

pMT-DEST48 Gateway® Vector

A destination vector for cloning and inducible expression in *Drosophila* cells

Catalog no. 12282-018

Version D 23 December 2008 25-0414

A Limited Label License covers this product (see Purchaser Notification). By use of this product, you accept the terms and conditions of the Limited Label License.

User Manual

Table of Contents

Important Information	v
Accessory Products	vi
Methods	1
Overview	1
Using pMT-DEST48	3
Transfection and Analysis	6
Appendix	8
Map and Features of pMT-DEST48	8
Technical Support	10
Purchaser Notification	12
References	14

Important Information

Shipping and Storage	pMT-DEST48 is shipped on wet ice. Upon receipt, store at -20°C. Product is guaranteed for six months from date of shipment when stored properly.
Contents	6 μg of pMT-DEST48 vector at 150 ng/μl in TE buffer, pH 8.0 (10 mM Tris-HCl, 1 mM EDTA, pH 8.0). Volume: 40 μl.
Quality Control	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 <u>www.invitrogen.com/support</u> and search for the Certificate of Analysis by product lot number, which is printed on the box.

Accessory Products

Additional Products

Additional products that may be used with pMT-DEST48 are available from Invitrogen. Ordering information is provided below.

Product	Amount	Catalog no.
DES® Inducible Kit with pCoHygro	1 kit	K4120-01
DES [®] Inducible Kit with pCoBlast	1 kit	K5120-01
DES [®] -Blasticidin Support Kit*	1 kit	K5150-01
Schneider (S2) Cells	1 ml vial, 1 x 10^7 cells/ml	R690-07
Schneider's Drosophila Medium	500 ml	11720-034
Fetal Bovine Serum	500 ml	16000-044
Penicillin-Streptomycin	100 ml	15070-063
Calcium Phosphate Transfection Kit	75 reactions	K2780-01
Hygromycin B	1 g	R220-05
Blasticidin S HCl	50 mg	R210-01
pENTR Directional TOPO [®] Cloning Kit	20 reactions	K2400-20
PureLink™ HiPure Plasmid Midiprep Kit	25 preps	K2100-04
One Shot [®] ccdB Survival [™] 2 T1 Phage-Resistant Cells	10 reactions	A10460
Gateway [®] LR Clonase [™] Enzyme Mix	20 reactions	11791-019
One Shot [®] TOP10 Chemically	10 reactions	C4040-10
Competent Cells	20 reactions	C4040-03
One Shot® TOP10 Electrocompetent	10 reactions	C4040-50
Cells	20 reactions	C4040-52

*The DES[®]-Blasticidin Support Kit contains pCoBlast, the Calcium Phosphate Transfection Kit, frozen S2 cells, DES[®] Expression Medium, and blasticidin (50 mg).

Accessory Products, continued

Detection of
RecombinantExpression of your recombinant fusion protein can be
detected using an antibody to the appropriate epitope. The
amount of antibody supplied is sufficient for 25 Westerns.

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

Purification of Recombinant Fusion Protein

If your gene of interest is in frame with the C-terminal peptide containing the V5 epitope and the polyhistidine (6xHis) tag, you may use Immobilized Metal Affinity Chromatography (IMAC) to purify your recombinant fusion protein. The ProBond[™] Purification System or bulk ProBond[™] resin are available separately from Invitrogen. See the table below for ordering information.

Product	Quantity	Catalog no.
ProBond [™] Nickel-chelating Resin	50 ml 150 ml	R801-01 R801-15
Purification Columns (10 ml polypropylene columns, empty)	50 each	R640-50
ProBond [™] Purification System	6 purifications	K850-01
Positope TM Control Protein	5 µg	R900-50

Accessory Products, continued

Gel Electrophoresis	A wide range of pre-cast NuPAGE [®] and Novex [®] Tris- Glycine polyacrylamide gels and electrophoresis apparatuses are available to facilitate separation and visualization of your recombinent fusion protein
	visualization of your recombinant fusion protein.

For more details on these products, visit our Web site at <u>www.invitrogen.com</u> or contact Technical Support (page 10).

