

Instruction Manual

Mammalian Lumio[™] Gateway[®] Vectors with Lumio[™] In-Cell Labeling Kits

Gateway[®] destination vectors for site-specific fluorescence labeling and detection of proteins in live mammalian cells

Catalog nos. 12589-016, 12589-024, and 12589-032

Version D 9 December 2010 *25-0699*

Table of Contents

Table of Contents .	iii
Kit Contents and Storage	v
Introduction	1
Overview.	1
Methods	4
Generating an Entry Clone .	4
Creating an Expression Clone	6
Performing the LR Recombination Reaction.	9
Transfecting Cells	
Creating Stable Cell Lines.	
Detecting Lumio [™] Fusion Proteins	
Appendix	
Recipes	
Blasticidin	
Map and Features of pcDNA [™] 6.2/cLumio [™] -DEST	20
Map and Features of pcDNA ^{TM} 6.2/nLumio ^{TM} -DEST	
Map of pcDNA [™] 6.2/nLumio [™] -GW/p64	24
Technical Service	
Purchaser Notification .	
Gateway [®] Clone Distribution Policy	
References	

Kit Contents and Storage

Types of Kits This manual is supplied with the following products. All products are also supplied with the Lumio[™] In-Cell Labeling Kits manual.

Product	Catalog no.
Mammalian Lumio [™] Gateway [®] Vectors with Lumio [™] Green In-Cell Labeling Kit	12589-016
Mammalian Lumio [™] Gateway [®] Vectors with Lumio [™] Red In-Cell Labeling Kit	12589-024
Mammalian Lumio [™] Gateway [®] Vectors with Lumio [™] Dual Green and Red In-Cell Labeling Kits	12589-032

Kit Components The Mammalian Lumio[™] Gateway[®] Vectors with Lumio[™] In-Cell Labeling Kits include the following components. For a detailed description of the contents of each component, see the next page.

Component	Catalog no.		
	<u>12589-016</u>	<u>12589-024</u>	<u>12589-032</u>
Mammalian Lumio [™] Gateway [®] Vectors	\checkmark	\checkmark	\checkmark
Lumio™ Green In-Cell Labeling Kit	\checkmark		\checkmark
Lumio [™] Red In-Cell Labeling Kit		\checkmark	\checkmark

Shipping and Storage

The Mammalian Lumio[™] Gateway[®] Vectors with Lumio[™] In-Cell Labeling Kits are shipped on blue ice. Upon receipt, store as detailed below.

Box	Storage
Mammalian Lumio [™] Gateway [®] Vectors	-20°C
Lumio [™] Green In-Cell Labeling Kit	-20°C, protected from light
Lumio [™] Red In-Cell Labeling Kit	-20°C, protected from light

Lumio[™] Gateway[®] The Lumio[™] Gateway[®] Vectors are listed below. Store vectors at -20°C. Vectors

Item	Concentration	Volume
pcDNA [™] 6.2/cLumio [™] -DEST	Lyophilized in TE, pH 8.0	6 µg
pcDNA [™] 6.2/nLumio [™] -DEST	Lyophilized in TE, pH 8.0	6 µg
pcDNA [™] 6.2/nLumio [™] -GW/p64 Control Plasmid	Lyophilized in TE, pH 8.0	10 µg

Kit Contents and Storage, continued

Blasticidin

Lumio [™] In-Cell Labeling Kits	The Mammalian Lumio [™] Gateway [®] Vectors are supplied with either the Lumio [™] Green, the Lumio [™] Red, or both the Lumio [™] Green and Lumio [™] Red In-Cell Labeling Kits. Refer to the Lumio [™] In-Cell Labeling Kits manual for detailed information pertaining to each kit and a description of the reagents provided in each kit.		
Product Qualification	pcDNA TM 6.2/cLumio TM -DEST, pcDNA TM 6.2/nLumio TM -DEST, and pcDNA TM 6.2/ nLumio TM -GW/p64 are qualified by restriction endonuclease digestion. The pcDNA TM 6.2/cLumio TM -DEST and pcDNA TM 6.2/nLumio TM -DEST vectors are further qualified in a recombination assay using Gateway [®] LR Clonase TM II enzyme mix. The <i>ccd</i> B gene is assayed by transformation using an appropriate <i>E. coli</i> strain.		
	For information about how the components o are qualified, refer to the Lumio [™] In-Cell Labo		ell Labeling Kits
Accessory	Additional products that may be used with the Mammalian Lumio [™] Gateway [®] Vectors are available from Invitrogen. Ordering information is provided below.		
Products		ng information is p	
		ng information is p Amount	
	Vectors are available from Invitrogen. Orderin		provided below.
	Vectors are available from Invitrogen. Orderin Product	Amount	Catalog no.
	Vectors are available from Invitrogen. Orderin Product Lumio [™] Green In-Cell Labeling Kit	Amount 1 kit	Catalog no.12589-057
	Vectors are available from Invitrogen. Orderin Product Lumio [™] Green In-Cell Labeling Kit Lumio [™] Red In-Cell Labeling Kit	Amount Amount 1 kit 1 kit	Catalog no. 12589-057 12589-040
	Vectors are available from Invitrogen. Orderin Product Lumio [™] Green In-Cell Labeling Kit Lumio [™] Red In-Cell Labeling Kit Lumio [™] Green Detection Kit	Amount1 kit1 kit100 reactions	Catalog no. 12589-057 12589-040 LC6090
	Vectors are available from Invitrogen. Orderin Product Lumio [™] Green In-Cell Labeling Kit Lumio [™] Red In-Cell Labeling Kit Lumio [™] Green Detection Kit BenchMark [™] Fluorescent Protein Standard	Amount1 kit1 kit100 reactions100 reactions	Catalog no. 12589-057 12589-040 LC6090 LC5928
	Vectors are available from Invitrogen. Orderin Product Lumio [™] Green In-Cell Labeling Kit Lumio [™] Red In-Cell Labeling Kit Lumio [™] Green Detection Kit BenchMark [™] Fluorescent Protein Standard	Amount1 kit1 kit100 reactions100 reactions20 reactions	Catalog no. 12589-057 12589-040 LC6090 LC5928 11791-020
	Vectors are available from Invitrogen. Orderin Product Lumio [™] Green In-Cell Labeling Kit Lumio [™] Red In-Cell Labeling Kit Lumio [™] Green Detection Kit BenchMark [™] Fluorescent Protein Standard Gateway [®] LR Clonase [™] II Enzyme Mix	Amount1 kit1 kit1 kit100 reactions100 reactions20 reactions100 reactions100 reactions	Catalog no. 12589-057 12589-040 LC6090 LC5928 11791-020 11791-100
	Vectors are available from Invitrogen. Orderin Product Lumio [™] Green In-Cell Labeling Kit Lumio [™] Red In-Cell Labeling Kit Lumio [™] Green Detection Kit BenchMark [™] Fluorescent Protein Standard Gateway [®] LR Clonase [™] II Enzyme Mix One Shot [®] TOP10 Chemically Competent	Amount1 kit1 kit1 kit100 reactions100 reactions20 reactions100 reactions100 reactions100 reactions10 reactions	Catalog no. 12589-057 12589-040 LC6090 LC5928 11791-020 11791-100 C4040-10
	Vectors are available from Invitrogen. Orderin Product Lumio [™] Green In-Cell Labeling Kit Lumio [™] Red In-Cell Labeling Kit Lumio [™] Green Detection Kit BenchMark [™] Fluorescent Protein Standard Gateway [®] LR Clonase [™] II Enzyme Mix One Shot [®] TOP10 Chemically Competent Cells	Amount1 kit1 kit1 kit100 reactions100 reactions20 reactions100 reactions100 reactions20 reactions20 reactions20 reactions	Catalog no. 12589-057 12589-040 LC6090 LC5928 11791-020 11791-100 C4040-10 C4040-03
	Vectors are available from Invitrogen. Orderin Product Lumio [™] Green In-Cell Labeling Kit Lumio [™] Red In-Cell Labeling Kit Lumio [™] Green Detection Kit BenchMark [™] Fluorescent Protein Standard Gateway [®] LR Clonase [™] II Enzyme Mix One Shot [®] TOP10 Chemically Competent Cells	Amount1 kit1 kit1 kit100 reactions100 reactions20 reactions100 reactions10 reactions20 reactions10 reactions10 reactions10 reactions10 reactions10 reactions	Catalog no. 12589-057 12589-040 LC6090 LC5928 11791-020 11791-100 C4040-10 C4040-03 C4040-50
	Vectors are available from Invitrogen. Orderin Product Lumio [™] Green In-Cell Labeling Kit Lumio [™] Red In-Cell Labeling Kit Lumio [™] Green Detection Kit BenchMark [™] Fluorescent Protein Standard Gateway [®] LR Clonase [™] II Enzyme Mix One Shot [®] TOP10 Chemically Competent Cells One Shot [®] TOP10 Electrocompetent Cells	Amount1 kit1 kit1 kit100 reactions100 reactions20 reactions100 reactions10 reactions20 reactions10 reactions20 reactions20 reactions20 reactions20 reactions20 reactions20 reactions20 reactions	Catalog no. 12589-057 12589-040 LC6090 LC5928 11791-020 11791-100 C4040-10 C4040-50 C4040-52

