

# BLOCK-iT<sup>™</sup> Lentiviral Pol II miR RNAi Expression System

### A Gateway<sup>®</sup>-adapted, lentiviral destination vector for high-level expression of microRNA in dividing and non-dividing mammalian cells

Cat. nos. K4937-00 and K4938-00

Rev. date: 15 September 2010 Manual part no. 25-0853 MAN0000523

**User Manual** 

### Contents

Kit Contents and Storage	iv
Introduction	1
System Summary	1
Using miRNA for RNAi Analysis	5
The BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Expression System	9
Gateway <sup>®</sup> Recombination Reactions	11
Green Fluorescent Protein	14
Biosafety Features of the System	16
Experimental Outline	
Methods	19
Cloning miRNA	
Creating Entry Clones for Use with pLenti6/V5-DEST	
Performing the Rapid BP/LR Recombination Reaction	
Transforming One Shot <sup>®</sup> Stbl3 <sup>™</sup> Competent <i>E. coli</i>	
Producing Lentivirus in 293FT Cells	
Titering Your Lentiviral Stock	
Transduction and Analysis	44
Detecting Fluorescence	
Examples of Expected Results	
Troubleshooting	
Appendix	57
Performing the Standard BP Recombination Reaction	
Performing the Standard LR Recombination Reaction	
Blasticidin	
Map and Features of pLenti6/V5-DEST	65
Map of pLenti6/V5-GW/lacZ	67
Map and Features of pDONR <sup>™</sup> 221	
Map and Features of pLP1	70
Map and Features of pLP2	72
Map and Features of pLP/VSVG	74
Map of pENTR <sup>™</sup> -gus	76
Additional Products	77
Technical Support	79
Purchaser Notification	
Gateway <sup>®</sup> Clone Distribution Policy	
References	

### Kit Contents and Storage

### **Types of Kits** This manual is supplied with the following products.

	Product		Cat. no.
	BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Expression Sys	stem	K4937-00
	BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Expression Sys EmGFP	stem with	K4938-00
Intended Use	<b>For research use only.</b> Not intended for human or anir therapeutic uses.	nal diagnosti	c or
Kit Components	The BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Kits include components. For a detailed description of the contents of following pages. For a detailed description of the BLOC Expression Vector Kit contents and how to use the reage BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kit man instructions to grow and maintain the 293FT Cell Line, s manual.	of each compo cK-iT <sup>™</sup> Pol II r ents supplied nual. For deta	onent, see the niR RNAi , see the iled
		Cat	. no.
	Components	K4937-00	K4938-00
	BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kit	✓	
	BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kit with EmGFP		~
	pLenti6/V5-DEST Gateway® Vector Kit	$\checkmark$	✓
	Gateway <sup>®</sup> BP Clonase <sup>®</sup> II Enzyme Mix	✓	✓
	Gateway <sup>®</sup> BP Clonase <sup>®</sup> II Enzyme Mix Gateway <sup>®</sup> LR Clonase <sup>®</sup> II Enzyme Mix	✓ ✓	✓ ✓
	Gateway <sup>®</sup> LR Clonase <sup>®</sup> II Enzyme Mix	* * *	* * *
	5	* * *	* * * *
	Gateway <sup>®</sup> LR Clonase <sup>®</sup> II Enzyme Mix One Shot <sup>®</sup> Stbl3 <sup>™</sup> Chemically Competent <i>E. coli</i>		
	Gateway <sup>®</sup> LR Clonase <sup>®</sup> II Enzyme Mix One Shot <sup>®</sup> Stbl3 <sup>™</sup> Chemically Competent <i>E. coli</i> One Shot <sup>®</sup> TOP10 Chemically Competent <i>E. coli</i>		

### Kit Contents and Storage, Continued

manual.

