

pcDNA[™]3.2/V5-DEST and pcDNA[™]6.2/V5-DEST Gateway[®] Vectors

Gateway[®]-adapted destination vectors for cloning and expression of C-terminal V5 fusion proteins in mammalian cells

Catalog nos. 12489-019 and 12489-027

Version E 27 October 2010 25-0500

User Manual

Table of Contents

Important Information	v
Accessory Products	vi
Methods	1
Overview	1
Generating an Entry Clone	3
Creating an Expression Clone	5
Performing the LR Recombination Reaction	8
Transfection	12
Expression and Analysis	14
Creating Stable Cell Lines	16
Appendix	19
Appendix	19 19
Appendix Recipes Blasticidin	19 19 21
Appendix Recipes Blasticidin Map of pcDNA™3.2/V5-DEST and pcDNA™6.2/V5-DEST	19 19 21 22
Appendix Recipes Blasticidin Map of pcDNA [™] 3.2/V5-DEST and pcDNA [™] 6.2/V5-DEST Features of pcDNA [™] 3.2/V5-DEST and pcDNA [™] 6.2/V5-DEST	19 19 21 22 23
Appendix Recipes Blasticidin Map of pcDNA [™] 3.2/V5-DEST and pcDNA [™] 6.2/V5-DEST Features of pcDNA [™] 3.2/V5-DEST and pcDNA [™] 6.2/V5-DEST Map of pcDNA [™] 3.2/V5/GW/CAT and pcDNA [™] 6.2/V5/GW/CA [™]	19 21 22 23 T24
Appendix Recipes Blasticidin Map of pcDNA [™] 3.2/V5-DEST and pcDNA [™] 6.2/V5-DEST Features of pcDNA [™] 3.2/V5-DEST and pcDNA [™] 6.2/V5-DEST Map of pcDNA [™] 3.2/V5/GW/CAT and pcDNA [™] 6.2/V5/GW/CA [™] Technical Service	19
Appendix Recipes Blasticidin Map of pcDNA [™] 3.2/V5-DEST and pcDNA [™] 6.2/V5-DEST Features of pcDNA [™] 3.2/V5-DEST and pcDNA [™] 6.2/V5-DEST Map of pcDNA [™] 3.2/V5/GW/CAT and pcDNA [™] 6.2/V5/GW/CA [™] Technical Service Purchaser Notification	19 19 22 23 T24 25 27
AppendixRecipesBlasticidinMap of $pcDNA^{TM}3.2/V5$ -DEST and $pcDNA^{TM}6.2/V5$ -DESTFeatures of $pcDNA^{TM}3.2/V5$ -DEST and $pcDNA^{TM}6.2/V5$ -DESTMap of $pcDNA^{TM}3.2/V5/GW/CAT$ and $pcDNA^{TM}6.2/V5/GW/CA'$ Technical ServicePurchaser NotificationGateway [®] Clone Distribution Policy	19

Important Information

Gateway [®] Vectors	This manual is supplied with the following products.									
	Product	Catalog no.								
	pcDNA [™] 3.2/V5-DEST Gateway [®] Vector	12489-019								
	pcDNA [™] 6.2/V5-DEST Gateway [®] Vector	12489-027								
Q Important	The pcDNA [™] 3.2/V5-DEST and pcDNA [™] 6.2/V5-DEST Gateway [®] Vectors have been renamed to be more descriptive and to better reflect the functionality of the vector.									
Shipping and Storage	The pcDNA [™] 3.2/V5-DEST and pcDNA [™] 6.2/ Gateway [®] Vectors are shipped on wet ice. Up store at –20°C.	'V5-DEST oon receipt,								
Contents	The pcDNA [™] 3.2/V5-DEST and pcDNA [™] 6.2/	'V5-DEST								

The pcDNA[™]3.2/V5-DEST and pcDNA[™]6.2/V5-DEST Gateway[®] Vector components are listed below.

Item	Concentration	Volume
Gateway [®] Destination Vector (pcDNA [™] 3.2/V5-DEST or pcDNA [™] 6.2/V5-DEST)	6 μg at 150 ng/μl, in TE buffer, pH 8.0 (10 mM Tris- HCl, 1 mM EDTA, pH 8.0)	40 µl
Control Plasmid (pcDNA [™] 3.2/V5/GW/CAT or pcDNA [™] 6.2/V5/GW/CAT)	10 μg at 0.5 μg/μl, in TE buffer, pH 8.0 (10 mM Tris- HCl, 1 mM EDTA, pH 8.0)	20 µl

Product	The Certificate of Analysis provides detailed quality control
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.

Accessory Products

Additional Products

Additional products that may be used with the pcDNA[™]3.2/V5-DEST and pcDNA[™]6.2/V5-DEST vectors are available from Invitrogen. Ordering information is provided below.

Product	Amount	Catalog no.			
Gateway [®] LR Clonase [™] II Enzyme Mix	20 reactions	11791-020			
	100 reactions	11791-100			
Tag-On-Demand [™] Suppressor	200 µl	K400-01			
Supernatant	5 x 200 μl	K405-01			
One Shot [®] TOP10 Chemically	10 reactions	C4040-10			
Competent Cells	20 reactions	C4040-03			
PureLink [™] HQ Mini Plasmid Purification Kit	100 preps	K2100-01			
PureLink [™] HiPure Plasmid Midiprep Kit	25 preps	K2100-04			
Lipofectamine [™] 2000	1.5 ml	11668-019			
	0.75 ml	11668-027			
Geneticin®	1 g	11811-023			
	5 g	11811-031			
Blasticidin	50 mg	R210-01			

Detection of Recombinant Proteins

You can detect expression of your recombinant fusion protein using the Anti-V5 antibodies available from Invitrogen. The amount of antibody supplied is sufficient for 25 Western blots or 25 immunostaining reactions (FITCconjugated antibody only).