Methods

Overview

Description	pMT-DEST48 is derived from pMT/V5-His and adapted for use with the Gateway [®] Technology. It is designed for use with the <i>Drosophila</i> Inducible Expression System (DES [®] ; Catalog no. K4120-01) available from Invitrogen. Upon transfection, the vector allows transient, inducible expression of your protein of interest in <i>Drosophila</i> cells. Co-transfecting pMT-DEST48 with the selection vector, pCoHygro or pCoBlast, allows selection of stable cell lines exhibiting inducible expression of the protein of interest.
	Note : The DES [®] Inducible Kit is available with either pCoHygro or pCoBlast. pCoBlast is also available in the DES [®] -Blasticidin Support Kit. For more information on these products, see page vi.
Features	 pMT-DEST48 contains the following elements: <i>Drosophila</i> metallothionein (MT) promoter for high-level, metal-inducible expression of the gene of interest in S2 cells (see next page for more information) Two recombination sites, <i>att</i>R1 and <i>att</i>R2, downstream of the MT promoter for recombinational cloning of the gene of interest from an entry clone Chloramphenicol resistance gene located between the two <i>att</i>R sites for counterselection The <i>ccd</i>B gene located between the two <i>att</i>R sites for negative selection The V5 epitope and 6xHis tag for detection and purification (optional) SV40 late polyadenylation sequence for proper termination and processing of the recombinant transcript The pUC origin for replication and maintenance of the plasmid in <i>E. coli</i>
	• The ampicillin (<i>bla</i>) resistance gene for selection in <i>E. coli</i> For a map of pMT-DEST48, see page 8.

Overview, continued

Gateway [®] Technology	Gateway [®] is a universal cloning technology that takes advantage of the site-specific recombination properties of bacteriophage lambda (Landy, 1989) to provide a rapid and highly efficient way to move your gene of interest into multiple vector systems. To express your gene of interest using Gateway [®] Technology, simply:
	1. Clone your gene of interest into a Gateway [®] entry vector to create an entry clone.
	2. Generate an expression clone by performing an LR recombination reaction between the entry clone and a Gateway [®] destination vector (<i>e.g.</i> pMT-DEST48).
	3. Introduce your expression clone into <i>Drosophila</i> S2 cells.
	For more information on the Gateway [®] Technology, refer to the Gateway [®] Technology Manual. This manual is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Support (page 10).
Description of MT Promoter	The <i>Drosophila</i> MT promoter allows high-level, inducible expression of the gene of interest in <i>Drosophila</i> S2 cells. When used to express heterologous proteins, the promoter is extremely efficient and tightly regulated, even at high copy number (Johansen <i>et al.</i> , 1989). The MT promoter is well characterized (Angelichio <i>et al.</i> , 1991; Bunch <i>et al.</i> , 1988; Maroni <i>et al.</i> , 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 <i>et al.,</i> 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.

Using pMT-DEST48

Q Important	The pMT-DEST48 vector is supplied as a supercoiled plasmid. Although Invitrogen has previously recommended using a linearized destination vector for more efficient recombination, further testing has found that linearization of this vector is NOT required to obtain optimal results for any downstream application.
Propagating pMT-DEST48	If you wish to propagate and maintain pMT-DEST48, we recommend using One Shot [®] <i>ccdB</i> Survival TM 2 T1 Phage-Resistant Cells for transformation (page vi). The <i>ccdB</i> Survival TM 2 T1 Phage-Resistant <i>E. coli</i> strain is resistant to CcdB effects and can support the propagation of plasmids containing the <i>ccdB</i> gene.
	Note: Do not use general <i>E. coli</i> cloning strains including TOP10 or DH5 α for propagation and maintenance as these strains are sensitive to CcdB effects.
Entry Clone	To recombine your gene of interest into pMT-DEST48, you should have an entry clone containing your gene of interest. For your convenience, Invitrogen offers the pENTR Directional TOPO [®] Cloning Kit (Catalog no. K2400-20) for 5 minute cloning of your gene of interest into an entry vector. For more information on entry vectors available from Invitrogen, refer to our Web site (<u>www.invitrogen.com</u>) or contact Technical Support (page 10).
	For detailed information on constructing an entry clone, refer to the specific entry vector manual. For detailed information on performing the LR recombination reaction, refer to the Gateway [®] Technology Manual.
	continued on next page

Using pMT-DEST48, continued

Points to Consider Before Recombining	If you wish to include the V5 epitope and 6xHis tag, your gene in the entry clone should not contain a stop codon. The gene should contain an ATG initiation codon and should also be designed to be in frame with the C-terminal epitope tag after recombination. Refer to the Recombination Region on the next page. If you DO NOT wish to include the V5 epitope and 6xHis tag, your gene should contain both an ATG initiation codon and a stop codon in the entry clone.
Recombining Your Gene of Interest	Each entry clone contains <i>att</i> L sites flanking the gene of interest. Genes in an entry clone are transferred to the destination vector backbone by mixing the DNAs with the Gateway [®] LR Clonase TM enzyme mix (see page vi for ordering information). The resulting recombination reaction is then transformed into <i>E. coli</i> and the expression clone selected. Recombination between the <i>att</i> R sites on the destination vector and the <i>att</i> L sites on the entry clone replaces the <i>ccd</i> B gene and the chloramphenicol (Cm ^R) gene with the gene of interest and results in the formation of <i>att</i> B sites in the expression clone. Follow the instructions in the Gateway [®] Technology Manual to set up the LR Clonase TM reaction, transform a <i>recA endA E</i> .
Confirming the Expression Clone	The <i>ccd</i> B gene mutates at a very low frequency, resulting in a very low number of false positives. True expression clones will be ampicillin-resistant and chloramphenicol-sensitive. Transformants containing a plasmid with a mutated <i>ccd</i> B gene will be both ampicillin- and chloramphenicol-resistant. To check your putative expression clone, test for growth on LB plates containing 30 µg/ml chloramphenicol. A true expression clone will not grow in the presence of chloramphenicol.