R210-01

50 mg

Introduction

Overview	
Description	The Mammalian Lumio [™] Gateway [®] Vector Kits contain Gateway [®] -adapted destination vectors designed for use with the Lumio [™] Technology. The pcDNA [™] 6.2/Lumio [™] -DEST vectors supplied with each kit facilitate <i>in vivo</i> fluorescence labeling and detection of recombinant proteins when used with a Lumio [™] In-Cell Labeling Kit.
Features	The pcDNA [™] 6.2/cLumio [™] -DEST and pcDNA [™] 6.2/nLumio [™] -DEST vectors contain the following elements:
	 Human cytomegalovirus immediate-early (CMV) promoter/enhancer for high-level expression in a wide range of mammalian cells
	 Lumio[™] tag for C-terminal (pcDNA[™]6.2/cLumio[™]-DEST) or N-terminal (pcDNA[™]6.2/nLumio[™]-DEST) fusion to the gene of interest for fluorescence detection
	• Two recombination sites, <i>att</i> R1 and <i>att</i> R2, downstream of the CMV 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 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 <i>E. coli</i>
	• SV40 early promoter and origin for expression of the Blasticidin resistance gene and stable propagation of the plasmid in mammalian hosts expressing the SV40 large T antigen
	Blasticidin resistance gene for selection of stable cell lines
	• The pUC origin for high copy replication and maintenance of the plasmid in <i>E. coli</i>
	• The ampicillin resistance gene for selection in <i>E. coli</i>
	For a map of pcDNA ^{TM} 6.2/cLumio ^{TM} -DEST and pcDNA ^{TM} 6.2/nLumio ^{TM} -DEST, refer to pages 20 and 22, respectively.
	continued on next page

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 in mammalian cells using Gateway [®] Technology, simply:
	 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> pcDNA [™] 6.2/ cLumio [™] -DEST or pcDNA [™] 6.2/nLumio [™] -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 Gateway [®] , 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).
Advantages of the Lumio [™] Technology	The Lumio [™] System is based on the FlAsH (<u>Fl</u> uorescein <u>Ars</u> enical <u>H</u> airpin) technology which uses biarsenical labeling reagents to bind and detect proteins containing a tetracysteine motif (<i>i.e.</i> Lumio [™] tag) (Griffin <i>et al.</i> , 1998). Using the Lumio [™] Technology and the Lumio [™] In-Cell Labeling Kits for fluorescence labeling of recombinant proteins provides the following advantages:
	• Small size of the Lumio [™] tag (6 amino acids, 585 Da) is less likely to interfere with the structure or biological activity of the protein of interest
	• Lumio [™] Labeling Reagents are membrane-permeable and readily cross the cell membrane, allowing labeling and detection of recombinant proteins in live mammalian cells
	• Lumio [™] Labeling Reagents bind the Lumio [™] tag with high specificity and high affinity (nanomolar or lower dissociation constant), allowing targeted labeling of the protein of interest
	• Lumio [™] Labeling Reagents become strongly fluorescent only upon binding the Lumio [™] tag, allowing specific detection of Lumio [™] -tagged proteins