Product	Epitope	Catalog no.
Anti-V5 Antibody	Detects 14 amino acid epitope	R960-25
Anti-V5-HRP Antibody	derived from the P and V	R961-25
Anti-V5-AP Antibody	SV5 (Southern <i>et al.</i> , 1991).	R962-25
Anti-V5-FITC Antibody	GKPIPNPLLGLDST	R963-25

Methods

Overview pcDNA[™]3.2/V5-DEST and pcDNA[™]6.2/V5-DEST are 7.7 kb Description and 7.3 kb vectors, respectively, that are adapted with the Gateway[®] Technology, and allow high-level, constitutive expression of the gene of interest in a variety of mammalian hosts. For more information on the Gateway® Technology, see the next page. pcDNA[™]3.2/V5-DEST and pcDNA[™]6.2/V5-DEST contain the Features following elements: Human cytomegalovirus immediate-early (CMV) promoter/enhancer for high-level expression in a wide range of mammalian cells Two recombination sites, attR1 and attR2, downstream of the CMV promoter for recombinational cloning of the gene of interest from an entry clone The *ccd*B gene located between the two *att*R sites for negative selection Chloramphenicol resistance gene located between the two attR sites for counterselection The V5 epitope tag for detection using Anti-V5 antibodies The Herpes Simplex Virus thymidine kinase polyadenylation signal for proper termination and processing of the recombinant transcript f1 intergenic region for production of single-strand DNA in F plasmid-containing E. coli SV40 early promoter and origin for expression of the neomycin (pcDNA[™]3.2/V5-DEST) or Blasticidin (pcDNA[™]6.2/V5-DEST) resistance gene and stable propagation of the plasmid in mammalian hosts expressing the SV40 large T antigen Neomycin (pcDNA[™]3.2/V5-DEST) or Blasticidin (pcDNA[™]6.2/V5-DEST) resistance gene for selection of stable cell lines The pUC origin for high copy replication and maintenance of the plasmid in E. coli The ampicillin (bla) resistance gene for selection in E. coli

Overview, continued

The Gateway[®] Technology

The Gateway[®] Technology is a universal cloning method 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.
- Generate an expression clone by performing an LR recombination reaction between the entry clone and a Gateway[®] destination vector (*e.g.* pcDNA[™]3.2/V5-DEST or pcDNA[™]6.2/V5-DEST).
- 3. Transfect your expression clone into the cell line of choice for transient or stable expression of your gene of interest.

For more information on the Gateway[®] Technology, refer to the Gateway[®] Technology with Clonase[™] II manual. This manual is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (page 25).

Generating an Entry Clone

Introduction	To recombine your gene of interest into pcDNA [™] 3.2/V5- DEST or pcDNA [™] 6.2/V5-DEST, you will need an entry clone containing the gene of interest (see below and the next page for recommendations). Many entry vectors including pENTR/D-TOPO [®] are available from Invitrogen to facilitate generation of entry clones. For more information, refer to our Web site (www.invitrogen.com) or contact Technical Service (page 25). Refer to the manual for the specific entry vector you are using for detailed instructions to construct an entry clone.
Tag-On- Demand [™] System	The pcDNA [™] 3.2/V5-DEST and pcDNA [™] 6.2/V5-DEST vectors are compatible with the Tag-On-Demand [™] System which allows expression of both native and C-terminally-tagged recombinant protein from the same expression construct.
	The System is based on stop suppression technology originally developed by RajBhandary and colleagues (Capone <i>et al.</i> , 1985) and consists of a recombinant adenovirus expressing a tRNA ^{ser} suppressor. When an expression vector encoding a gene of interest with the TAG (amber stop) codon is transfected into mammalian cells, the stop codon will be translated as serine, allowing translation to continue and resulting in production of a C-terminally-tagged fusion protein.
	For more information, refer to the Tag-On-Demand [™] Suppressor Supernatant manual. This manual is available for downloading from our Web site (www.invitrogen.com) or contact Technical Service (page 25).
Note	If you wish to express a human gene of interest from pcDNA [™] 3.2/V5-DEST or pcDNA [™] 6.2/V5-DEST, we recommend using an Ultimate [™] Human ORF (hORF) Clone available from Invitrogen. Each Ultimate [™] hORF Clone is a fully sequenced clone provided in a Gateway [®] entry vector that is ready-to-use in an LR recombination reaction with

pcDNA[™]3.2/V5-DEST or pcDNA[™]6.2/V5-DEST. In addition, each Ultimate[™] hORF Clone contains a **TAG** stop codon, making it fully compatible for use in the Tag-On-Demand[™] System. For more information about the Ultimate[™] hORF Clones available, refer to our Web site (www.invitrogen.com) or contact Technical Service (page 25).

Generating an Entry Clone, continued

Kozak Consensus Sequence	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. The ATG initiation codon is shown underlined.
	(G/A)NN <u>ATG</u> G
	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 consensus sequence.
Points to Consider Before Recombining	pcDNA [™] 3.2/V5-DEST and pcDNA [™] 6.2/V5-DEST are C-terminal fusion vectors; however, you may use these vectors to express native proteins or C-terminal fusion proteins. You may also use these vectors in the Tag-On- Demand [™] System (see previous page). Consider the following when generating your entry clone.