# Shipping and<br/>StorageThe BLOCK-iT<sup>™</sup> Lentiviral Pol II miR RNAi Kits are shipped as described below.<br/>Upon receipt, store each item as detailed below. For more detailed information<br/>about the reagents supplied in the BLOCK-iT<sup>™</sup> Pol II miR RNAi Expression<br/>Vector Kits, refer to the BLOCK-iT<sup>™</sup> Pol II miR RNAi Expression Vector Kit

Box	Component	Shipping	Storage
1	BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kit	Dry ice	-20°C
	or		
	BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kit with EmGFP		
2	One Shot® TOP10 Chemically Competent E. coli	Dry ice	-80°C
3	pLenti6/V5-DEST Gateway® Vector Kit	Blue ice	-20°C
4	pDONR <sup>™</sup> 221 Vector	Room	-20°C
		temperature	
5–7	ViraPower <sup>™</sup> Bsd Lentiviral Support Kit:		
	ViraPower <sup>™</sup> Packaging Mix	Blue ice	-20°C
	Lipofectamine <sup>®</sup> 2000	Blue ice	4°C (do not freeze)
	Blasticidin	Room	-20°C
		temperature	
8	One Shot <sup>®</sup> Stbl3 <sup>™</sup> Chemically Competent <i>E. coli</i>	Dry ice	-80°C
9–10	Gateway <sup>®</sup> BP Clonase <sup>®</sup> II Enzyme Mix	Dry ice	-20°C
	Gateway <sup>®</sup> LR Clonase <sup>®</sup> II Enzyme Mix	Dry ice	-20°C
11	293FT Cell Line	Dry ice	Liquid nitrogen

### BLOCK-iT<sup>™</sup> Pol II miR RNAi Expression Vector Kit

The BLOCK-iT<sup>™</sup> Lentiviral Pol II miR RNAi Expression Kits include an expression vector for expression of your microRNA (miRNA) sequence. The BLOCK-iT<sup>™</sup> Pol II miR RNAi Expression Vector Kits contain:

- BLOCK-iT<sup>™</sup> Pol II miR RNAi Expression Vector **or** BLOCK-iT<sup>™</sup> Pol II miR RNAi Expression Vector with EmGFP Reagents (Box 1)
- One Shot<sup>®</sup> TOP10 Chemically Competent *E. coli* (Box 2)

Refer to the BLOCK-iT<sup>™</sup> Pol II miR RNAi Expression Vector Kit manual for a detailed description of the reagents provided with the kit and instructions to produce the Gateway<sup>®</sup> expression vector construct.

### Kit Contents and Storage, Continued

#### Vectors

The following vectors are included with the BLOCK-iT<sup>™</sup> Lentiviral Pol II miR RNAi Expression Kits (Boxes 3 and 4). **Store the vectors at –20°C.** 

Vector	Composition	Amount
pLenti6/V5-DEST Gateway® Vector	40 μL of vector at 150 ng/μL in 10 mM Tris- HCl, 1 mM EDTA, pH 8.0	6 µg
pLenti6/V5-GW/lacZ control	20 μL of vector at 500 ng/μL in 10 mM Tris- HCl, 1 mM EDTA, pH 8.0	10 µg
pDONR <sup>™</sup> 221 Vector	40 μL of vector at 150 ng/μL in 10 mM Tris- HCl, 1 mM EDTA, pH 8.0	6 µg

### ViraPower<sup>™</sup> Bsd Lentiviral Support Kit Contents

The following reagents are included with the ViraPower<sup>™</sup> Bsd Lentiviral Support Kit (Boxes 5–7). Store the ViraPower<sup>™</sup> Packaging Mix and Blasticidin at –20°C. Store Lipofectamine<sup>®</sup> 2000 Reagent at 4°C.

Important: Do not freeze Lipofectamine<sup>®</sup> 2000 Reagent.