Overview, continued

Components of	The Lumio [™] System consists of two major components:
the Lumio [™] System	• The tetracysteine Lumio [™] tag (Cys-Cys-Pro-Gly-Cys-Cys) in the pcDNA [™] 6.2/ Lumio [™] -DEST vector. When fused to a gene of interest, the Lumio [™] tag allows the expressed fusion protein to be specifically recognized by a biarsenical labeling reagent. For more information on the tetracysteine motif, see below.
	• A biarsenical labeling reagent, Lumio [™] Green or Lumio [™] Red, which becomes fluorescent upon binding to recombinant proteins containing the Lumio [™] tag. The Lumio [™] Green and Lumio [™] Red Labeling Reagents are supplied precomplexed to the dithiol EDT (1,2-ethanedithiol) which stabilizes and solubilizes the biarsenic reagents.
Tetracysteine Motif	The Lumio [™] Reagents bind a tetracysteine motif consisting of Cys-Cys-Xaa-Xaa-Cys-Cys where Cys equals cysteine and Xaa equals any amino acid other than cysteine. This motif is rarely seen in naturally occurring proteins allowing specific fluorescence labeling of recombinant proteins fused to the Lumio [™] tag. In the Lumio [™] System, the optimized Cys-Cys-Pro-Gly-Cys-Cys tetracysteine motif is used as this motif has been shown to have a higher affinity for and more rapid binding to biarsenic compounds as well as enhanced stability compared to other characterized motifs (Adams <i>et al.</i> , 2002).
Lumio [™] Green Detection Kit	For sensitive and specific in-gel detection of Lumio [™] -tagged fusion proteins, we recommend the Lumio [™] Green Detection Kit available from Invitrogen (Catalog no. LC6090). The Lumio [™] Green Detection Kit enables immediate visualization of Lumio [™] -tagged proteins in polyacrylamide gels using a UV transilluminator or a visible light laser-based scanner and without the need for staining or western blotting. In addition, the BenchMark [™] Fluorescent Protein Standard (Catalog no. LC5928) allows you to easily visualize molecular weight ranges of proteins labeled with Lumio [™] Green Detection Reagent. For more information on these products and other products that may be used with the Lumio [™] Technology, visit our Web site (www.invitrogen.com) or contact Technical Service (page 25).

Methods

Generating an Entry Clone

Introduction	To recombine your gene of interest into pcDNA [™] 6.2/cLumio [™] -DEST or pcDNA [™] 6.2/nLumio [™] -DEST, you will need an entry clone containing the gene of interest. Many entry vectors including pENTR/D-TOPO [®] (Catalog no. K2400-20) 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 [™] 6.2/cLumio [™] -DEST vector is 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 or mouse gene of interest, we recommend using an Ultimate [™] Human ORF (hORF) or Ultimate [™] Mouse ORF (mORF) Clone available from Invitrogen. Each Ultimate [™] ORF Clone is a fully sequenced clone provided in a Gateway [®] entry vector that is ready-to-use in an LR recombination reaction with pcDNA [™] 6.2/cLumio [™] -DEST. In addition, each clone contains a TAG stop codon, making it fully compatible for use in the Tag-On-Demand [™] System. For more information about the Ultimate [™] ORF Clones available, refer to our Web site (www.invitrogen.com) or contact Technical Service (page 25).
Kozak Consensus Sequence	If you will be expressing your protein from pcDNA [™] 6.2/cLumio [™] -DEST, your insert in the entry clone 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 are the most critical for function (shown in bold).

Generating an Entry Clone, continued

Points to Consider for pcDNA[™]6.2/ cLumio[™]-DEST

pcDNA[™]6.2/cLumio[™]-DEST allows expression of recombinant proteins with a C-terminal peptide containing the V5 epitope and the Lumio[™] tag; however, you may use this vector to express a native protein, if desired. You may also use this vector 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 and Lumio [™] tag	 should contain a Kozak initiation sequence (see previous page)
	• should not contain a stop codon
	 should be in frame with the V5 epitope and Lumio[™] tag after recombination (see page 7 for a diagram)
include the V5 epitope and Lumio™ tag for use in the Tag-	 should contain a Kozak initiation sequence (see previous page)
On-Demand [™] System	• should contain a TAG stop codon
	 should be in frame with the V5 epitope and Lumio[™] tag after recombination (see page 7 for a diagram)
not include the V5 epitope and Lumio [™] tag	 should contain a Kozak initiation sequence (see previous page)
	should contain a stop codon

Points to Consider for pcDNA[™]6.2/ nLumio[™]-DEST

pcDNA[™]6.2/nLumio[™]-DEST allows expression of recombinant proteins with an N-terminal peptide containing the Lumio[™] and V5 epitope tags and contains an ATG initiation codon within the context of a Kozak consensus sequence (see previous page). To include the Lumio[™] and V5 epitope tags, your insert in the entry clone should:

- not contain a Kozak initiation sequence
- be in frame with the Lumio[™] and V5 epitope tags after recombination (see page 8 for a diagram)
- 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 your pcDNA [™] 6.2/Lumio [™] -DEST vector to create your expression clone. To ensure that you obtain the best possible results, we recommend that you read this section and the next section entitled Performing the LR Recombination Reaction (pages 9-11) 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 [™] 6.2/Lumio [™] -DEST vector.
	2. Transform the reaction mixture into a suitable <i>E. coli</i> host.
	3. Select for expression clones (refer to pages 7-8 for diagrams of the recombination region of the resulting expression clones).
Resuspending the Vectors	The pcDNA ^{TM} 6.2/Lumio ^{TM} -DEST vectors are supplied as 6 µg of plasmid, lyophilized in TE, pH 8.0. To use, simply resuspend the destination vector in 40 µl of sterile water to a final concentration of 150 ng/µl.
Propagating the Vectors	If you wish to propagate and maintain the pcDNA TM 6.2/Lumio TM -DEST vectors, we recommend using One Shot [®] <i>ccd</i> B Survival T1 ^R Chemically Competent <i>E. coli</i> (Catalog no. C7510-03) from Invitrogen for transformation. The <i>ccd</i> B Survival T1 ^R <i>E. coli</i> strain is resistant to CcdB effects and can support the propagation of plasmids containing the <i>ccd</i> B 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 DH5α for propagation and maintenance of the pcDNA [™] 6.2/Lumio [™] -DEST vectors as these strains are sensitive to CcdB effects.