If you wish to	Then your insert						
include the V5 epitope tag and NOT use the Tag-On-Demand [™] System	 should NOT contain a stop codon should be in frame with the V5 epitope tag after recombination (see page 6 for a diagram) 						
include the V5 epitope tag for use in the Tag-On-Demand [™] System	 should contain a TAG stop codon should be in frame with the V5 epitope tag after recombination (see page 7 for a diagram) 						
not include the V5 epitope tag	should contain a stop codon						

Creating an Expression Clone

Introduction	After you have generated an entry clone, you will perform the LR recombination reaction to transfer the gene of interest into the pcDNA [™] 3.2/V5-DEST or pcDNA [™] 6.2/V5- DEST vector to create your expression clone. To ensure that you obtain the best results, we recommend that you read this section and the next section entitled Performing the LR Recombination Reaction (pages 8-11) before beginning.							
Experimental	To generate an expression clone, you will:							
Outline	1. Perform an LR recombination reaction using the <i>att</i> L- containing entry clone and the <i>att</i> R-containing pcDNA [™] 3.2/V5-DEST or pcDNA [™] 6.2/V5-DEST vector.							
	2. Transform the reaction mixture into a suitable <i>E. coli</i> host.							
	3. Select for expression clones (refer to pages 6-7 for a diagram of the recombination region of the resulting expression clones).							
Propagating the Vectors	If you wish to propagate and maintain pcDNA TM 3.2/V5- DEST or pcDNA TM 6.2/V5-DEST, we recommend using One Shot [®] ccdB Survival TM 2 T1 Phage-Resistant Cells (Catalog no. A10460) from Invitrogen for transformation. The ccdB 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 ccdB gene. To maintain the integrity of the vector, select for transformants in media containing 50– 100 µg/ml ampicillin and 15–30 µg/ml chloramphenicol. 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.							

Creating an Expression Clone, continued

Rec Rec	comb gion	inatio	on	Th fro pcl	e reo m p DNA	coml cDN A™6.	oinat IA™3 2/V§	tion 1 3.2/V 5-DE	n region of the expression clone resulting /V5-DEST × entry clone or DEST × entry clone is shown below.								
				Fea	atur	es of	f the	Reco	ombi	nati	on R	egio	n:				
	 Shaded regions correspond to DNA sequences transferred from the entry clone into pcDNA[™]3.2/V5- DEST or pcDNA[™]6.2/V5-DEST by recombination. Non- shaded regions are derived from the pcDNA[™]3.2/V5- DEST or pcDNA[™]6.2/V5-DEST vector. 														- on- -		
	• The underlined nucleotides flanking the shaded region correspond to bases 918 and 3161 of the pcDNA [™] 3.2/V5 DEST or pcDNA [™] 6.2/V5-DEST vector sequence.													on V5-			
771	CAAT											start T					
0.21	70707	ACCCI		مەرىپەر	T A C T		· ۳ ۳ ۳ ۸	TCCA		ר א א שי	T7 pror	noter/p		ite		VCCCZ	. 7.
031	AGAGA	ACCCF	1 010	JC I .	IACI	G G(JIIA	918	attB	1	ACGA	CICI	AC I A I	AG	GGAGA	ICCCF	171
891	GCTGG	CTAG	TAZ	AGC	ΓΑΤΟ	AA	CAAG	TT <u>T</u> G	F ACA		AAGC	AGG	CTN				NAC
		3	161	attB	2	ΤC	5110	AACI	A IG.		LICG	ICC	GAN		OFILE		NIG
3153	CCA G GGT C Pro A	CT TI CGA AA Ala Pi	AG AZ	IG I AC J Bu '	TAC ATG Tyr	AAA TTT Lys	GTG CAC Val	GTT CAA Val	GAT CTA Asp	CTA GAT Leu	GAG CTC Glu	GGC CCG Gly	CCG GGC Pro	CGG GCC Arg	TTC AAG Phe	GAA CTT Glu	
	V5	epitope	_			V5 re	verse p	riming s	site		-						
3201	GGT A Gly I	AG CO Jys Pi	CT A	rc (le 1	CCT Pro	AAC Asn	CCT Pro	CTC Leu	CTC Leu	GGT Gly	CTC Leu	GAT Asp	TCT Ser	ACG Thr	CGT Arg	ACC Thr	
3249	GGT T Gly *	'AG T#	AA T(GA (* *	GTTI	AAA	CGG (GGGA	GGCT2	AA CI	[GAA	ACAC	g ga <i>i</i>	AGGA	GACA		

Creating an Expression Clone, continued

Recombination Region for Use in the Tag-On-Demand[™] System The recombination region of the expression clone resulting from pcDNA[™]3.2/V5-DEST × entry clone or pcDNA[™]6.2/V5-DEST × entry clone is shown below.

Note: The gene of interest must contain a **TAG** stop codon for use in the Tag-On-Demand[™] System (see page 3 for more information).

Features of the Recombination Region:

- Shaded regions correspond to DNA sequences transferred from the entry clone into pcDNA[™]3.2/V5-DEST or pcDNA[™]6.2/V5-DEST by recombination. Nonshaded regions are derived from the pcDNA[™]3.2/V5-DEST or pcDNA[™]6.2/V5-DEST vector.
- The underlined nucleotides flanking the shaded region correspond to bases 918 and 3161 of the pcDNA[™]3.2/V5-DEST or pcDNA[™]6.2/V5-DEST vector sequence.

	CAAT										TATA 3'end of CMV promoter					Putative transcriptional start			
771	CAAATGGGCG GTAGGCGTGT					GT AC	CGGT	GGGA	G GT(TAT	FATATAA GCAGAGCTC					CTGGCTAACT			
831	AGAG	GAACO	CCA (CTGCI	TACI	rg go	T7 promoter/priming						iming s	site					
	918 attB 1																		
891	GCTGGCTAGT TAAGCTATC				CA AC	CAAG: GTTC2	rt <u>t</u> gi Aaacz	r acz	AAAA TTTTT	AAGC FTCG	AGG TCC	CTN GAN	GE	NE_	TAG ATC	NAC NTG			
			316 ⁻	1 _{at}	fB 2														
3153	CCA GGT Pro	GCT CGA Ala	TT <u>C</u> AAG Phe	TTG AAC Leu	TAC ATG Tyr	AAA TTT Lys	GTG CAC Val	GTT CAA Val	GAT CTA Asp	CTA GAT Leu	GAG CTC Glu	GGC CCG Gly	CCG GGC Pro	CGG GCC Arg	TTC AAG Phe	GAA CTT Glu			
	,	V5 epito	pe			V5 rev	verse p	riming s	ite		-								
3201	GGT Gly	AAG Lys	CCT Pro	ATC Ile	CCT Pro	AAC Asn	CCT Pro	CTC Leu	CTC Leu	GGT Gly	CTC Leu	GAT Asp	TCT Ser	ACG Thr	CGT Arg	ACC Thr			
3249	GGT Glv	TAG * * *	TAA * * *	TGA ***	GTTI	TAAA	CGG (GGGAC	GGCTI	AA CI	[GAA]	ACACO	g ga <i>i</i>	AGGA	GACA				