Reagent	Composition	Amount
ViraPower <sup>™</sup> Packaging Mix	Contains a mixture of the pLP1, pLP2, and pLP/VSVG plasmids, 1 µg/µL in TE, pH 8.0	195 µg
Lipofectamine <sup>®</sup> 2000 Reagent	Proprietary	0.75 mL
Blasticidin	Powder	50 mg

### One Shot<sup>®</sup> Stbl3<sup>™</sup> Chemically Competent *E. coli*

The following reagents are included with the One Shot<sup>®</sup> Stbl3<sup>TM</sup> Chemically Competent *E. coli* kit (Box 8). Transformation efficiency is  $\geq 1 \times 10^8$  cfu/µg plasmid DNA. Store Box 8 at -80°C.

Reagent	Composition	Amount
S.O.C. Medium	2% Tryptone	6 mL
	0.5% Yeast Extract	
	10 mM NaCl	
	2.5 mM KCl	
	10 mM MgCl <sub>2</sub>	
	10 mM MgSO <sub>4</sub>	
	20 mM glucose	
Stbl3 <sup>™</sup> Cells		$21 \times 50 \ \mu L$
pUC19 Control DNA	10 pg/ $\mu$ L in 5 mM Tris-HCl, 0.5 mM EDTA, pH 8	50 µL

### Kit Contents and Storage, Continued

Genotype of Stbl3 <sup>™</sup> Cells	F <sup>-</sup> mcrB mrr hsdS20( $r_B^-$ , $m_B^-$ ) recA13 supE44 ara-14 galK2 lacY1 proA2 rpsL20(Str <sup>R</sup> ) xyl-5 $\lambda^-$ leu mtl-1
Gateway <sup>®</sup> LR Clonase <sup>®</sup> II	The following reagents are included with the Gateway <sup>®</sup> Clonase <sup>®</sup> II Enzyme Mixes (Boxes 9–10).
Enzyme Mixes	<b>Store Box 9 and 10 at –20°C for up to 6 months.</b> For long-term storage, store at –80°C.
	Gateway <sup>®</sup> BP Clonase <sup>®</sup> II

Reagent	Composition	Amount
Gateway <sup>®</sup> BP Clonase <sup>®</sup> II Enzyme Mix	Proprietary	40 µL
Proteinase K Solution	2 μg/μL in: 10 mM Tris-HCl, pH 7.5 20 mM CaCl <sub>2</sub> 50% glycerol	40 µL
PEG Solution	30% PEG 8000 30 mM MgCl <sub>2</sub>	1 mL
pEXP7-tet Positive Control	50 ng/μL in TE Buffer, pH 8.0	20 µL

### Gateway® LR Clonase® II

Reagent	Composition	Amount
Gateway <sup>®</sup> LR Clonase <sup>®</sup> II Enzyme Mix	Proprietary	40 µL
Proteinase K Solution	2 μg/μL in: 10 mM Tris-HCl, pH 7.5 20 mM CaCl <sub>2</sub> 50% glycerol	40 µL
pENTR <sup>™</sup> -gus Positive Control	50 ng/μL in TE Buffer, pH 8.0	20 µL

**Note:** The pENTR<sup>™</sup>-gus control included with the Gateway<sup>®</sup> LR Clonase<sup>®</sup> II Enzyme Mix may be used as a positive control for the LR recombination reaction **only** (see page 61).

#### 293FT Cell Line

The 293FT Cell Line (Box 11) is used for the production of lentiviral stocks. The 293FT Cell Line is supplied as one vial containing  $3 \times 10^6$  frozen cells in 1 mL of Freezing Medium. **Upon receipt, store in liquid nitrogen.** 

For instructions to thaw, culture, and maintain the 293FT Cell Line, see the 293FT Cell Line manual.