Creating an Expression Clone, continued

Recor Regio pcDN cLum	n for A [™] 6.2	2/		 The recombination region of the expression clone resulting from pcDNA[™]6.2/ cLumio[™]-DEST × entry clone is shown below. Note: If you are using pcDNA[™]6.2/cLumio[™]-DEST in the Tag-On-Demand[™] System, your gene of interest must contain a TAG stop codon (see page 4 for more information). Features of the Recombination Region: Shaded regions correspond to DNA sequences transferred from the entry clone into pcDNA[™]6.2/cLumio[™]-DEST by recombination. Non-shaded regions are derived from the pcDNA[™]6.2/cLumio[™]-DEST vector. Bases 922 and 2605 of the pcDNA[™]6.2/cLumio[™]-DEST vector sequence are marked. 						ır							
771	CAA			CTT A CI	GCGTO	ርጥ ልር	CCT	CCA	2 CTC	TATA	7			¥		e ptional sta	>
,,1	Chhi	11000	300 (JIAG	30310		.9910	JADDU	3 010						1990	, I AAC I	
831	AGA	GAAC	CCA (CTGC	FTACI	rg go	CTTAI	CGA	A ATT		T7 pron ACGA		-		GAGA	CCCAA	7
									922	ć	attB1						
891	GCT	GGCTZ	AGT 1	TAAG(CTGA				г т <u>т</u> с А ААС		AAA				GEN	E NA	
			260	5 at	ťB2	μŗ	49110	51101	A AAG	JAIGI		TICC	JICCC	- 1171			G
2597	CCA	GCT	TTC	TTG	TAC	AAA	GTG	GTT	GAT	GCT	GTT	AAC	GGG	AAG	CCT	ATC	
		CGA	_														
									СТА		CAA	TTG	CCC	TTC	GGA	-	
	Pro	Ala			Tyr						CAA	TTG	CCC	TTC		-	
	Pro	Ala			Tyr						CAA	TTG	CCC	TTC		-	
2645	CCT	AAC	Phe CCT	Leu V5 epi CTC	Tyr cope CTC	Lys GGT	Val CTC	Val GAT	Asp TCT	Ala ACG	CAA Val	TTG Asn ACC	CCC Gly GGT	TTC Lys GCT	Pro GGT	Ile GGC	
2645	CCT GGA	AAC TTG	Phe CCT GGA	Leu V5 epit CTC GAG	Tyr cope	Lys GGT CCA	Val CTC GAG	Val GAT CTA	Asp TCT AGA	Ala ACG TGC	CAA Val CGT GCA	TTG Asn ACC TGG	CCC Gly GGT CCA	TTC Lys GCT CGA	Pro GGT CCA	Ile GGC CCG	
2645	CCT GGA	AAC TTG	Phe CCT GGA Pro	Leu V5 epit CTC GAG Leu	Tyr cope CTC GAG	Lys GGT CCA	Val CTC GAG	Val GAT CTA	Asp TCT AGA	Ala ACG TGC	CAA Val CGT GCA	TTG Asn ACC TGG	CCC Gly GGT CCA	TTC Lys GCT CGA	Pro GGT CCA	Ile GGC CCG	
	CCT GGA Pro	AAC TTG Asn	Phe CCT GGA Pro Lumio	Leu V5 epit CTC GAG Leu ™ tag	Tyr cope CTC GAG Leu	Lys GGT CCA Gly	Val CTC GAG Leu	Val GAT CTA Asp	Asp TCT AGA Ser	Ala ACG TGC Thr	CAA Val CGT GCA Arg	TTG Asn ACC TGG Thr	CCC Gly GGT CCA Gly	TTC Lys GCT CGA Ala	Pro GGT CCA Gly	Ile GGC CCG	
2645 2693	CCT GGA Pro TGT	AAC TTG Asn TGT	Phe CCT GGA Pro Lumio	Leu V5 epit GAG Leu ™ tag GGC	Tyr cope CTC GAG	Lys GGT CCA Gly TGC	Val CTC GAG Leu GGT	Val GAT CTA Asp GGC	Asp TCT AGA Ser GGC	Ala ACG TGC	CAA Val CGT GCA	TTG Asn ACC TGG	CCC Gly GGT CCA Gly	TTC Lys GCT CGA	Pro GGT CCA Gly	Ile GGC CCG	

TK polyA Reverse priming site

2739 GGGAGGCTAA CTGAAACACG GAAGGAGACA ATACCGGAAG GAACCCGCGC TATGACGGCA

Creating an Expression Clone, continued

Regi pcDN	ombination on for NA [™] 6.2/ nio [™] -DEST	nLumio [™] - Features o • Shade clone i region	DEST × entry cl of the Recombin d regions correst into pcDNA [™] 6.2 as are derived fr 1018 and 2680 c	one is shown b nation Region: spond to DNA 2/nLumio [™] -DH om the pcDNA		l from the entry . Non-shaded ? vector.
	CAAT			TATA	A 3'end of CMV promoter	Putative transcriptional start
771	CAAATGGGCG	GTAGGCGT	GT ACGGTGG	GAG GTCTAT.	ATAA GCAGAGCTCI	CTGGCTAACT
					T7 promoter/priming site	
831	AGAGAACCCA	CTGCTTAC	TG GCTTATC	GAA ATTAAT.	ACGA CTCACTATAG	GGAGACCCAA
891	GCTGGCTAGT	TAAGCTGC	TAC CG.	A CCA CCG .	Lumio [™] tag TGT TGT CCT GGC ACA ACA GGA CCG Cys Cys Pro Gly	G ACA ACG CCA
				V5	o epitope	
945	CCG CCG TTO	C GAC CCC	TTC GGA T	AG GGA TTG	CCT CTC CTC GG GGA GAG GAG CC Pro Leu Leu Gl	CA GAG CTA AGA
			10	0 18 attB1		
996		A CAA TAG	TGT TCA A	AC ATG TTT	AAA GCA GGC TN TTT CGT CCG AN Lys Ala Gly	
	:	2680 attB2				
2670	ACCCAGCTTT TGGGTCGAAA				TTAG TAATGAGTTI	AAACGGGGGA

TK polyA Reverse priming site

2730 GGCTAACTGA AACACGGAAG GAGACAATAC CGGAAGGAAC CCGCGCTATG ACGGCAATAA

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 your pcDNA [™] 6.2/Lumio [™] -DEST vector, and transform the reaction mixture into a suitable <i>E. coli</i> host (see below) to select for an expression clone. We recommend including the pENTR [™] -gus positive control supplied with the LR Clonase [™] II enzyme mix in your experiments 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 α^{TM} , or equivalent for transformation. 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>ccd</i> A gene and will prevent negative selection with the <i>ccd</i> B gene.
Note	The presence of the EM7 promoter and the Blasticidin resistance gene in the pcDNA ^{TM} 6.2/Lumio ^{TM} -DEST vectors 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 the Appendix , page 18 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 the Appendix , page 19 for instructions on how to prepare and store Blasticidin.
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 provided in this manual as reaction conditions differ.

