



## pSecTag2/Hygro A, B, and C

Catalog number V910-20

Revision date 19 January 2012

Publication Part number 28-0171

MAN0000655



## **Contents**

Kit Contents and Storage	iv
Introduction	
Overview	1
pSecTag2/Hygro Vector	
Methods	
Cloning into pSecTag2/Hygro	
Transformation into <i>E. coli</i>	
Transfection into Mammalian Cells	9
Expression and Purification	10
Appendix	11
pSecTag2/Hygro/PSA	11
Additional Products	
Technical Support	
Purchaser Notification	14
References	15

## **Kit Contents and Storage**

Kit Contents	Each kit contains 20 $\mu g$ each of pSecTag2/Hygro A, B, and C and pSecTag2/Hygro/PSA. Each vector is supplied at a concentration of 0.5 $\mu g/\mu L$ in 40 $\mu L$ of 10 mM Tris-HCl, 1 mM EDTA, pH 8.0.
Shipping and Storage	Plasmids are shipped on ice and should be stored at -20°C.
Intended Use	For research use only. Not intended for any human or animal diagnostic or therapeutic uses.

#### Introduction

#### **Overview**

#### Introduction

pSecTag2/Hygro A, B, and C are 5.7 kb expression vectors designed for high-level expression and secretion in mammalian hosts. The pSecTag2/Hygro vectors are identical to the pSecTag2 vectors except that the Zeocin<sup>™</sup> resistance gene is replaced with the Hygromycin B resistance gene (Gritz and Davies, 1983) for selection in mammalian cells. Like pSecTag2, pSecTag2/Hygro contains the gene encoding β-lactamase for bacterial selection on ampicillin.

Proteins expressed from pSecTag2/Hygro are fused at the N-terminus to the murine Ig kappa chain leader sequence for protein secretion and at the C-terminus to a peptide containing the *c-myc* epitope and six tandem histidine residues for detection and purification.

The pSecTag2/Hygro vector is supplied in three different versions (A, B, and C) to facilitate correct in-frame fusion with the Ig kappa chain leader sequence.

For more information on the pSecTag2/Hygro vector, see page 2.

To get started with cloning into pSecTag2/Hygro, see page 4.

#### E. coli Strain

We recommend that you propagate pSecTag2/Hygro A, B and C in *E. coli* strains that are recombination deficient (recA) and endonuclease A deficient (endA) such as TOP10F′, DH5 $\alpha$ F′, and INV $\alpha$ F′.

For your convenience, TOP10F´ is available as chemically competent or electrocompetent cells from Life Technologies.

Item	Quantity	Catalog no.
Electrocomp <sup>™</sup> TOP10F′	$5 \times 80 \mu L$	C665-55
One Shot® TOP10F' (chemically competent cells)	21 × 50 μL	C3030-03

## pSecTag2/Hygro Vector

## Features of pSecTag2/Hygro

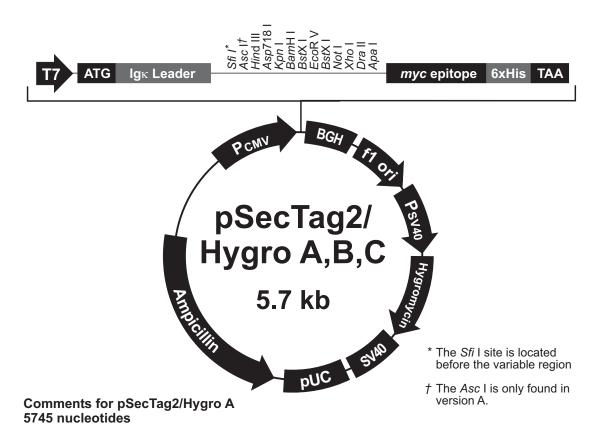
pSecTag2/Hygro A (5745 bp), pSecTag2/Hygro B (5749 bp), and pSecTag2/Hygro C (5753 bp) contain the following elements. All features have been functionally tested.