### Introduction

# System Summary

Description of the System	The BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Expression System combines BLOCK-iT <sup>™</sup> Pol II miR RNAi and ViraPower <sup>™</sup> Lentiviral technologies to facilitate creation of a replication-incompetent lentivirus that delivers a microRNA (miRNA) sequence of interest to dividing or non-dividing mammalian cells for RNA interference (RNAi) analysis.
System Components	<ul> <li>The BLOCK-iT<sup>™</sup> Lentiviral Pol II miR RNAi Expression System includes:</li> <li>A BLOCK-iT<sup>™</sup> Pol II miR RNAi Expression Vector Kit for producing an expression clone containing a double-stranded oligonucleotide (ds oligo) encoding a pre-miRNA sequence (see page 8) for expression in mammalian cells using an RNA Polymerase II (Pol II) promoter, the human cytomegalovirus (CMV) immediate early promoter. The BLOCK-iT<sup>™</sup> Pol II miR RNAi Expression Vector Kits are supplied with a choice of pcDNA<sup>™</sup>6.2-GW/miR or pcDNA<sup>™</sup>6.2-GW/EmGFP-miR expression vectors (collectively referred to as pcDNA<sup>™</sup>6.2-GW/± EmGFP-miR).</li> <li>The pDONR<sup>™</sup>221 vector is used as an intermediate to transfer the pre-miRNA expression cassette (see page 8 for details) into the lentiviral expression plasmid (see below) using Gateway<sup>®</sup> Technology.</li> <li>A pLenti6/V5-DEST destination vector into which the pre-miRNA cassette from the expression clone is transferred using Gateway<sup>®</sup> Technology (see below). This expression plasmid contains elements that allow packaging of the construct into virions and the Blasticidin resistance marker for selection of stably transduced cell lines.</li> <li>Gateway<sup>®</sup> BP and LR Clonase<sup>®</sup> II Enzyme Mixes that facilitate the transfer of the pre-miRNA expression cassette from the expression vector.</li> <li>Components of the ViraPower<sup>™</sup> Lentiviral System for production of a replication-incompetent lentivirus that stably expresses the miRNA of interest in both dividing and non-dividing mammalian cells.</li> </ul>
	ViraPower <sup>™</sup> Lentiviral Technology, and Gateway <sup>®</sup> Technology, see pages 2–3.

# System Summary, Continued

Advantages of the BLOCK-iT <sup>™</sup> Lentiviral Pol II	Using BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Expression Kits to facilitate lentiviral-based delivery of miRNA to mammalian cells provides the following advantages:
miR RNAi Expression System	<ul> <li>The BLOCK-iT<sup>™</sup> Pol II miR RNAi Expression Vectors provide a rapid and efficient way to clone ds oligo duplexes encoding a desired miRNA target sequence into a vector containing a Pol II promoter for use in RNAi analysis.</li> </ul>
	<ul> <li>Gateway<sup>®</sup>-adapted vectors for easy transfer of the miRNA of interest from one expression vector (pcDNA<sup>™</sup>6.2-GW/<u>+</u>EmGFPmiR) into another (pLenti6/V5-DEST vector).</li> </ul>
	• Generates a replication-incompetent lentivirus that effectively transduces both dividing and non-dividing mammalian cells, thus broadening the potential RNAi applications beyond those of other traditional retroviral systems (Naldini, 1998).
	• Efficiently delivers the miRNA of interest to mammalian cells in culture or <i>in vivo</i> .
	• Provides stable, long-term expression of the miRNA of interest beyond that offered by traditional adenoviral-based systems.
	• Produces a pseudotyped virus with a broadened host range (Yee, 1999).
	• Includes multiple features designed to enhance the biosafety of the system.
The BLOCK-iT <sup>™</sup> Pol II miR RNAi Technology	The BLOCK-iT <sup>™</sup> Pol II miR RNAi Technology is a next generation RNAi technology employing miRNA expression vectors that allow flexible expression of miRNA-based knockdown cassettes driven by RNA Polymerase II (Pol II) promoters in mammalian cells.
	The BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vectors are specifically designed to allow expression of miRNA sequences and contain specific miR flanking sequences that allow proper processing of the miRNA. The expression vector design is based on the miRNA vector system developed in the laboratory of David Turner (U.S. Patent Publication No. 2004/0053876) and includes the use of endogenous murine miR-155 flanking sequences (see page 8 for details).
	A variety of BLOCK-iT <sup>™</sup> RNAi products are available to facilitate RNAi analysis in mammalian and invertebrate systems. For more information about any of the BLOCK-iT <sup>™</sup> RNAi products, see the RNAi Central application portal at <u>www.invitrogen.com/rnai</u> or contact Technical Support (see page 79).
	Continued on next page