Performing the LR Recombination Reaction, continued

Materials Needed	Yo	u should have the following materials	on hand before begir	nning:			
	٠	Purified plasmid DNA of your entry clone (50-150 ng/ μ l in TE, pH 8.0)					
	•	 pcDNA[™]6.2/cLumio[™]-DEST or pcDNA[™]6.2/nLumio[™]-DEST vector (150 ng/µl in TE, pH 8.0) 					
	•	LR Clonase [™] II enzyme mix (Invitrog until immediately before use)	gen, Catalog no. 11791	1-020; keep at -20°C			
	•	pENTR [™] -gus positive control, option the LR Clonase [™] II enzyme mix)	nal (50 ng/μl in TE, pl	H 8.0; supplied with			
	٠	TE Buffer, pH 8.0 (10 mM Tris-HCl,	pH 8.0, 1 mM EDTA)				
	•	$2 \ \mu g/\mu l$ Proteinase K solution (supplet thaw and keep on ice until use)	lied with the LR Clona	ase™ II enzyme mix;			
	٠	Appropriate competent <i>E. coli</i> host and growth media for expression					
	•	S.O.C. Medium					
	•	LB agar plates with the appropriate antibiotic to select for expression clones					
Setting Up the LR Recombination	1.	Add the following components to 1.5 ml microcentrifuge tubes at room temperature and mix.					
Reaction		Note: To include a negative control, set u TE Buffer, pH 8.0 for the LR Clonase™ II					
		Component	Sample	Positive Control			
		Entry clone (50-150 ng/reaction)	1-7 µl				
		Destination vector (150 ng/ μ l)	1 µl	1 µl			
		pENTR [™] -gus (50 ng/µl)		2 µl			
		TE Buffer, pH 8.0	to 8 μl	5 μl			
	2	Remove the LR Clonase [™] II enzvme	mix from -20°C and th	haw on ice			

- Remove the LR Clonase[™] II enzyme mix from -20°C and thaw on ice (~ 2 minutes).
- 3. Vortex the LR Clonase[™] II enzyme mix briefly twice (2 seconds each time).
- 4. 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.

- Incubate reactions at 25°C for 1 hour.
 Note: Extending the incubation time to 18 hours typically yields more colonies.
- 6. Add 1 μl of the Proteinase K solution to each reaction. Incubate for 10 minutes at 37°C.
- 7. 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 $^\circ C$ for up to 1 week before transformation.

Performing the LR Recombination Reaction, 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 I reaction should give > 5000 colonies if the entire reaction is transformed and plated.				
Confirming the Expression Clone	• • • • • •				
Sequencing	appropriate N-terminal or C-ter T7 Promoter and TK polyA Rev	uct to confirm that your gene is in frame with the rminal fusion tag, if desired. We suggest using the rerse primer sequences (see below). Refer to the r the location of the primer binding sites.			
	Primer	Sequence			
	T7 Promoter Primer	5′-TAATACGACTCACTATAGGG-3′			
	TK polyA Reverse Primer	5'-CTTCCGTGTTTCAGTTAGC-3'			

Transfecting Cells

Introduction	This section provides general information for transfecting your expression clone into the mammalian cell line of choice. We recommend that you include the pcDNA [™] 6.2/nLumio [™] -GW/p64 positive control vector and a non-Lumio [™] control in your transfection experiment to help you evaluate your results. The non-Lumio [™] control should contain cells transfected with a vector that does not encode the Lumio [™] tag. Using untransfected cells (no transfection reagent, no DNA) as a control is not optimal since transfection may introduce fluorescent artifacts or alter cell morphology that could affect the labeling and detection of your protein.
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 from contamination with 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). Other plasmid purification kits or methods are suitable.
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 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 the other transfection reagents available from Invitrogen, refer to our Web site (www.invitrogen.com) or contact Technical Service (page 25).
Cell Density	If you plan to label your recombinant protein using the Lumio TM In-Cell Labeling Kit, we recommend designing your transfection experiment so that cells will be at optimal density at the time of labeling. Suspension cells typically label most efficiently at a density of 1-2 x 10 ⁶ cells/ml. Adherent cells label most efficiently when they are 60-90% confluent at the time of labeling.

Transfecting Cells, continued

Positive Control pcDNA[™]6.2/nLumio[™]-GW/p64 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. This vector allows expression of the human c-myc (p64) protein with an N-terminal fusion to the Lumio[™] and V5 epitope tags.

To propagate and maintain the plasmid:

- Resuspend the vector in 10 µl sterile water to prepare a 1 µg/µl stock solution. Use the stock solution to transform a *recA*, *endA E*. *coli* strain like TOP10, DH5α, or equivalent.
- 2. Select transformants on LB agar plates containing $50-100 \mu g/ml$ ampicillin.
- 3. Prepare a glycerol stock of a positive transformant for long-term storage.

Creating Stable Cell Lines

Introduction

The pcDNA[™]6.2/Lumio[™]-DEST vectors contain the Blasticidin resistance gene 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 Blasticidin. General information and guidelines are provided below.



The Lumio[™] Technology works best for labeling proteins that are expressed at high levels or are concentrated in a subcellular region. If you are expressing proteins at low levels (*e.g.* from a weak promoter) or if you are expressing a cytoplasmic protein, we recommend conducting initial labeling and detection studies using the Lumio[™] In-Cell Labeling Kits in transiently transfected cells where protein expression levels are higher. Once labeling conditions are optimized, you may analyze your protein in stable cell lines.



To obtain stable transfectants, we recommend that you linearize your pcDNA[™]6.2/Lumio[™]-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.

Determining Blasticidin Sensitivity

To successfully generate a stable cell line expressing your protein of interest, you need to determine the minimum concentration of Blasticidin required to kill your untransfected host cell line by performing a kill curve experiment (see below). Typically, concentrations ranging from 2.5 to $10 \,\mu$ g/ml Blasticidin are sufficient to kill most untransfected mammalian cell lines. Blasticidin is available separately from Invitrogen (Catalog no. R210-01). Refer to the **Appendix**, page 19 for instructions on how to prepare and store Blasticidin.

- 1. Plate cells at approximately 25% confluence. Prepare a set of 6 plates.
- 2. On the following day, replace the growth medium with fresh growth medium containing varying concentrations of Blasticidin (*e.g.* 0, 1, 3, 5, 7.5, and 10μ g/ml Blasticidin).
- 3. Replenish the selective media every 3-4 days, and observe the percentage of surviving cells.
- 4. Count the number of viable cells at regular intervals to determine the appropriate concentration of Blasticidin that prevents growth within 10-14 days after addition of Blasticidin.

