V5 epitope and the polyhistidine (6xHis) tag for detection and purification of your recombinant protein, you must clone your gene in frame with the C-terminal peptide. The vector is supplied in three reading frames to facilitate cloning. Refer to the diagrams on pages 5-7 to develop a cloning strategy. **Be sure that your gene does not contain a stop codon upstream of the C-terminal peptide.**
- If you do not wish to include the C-terminal peptide, include the native stop codon for your gene of interest.

Continued on next page

Cloning into pMT/V5-His A, B, and C, Continued

Multiple Cloning Site of pMT/V5-His A

Below is the multiple cloning site for pMT/V5-His A. The metal regulatory regions are marked as per Maroni, *et al.*, 1986. The start of transcription is at nucleotide 778. Restriction sites are labeled to indicate the actual cleavage site. The boxed nucleotides indicate the variable region. The multiple cloning site has been confirmed by sequencing and functional testing. The nucleotide sequence of pMT/V5-His A is available from www.lifetechnologies.com or from Technical Support (see page 15). For a map and a description of the features of pMT/V5-His A, refer to pages 10–11.

```

411  ┌ 5' end of metallothionein promoter
      CGTTGCAGGA CAGGATGTGG TGCCCGATGT GACTAGCTCT TTGCTGCAGG CCGTCCTATC
      └──────────────────────────────────────────────────────────────────────────────────┘

      Metal regulatory region
471  ┌──────────────────────────────────────────────────────────────────────────────────┘
      CTCTGGTTCC GATAAGAGAC CCAGAACTCC GGCCCCCCAC CGCCCACCGC CACCCCCATA

831  ┌──────────────────────────────────────────────────────────────────────────────────┘
      CATATGTGGT ACGCAAGTAA GAGTGCCTGC GCATGCCCCA TGTGCCCCAC CAAGAGTTTT
      ┌──────────────────────────────────────────────────────────────────────────────────┘
      region
591  GCATCCCATA CAAGTCCCCA AAGTGGAGAA CCGAACCAAT TCTTCGCGGG CAGAACAAAA

      Metal regulatory regions
651  ┌──────────────────────────────────────────────────────────────────────────────────┘
      GCTTCTGCAC ACGTCTCCAC TCGAATTTGG AGCCGGCCGG CGTGTGCAAA AGAGGTGAAT
      ┌──────────────────────────────────────────────────────────────────────────────────┘
      Metal regulatory region
      TATA box
711  CGAACGAAAG ACCCGTGTGT AAAGCCGCGT TTCCAAAATG TATAAAACCG AGAGCATCTG
      ┌──────────────────────────────────────────────────────────────────────────────────┘
      Metal regulatory regions
      ┌──────────────────────────────────────────────────────────────────────────────────┘
      Start of transcription
771  GCCAATGTGC ATCAGTTGTG GTCAGCAGCA AAATCAAGTG AATCATCTCA GTGCAACTAA

831  Xba I#           Kpn I Spe I           BstX I* EcoR I           EcoR V
      A G G G G G G G A T C TAGATCGGGGT ACC TAC TAG TCC AGT GTG GTG GAA TTC TGC AGA TAT
      Thr Tyr *** Ser Ser Val Val Glu Phe Cys Arg Tyr

888  BstX I* Not I   Xho I   Xba I           Apa I BstB I           V5 epitope
      CCA GCA CAG TGG CGG CCG CTC GAG TCT AGA GGG CCC TTC GAA GGT AAG CCT ATC
      Pro Ala Gln Trp Arg Pro Leu Glu Ser Arg Gly Pro Phe Glu Gly Lys Pro Ile

942  Age I           Polyhistidine region
      CCT AAC CCT CTC CTC GGT CTC GAT TCT ACG CGT ACC GGT CAT CAT CAC CAT CAC
      Pro Asn Pro Leu Leu Gly Leu Asp Ser Thr Arg Thr Gly His His His His His

996  Pme I           BGH Reverse priming site
      CAT TGA GTTTAAACC CGCTGATCAG CCTCGACTGT GCCTTCTAAG ATCCAGACAT GATAAGATAC
      His ***

1061  ATTGATGAGT TTGACAAAC CACAAC TAGA ATGCAGTGAA AAAAATGCTT TATTTGTGAA

1121  SV40 late polyadenylation signal
      A T T T G T G A T G C T A T T G C T T T A T T T G T A A C C A T T A T A A G C T G C A A T A A A C A A G T T A A C A A C
  
```

This Xba I site is *dam* methylated in certain strains of *E. coli*.