# System Summary, Continued

ViraPower <sup>™</sup> Lentiviral Technology	ViraPower <sup>™</sup> Lentiviral Technology facilitates highly efficient, <i>in vitro</i> or <i>in vivo</i> delivery of a target gene or RNA to dividing and non-dividing mammalian cells using a replication-incompetent lentivirus. Based on the lentikat <sup>™</sup> system developed by Cell Genesys (Dull <i>et al.</i> , 1998), the ViraPower <sup>™</sup> Lentiviral Technology possesses features which enhance its biosafety while allowing high-level expression in a wider range of cell types than traditional retroviral systems. The main components of the ViraPower <sup>™</sup> Lentiviral Expression System include:
	• A pLenti-based expression vector ( <i>e.g.</i> , pLenti6/V5-DEST) into which the DNA sequence of interest will be cloned. This vector contains elements required to allow packaging of the expression construct into virions and an antibiotic resistance marker to allow selection of stably transduced cell lines. For more information, see page 9.
	<ul> <li>The ViraPower<sup>™</sup> Packaging Mix, an optimized mixture of the three packaging plasmids required for production of the lentivirus.</li> </ul>
	• An optimized 293FT producer cell line to facilitate optimal production of virus.
	For more information about these components, see page 9. For more information about the biosafety features of the system, see page 16.
Gateway <sup>®</sup> Technology	Gateway <sup>®</sup> 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 DNA sequence of interest into multiple vector systems.
	To express your miRNA of interest in mammalian cells using the BLOCK-iT™ Lentiviral Pol II miR RNAi Expression System and Gateway® Technology, simply:
	<ol> <li>Clone a double-stranded oligonucleotide encoding your miRNA sequence of interest into the pcDNA<sup>™</sup>6.2-GW/miR or pcDNA<sup>™</sup>6.2-GW/EmGFP-miR expression vector to create an expression clone. Transfect this expression clone directly into mammalian cells for initial screening, if desired.</li> </ol>
	<ol> <li>To transfer your pre-miRNA expression cassette into pLenti6/V5-DEST vector, generate an entry clone by performing a BP recombination reaction between the pcDNA<sup>™</sup>6.2-GW/miR or pcDNA<sup>™</sup>6.2-GW/EmGFP-miR expression clone and pDONR<sup>™</sup>221 donor vector, then perform a LR recombination reaction between the resulting entry clone (pENTR<sup>™</sup>221/miR) and pLenti6/V5-DEST. See page 25 for more details.</li> </ol>
	<ol><li>Use your lentiviral expression clone and the reagents supplied in the kit to produce a lentiviral construct.</li></ol>
	<ol> <li>Transduce the lentiviral construct into mammalian cells to express the miRNA. Select for stably transduced cells, if desired.</li> </ol>
	For detailed information about the Gateway <sup>®</sup> Technology, refer to the Gateway <sup>®</sup> Technology with Clonase <sup>®</sup> II manual which is available at <u>www.invitrogen.com</u> or by contacting Technical Support (see 79).
	Continued on next page