Feature	Benefit
Human cytomegalovirus (CMV) immediate-early promoter/enhancer	Allows efficient, high-level expression of your recombinant protein (Andersson <i>et al.</i> , 1989; Boshart <i>et al.</i> , 1985; Nelson <i>et al.</i> , 1987)
T7 promoter/priming site	Allows for <i>in vitro</i> transcription in the sense orientation and sequencing through the insert
ATG initiation codon	Allows initiation of translation of the pSecTag2/Hygro fusion protein
Murine Ig κ-chain leader sequence	Allows secretion of the fusion protein (Coloma <i>et al.</i> , 1992)
Multiple cloning site	Allows insertion of your gene and facilitates cloning
<i>c-myc</i> epitope (Glu-Gln-Lys-Leu-Ile-Ser-Glu-Glu-Asp-Leu)	Allows detection of pSecTag2/Hygro fusion protein with the Anti- <i>myc</i> Antibody (Catalog no. R950-25) (Evans <i>et al.</i> , 1985)
Polyhistidine tag	For high affinity binding to Ni <sup>2+</sup> -chelating resin (i.e. ProBond <sup>™</sup> ) and easy purification
	In addition, it allows detection of pSecTag2/Hygro fusion proteins with the Anti-His(C-term) Antibody (Catalog no. R930-25) (Lindner <i>et al.</i> , 1997)
BGH reverse priming site	Allows sequencing through the insert
Bovine growth hormone (BGH) polyadenylation signal	Efficient transcription termination and polyadenylation of mRNA (Goodwin and Rottman, 1992)
f1 origin	Allows rescue of single-stranded DNA
SV40 early promoter and origin	Allows efficient, high-level expression of the Hygromycin resistance gene and episomal replication in cells expressing the SV40 large T antigen
Hygromycin resistance gene (Hygromycin B- phosphotransferase)	Selection of stable transfectants in mammalian cells (Gritz and Davies, 1983; Palmer <i>et al.</i> , 1987)
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 in E. coli

### pSecTag2/Hygro Vector, continued

## Map of pSecTag2/Hygro

The figure below shows where the features of pSecTag2/Hygro are located in the vector. The sequences for pSecTag2/Hygro A, B, and C are available for downloading from our website (www.lifetechnologies.com) or from Technical Support (page 13). Details of the multiple cloning sites for all three vectors, including the variable region that determines the reading frame, are shown on pages 5–7.



CMV promoter: bases 209-863

T7 promoter/priming site: bases 863-882

Murine Ig kappa-chain V-J2-C signal peptide: bases 905-967

Multiple cloning site: bases 970-1081 *c-myc* epitope: bases 1082-1111 Polyhistidine tag: bases 1127-1144

BGH reverse priming site: bases 1167-1184 BGH polyadenylation site: bases 1166-1380

f1 origin: bases 1443-1856

SV40 promoter and origin: bases 1924-2245

Hygromycin B phosphotransferase ORF (HygR): bases 2263-3288

SV40 polyadenylation site: bases 3418-3547

pUC origin: bases 3931-4604

β-lactamase ORF (Amp<sup>R</sup>): bases 4749-5609

#### **Methods**

## Cloning into pSecTag2/Hygro

#### Introduction

This section contains information on cloning your insert into the pSecTag2/Hygro vectors. Details of the multiple cloning sites are found on pages 5–7. Brief information on transforming into *E. coli* is located 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, see *Molecular Cloning: A Laboratory Manual* (Sambrook *et al.*, 1989) or *Current Protocols in Molecular Biology* (Ausubel *et al.*, 1994)

## Maintenance of pSecTag2/Hygro

To propagate and maintain pSecTag2/Hygro A, B, and C, we recommend that you transform the plasmids into *E. coli* and prepare glycerol stocks for long-term storage as described below:

- 1. Use the supplied  $0.5 \,\mu\text{g}/\mu\text{L}$  stock solution (in TE, pH 8.0) to transform a recA, endA E. coli strain like INV $\alpha$ F′, TOP10F′, DH5 $\alpha$ F′, or equivalent.
- 2. Select transformants on LB plates containing 50–100 µg/mL ampicillin.
- 3. Analyze transformants for the appropriate plasmid and prepare glycerol stocks by mixing 0.85 mL of an overnight culture with 0.15 mL of sterile glycerol. Transfer the resulting solution to a cryovial and store at -80°C.