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

Continued on next page

Cloning into pMT/V5-His A, B, and C, Continued

Multiple Cloning Site of pMT/V5-His B

Below is the multiple cloning site for pMT/V5-His B. The metal regulatory regions are marked as per Maroni, *et al.*, 1986. The start of transcription is at nucleotide 778. Restriction sites are labeled to indicate the actual cleavage site. The boxed nucleotides indicate the variable region. The multiple cloning site has been confirmed by sequencing and functional testing. The nucleotide sequence of pMT/V5-His B is available from www.lifetechnologies.com or from Technical Support (see page 15). For a map and a description of the features of pMT/V5-His B, refer to pages 10–11.

```

411      ┌ 5' end of metallothionein promoter
CGTTGCAGGA CAGGATGTGG TGCCCCGATGT GACTAGCTCT TTGCTGCAGG CCGTCCTATC

Metal regulatory region
471      └
CTCTGGTTCC GATAAGAGAC CCAGAACTCC GGCCCCCCAC CGCCCACCGC CACCCCCATA

531      ┌ Metal regulatory
CATATGTGGT ACGCAAGTAA GAGTGCCTGC GCATGCCCCA TGTGCCCCAC CAAGAGTTTT
region
591      └
GCATCCATA CAAGTCCCA AAGTGGAGAA CCGAACCAAT TCTTCGCGGG CAGAACAAAA

Metal regulatory regions
651      └
GCTTCTGCAC ACGTCTCCAC TCGAATTGG AGCCGGCCGG CGTGTGCAAA AGAGGTGAAT
Metal regulatory region
711      ┌ TATA box
CGAACGAAAG ACCCGTGTGT AAAGCCGCGT TTCCAAAATG TATAAAACCG AGAGCATCTG
Metal regulatory regions
771      ┌ Start of transcription
GCCAATGTGC ATCAGTTGTG GTCAGCAGCA AAATCAAGTG AATCATCTCA GTGCAACTAA
MT Forward priming site

831      ┌ Xba I#
AGGGGGGATC TAGATCGGGG TA CCT ACT AGT CCA GTG TGG TGG AAT TCT GCA GAT ATC
Pro Thr Ser Pro Val Trp Trp Asn Ser Ala Asp Ile

889      ┌ BstX I* Not I Xho I Xba I Apa I Sac II BstB I
CAG CAC AGT GGC GGC CGC TCG AGT CTA GAG GGC CCG CGG TTC GAA GGT AAG CCT
Gln His Ser Gly Gly Arg Ser Ser Leu Glu Gly Pro Arg Phe Glu Gly Lys Pro
V5 epitope
943      ┌ Age I Polyhistidine region
ATC CCT AAC CCT CTC CTC GGT CTC GAT TCT ACG CGT ACC GGT CAT CAT CAC CAT
Ile Pro Asn Pro Leu Leu Gly Leu Asp Ser Thr Arg Thr Gly His His His His

997      ┌ Pme I BGH Reverse priming site
CAC CAT TGA GTTTA AACCCGCTGA TCAGCCTCGA CTGTGCCTTC TAAGATCCAG ACATGATAAG
His His ***

1061      ATACATTGAT GAGTTTGGAC AAACCACAAC TAGAATGCAG TGAAAAAAT GCTTTATTG

1121      ┌ SV40 late polyadenylation signal
TGAAATTTGT GATGCTATTG CTTTATTTGT AACCATTATA AGCTGCAATA AACAAGTTAA

```

This *Xba* I site is *dam* methylated in certain strains of *E. coli*.