Using pMT-DEST48, continued

Reco Regio	mbination on	from pMT	-DEST48 ×	0	is shown bel	lone resulting low.
		Shaded transfe recomb	d regions co rred from t	orrespond to he entry clor on-shaded re	those DNA ne into pMT	
	• The underlined nucleotides flanking the shaded region correspond to bases 866 and 2717, respectively, of the pMT-DEST48 vector sequence.					
	Metallothionein pr	omoter	TATA box		→	
725	AAAGCCGCGT	TTCCAAAATG			GCCAATGTGC CGGTTACACG	
785					AGGGGGGGATC TCCCCCCTAG	
845	TACATCAAAC ATGTAGTT <mark>TG</mark>	_	TTTTTTCGAC	CEN	IF NAC CCA	2543 ala phe GCT TTC CGA AAG attB2
					V5 epitope	
2544	TTG TAC AAA	GTG GTT CG	GA TGG TTC	GAA GGT AÂG	pro ile pro CCT ATC CCT GGA TAG GGA	AAC CCT
					6xHis tag	5
2592	CTC CTC GGT	CTC GAT TC	CT ACG CGT	ACC GGT CAT	his his his CAT CAC CAT GTA GTG GTA	

Transfection and Analysis

Introduction

The section provides general guidelines for transfecting *Drosophila* S2 cells. If you are assaying for transient, inducible expression of your gene of interest, you may transfect your pMT-DEST48 construct alone into S2 cells. If you wish to generate stable cell lines, you **must** cotransfect your pMT-DEST48 construct with pCoHygro or pCoBlast into S2 cells. We recommend that you include a mock transfection (negative control) in your experiments to evaluate your results.

For specific guidelines on culturing S2 cells, performing transient transfection, and generating stable cell lines, refer to the *Drosophila* Expression System (DES[®]) manual. This manual also contains detailed information about pCoHygro and pCoBlast and is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Support (page 10).



Note that pMT-DEST48 does not contain a resistance marker for selection in *Drosophila* cells.

Plasmid Preparation

Plasmid DNA for transfection into S2 cells must be very clean and free from phenol and sodium chloride. Contaminants will kill the cells, and salt will interfere with lipid complexing, decreasing transfection efficiency. We recommend isolating plasmid DNA using the S.N.A.P.[™] MiniPrep Kit (10-15 µg DNA, Catalog no. K1900-01), the S.N.A.P.[™] MidiPrep Kit (10-200 µg DNA, Catalog no. K1910-01), or CsCl gradient centrifugation.

Transfection and Analysis, continued

Induction of Recombinant Protein Expression	Once you have transfected your pMT-DEST48 construct into S2 cells, you will use copper sulfate to induce expression of your recombinant protein. In general, we recommend that you add copper sulfate directly to the culture medium to a final concentration of 500 μ M and incubate the cells for 24 hours to obtain maximal induction of your recombinant protein of interest. Expression conditions may vary depending on the nature of the recombinant protein of interest. You may want to perform a time course to optimize expression of the recombinant protein. For more details, refer to the DES [®] manual. Copper sulfate is provided with the DES [®] Inducible Kit available from Invitrogen (see page vi for ordering information).
Detecting Recombinant Fusion Proteins	To detect expression of your recombinant fusion protein by Western blot analysis, you may use the Anti-V5 antibodies or the Anti-His(C-term) antibodies available from Invitrogen (see page vii for ordering information) or an antibody to your protein of interest. In addition, the Positope [™] Control Protein (Catalog no. R900-50) is available from Invitrogen for use as a positive control for detection of fusion proteins containing a V5 epitope or a polyhistidine (6xHis) tag.
Note	The C-terminal peptide containing the V5 epitope and the 6xHis tag will add approximately 4 kDa to your protein.
Purification of Recombinant Fusion Proteins	The presence of the C-terminal polyhistidine (6xHis) tag in your recombinant fusion protein allows use of a metal- chelating resin such as ProBond [™] to purify your fusion protein. The ProBond [™] Purification System and bulk ProBond [™] resin are available from Invitrogen (see page vii for ordering information). Refer to the ProBond [™] Purification System manual for protocols to purify your fusion protein. Invitrogen also offers Ni-NTA Agarose (Catalog no. R901-01) for purification of proteins containing a polyhistidine (6xHis) tag. Note: Other metal-chelating resins and purification methods are suitable.