Creating Stable Cell Lines, continued

Generating Stable Cell Lines	sele	ce you have determined the appropriate Blasticidin concentration to use for ection, you can generate a stable cell line expressing your pcDNA™6.2/Lumio™- ST construct.
	1.	Transfect the mammalian cell line of interest with the pcDNA [™] 6.2/ cLumio [™] -DEST or pcDNA [™] 6.2/nLumio [™] -DEST expression construct using your transfection method of choice.
	2.	24 hours after transfection, wash the cells and add fresh growth medium.
	3.	48 hours after transfection, split the cells into fresh growth medium such that they are no more than 25% confluent. If the cells are too dense, the antibiotic will not kill the cells. Antibiotics work best on actively dividing cells.
	4.	Incubate the cells at 37°C for 2-3 hours until they have attached to the culture dish.
	5.	Remove the growth medium and replace with fresh growth medium containing Blasticidin at the predetermined concentration required for your cell line.
	6.	Feed the cells with selective media every 3-4 days until Blasticidin-resistant colonies can be identified.
	7.	Pick at least 10 Blasticidin-resistant colonies and expand them to assay for recombinant protein expression.

Detecting Lumio[™] Fusion Proteins

Introduction	Once you have transfected your expression clone into mammalian cells, you may:
	 Detect protein expression and localization in live cells by fluorescence microscopy using the Lumio[™] In-Cell Labeling Kit. For detailed guidelines and protocols, refer to the Lumio[™] In-Cell Labeling Kits manual.
	 Detect protein expression directly in polyacrylamide gels using the Lumio[™] Green Detection Kit (see below).
	• Detect protein expression by Western blot analysis using Anti-V5 Antibodies available from Invitrogen (see below).
Lumio [™] Green Detection Kit	The Lumio [™] Green Detection Kit (Catalog no. LC6090) enables immediate and specific visualization of Lumio [™] -tagged proteins in polyacrylamide gels using a UV transilluminator or a visible light laser-based scanner and without the need for staining or western blotting. Refer to the Lumio [™] Green Detection Kit manual for detailed protocols to prepare lysate samples specifically for detection with the Lumio [™] Green Detection Reagent. This manual is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (page 25).
Western Blotting	You may detect expression of your recombinant fusion protein using the Anti-V5 Antibody (Catalog no. R960-25), Anti-V5-HRP Antibody (Catalog no. R961-25), or Anti-V5-AP Antibody (Catalog no. R962-25) available from Invitrogen. You may use any method of choice to prepare your mammalian cell lysates for Western blot analysis. We recommend the following guidelines:
	 If you plan to analyze your samples using the Lumio[™] Green Detection Kit (see above) in addition to Western blotting, you will need to prepare your samples using lysis buffer. Lysates containing standard Laemmli SDS-PAGE sample buffer will not be suitable for in-gel detection with the Lumio[™] Green Detection Kit. Refer to the Lumio[™] Green Detection Kit manual for a protocol to prepare cell lysates that are compatible with both in-gel detection and Western blot analysis.
	• For cells transfected with the pcDNA [™] 6.2/nLumio [™] -GW/p64 positive control vector, you will need to prepare lysates using RIPA or SDS-PAGE sample buffer to adequately release p64 from the nucleoli. If you are preparing samples using lysis buffer, you may sonicate your samples to release p64.
	• To detect p64 (human c-myc) expression, you may use any of the Anti-V5 Antibodies or the Anti- <i>myc</i> Antibodies available from Invitrogen.
	Note: The <i>c-myc</i> gene encodes a protein with an expected molecular weight of 48 kDa,

Detecting Lumio[™] Fusion Proteins, continued

Polyacrylamide Gels Available from Invitrogen

To facilitate separation and visualization of your recombinant 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. In addition, Invitrogen carries a large selection of molecular weight protein standards and staining kits. For more information, refer to our Web site (www.invitrogen.com) or contact Technical Service (page 25).



Expression of your protein fused to the Lumio[™] and V5 epitope tags will increase the size of your recombinant protein. The table below lists the expected size increase in molecular weight for your recombinant protein. Note that the expected sizes take into account any additional amino acids between the gene of interest and the fusion peptide (see pages 7 and 8 for a diagram).

Vector	Fusion	Expected Size Increase
pcDNA [™] 6.2/cLumio [™] -DEST	C-terminal	4 kDa
pcDNA [™] 6.2/nLumio [™] -DEST	N-terminal	4 kDa

Appendix

Recipes

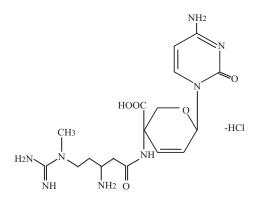
LB (Luria-Bertani) Medium and Plates	Composition: 1.0% Tryptone 0.5% Yeast Extract 1.0% NaCl pH 7.0					
	1.	For 1 liter, dissolve 10 g tryptone, 5 g yeast extract, and 10 g NaCl in 950 ml deionized water.				
	2.	Adjust the pH of the solution to 7.0 with NaOH and bring the volume up to 1 liter.				
	3.	Autoclave on liquid cycle for 20 minutes at 15 psi. Allow solution to cool to 55° C and add antibiotic (100 µg/ml ampicillin) if needed.				
	4.	Store at room temperature or at +4°C.				
	LB	agar plates				
	1.	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 (100 μ g/ml of ampicillin), and pour into 10 cm plates.				
	4.	Let harden, then invert and store at +4°C.				
Low Salt LB Medium with Blasticidin	10 5 ք	w Salt LB Medium: g Tryptone g NaCl g Yeast Extract				
	1.	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.				
	2.	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.				
	4.	Store plates at +4°C in the dark. Plates containing Blasticidin are stable for up to 2 weeks.				

Blasticidin

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

Molecular Weight, Formula, and Structure

The formula for Blasticidin 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.