# Cloning into the pSecTag2/Hygro Vectors

pSecTag2/Hygro A, B, and C vectors are fusion vectors requiring that you clone your gene of interest in frame with the initiation ATG of the N-terminal Ig kappa chain leader sequence and/or the C-terminal *myc* epitope/polyhistidine tag. Three versions of this vector are provided to facilitate cloning. For proper expression, first determine which restriction sites are appropriate for ligation and then which vector will preserve the reading frame at **BOTH** the 5´ and the 3´ ends. It may be necessary to PCR your gene product to create a fragment with the appropriate restriction sites to clone in frame at both ends. Carefully inspect your gene and the multiple cloning site of each vector before cloning your gene of interest. Be sure to remove the stop codon in your gene if you wish to express your protein with the C-terminal tag. See pages 5–7 for details of the multiple cloning sites.

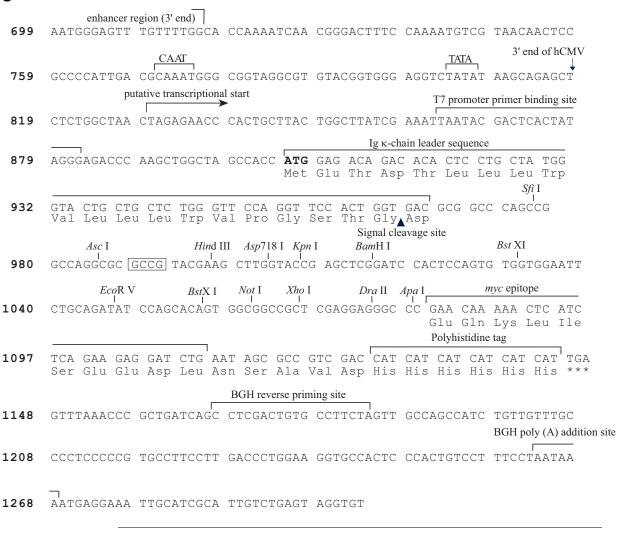
#### **Note**

If you wish to express and secrete your protein without the C-terminal tag, include a termination codon in your gene of interest. Note that you will be unable to detect the fusion protein with the Anti-myc Antibody or Anti-His (C-term) Antibody or purify it using nickel-chelating resin (i.e. ProBond<sup>TM</sup>).

### Cloning into pSecTag2/Hygro, continued

#### Multiple Cloning Site of pSecTag2/ Hygro A

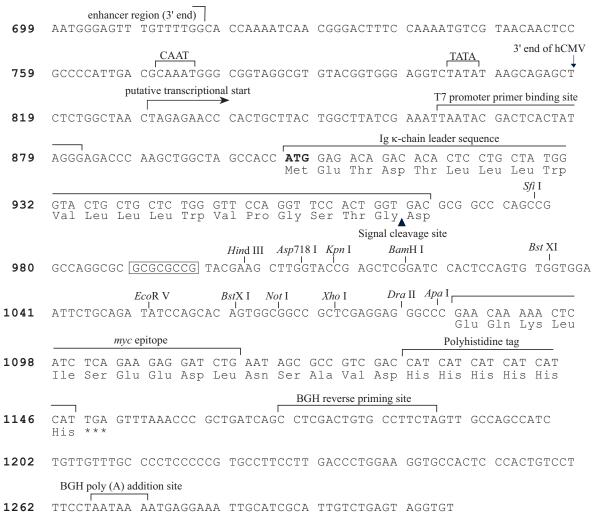
Below is the multiple cloning site for pSecTag2/Hygro A. Restriction sites are labeled to indicate the cleavage site. The variable region is the boxed region located after the Ig kappa chain leader sequence. The multiple cloning site has been confirmed by sequencing and functional testing.