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

Continued on next page

Cloning into pMT/V5-His A, B, and C, Continued

Multiple Cloning Site of pMT/V5-His C

Below is the multiple cloning site for pMT/V5-His C. The metal regulatory regions are marked as per Maroni, *et al.*, 1986. The start of transcription is at nucleotide 778. Restriction sites are labeled to indicate the actual cleavage site. The boxed nucleotides indicate the variable region. The multiple cloning site has been confirmed by sequencing and functional testing. The nucleotide sequence of pMT/V5-His C is available from www.lifetechnologies.com or from Technical Support (see page 15). For a map and a description of the features of pMT/V5-His C, refer to pages 10–11.

```

┌ 5' end of metallothionein promoter
411  CGTTGCAGGA CAGGATGTGG TGCCCGATGT GACTAGCTCT TTGCTGCAGG CCGTCCTATC

Metal regulatory region
471  CTCTGGTTCC GATAAGAGAC CCAGAACTCC GGCCCCCACC CGCCCACCGC CACCCCCATA

Metal regulatory
531  CATATGTGGT ACGCAAGTAA GAGTGCCTGC GCATGCCCCA TGTGCCCCAC CAAGAGTTTT
region
591  GCATCCCATA CAAGTCCCCA AAGTGGAGAA CCGAACCAAT TCTTCGCGGG CAGAACAAAA

Metal regulatory regions
651  GCTTCTGCAC ACGTCTCCAC TCGAATTTGG AGCCGGCCGG CGTGTGCAAA AGAGGTGAAT
Metal regulatory region
711  CGAACGAAAG ACCCGTGTGT AAAGCCGCGT TTCCAAAATG TATAAAACCG AGAGCATCTG
TATA box
Metal regulatory regions
771  GCCAATGTGC ATCAGTTGTG GTCAGCAGCA AAATCAAGTG AATCATCTCA GTGCAACTAA
└ Start of transcription
MT Forward priming site

Xba I# Kpn I Spe I BstX I* EcoR I EcoR V
831  AGGGGGGATC TAGATCGGGG TAC CTA CTA GTC CAG TGT GGT GGA ATT CTG CAG ATA
Tyr Leu Leu Val Gln Cys Gly Gly Ile Leu Gln Ile

BstX I* Not I Xho I BstE II BstB I V5 epitope
887  TCC AGC ACA GTG GCG GCC GCT CGA GGT CAC CCA TTC GAA GGT AAG CCT ATC CCT
Ser Ser Thr Val Ala Ala Ala Arg Gly His Pro Phe Glu Gly Lys Pro Ile Pro

Age I Polyhistidine region
941  AAC CCT CTC CTC GGT CTC GAT TCT ACG CGT ACC GGT CAT CAT CAC CAT CAC CAT
Asn Pro Leu Leu Gly Leu Asp Ser Thr Arg Thr Gly His His His His His His

Pme I BGH Reverse priming site
995  TGA GTT TAAACCCGCT GATCAGCCTC GACTGTGCCT TCTAAGATCC AGACATGATA AGATACATTG
***

1061  ATGAGTTTGG ACAAACCACA ACTAGAATGC AGTGAAAAAA ATGCTTTATT TGTGAAATTT

SV40 late polyadenylation signal
1121  GTGATGCTAT TGCTTTATTT GTAACCATTA TAAGCTGCAA TAAACAAGTT AACAACAACAA

```

This *Xba* I site is *dam* methylated in certain strains of *E. coli*.

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

Continued on next page

Cloning into pMT/V5-His A, B, and C, Continued

E. coli **Transformation**

Transform your ligation mixtures into a competent *recA*, *endA* *E. coli* strain (e.g. TOP10, DH5 α [™]-T1[®]) and select on LB agar plates containing 50–100 $\mu\text{g}/\text{mL}$ ampicillin. Select 10–20 clones and analyze for the presence and orientation of your insert.