Performing the LR Recombination Reaction

Introduction	Once you have obtained an entry clone containing your gene of interest, you may perform an LR recombination reaction between the entry clone and pcDNA [™] 3.2/V5-DEST or pcDNA [™] 6.2/V5-DEST, and transform the reaction mixture into a suitable <i>E. coli</i> host (see below) to select for an expression clone. We recommend including a negative control (no LR Clonase [™] II) in your experiment to help you evaluate your results.
<i>E. coli</i> Host	You may use any <i>recA</i> , <i>endA E</i> . <i>coli</i> strain including TOP10, $DH5\alpha^{TM}$, or equivalent for transformation (see page vi for ordering information). Do not transform the LR reaction mixture into <i>E</i> . <i>coli</i> strains that contain the F' episome (<i>e.g.</i> TOP10F'). These strains contain the <i>ccdA</i> gene and will prevent negative selection with the <i>ccdB</i> gene.
Note	The presence of the EM7 promoter and the Blasticidin resistance gene in pcDNA [™] 6.2/V5-DEST allows for selection of <i>E. coli</i> transformants using Blasticidin. For selection, use Low Salt LB agar plates containing 100 µg/ml Blasticidin (see page 19 for a recipe). For Blasticidin to be active, the salt concentration of the medium must remain low (< 90 mM) and the pH must be 7.0. Blasticidin is available separately from Invitrogen (see page vi for ordering information). Refer to page 21 for instructions on how to prepare and store Blasticidin.

Performing the LR Recombination Reaction, continued

LR Clonase [™] II Enzyme Mix	LR Clonase [™] II enzyme mix is available separately from Invitrogen (Catalog no. 11791-020) to catalyze the LR recombination reaction. The LR Clonase [™] II enzyme mix combines the proprietary enzyme formulation and 5X LR Clonase [™] Reaction Buffer previously supplied as separate components in LR Clonase [™] enzyme mix into an optimized single-tube format for easier set-up of the LR recombination reaction. Use the protocol provided on page 10 to perform the LR recombination reaction using LR Clonase [™] II enzyme mix. Note: You may perform the LR recombination reaction using LR Clonase [™] enzyme mix, if desired. To use LR Clonase [™] enzyme mix, follow the protocol provided with the product. Do not use the protocol for LR Clonase [™] II enzyme mix as reaction conditions differ.	
Materials Needed	You should have the following materials on hand before beginning:	
	 Purified plasmid DNA of your entry clone (50– 150 ng/μl in TE, pH 8.0) 	
	 pcDNA[™]3.2/V5-DEST or pcDNA[™]6.2/V5-DEST (150 ng/µl in TE, pH 8.0) 	
	 LR Clonase[™] II enzyme mix (Invitrogen, Catalog no. 11791-020; keep at -20°C until immediately before use) 	
	• TE Buffer, pH 8.0 (10 mM Tris-HCl, pH 8.0, 1 mM EDTA)	
	 2 µg/µl Proteinase K solution (supplied with LR Clonase[™] II enzyme mix; thaw and keep on ice until use) 	
	 pENTR[™]-gus (supplied with LR Clonase[™] II enzyme mix; use as a control for the LR reaction; 50 ng/µl) 	
	• Appropriate competent <i>E. coli</i> host and growth media for expression	
	• S.O.C. Medium	
	• LB agar plates containing 100 μ g/ml ampicillin or Low Salt LB plates containing 100 μ g/ml Blasticidin)	

Performing the LR Recombination Reaction, continued

Setting Up the LR Reaction

Follow this procedure to perform the LR reaction between your entry clone and a destination vector. To include a negative control, set up a second sample reaction, but omit the LR Clonase[™] II enzyme mix.

 Add the following components to 1.5 ml microcentrifuge tubes at room temperature and mix.

Component	Sample	Positive Control
Entry clone (50–150 ng/rxn)	1–7 µl	
Destination vector (150 ng/µl)	1 µl	1 μl
pENTR [™] -gus (50 ng/µl)		2 µl
TE Buffer, pH 8.0	to 8 µl	5 µl

- Remove the LR Clonase[™] II enzyme mix from -20°C and thaw on ice (~ 2 minutes).
- Vortex the LR Clonase[™] II enzyme mix briefly twice (2 seconds each time).
- To each sample above, add 2 µl of LR Clonase[™] II enzyme mix. Mix well by pipetting up and down.
 Reminder: Return LR Clonase[™] II enzyme mix to -20°C immediately after use.
- 5. Incubate reactions at 25°C for 1 hour.

Note: Extending the incubation time to 18 hours typically yields more colonies.

- Add 1 μl of the Proteinase K solution to each reaction. Incubate for 10 minutes at 37°C.
- Transform 1 μl of the LR recombination reaction into a suitable *E. coli* host (follow the manufacturer's instructions) and select for expression clones.

Note: You may store the LR reaction at -20°C for up to 1 week before transformation, if desired.