### Cloning into pSecTag2/Hygro, continued

#### Multiple Cloning Site of pSecTag2/ Hygro B

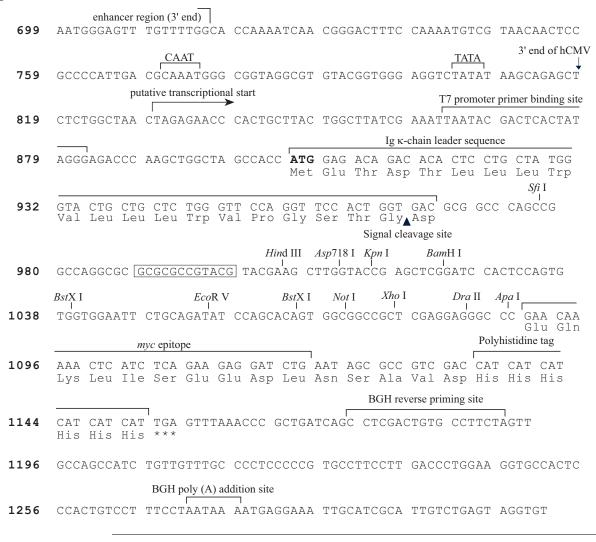
Below is the multiple cloning site for pSecTag2/Hygro B. Restriction sites are labeled to indicate the cleavage site. The variable region is the boxed region located after the Ig kappa chain leader sequence. The multiple cloning site has been confirmed by sequencing and functional testing.



### Cloning into pSecTag2/Hygro, continued

#### Multiple Cloning Site of pSecTag2/ Hygro C

Below is the multiple cloning site for pSecTag2/Hygro C. Restriction sites are labeled to indicate the cleavage site. The variable region is the boxed region located after the Ig kappa chain leader sequence. The multiple cloning site has been confirmed by sequencing and functional testing.



#### Transformation into E. coli

#### Introduction

At this point you should have ligation mixtures that are ready to be transformed into competent *E. coli*. The following guidelines and recommendations are provided for your convenience. If you need more details about the techniques discussed, refer to the general molecular biology references in the **Reference** section.

## E. coli Transformation

Transform your ligation mixtures into a competent recA, endA E. coli strain (e.g. INV $\alpha$ F´, TOP10F´, DH5 $\alpha$ <sup>TM</sup>) and select on LB plates containing 50–100  $\mu$ g/mL ampicillin. Select 10–20 clones and analyze for the presence and orientation of your insert.



We recommend that you sequence your construct with the T7 forward and BGH reverse primer binding sites to confirm that your gene is correctly fused to the Ig kappa chain leader sequence at the N-terminal and the C-terminal tag. Refer to diagrams on pages 5–7 for sequence and location of primer binding sites.

For your convenience, Life Technologies offers a custom primer synthesis service. For more information, refer to our website (www.lifetechnologies.com) or contact Technical Support (page 13).

## Plasmid Preparation

Once you have confirmed that your gene is in the correct reading frame, prepare plasmid DNA for transfection. 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 Plasmid MiniPrep or Plasmid MidiPrep kits, or the ChargeSwitch®-Pro Plasmid Miniprep Kit (see page 12, Additional Products).

#### **Transfection into Mammalian Cells**

#### Methods of Transfection

For established cell lines (e.g., HeLa), consult original references or the supplier of your cell line for the optimal method of transfection. It is recommended that you follow exactly the protocol for your cell line. Pay particular attention to medium requirements, when to pass the cells, and at what dilution to split the cells. Further information is provided in *Current Protocols in Molecular Biology*.

#### Hygromycin B Selection Guidelines

If you wish to create stable cell lines expressing your gene of interest you will need to select using Hygromycin B. Use Hygromycin B (527.5 MW) as follows:

- Test varying concentrations of Hygromycin on your cell line to determine the concentration that kills your cells (kill curve). Cells differ in their susceptibility to Hygromycin.
- Prepare complete medium supplemented with 100–1000 μg/mL
   Hygromycin B. Generally, concentrations that kill mammalian cells are in the 150–400 μg/mL range.
- Calculate concentration based on the amount of active drug (check the lot label).

Cells will divide once or twice in the presence of lethal doses of Hygromycin B, so the effects of the drug take several days to become apparent. Complete inhibition of cell growth can take 2–3 weeks of growth in selective medium. Selection and expansion of clones will take additional time.

## Hygromycin B Activity

Hygromycin B is an aminocyclitol that inhibits protein synthesis by disrupting translocation and promoting mistranslation. It is used for selection of stable mammalian cell lines. The resistance gene encodes Hygromycin B-phosphotransferase, which detoxifies Hygromycin B by phosphorylation.

### **Expression and Purification**

#### Introduction

Expression of your recombinant protein can be detected using an antibody to the c-myc epitope encoded in the C-terminal fusion peptide. In addition, the metal binding domain allows simple, one-step purification of your recombinant protein by Immobilized Metal Affinity Chromatography (IMAC) using a nickel-chelating resin. (i.e., ProBond $^{\text{TM}}$ ). See **Additional Products** on page 12 for detection and purification products from Life Technologies.