We recommend that you sequence your construct with the MT Forward and BGH Reverse primers to confirm that your gene is in the correct orientation for expression and is cloned in frame with the C-terminal peptide. The MT Forward and BGH Reverse primers are included in each DES[®] Inducible Kit. Refer to the diagrams on pages 5–7 for the sequences and location of the priming sites.

Note: For your convenience, we offer a custom primer synthesis service. For more information, see www.lifetechnologies.com or call Technical Support (see page 15).

Preparing a **Glycerol Stock**

Once you have identified the correct clone, purify the colony and make a glycerol stock for long-term storage. You should keep a DNA stock of your plasmid at -20°C .

1. Streak the original colony out on an LB plate containing 50 $\mu\text{g}/\text{mL}$ ampicillin. Incubate the plate at 37°C overnight.
 2. Isolate a single colony and inoculate into 1–2 mL of LB containing 50 $\mu\text{g}/\text{mL}$ ampicillin.
 3. Grow the culture to mid-log phase ($\text{OD}_{600} = 0.5\text{--}0.7$).
 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 cloned your gene of interest into pMT/V5-His and have prepared purified plasmid DNA, you are ready to transfect your construct into S2 cells. If you are assaying for transient, inducible expression of your gene of interest, you may transfect your pMT/V5-His construct alone into S2 cells. If you wish to generate stable cell lines, you **must** cotransfect your pMT/V5-His construct with pCoHygro or pCoBlast into S2 cells. Note that the pMT/V5-His vector does not contain a resistance marker for selection in *Drosophila* cells. We recommend that you include the pMT/V5-His/*lacZ* positive control vector and a mock transfection (negative control) in your experiments to evaluate your results. Specific guidelines and protocols for transient transfection and generation of stable cell lines can be found in the DES[®] manual.

Note: The pCoHygro or pCoBlast selection vector is supplied with the appropriate DES[®] Inducible Kit. For more information about each vector, refer to the DES[®] manual.

Plasmid Preparation

Plasmid DNA for transfection into eukaryotic cells must be pure and free from phenol and sodium chloride. Contaminants will kill the cells, decreasing transfection efficiency. We recommend isolating plasmid DNA using the PureLink[®] HiPure MiniPrep Kit (up to 30 µg DNA), the PureLink[®] HiPure MidiPrep Kit (up to 150 µg DNA) (see page 13 for ordering), or CsCl gradient centrifugation.

Positive Control

pMT/V5-His/*lacZ* is provided as a positive control vector for *Drosophila* cell transfection and expression (see page 12 for a map) and may be used to optimize transfection conditions for S2 cells. Transfection of pMT/V5-His/*lacZ* results in induction of β-galactosidase expression upon addition of copper sulfate. A successful transfection will result in β-galactosidase expression that can be easily assayed by staining with X-gal.

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. The β-Gal Assay Kit and the β-Gal Staining Kit are available for fast and easy detection of β-galactosidase expression (see page 13 for ordering information).

Induction of Recombinant Protein Expression

Once you have transfected your pMT/V5-His construct into S2 cells, you will induce expression of recombinant protein using copper sulfate. In general, we recommend that you add copper sulfate directly to the culture medium to a final concentration of 500 µM and incubate the cells for 24 hours to obtain maximal induction of your protein of interest. Refer to the DES[®] manual for more details. Copper sulfate is provided in each DES[®] Inducible Kit.