Creating an Expression Clone, continued

What You Should See	If you use <i>E. coli</i> cells with a transformation efficiency of $\ge 1 \times 10^8$ cfu/µg, the LR reaction should give > 5,000 colonies if the entire reaction is transformed and plated.		
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.		
Sequencing	To confirm that your gene of interest is in frame with the C-terminal V5 epitope, you may sequence your expression construct, if desired. We suggest using the following primer sequences. Refer to the diagram on page 6 for the location of the primer binding sites. For your convenience, Invitrogen offers a custom primer synthesis service. For more information, refer to our Web		
	site (www.invitrogen.com) or contact Technical Service (page 25).		
	Primer	Sequence	
	T7 Promoter	5'-TAATACGACTCACTATAGGG-3'	
	V5 Reverse	5'-ACCGAGGAGAGGGTTAGGGAT-3'	

Transfection

Introduction	This section provides general information for transfecting your expression clone into the mammalian cell line of choice. We recommend that you include a positive control vector (pcDNA [™] 3.2/V5/GW/CAT or pcDNA [™] 6.2/V5/GW/CAT) and a mock transfection (negative control) in your experiments to evaluate your results.
Plasmid Preparation	Once you have generated your expression clone, you must isolate plasmid DNA for transfection. Plasmid DNA for transfection into eukaryotic cells must be clean and free contamination with 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 PureLink [™] HQ Mini Plasmid Purification Kit (Catalog no. K2100-01), the PureLink [™] HiPure Plasmid Midiprep Kit (Catalog no. K2100-04), or CsCl gradient centrifugation.
Methods of Transfection	For established cell lines (<i>e.g.</i> HeLa), consult original references or the supplier of your cell line for the optimal method of transfection. We recommend that you follow exactly the protocol for your cell line. Pay particular attention to medium requirements, when to pass the cells, and at what dilution to split the cells. Further information is provided in <i>Current Protocols in Molecular Biology</i> (Ausubel <i>et al.</i> , 1994).
	Methods for transfection include calcium phosphate (Chen and Okayama, 1987; Wigler <i>et al.</i> , 1977), lipid-mediated (Felgner <i>et al.</i> , 1989; Felgner and Ringold, 1989) and electroporation (Chu <i>et al.</i> , 1987; Shigekawa and Dower, 1988). For high efficiency transfection in a broad range of mammalian cell lines, we recommend using Lipofectamine [™] 2000 Reagent (Catalog no. 11668-027) available from Invitrogen. For more information about Lipofectamine [™] 2000 and other transfection reagents, refer to our Web site (www.invitrogen.com) or contact Technical Service (page 25).

Transfection, continued

Positive Control pcDNA[™]3.2/V5/GW/CAT or pcDNA[™]6.2/V5/GW/CAT is provided as a positive control vector for mammalian cell transfection and expression (see page 24 for a map) and may be used to optimize recombinant protein expression levels in your cell line. These vectors allow expression of a C-terminally tagged chloramphenicol acetyl transferase (CAT) fusion protein that may be detected by Western blot or functional assay.

To propagate and maintain the plasmid:

- Prepare a 1:50 dilution of the positive control vector in sterile water (i.e. 1 µl vector + 49 µl water) for a 10 ng/ul stock solution. Use 10 ng of the stock solution to transform a *recA*, *endA E. coli* strain like TOP10, DH5α, JM109, or equivalent.
- Select transformants on LB agar plates containing 50– 100 μg/ml ampicillin.
- 3. Prepare a glycerol stock of a transformant containing plasmid for long-term storage.

Expression and Analysis

Introduction	Expression of your gene of interest from the expression clone can be performed in either transiently transfected cells or stable cell lines (see page 16 for guidelines to create stable cell lines). You may use a functional assay or a Western blot analysis to detect your recombinant protein (see below).		
Preparing Cell Lysates	To detect your fusion protein by Western blot, you to prepare a cell lysate from transfected cells. A san protocol is provided below. Other protocols are sui lyse cells:		
	1.	Wash cell monolayer (\sim 5 x 10 ⁵ to 1 x 10 ⁶ cells) once with phosphate-buffered saline (PBS; Invitrogen Catalog no. 10010-023).	
	2.	Scrape cells into 1 ml PBS and pellet the cells at $1500 \times g$ for 5 minutes.	
	3.	Resuspend in 50 µl Cell Lysis Buffer (see page 20 for a recipe). Other cell lysis buffers are suitable. Vortex.	
	4.	Incubate cell suspension at 37°C for 10 minutes to lyse the cells. Note: You may prefer to lyse the cells at room temperature or on ice if degradation of your protein is a potential problem.	
	5.	Centrifuge the cell lysate at 10,000 x g for 10 minutes at +4°C to pellet nuclei and transfer the supernatant to a fresh tube. Assay the lysate for protein concentration. Note: Do not use protein assays utilizing Coomassie Blue or other dyes. NP-40 interferes with the binding of the dye with the protein.	
	6.	Add SDS-PAGE sample buffer (see page 20 for a recipe) to a final concentration of 1X and boil the sample for 5 minutes.	
	7.	Load 20 μ g of lysate onto an SDS-PAGE gel and electrophorese. Use the appropriate percentage of acrylamide to resolve your fusion protein.	

Expression and Analysis, continued

Polyacrylamide Gel Electrophoresis	To facilitate separation and visualization of your recombinar fusion protein by polyacrylamide gel electrophoresis, a wide range of pre-cast NuPAGE [®] and Novex [®] Tris-Glycine polyacrylamide gels and electrophoresis apparatus are available from Invitrogen. For more information, refer to our Web site (www.invitrogen.com) or contact Technical Service (page 25).	
Detecting Recombinant Fusion Proteins	To detect expression of your recombinant fusion protein by Western blot analysis, you may use the Anti-V5 antibodies available from Invitrogen (see page vi 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. The ready-to-use WesternBreeze [®] Chromogenic Kits and WesternBreeze [®] Chemiluminescent Kits are available from Invitrogen to facilitate detection of antibodies by colorimetric or chemiluminescent methods. For more information, refer to our Web site (www.invitrogen.com) or contact Technical Service (page 25).	
Note	The C-terminal peptide containing the V5 epitope will add approximately 4 kDa to your protein.	
Detecting CAT Protein	If you use the provided positive control vector in your experiment, you may assay for CAT expression using your method of choice. Note that CAT is fused to the C-terminal V5 epitope tag so you can use Western blot analysis and an Anti-V5 antibody to detect expression of CAT. Other commercial kits are available for assaying CAT expression. The molecular weight of the CAT fusion protein is approximately 30 kDa.	