Appendix

Map and Features of pMT-DEST48

Map of pMT-DEST48

The map below shows the elements of pMT-DEST48. DNA from the entry clone replaces the region between bases 860 and 2543. The complete sequence of pMT-DEST48 is available from our Web site (www.invitrogen.com) or by contacting Technical Support (page 10).



Map and Features of pMT-DEST48, continued

Features of
pMT-DEST48pMT-DEST48 (5179 bp) contains the following elements. All
features have been functionally tested.

Feature	Benefit
<i>Drosophila</i> metallothionein (MT) promoter	Allows strong, inducible expression of your recombinant protein (Angelichio <i>et al.</i> , 1991; Bunch <i>et al.</i> , 1988; Maroni <i>et al.</i> , 1986; Olsen, 1992)
attR1 and attR2 sites	Allows recombinational cloning of the gene of interest from an entry clone
Chloramphenicol resistance gene	Allows counterselection of expression clones
<i>ccd</i> B gene	Allows negative selection of expression clones
V5 epitope	Allows detection of recombinant fusion proteins by the Anti-V5 antibodies (Southern <i>et al.</i> , 1991)
C-terminal polyhistidine tag	Allows purification of recombinant fusion proteins on metal-chelating resin (<i>e.g.</i> ProBond [™])
	Allows detection of the recombinant protein by the Anti-His (C-term) antibodies (Lindner <i>et al.,</i> 1997)
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 resistance gene (β-lactamase)	Allows selection of transformants in <i>E. coli</i>

Technical Support

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

Web

World Wide

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

Contact Us

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

Corporate Headquarters:

Invitrogen Corporation 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:

Invitrogen Japan 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:

Invitrogen Ltd 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 <u>www.invitrogen.com/msds</u>.

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 <u>www.invitrogen.com/support</u> and search for the Certificate of Analysis by product lot number, which is printed on the box.

Technical Support, continued

Limited Invitrogen is committed to providing our customers with Warranty 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. Invitrogen warrants that all of its products will 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 Invitrogen Corporation's liability only to the cost 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. Invitrogen reserves the right to select the method(s) used to analyze a product unless Invitrogen 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 Invitrogen 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. Invitrogen assumes 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. 19: Gateway[®] Cloning Products

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 purchase of this product does not convey a license under any method claims in the foregoing patents or patent applications, or to use this product with any recombination sites other than those purchased from Invitrogen Corporation or its authorized distributor. The right to use methods claimed in the foregoing patents or patent applications with this product for research purposes only can only be acquired by the use of Clonase[™] purchased from Invitrogen Corporation or its authorized distributors. The buyer cannot modify the recombination sequence(s) contained in this product for any purpose. The buyer cannot sell or otherwise transfer (a) this product, (b) its components, or (c) materials made by the employment of this product or its components to a third party or otherwise use this product or its components or materials made by the employment of this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the employment 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. Notwithstanding the preceding, any buyer who is employed in an academic or government institution may transfer materials made with this product to a third party who has a license from Invitrogen under the patents identified above to distribute such materials. Transfer of such materials and/or information to collaborators does not convey rights to practice any methods claimed in the foregoing patents or patent applications. 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. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents 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 none of (i) this product, (ii) any of its components, or (iii) a method claim of the foregoing patents, was used in the manufacture of such product. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents based upon the use of this product to manufacture a protein for sale, provided that no method claim in the above patents was used in the manufacture of such protein. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to use this product for purposes other than those permitted above, contact Licensing Department, Invitrogen Corporation, 5791 Van Allen Way, Carlsbad, California 92008. Phone (760) 603-7200.

Purchaser Notification, continued

Limited Use Label License No. 56: Drosophila Expression System (DES[®]) This product is licensed to you under patent license from SmithKline Beecham Corporation for your research use only. This product may not be sold or transferred to any other person. In using this product, you agree to make no commercial use of this product. Requests for licenses for commercial manufacture or use may be made directly to the Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, CA. Tel: 760-603-7200; Fax: 760-602-6500.

Please note that you need a license to practice this technology. If you do not have a signed agreement on file, please use the contact information above to obtain a license agreement.

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.

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.

Landy, A. (1989). Dynamic, Structural, and Regulatory Aspects of Lambda Sitespecific Recombination. Annu. Rev. Biochem. 58, 913-949.

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.

Olsen, M. K. (1992). Stable Production of an Analog of Human Tissue Plasminogen Activator from Cultured *Drosophila* Cells. Cytotechnology 10, 157-167.

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.

©2002–2008 Invitrogen Corporation. All rights reserved.

invitrogen

Corporate Headquarters Invitrogen Corporation 5791 Van Allen Way Carlsbad, CA 92008 T: 1 760 603 7200 F: 1 760 602 6500 E: tech_support@invitrogen.com

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