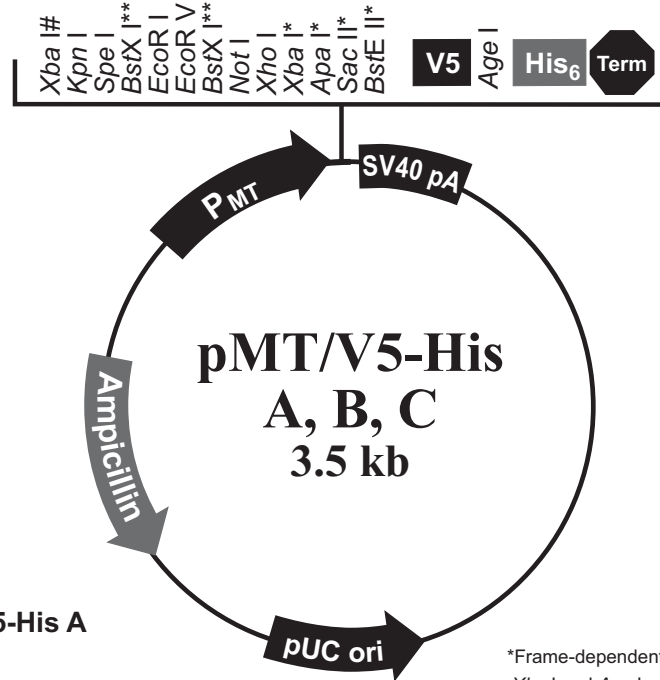
Detection and Purification of Recombinant Fusion Proteins

If you have cloned your gene of interest in frame with the C-terminal peptide containing the V5 epitope and the polyhistidine (6xHis) tag, you may use the Anti-V5 antibodies or Anti-His(C-term) antibodies to detect expression of your recombinant fusion protein by western blot analysis (see page 13 for ordering). The 6xHis tag also allows purification of recombinant protein using metal-chelating resins including ProBond[™]. Refer to the DES[®] manual for more detailed guidelines and instructions to detect and purify your recombinant fusion protein.

Appendix

pMT/V5-His Vector

Map of pMT/V5-His The figure below summarizes the features of the pMT/V5-His A, B, and C vectors. For a more detailed description of each feature, see page 11. The nucleotide sequences of pMT/V5-His A, B, and C are available from www.lifetechnologies.com or from Technical Support (see page 15).



Comments for pMT/V5-His A 3538 nucleotides

Metallothionein promoter: bases 412-778
 Start of transcription: base 778
 MT Forward priming site: bases 814-831
 Multiple cloning site: bases 854-923
 V5 epitope tag: bases 930-971
 Polyhistidine region: bases 981-1001
 BGH Reverse priming site: bases 1021-1038
 SV40 late polyadenylation signal: bases 1163-1168
 pUC origin: bases 1601-2334 (complementary strand)
bla promoter: bases 3340-3438 (complementary strand)
 Ampicillin (*bla*) resistance gene ORF: bases 2479-3339 (complementary strand)

*Frame-dependent variations
Xba I and *Apa* I are found only in versions A and B.
Sac II is found only in version B.
BstE II is found only in version C.
 #This *Xba* I site is *dam* methylated.
 **There are two *BstX* I sites in the polylinker.

Continued on next page

pMT/V5-His Vector, Continued

Features of pMT/V5-His

The features of pMT/V5-His A (3538 bp), pMT/V5-His B (3542 bp), and pMT/V5-His C (3534 bp) are described below. All features have been functionally tested. The multiple cloning site has been tested by restriction enzyme analysis.

Feature	Benefit
<i>Drosophila</i> metallothionein (MT) promoter	Permits high-level, inducible expression of heterologous proteins (Bunch <i>et al.</i> , 1988; Maroni <i>et al.</i> , 1986)
MT Forward priming site	Allows sequencing in the sense orientation
Multiple cloning site	Allows insertion of your gene of interest
V5 epitope (Gly-Lys-Pro-Ile-Pro-Asn-Pro-Leu-Leu-Gly-Leu-Asp-Ser-Thr)	Allows detection of your recombinant protein with the Anti-V5, Anti-V5-HRP, or Anti-V5-AP antibodies (Southern <i>et al.</i> , 1991)
Polyhistidine (6xHis) tag	Permits purification of your recombinant protein on metal-chelating resin such as ProBond™. In addition, the C-terminal 6xHis tag is the epitope for the Anti-His(C-term) Antibody and the Anti-His(C-term)-HRP Antibody (Lindner <i>et al.</i> , 1997) (see page 13 for ordering)
BGH Reverse priming site	Permits sequencing of the non-coding strand
SV40 late polyadenylation signal	Allows efficient transcription termination and polyadenylation of mRNA (Angelichio <i>et al.</i> , 1991)
pUC origin	Allows high-copy number replication and growth in <i>E. coli</i>
<i>bla</i> promoter	Allows expression of the ampicillin (<i>bla</i>) resistance gene
Ampicillin (<i>bla</i>) resistance gene (β-lactamase)	Permits selection of transformants in <i>E. coli</i>