Creating Stable Cell Lines

Introduction

The pcDNA[™]3.2/V5-DEST and pcDNA[™]6.2/V5-DEST vectors contain the neomycin and Blasticidin resistance genes, respectively, to allow selection of stable cell lines. If you wish to create stable cell lines, transfect your construct into the mammalian cell line of choice and select for foci using Geneticin[®] (pcDNA[™]3.2/V5-DEST only) or Blasticidin (pcDNA[™]6.2/V5-DEST only). General information and guidelines are provided below.



To obtain stable transfectants, we recommend that you linearize your pcDNA[™]3.2/V5-DEST or pcDNA[™]6.2/V5-DEST construct before transfection. While linearizing the vector may not improve the efficiency of transfection, it increases the chances that the vector does not integrate in a way that disrupts elements necessary for expression in mammalian cells. To linearize your construct, cut at a unique site that is not located within a critical element or within your gene of interest.

Geneticin[®]

Geneticin[®] blocks protein synthesis in mammalian cells by interfering with ribosomal function. It is an aminoglycoside, similar in structure to neomycin, gentamycin, and kanamycin. Expression in mammalian cells of the bacterial aminoglycoside phosphotransferase gene (APH), derived from Tn*5*, results in detoxification of Geneticin[®] (Southern and Berg, 1982).

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 nontoxic deaminohydroxy derivative (Izumi *et al.*, 1991).

Creating Stable Cell Lines, continued

Determining Antibiotic Sensitivity

To successfully generate a stable cell line expressing your protein of interest, you need to determine the minimum concentration of antibiotic (Geneticin[®] or Blasticidin) required to kill your untransfected host cell line. Test a range of concentrations (see protocol below) to ensure that you determine the minimum concentration necessary for your cell line. Refer to page 21 for instructions on how to prepare and store Blasticidin.

- Plate or split a confluent plate so the cells will be approximately 25% confluent. For each antibiotic, prepare a set of 6–7 plates. Add the following concentrations of antibiotic to each plate:
 - For Blasticidin selection, test 0, 1, 3, 5, 7.5, and 10 $\mu g/ml$ Blasticidin
 - For Geneticin[®] selection, test 0, 50, 125, 250, 500, 750, and 1000 μg/ml Geneticin[®].
- 2. Replenish the selective media every 3-4 days, and observe the percentage of surviving cells.
- 3. Count the number of viable cells at regular intervals to determine the appropriate concentration of antibiotic that prevents growth within 1–3 weeks after addition of the antibiotic.

Geneticin[®] Selection Guidelines

Once you have determined the appropriate Geneticin[®] concentration to use for selection, you can generate a stable cell line expressing your **pcDNA**[™]**3.2/V5-DEST** construct. Geneticin[®] is available separately from Invitrogen (see page vi for ordering information). Use as follows:

- 1. Prepare Geneticin[®] in a buffered solution (*e.g.* 100 mM HEPES, pH 7.3).
- 2. Use the predetermined concentration of Geneticin[®] in complete medium.
- 3. Calculate concentration based on the amount of active drug.
- 4. Cells will divide once or twice in the presence of lethal doses of Geneticin[®], so the effects of the drug take several days to become apparent. Complete selection can take from 2 to 3 weeks of growth in selective medium.

Creating Stable Cell Lines, continued

Blasticidin Selection Guidelines

Once you have determined the appropriate Blasticidin concentration to use for selection, you can generate a stable cell line expressing your **pcDNA**[™]**6.2/V5-DEST** construct. Blasticidin is available separately from Invitrogen (see page vi for ordering information). Use as follows:

- 1. Prepare a stock solution of 5–10 mg/ml of Blasticidin in sterile water. Filter-sterilize the solution.
- 2. Use the predetermined concentration of Blasticidin in complete medium.
- 3. Cells differ in their susceptibility to Blasticidin. Complete selection can take up to 10 days of growth in selective medium.

Refer to page 21 for instructions on how to prepare and store Blasticidin.

Appendix

Recipes

LB (Luria- Bertani) Medium and Plates	Composition: 1.0% Tryptone 0.5% Yeast Extract 1.0% NaCl pH 7.0	
	 For 1 liter, dissolve 10 g tryptone, 5 g yeast extract, and 10 g NaCl in 950 ml deionized water. 	
	2. Adjust the pH of the solution to 7.0 with NaOH and bring the volume up to 1 liter.	
	3. Autoclave on liquid cycle for 20 minutes at 15 psi. Allow solution to cool to 55°C and add antibiotic if needed.	
	4. Store at room temperature or at $+4^{\circ}$ C.	
	LB agar plates1. Prepare LB medium as above, but add 15 g/L agar before autoclaving.	
	2. Autoclave on liquid cycle for 20 minutes at 15 psi.	
	3. After autoclaving, cool to ~55°C, add antibiotic if needed, and pour into 10 cm plates.	
	4. Let harden, then invert and store at +4°C.	
Low Salt LB Medium with Blasticidin	Low Salt LB Medium: 10 g Tryptone 5 g NaCl 5 g Yeast Extract	
	 Combine the dry reagents above and add deionized, distilled water to 950 ml. Adjust pH to 7.0 with 1 N NaOH. Bring the volume up to 1 liter. For plates, add 15 g/L agar before autoclaving. 	
	 Autoclave on liquid cycle at 15 psi and 121°C for 20 minutes. 	
	3. Allow the medium to cool to at least 55°C before adding the Blasticidin to $100 \mu g/ml$ final concentration.	
	 Store plates at +4°C in the dark. Plates containing Blasticidin are stable for up to 2 weeks. 	