#### Detection of Secreted Protein from Medium

The medium in which transfected cells are grown can be analyzed for secreted, recombinant protein by functional assay or western blot analysis. It may be necessary to perform a time course analysis to determine the optimal time for expression of your recombinant protein. If you do not have an antibody to your particular protein, you can use the Anti-*myc* Antibody to detect your protein.

A positive control vector, pSecTag2/Hygro/PSA, is included to test for expression and secretion in your particular cell line. Prostate-specific antigen (PSA) is fused to the *c-myc* epitope and the polyhistidine tag. The resulting fusion protein is ~33 kDa which includes the N-terminal secretion signal. When the secretion signal is cleaved off, the size of the fusion protein should decrease by 2.4 kDa. **Note**: There are glycosylation sites in PSA. The secreted fusion protein expressed in human breast carcinoma cells migrates at ~45 kDa.

#### Purification of Secreted Recombinant Protein

To purify secreted, recombinant protein from the medium, follow the manufacturer's instructions for the nickel-chelating resin that you are using. Start with about 3–5 mL of medium and load onto 1–2 mL of resin. Scale up or down depending on the level of expression.

# Analyzing Cells for Recombinant Protein

If you do not detect any secreted protein in the medium, use the procedure below to check cells for production of recombinant protein. You will need  $5 \times 10^6$  to  $1 \times 10^7$  cells for purification on a 2 mL ProBond<sup>TM</sup> column (see ProBond<sup>TM</sup> Protein Purification manual).

- 1. Seed cells in either five T-75 flasks or 2–3 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 2–5 minutes or by scraping the cells in PBS.
- 4. Inactivate the trypsin by diluting with complete medium. Transfer the cells to a sterile conical tube.
- 5. Centrifuge the cells at 1500 rpm for 5 minutes. You may wish to wash the cells one time with PBS prior to lysis. The cells may be lysed immediately or frozen in liquid nitrogen and stored at –80°C until needed.

#### **Lysis of Cells**

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

If you are using other resin, refer to the manufacturer's instruction for recommendations on sample preparation.

## **Appendix**

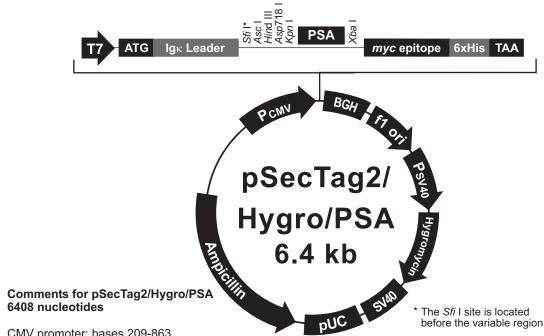
## pSecTag2/Hygro/PSA

#### **Description**

pSecTag2/Hygro/PSA is a 6408 bp positive control vector expressing and secreting the prostate-specific antigen (PSA) fused to the *c-myc* epitope and the polyhistidine tag. The vector was constructed by amplifying the PSA gene and cloning it into pCR<sup>™</sup>II. The fragment of DNA containing the PSA gene was excised using Kpn I and Apa I and cloned into Kpn I/Apa I digested pSecTag2/Hygro A.

#### Map

The figure below shows the features of pSecTag2/Hygro/PSA. The complete nucleotide sequence for pSecTag2/Hygro/PSA is available for downloading from our website (www.lifetechnologies.com) or from Technical Support (page 13).



CMV promoter: bases 209-863

T7 promoter/priming site: bases 863-882

Murine Ig kappa-chain V-J2-C signal peptide: bases 905-967

PSA gene: bases 1042-1732 c-myc epitope: bases 1745-1777 Polyhistidine tag: bases 1790-1807

BGH reverse priming site: bases 1830-1847 BGH polyadenylation sequence: bases 1829-2043

f1 origin: bases 1954-2367

SV40 promoter and origin: bases 2587-2908

Hygromycin B phosphotransferase ORF (HygR): bases 2926-3951

SV40 polyadenylation site: bases 4144-4210

pUC origin: bases 4594-5267

β-lactamase ORF (Amp<sup>R</sup>): bases 5412-6272

## **Additional Products**

#### **Kit Contents**

The following products are available separately from Life Technologies. To order, visit our website at **www.lifetechnologies.com** or contact Technical Support (see page 13).