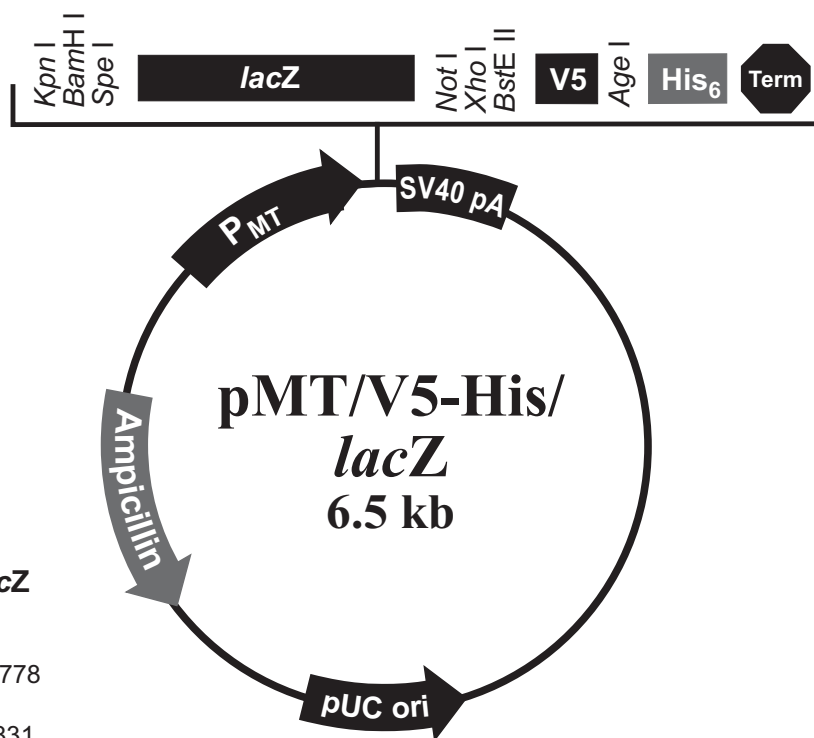
pMT/V5-His/lacZ Vector

Description

pMT/V5-His/lacZ is a 6596 bp control vector expressing β -galactosidase. The plasmid was constructed by digesting pMT/V5-His B with *Kpn* I and *Age* I and ligating a 3.2 kb *Kpn* I-*Age* I fragment containing the *lacZ* gene and the V5 epitope in frame with the polyhistidine tag.

Map of pMT/V5-His/lacZ

The figure below summarizes the features of the pMT/V5-His/lacZ vector. The nucleotide sequence for pMT/V5-His/lacZ is available for downloading from www.lifetechnologies.com or by contacting Technical Support (see page 15).



Comments for pMT/V5-His/lacZ

6596 nucleotides

Metallothionein promoter: bases 412-778

Start of transcription: base 778

MT Forward priming site: bases 814-831

LacZ portion of fusion: bases 904-3987

V5 epitope tag: bases 3988-4029

Polyhistidine region: bases 4039-4059

BGH Reverse priming site: bases 4079-4096

SV40 late polyadenylation signal: bases 4221-4226

pUC origin: bases 4719-5392 (complementary strand)

bla promoter: bases 6398-6496 (complementary strand)

Ampicillin (*bla*) resistance gene ORF: bases 5537-6397 (complementary strand)

Accessory Products

Additional Products

Many of the reagents supplied with the pMT/V5-His vector and the DES[®] Inducible Kits, and other reagents suitable for use with the kits are available separately. Ordering information for these reagents is provided below. For more information, refer to www.lifetechnologies.com or call Technical Support (see page 15).