Recipes, continued

Cell Lysis Buffer	50 mM Tris, pH 7.8 150 mM NaCl 1% Nonidet P-40				
	1. This stoc	. This solution can be prepared from the following c stock solutions. For 100 ml, combine			
	1 M 5 M Nor	Tris base NaCl iidet P-40	5 ml 3 ml 1 ml		
	2. Brin adju	ig the volume up to ist the pH to 7.8 with	90 ml with deionized water and h HCl.		
	3. Brin	g the volume up to	100 ml. Store at room temperature.		
	To prevent proteolysis, you may add 1 mM PMSF, 1 μM leupeptin, or 0.1 μM aprotinin before use.				
4X SDS-PAGE	1. Con	nbine the following	reagents:		
Sample Buffer	0.5 M Glyo β-m Bron SDS	M Tris-HCl, pH 6.8 cerol (100%) ercaptoethanol mophenol Blue	5 ml 4 ml 0.8 ml 0.04 g 0.8 g		
	2. Brin	ig the volume to 10 i	nl with sterile water.		
	3. Alic	µuot and freeze at −2	0°C until needed.		

Blasticidin

Molecular Weight, Formula, and Structure The formula for Blasticidin S is $C_{17}H_{26}N_8O_5$ -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 (<i>e.g.</i> a laboratory coat) when handling Blasticidin. Weigh out Blasticidin and prepare solutions in a hood.		
Preparing and Storing Stock Solutions	Blasticidin may be obtained separately from Invitrogen (Catalog no. R210-01) in 50 mg aliquots. Blasticidin is soluble in water. Sterile water is generally used to prepare stock solutions of 5 to 10 mg/ml.		
	• Dissolve Blasticidin in sterile water and filter-sterilize the solution.		
	• Aliquot in small volumes suitable for one time use (see next to 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.		
	• pH of the aqueous solution should be 7.0 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 store the thawed stock solution at +4°C for up to 2 weeks. 		
	• Medium containing Blasticidin may be stored at +4°C for up to 2 weeks.		

Map of pcDNA[™]3.2/V5-DEST and pcDNA[™]6.2/V5-DEST

Map

The map below shows the elements of pcDNA[™]3.2/V5-DEST and pcDNA[™]6.2/V5-DEST. DNA from the entry clone replaces the region between bases 918 and 3161. The complete sequences of these vectors are available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (page 25).



continued on next page

f1 origin:

Features of pcDNA[™]3.2/V5-DEST and pcDNA[™]6.2/V5-DEST

FeaturespcDNA[™]3.2/V5-DEST (7711 bp) and pcDNA[™]6.2/V5-DEST
(7341 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 <i>in vitro</i> transcription in the sense orientation and sequencing through the insert
attR1 and attR2 sites	Allows recombinational cloning of the gene of interest from an entry clone
<i>ccd</i> B gene	Allows negative selection of plasmid
Chloramphenicol resistance gene	Allows counterselection of plasmid
V5 epitope	Allows detection of the recombinant fusion protein by the Anti-V5 antibodies (Southern <i>et al.</i> , 1991).
V5 reverse priming site	Allows sequencing of the insert
Herpes Simplex Virus Thymidine Kinase (TK) polyadenylation signal	Allows efficient transcription termination and polyadenylation of mRNA (Cole and Stacy, 1985)
f1 origin	Allows rescue of single-stranded DNA
SV40 early promoter and origin	Allows efficient, high-level expression of the neomycin or Blasticidin resistance gene and episomal replication in cells expressing the SV40 large T antigen
Neomycin resistance gene (pcDNA [™] 3.2/V5-DEST only)	Allows selection of stable transfectants in mammalian cells (Southern and Berg, 1982)
EM7 promoter (pcDNA [™] 6.2/V5-DEST only)	Allows expression of the Blasticidin resistance gene in <i>E. coli</i>
Blasticidin (<i>bsd</i>) resistance gene (pcDNA [™] 6.2/V5-DEST only)	Allows selection of stable transfectants in mammalian cells (Kimura <i>et al.,</i> 1994)
SV40 early polyadenylation signal	Allows efficient transcription termination and polyadenylation of mRNA
pUC origin	Allows high-copy number replication and growth in <i>E. coli</i>
Ampicillin (<i>bla</i>) resistance gene (β-lactamase)	Allows selection of transformants in <i>E. coli</i>

Map of pcDNA[™]3.2/V5/GW/CAT and pcDNA[™]6.2/V5/GW/CAT

Description

pcDNA[™]3.2/V5/GW/CAT (6188 bp) and pcDNA[™]6.2/V5/ GW/CAT (5818 bp) are control vectors expressing chloramphenicol acetyltransferase (CAT). Each vector was constructed using the LR recombination reaction between an entry clone containing the CAT gene and the respective destination vector. **Note:** The CAT gene is in frame with the C-terminal V5 epitope and does not contain a stop codon. The molecular weight of the CAT fusion protein is ~30 kDa.

Мар

The map below shows the elements of pcDNA[™]3.2/V5/GW/ CAT and pcDNA[™]6.2/V5/GW/CAT. The complete sequences of these vectors are available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (page 25).



Technical Service

World Wide Web	 Visit the Invitrogen website at <u>www.invitrogen.com</u> for: Technical resources, including manuals, vector maps and sequences, application notes, MSDSs, FAQs, formulations, citations, handbooks, etc. Complete technical support contact information Access to the Invitrogen Online Catalog Additional product information and special offers 		
Contact Us	For more information or technical assistance, call, write, for email. Additional international offices are listed on our Web page (www.invitrogen.com).Corporate Headquarters:European Headquarters:Invitrogen CorporationInvitrogen Ltd5791 Van Allen WayInchinnan Business ParkCarlsbad, CA 92008 USA3 Fountain DriveTel: 1 760 603 7200Paisley PA4 9RF, UKTel (Toll Free): 1 800 955 6288Tel: +44 (0) 141 814 6100Fax: 1 760 602 6500Tech Fax: +44 (0) 141 814 6100E-mail:E-mail:techsupport@invitrogen.comE-mail:		
MSDS	Material Safety Data Sheets (M website at <u>www.invitrogen.co</u>	ISDSs) are available on our m/msds.	