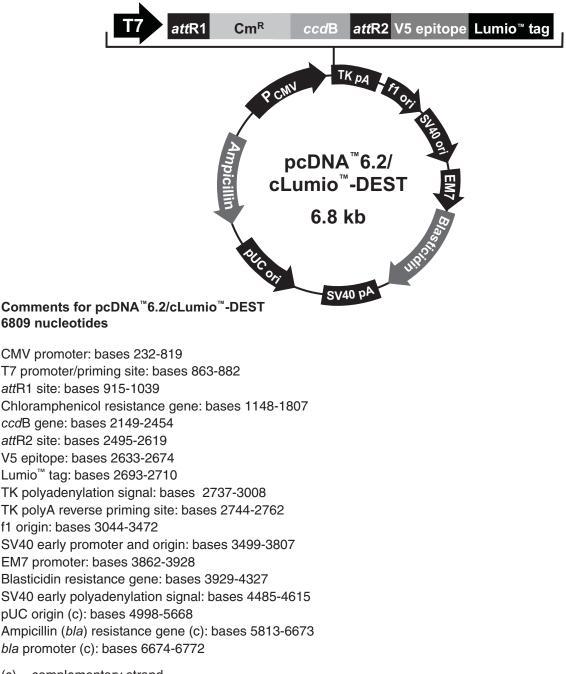


Handling Blasticidin	Always wear gloves, mask, goggles, and protective clothing (<i>e.g.</i> a laboratory coat) when handling Blasticidin. Weigh out Blasticidin and prepare solutions in a hood.
Preparing and Storing Stock Solutions	Blasticidin may be obtained separately from Invitrogen (Catalog no. R210-01) 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 and Features of pcDNA[™]6.2/cLumio[™]-DEST

Мар

The map below shows the elements of pcDNA[™]6.2/cLumio[™]-DEST. DNA from the entry clone replaces the region between bases 922 and 2605. **The complete sequence of this vector is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (page 25).**



(c) = complementary strand

Map and Features of pcDNA[™]6.2/cLumio[™]-DEST, continued

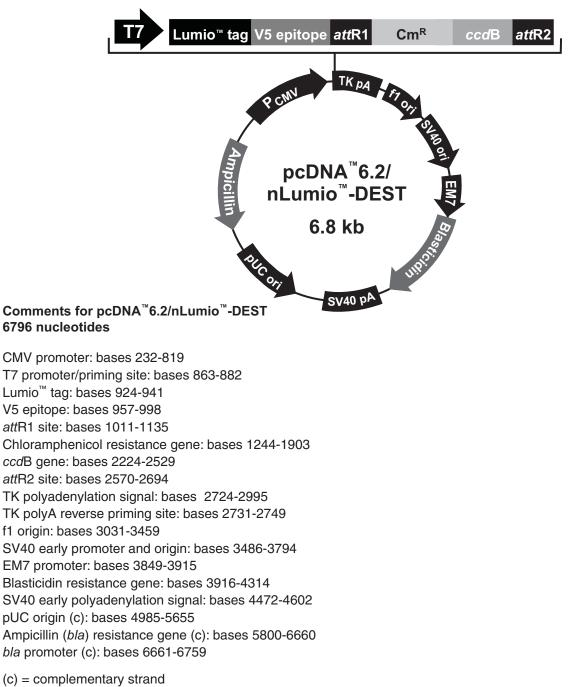
Features

pcDNA[™]6.2/cLumio[™]-DEST (6809 bp) contains 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
Chloramphenicol resistance gene	Allows counterselection of plasmid
ccdB gene	Allows negative selection of plasmid
V5 epitope	Allows detection of the recombinant fusion protein by the Anti-V5 antibodies (Southern <i>et al.</i> , 1991)
Lumio [™] tag	Allows binding of the Lumio [™] Labeling Reagents to facilitate <i>in vivo</i> fluorescence detection of the recombinant fusion protein (Adams <i>et al.,</i> 2002)
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 Blasticidin resistance gene and episomal replication in cells expressing the SV40 large T antigen
EM7 promoter	Allows expression of the Blasticidin resistance gene in <i>E. coli</i>
Blasticidin (bsd) resistance gene	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 resistance gene	Allows selection of transformants in E. coli

Мар

The map below shows the elements of pcDNA[™]6.2/nLumio[™]-DEST. DNA from the entry clone replaces the region between bases 1018 and 2680. **The complete sequence of this vector is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (page 25).**



Map and Features of pcDNA[™]6.2/nLumio[™]-DEST, continued

Features

pcDNA[™]6.2/nLumio[™]-DEST (6796 bp) contain the following elements. All features have been functionally tested.

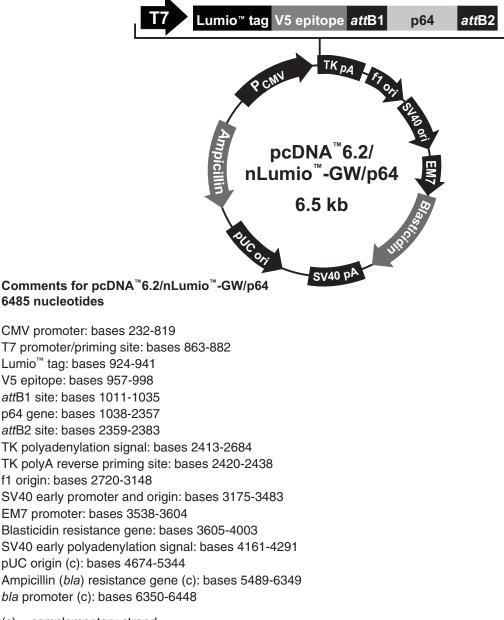
Feature	Benefit
Human cytomegalovirus (CMV) immediate-early promoter/enhancer	Allows efficient, high-level expression of your recombinant protein (Andersson <i>et al.,</i> 1989; Boshart <i>et al.,</i> 1985; Nelson <i>et al.,</i> 1987)
T7 promoter	Allows <i>in vitro</i> transcription in the sense orientation
Lumio [™] tag	Allows binding of the Lumio [™] Labeling Reagents to facilitate <i>in vivo</i> fluorescence detection of the recombinant fusion protein (Adams <i>et al.</i> , 2002)
V5 epitope	Allows detection of the recombinant fusion protein by the Anti-V5 antibodies (Southern <i>et al.,</i> 1991)
attR1 and attR2 sites	Allows recombinational cloning of the gene of interest from an entry clone
Chloramphenicol resistance gene	Allows counterselection of plasmid
<i>ccd</i> B gene	Allows negative selection of plasmid
Herpes Simplex Virus Thymidine Kinase (TK) polyadenylation signal	Allows efficient transcription termination and polyadenylation of mRNA (Cole and Stacy, 1985)
TK polyA reverse priming site	Allows sequencing through the insert
f1 origin	Allows rescue of single-stranded DNA
SV40 early promoter and origin	Allows efficient, high-level expression of the Blasticidin resistance gene and episomal replication in cells expressing the SV40 large T antigen
EM7 promoter	Allows expression of the Blasticidin resistance gene in <i>E. coli</i>
Blasticidin (bsd) resistance gene	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 resistance gene	Allows selection of transformants in E. coli

Map of pcDNA[™]6.2/nLumio[™]-GW/p64

Description

pcDNA[™]6.2/nLumio[™]-GW/p64 (6485 bp) is a control vector expressing the p64 gene, and was generated using the Gateway[®] LR recombination reaction between an entry clone containing the human *c-myc* gene (Bernard *et al.*, 1983; Colby *et al.*, 1983; Watt *et al.*, 1983) and pcDNA[™]6.2/nLumio[™]-DEST. **The complete sequence of this vector is available for downloading from our Web site** (www.invitrogen.com) or by contacting Technical Service (page 25).