Item	Quantity	Catalog no.
<i>Drosophila</i> Inducible Expression System (DES)	1 kit	K4120-01
	1 kit	K5120-01
One Shot [®] TOP10 (chemically competent cells)	20 reactions	C4040-03
One Shot [®] TOP10 Electrocomp [™] (electrocompetent cells)	20 reactions	C4040-52
One Shot [®] DH5 α [™] -T1 ^R Max Efficiency [®] Chemically Competent <i>E. coli</i>	20 × 50 μ L	12297-016
β -Gal Assay Kit	80 mL	K1455-01
β -Gal Staining Kit	1 kit	K1465-01
PureLink [®] HiPure MiniPrep Kit	25 preps	K2100-02
PureLink [®] HiPure MidiPrep Kit	25 preps	K2100-04
BGH Reverse Primer	2 μ g, lyophilized in TE	N575-02
Hygromycin B	20 mL	10687-010
Blasticidin S HCl	50 mg	R210-01
Schneider's <i>Drosophila</i> Medium	500 mL	21720-024
Calcium Phosphate Transfection Kit	75 reactions	K2780-01

Continued on next page

Accessory Products, Continued

Detection of Recombinant Proteins

Expression of your recombinant fusion protein can be detected using an antibody to the appropriate epitope. The table below describes the antibodies available for detection of C-terminal fusion proteins expressed using pMT/V5-His. Horseradish peroxidase (HRP)- and alkaline phosphatase (AP)-conjugated antibodies allow one-step detection using colorimetric or chemiluminescent detection methods.

The amount of antibody supplied is sufficient for 25 western blots.

Product	Epitope	Catalog no.
Anti-V5 Antibody	Detects 14 amino acid epitope derived from the P and V proteins of the paramyxovirus, SV5 (Southern <i>et al.</i> , 1991) GKPIP NPLLGLDST	R960-25
Anti-V5-HRP Antibody		R961-25
Anti-V5-AP Antibody		R962-25
Anti-His (C-term) Antibody	Detects the C-terminal polyhistidine (6xHis) tag (requires the free carboxyl group for detection (Lindner <i>et al.</i> , 1997) HHHHHH-COOH	R930-25
Anti-His(C-term)-HRP Antibody		R931-25
Anti-His(C-term)-AP Antibody		R932-25

Purification of Recombinant Protein

The metal binding domain encoded by the polyhistidine tag allows simple, easy purification of your recombinant protein by Immobilized Metal Affinity Chromatography (IMAC) using the ProBond™ Resin. To purify proteins expressed from pMT/V5-His, the Xpress™ Purification System or the ProBond™ resin in bulk are available separately. See the table below for ordering.

Product	Quantity	Catalog no.
ProBond™ Metal-Binding Resin (precharged resin provided as a 50% slurry in 20% ethanol)	50 mL	R801-01
	150 mL	R801-15
Xpress™ Purification System (includes six 2 mL precharged, prepacked ProBond™ resin columns and buffers for native and denaturing purification)	6 purifications	K850-01
Xpress™ Purification System with Anti-V5-HRP Antibody	1 kit	K854-01
Xpress™ Purification System with Anti-His(C-term)-HRP Antibody	1 kit	K853-01
Purification Columns (10 mL polypropylene columns)	50	R640-50

Technical Support

Obtaining support For the latest services and support information for all locations, go to www.lifetechnologies.com/support.
At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support (techsupport@lifetech.com)
- Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents
- Obtain information about customer training
- Download software updates and patches

Safety Data Sheets (SDS) Safety Data Sheets (SDSs) are available at www.lifetechnologies.com/support.

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.lifetechnologies.com/support and search for the Certificate of Analysis by product lot number, which is printed on the box.