Technical Service, continued

Certificate of Analysis	The Certificate of Analysis (CofA) provides detailed quality control information for each product. CofAs are available on our website at <u>www.invitrogen.com/support</u> , and are searchable by product lot number, which is printed on each box.
Limited Warranty	Invitrogen 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 Service 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. <u>This warranty limits Invitrogen Corporation's liability only to the cost of the product</u> . 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 Service 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

Introduction

Use of the pcDNA[™]3.2/V5-DEST and pcDNA[™]6.2/V5-DEST Gateway[®] Vectors is covered under the licenses detailed below.

Limited Use Label License No. 19: Gateway[®] Cloning Products The purchase of this product conveys to the buyer the nontransferable 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 Life Technologies 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 ClonaseTM purchased from Life Technologies 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 Life Technologies 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.

Purchaser Notification, continued

Limited Use Label License No. 19: Gateway [®] Cloning Products, continued	Life Technologies 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 de- veloped 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. Life Technologies 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, Life Technologies is willing to accept return of the product with a full refund. For information on purchasing a license to
Gateway [®] Clone Distribution Policy	For additional information about Invitrogen's policy for the use and distribution of Gateway [®] clones, see the section entitled Gateway[®] Clone Distribution Policy , page 29.
Limited Use Label License No. 51: Blastici- din & the Blasti- cidin 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.

Gateway[®] Clone Distribution Policy

Introduction	The information supplied in this section is intended to provide clarity concerning Invitrogen's policy for the use and distribution of cloned nucleic acid fragments, including open reading frames, created using Invitrogen's commercially available Gateway [®] Technology.
Gateway [®] Entry Clones	Invitrogen understands that Gateway [®] entry clones, containing <i>att</i> L1 and <i>att</i> L2 sites, may be generated by academic and government researchers for the purpose of scientific research. Invitrogen agrees that such clones may be distributed for scientific research by non-profit organizations and by for-profit organizations without royalty payment to Invitrogen.
Gateway [®] Expression Clones	Invitrogen also understands that Gateway [®] expression clones, containing <i>att</i> B1 and <i>att</i> B2 sites, may be generated by academic and government researchers for the purpose of scientific research. Invitrogen agrees that such clones may be distributed for scientific research by academic and government organizations without royalty payment to Invitrogen. Organizations other than academia and government may also distribute such Gateway [®] expression clones for a nominal fee (\$10 per clone) payable to Invitrogen.
Additional Terms and Conditions	We would ask that such distributors of Gateway [®] entry and expression clones indicate that such clones may be used only for research purposes, that such clones incorporate the Gateway [®] Technology, and that the purchase of Gateway [®] Clonase [™] from Invitrogen is required for carrying out the Gateway [®] recombinational cloning reaction. This should allow researchers to readily identify Gateway [®] containing clones and facilitate their use of this powerful technology in their research. Use of Invitrogen's Gateway [®] Technology, including Gateway [®] clones, for purposes other than scientific research may require a license and questions concerning such commercial use should be directed to Invitrogen's licensing department at 760-603-7200.

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.

Capone, J. P., Sharp, P. A., and RajBhandary, U. L. (1985). Amber, Ochre and Opal Suppressor tRNA Genes Derived from a Human Serine tRNA Gene. EMBO J. 4, 213-221.

Chen, C., and Okayama, H. (1987). High-Efficiency Transformation of Mammalian Cells by Plasmid DNA. Mol. 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.

Cole, C. N., and Stacy, T. P. (1985). Identification of Sequences in the Herpes Simplex Virus Thymidine Kinase Gene Required for Efficient Processing and Polyadenylation. Mol. Cell. Biol. *5*, 2104-2113.

Felgner, P. L., Holm, M., and Chan, H. (1989). Cationic Liposome Mediated Transfection. Proc. West. Pharmacol. Soc. 32, 115-121.

Felgner, P. L. a., and Ringold, G. M. (1989). Cationic Liposome-Mediated Transfection. Nature 337, 387-388.

Izumi, M., Miyazawa, H., Kamakura, T., Yamaguchi, I., Endo, T., and Hanaoka, F. (1991). Blasticidin 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). Blasticidin S Deaminase Gene from *Aspergillus terreus* (*BSD*): A New Drug Resistance Gene for Transfection of Mammalian Cells. Biochim. Biophys. ACTA *1219*, 653-659.

References, continued

Kozak, M. (1987). An Analysis of 5´-Noncoding Sequences from 699 Vertebrate Messenger RNAs. Nucleic Acids Res. *15*, 8125-8148.

Kozak, M. (1991). An Analysis of Vertebrate mRNA Sequences: Intimations of Translational Control. J. Cell Biology 115, 887-903.

Kozak, M. (1990). Downstream Secondary Structure Facilitates Recognition of Initiator Codons by Eukaryotic Ribosomes. Proc. Natl. Acad. Sci. USA *87*, 8301-8305.

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

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.

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.

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.

Southern, P. J., and Berg, P. (1982). Transformation of Mammalian Cells to Antibiotic Resistance with a Bacterial Gene Under Control of the SV40 Early Region Promoter. J. Molec. Appl. Gen. 1, 327-339.

Takeuchi, S., Hirayama, K., Ueda, K., Sakai, H., and Yonehara, H. (1958). Blasticidin S, A New Antibiotic. The Journal of Antibiotics, Series A 11, 1-5.

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.

Yamaguchi, H., Yamamoto, C., and Tanaka, N. (1965). Inhibition of Protein Synthesis by Blasticidin S. I. Studies with Cell-free Systems from Bacterial and Mammalian Cells. J. Biochem (Tokyo) 57, 667-677.

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

^{©2002–2008, 2010} 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