Note: The *c-myc* gene encodes a protein with an expected molecular weight of 48 kDa, however, the native protein actually runs at a range of 55-64 kDa on an SDS-PAGE gel.



(c) = complementary strand

Technical Service

World Wide Web



Visit the Invitrogen Web Resource using your World Wide Web browser. At the site, you can:

- Get the scoop on our hot new products and special product offers
- View and download vector maps and sequences
- Download manuals in Adobe® Acrobat® (PDF) format
- Explore our catalog with full color graphics
- Obtain citations for Invitrogen products
- Request catalog and product literature

Once connected to the Internet, launch your Web browser (Internet Explorer 5.0 or newer or Netscape 4.0 or newer), then enter the following location (or URL):

http://www.invitrogen.com

...and the program will connect directly. Click on underlined text or outlined graphics to explore. Don't forget to put a bookmark at our site for easy reference!

Contact Us For more information or technical assistance, call, write, fax, or email. Additional international offices are listed on our Web page (www.invitrogen.com).

Corporate Headquarters:	Japanese Headquarters:	European Headquarters:
Invitrogen Corporation 1600 Faraday Avenue Carlsbad, CA 92008 USA Tel: 1 760 603 7200 Tel (Toll Free): 1 800 955 6288	Invitrogen Japan K.K. Nihonbashi Hama-Cho Park Bldg. 4F 2-35-4, Hama-Cho, Nihonbashi Tel: 81 3 3663 7972	Invitrogen Ltd Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF, UK Tel: +44 (0) 141 814 6100
Fax: 1 760 602 6500 E-mail: tech_service@invitrogen.com	Fax: 81 3 3663 8242 E-mail: jpinfo@invitrogen.com	Tech Fax: +44 (0) 141 814 6117 E-mail: eurotech@invitrogen.com

MSDS Requests

To request an MSDS, visit our Web site at www.invitrogen.com. On the home page, go to 'Technical Resources', select 'MSDS', and follow instructions on the page.

Technical Service, continued

Invitrogen is committed to providing our customers with high-quality goods and Limited Warranty 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, please contact our Technical Service Representatives. Invitrogen warrants that all of its products will perform according to the 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 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

Limited Use Label

Gateway[®] Cloning

License No. 19:

Products

Use of the Mammalian Lumio[™] Gateway[®] Vector Kits with Lumio[™] Technology are covered under the licenses detailed below.

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 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. 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 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. 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 use this product for purposes other than those permitted above, contact Licensing Department, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008. Phone (760) 603-7200.

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.

Purchaser Notification, continued

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

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

Limited Use Label License No. 167: Target Sequences for Synthetic Molecules

6,008,378, 6,054,271, and 6,451,569 and foreign equivalents owned by and/or licensed to Life Technologies Corporation.

This product and/or its use is the subject of one or more of U.S. Patent Nos. 5,932,474,

Limited Use Label License No. 182: Cloning Vectors

The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of patents owned by Invitrogen and claiming this product based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, 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

Adams, S. R., Campbell, R. E., Gross, L. A., Martin, B. R., Walkup, G. K., Yao, Y., Llopis, J., and Tsien, R. Y. (2002). New Biarsenical Ligands and Tetracysteine Motifs for Protein Labeling in Vitro and in Vivo: Synthesis and Biological Applications. J. Am. Chem. Soc. *124*, 6063-6076.

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).

Bernard, O., Cory, S., Gerondakis, S., Webb, E., and Adams, J. M. (1983). Sequence of the Murine and Human Cellular myc Oncogenes and Two Modes of myc Transcription Resulting from Chromosome Translocation in B Lymphoid Tumors. EMBO J. 2, 2375-2383.

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.

Colby, W. W., Chen, E. Y., Smith, D. H., and Levinson, A. D. (1983). Identification and Nucleotide Sequence of a Human Locus Homologous to the v-myc Oncogene of Avian Myelocytomatosis Virus MC29. Nature *301*, 722-725.

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.

Griffin, B. A., Adams, S. R., and Tsien, R. Y. (1998). Specific Covalent Labeling of Recombinant Protein Molecules Inside Live Cells. Science 281, 269-272.

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.

References, continued

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.

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 Site-specific 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.

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.

Watt, R., Stanton, L. W., Marcu, K. B., Gallo, R. C., Croce, C. M., and Rovera, G. (1983). Nucleotide Sequence of Cloned cDNA of Human c-myc Oncogene. Nature *303*, 725-728.

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 or human therapeutic or diagnostic use.

^{©2003-2004, 2010} Invitrogen Corporation. All rights reserved.



Corporate Headquarters:

Invitrogen Corporation 1600 Faraday Avenue Carlsbad, California 92008 Tel: 1 760 603 7200 Tel (Toll Free): 1 800 955 6288 Fax: 1 760 602 6500 Email: tech_service@invitrogen.com

European Headquarters:

Invitrogen Ltd Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF, UK Tel (Free Phone Orders): 0800 269 210 Tel (General Enquiries): +44 (0) 141 814 6100 Fax: +44 (0) 141 814 6260 Email: eurotech@invitrogen.com

International Offices:

Argentina 5411 4556 0844 Australia 1 800 331 627 Austria 0800 20 1087 Belgium 0800 14894 Brazil 55 11 5051 7422 Canada 800 263 6236 China 10 6849 2578 Denmark 80 30 17 40

France 0800 23 20 79 Germany 0800 083 0902 Hong Kong 2407 8450 Italy 02 98 22 201 Japan 03 3663 7974 The Netherlands 0800 099 3310 New Zealand 0800 600 200 Norway 00800 5345 5345

Singapore 65 686 186 38 Spain & Portugal 900 181 461 Sweden 020 26 34 52 Switzerland 0800 848 800 Taiwan 2 2651 6156 UK 0800 269 010 For other countries see our website

www.invitrogen.com