Limited Product Warranty Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

Purchaser Notification

**Limited Use Label
License No: 358
Research Use
Only**

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

**Limited Use Label
License:
6X His Tag**

This product is licensed from Hoffmann-La Roche, 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:
Drosophila
Expression
System (DES)**

THIS PRODUCT IS LICENSED TO YOU UNDER PATENT LICENSE FROM SMITHKLINE BEECHAM CORPORATION (SB) FOR YOUR RESEARCH USE ONLY. THIS PRODUCT MAY NOT BE SOLD OR TRANSFERRED TO ANY OTHER PERSON. IN USING THIS PRODUCT, YOU AGREE TO (1) MAKE NO COMMERCIAL USE OF THIS PRODUCT; (2) INFORM INVITROGEN OF ANY INVENTIONS OR DISCOVERIES ARISING FROM YOUR USE OF THIS PRODUCT; AND (3) GRANT TO SB THE FIRST RIGHT TO NEGOTIATE A LICENSE TO SUCH INVENTIONS AND DISCOVERIES. YOU AGREE NOT TO GRANT A LICENSE TO ANY THIRD PARTY ON TERMS MORE FAVORABLE THAN THOSE OFFERED TO SB. REQUESTS FOR LICENSES FOR COMMERCIAL MANUFACTURE OR USE MAY BE OBTAINED DIRECTLY BY CONTACTING: OUTLICENSING@LIFETECH.COM OR OUT LICENSING, LIFE TECHNOLOGIES, 5791 VAN ALLEN WAY, CARLSBAD, CALIFORNIA 92008.

References

- Angelichio, P. L., Beck, J. A., Johansen, H., and Ivey-Hoyle, M. (1991). Comparison of Several Promoters and Polyadenylation Signals for Use in Heterologous Gene Expression in Cultured *Drosophila* Cells. *Nucleic Acids Res.* 19, 5037-5043.
- 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).
- Bunch, T. A., Grinblat, Y., and Goldstein, L. S. (1988). Characterization and Use of the *Drosophila* Metallothionein Promoter in Cultured *Drosophila melanogaster* Cells. *Nucleic Acids Res.* 16, 1043-1061.
- Johansen, H., van der Straten, A., Sweet, R., Otto, E., Maroni, G., and Rosenberg, M. (1989). Regulated Expression at High Copy Number Allows Production of a Growth Inhibitory Oncogene Product in *Drosophila* Schneider Cells. *Genes and Development* 3, 882-889.
- Kozak, M. (1987). An Analysis of 5'-Noncoding Sequences from 699 Vertebrate Messenger RNAs. *Nuc. Acids Res.* 15, 8125-8148.
- Kozak, M. (1991). An Analysis of Vertebrate mRNA Sequences: Intimations of Translational Control. *J. Cell Biol.* 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.
- Maroni, G., Otto, E., and Lastowski-Perry, D. (1986). Molecular and Cytogenetic Characterization of a Metallothionein Gene of *Drosophila*. *Genetics* 112, 493-504.
- Miller, J. H. (1972). *Experiments in Molecular Genetics* (Cold Spring Harbor, New York: Cold Spring Harbor Laboratory).
- Olsen, M. K. (1992). Stable Production of an Analog of Human Tissue Plasminogen Activator from Cultured *Drosophila* Cells. *Cytotechnology* 10, 157-167.
- Sambrook, J., Fritsch, E. F., and Maniatis, T. (1989). *Molecular Cloning: A Laboratory Manual*, Second Edition (Plainview, New York: Cold Spring Harbor Laboratory Press).
- Southern, J. A., Young, D. F., Heaney, F., Baumgartner, W., and Randall, R. E. (1991). Identification of an Epitope on the P and V Proteins of Simian Virus 5 That Distinguishes Between Two Isolates with Different Biological Characteristics. *J. Gen. Virol.* 72, 1551-1557.

©2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the properties of Life Technologies Corporation or their respective owners.

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

Notes



Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA | Phone +1 760 603 7200 | Toll Free in USA 800 955 6288

For support visit www.invitrogen.com/support or email techsupport@invitrogen.com

www.lifetechnologies.com