Product	Amount	Catalog number
PureLink® HiPure Plasmid MiniPrep Kits	25 preps 100 preps	K2100-02 K2100-03
PureLink® HiPure Plasmid MidiPrep Kits	25 preps 50 preps	K2100-04 K2100-05
ChargeSwitch®-Pro Plasmid Miniprep Kit	50 preps 250 preps	CS30050 CS30250
Anti-myc Antibody	25 westerns	R950-25
Anti-His (C-term) Antibody	25 westerns	R930-25
ProBond <sup>™</sup> Purification System	6 purifications	K850-01
ProBond™ Metal-Binding Resin	50 mL 150 mL	R801-01 R801-15
Purification Columns (10-mL polypropylene columns)	50 each	R640-50
Hygromycin B	20 mL	10687-010

### **Technical Support**

#### **Obtaining support**

For the latest services and support information for all locations, go to www.lifetechnologies.com.

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 Sata Sheets (SDS)

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

#### 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 <a href="https://www.lifetechnologies.com/support">www.lifetechnologies.com/support</a> and search for the Certificate of Analysis by product lot number, which is printed on the box.

#### **Limited warranty**

Life Technologies and/or its affiliate(s) warrant their products as set forth in the Life Technologies General Terms and Conditions of Sale found on the Life Technologies web site at <a href="www.lifetechnologies.com/termsandconditions">www.lifetechnologies.com/termsandconditions</a>. If you have any questions, please contact Life Technologies.

Life Technologies and/or its affiliate(s) disclaim all warranties with respect to this document, expressed or implied, including but not limited to those of merchantability or fitness for a particular purpose. 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.

### **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 <code>outlicensing@lifetech.com</code> Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

#### References

- Andersson, S., Davis, D. L., Dahlbäck, H., Jörnvall, H., and Russell, D. W. (1989). Cloning, Structure, and Expression of the Mitochondrial Cytochrome P-450 Sterol 26-Hydroxylase, a Bile Acid Biosynthetic Enzyme. J. Biol. Chem. 264, 8222-8229.
- Ausubel, F. M., Brent, R., Kingston, R. E., Moore, D. D., Seidman, J. G., Smith, J. A., and Struhl, K. (1994). Current Protocols in Molecular Biology (New York: Greene Publishing Associates and Wiley-Interscience).
- Boshart, M., Weber, F., Jahn, G., Dorsch-Häsler, K., Fleckenstein, B., and Schaffner, W. (1985). A Very Strong Enhancer is Located Upstream of an Immediate Early Gene of Human Cytomegalovirus. Cell 41, 521-530.
- Chen, C., and Okayama, H. (1987). High-Efficiency Transformation of Mammalian Cells by Plasmid DNA. Molec. Cell. Biol. 7, 2745-2752.
- Chu, G., Hayakawa, H., and Berg, P. (1987). Electroporation for the Efficient Transfection of Mammalian Cells with DNA. Nucleic Acids Res. *15*, 1311-1326.
- Coloma, M. J., Hastings, A., Wims, L. A., and Morrison, S. L. (1992). Novel Vectors for the Expression of Antibody Molecules Using Variable Regions Generated by Polymerase Chain Reaction. J. Imm. Methods 152, 89-104.
- 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.
- Gritz, L., and Davies, J. (1983). Plasmid-Encoded Hygromycin-B Resistance: The Sequence of Hygromycin-B-Phosphotransferase Gene and its Expression in *E. coli* and *S. Cerevisiae*. Gene 25, 179-188
- 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.
- 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.
- Palmer, T. D., Hock, R. A., Osborne, W. R. A., and Miller, A. D. (1987). Efficient Retrovirus-Mediated Transfer and Expression of a Human Adenosine Deaminase Gene in Diploid Skin Fibroblasts from an Adenosine-Deficient Human. Proc. Natl. Acad. Sci. U.S.A. *84*, 1055-1059.
- 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.
- ©2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners. Zeocin is a trademark of Cayla.

## Notes

## Notes

## Notes



#### Headquarters