### System Summary, Continued

Purpose of this Manual	This manual provides an overview of the BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Expression Kits and provides instructions and guidelines to:
	<ol> <li>Use the pLenti6/V5-DEST vector and expression clone (pcDNA<sup>™</sup>6.2- GW/<u>+</u>EmGFP-miR) in a Rapid BP/LR recombination reaction to generate a lentiviral expression clone containing the pre-miRNA sequence of interest.</li> </ol>
	<ol> <li>Co-transfect the pLenti6/V5-GW/<u>+</u>EmGFP-miR expression construct and the ViraPower<sup>™</sup> Packaging Mix into the 293FT Cell Line to produce a lentiviral stock.</li> </ol>
	3. Titer the lentiviral stock.
	<ol> <li>Transduce the lentiviral construct into mammalian cells and perform "transient" RNAi analysis</li> </ol>
	5. Generate a stably transduced cell line, if desired.
	For detailed instructions to generate a pcDNA6.2 <sup>™</sup> -GW/ <u>+</u> EmGFP-miR expression clone containing the pre-miRNA expression cassette, refer to the BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kit manual. For instructions to culture and maintain the 293FT producer cell line, refer to the 293FT Cell Line manual. These manuals are supplied with the BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Expression Kits, but are also available at <u>www.invitrogen.com</u> or by contacting Technical Support (see page 79).
Note	The One Shot <sup>®</sup> Stbl3 <sup>™</sup> Chemically Competent <i>E. coli</i> , Gateway <sup>®</sup> LR Clonase <sup>®</sup> II Enzyme Mix, and Lipofectamine <sup>®</sup> 2000 Reagent included in the BLOCK-iT <sup>™</sup> Lentiviral RNAi Expression System are available separately and are supplied with individual documentation detailing general use of the product. For instructions to use these products specifically with the BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Kits, follow the recommended protocols in this manual.
Important	The BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Expression Kit is designed to help you create a lentivirus to deliver and express an miRNA sequence in mammalian cells for RNAi analysis. Although the system is designed to help you express your miRNA sequence in the simplest, most direct fashion, use of the system is intended for users familiar with the principles of retrovirus biology and gene silencing. We highly recommend that users possess a working knowledge of viral and tissue culture techniques, lipid-mediated transfection, Gateway <sup>®</sup> Technology, and the RNAi pathway. For more information about the following topics, refer to these references:
	<ul> <li>Retrovirus biology and the retroviral replication cycle: see Buchschacher and Wong-Staal, 2000 and Luciw, 1996.</li> </ul>
	Retroviral and lentiviral vectors: see Naldini 1999 Naldini 1998 and Yee 1999

- Retroviral and lentiviral vectors: see Naldini, 1999, Naldini, 1998, and Yee, 1999.
- RNAi pathway and expression of miRNA in mammalian cells: see published references (Brummelkamp *et al.*, 2002; Cullen, 2004; Kim, 2005; McManus & Sharp, 2002; Sui *et al.*, 2002; Yu *et al.*, 2002; Zeng *et al.*, 2002)

# Using miRNA for RNAi Analysis

Introduction	RNA interference (RNAi) describes the phenomenon by which short, homologous RNA duplexes induce potent and specific inhibition of eukaryotic gene expression via the degradation of complementary messenger RNA (mRNA), and is functionally similar to the processes of post-transcriptional gene silencing (PTGS) or cosuppression in plants (Cogoni <i>et al.</i> , 1994; Napoli <i>et al.</i> , 1990; Smith <i>et al.</i> , 1990; van der Krol <i>et al.</i> , 1990) and quelling in fungi (Cogoni & Macino, 1997; Cogoni & Macino, 1999; Romano & Macino, 1992). In plants, the PTGS response is thought to occur as a natural defense against viral infection or transposon insertion (Anandalakshmi <i>et al.</i> , 1998; Jones <i>et al.</i> , 1998; Li & Ding, 2001; Voinnet <i>et al.</i> , 1999). In experimental settings, RNAi is widely used to silence genes through transfection of RNA duplexes or introduction of vector- expressed short hairpin RNA (shRNA).	
The RNAi Pathway	In eukaryotic organisms, dsRNA produced <i>in vivo</i> , introduced by pathogens, or through research, is processed into 21–23 nucleotide double-stranded short interfering RNA duplexes (siRNA) by an enzyme called Dicer, a member of the RNase III family of double-stranded RNA-specific endonucleases (Bernstein <i>et al.</i> , 2001; Ketting <i>et al.</i> , 2001). Each siRNA then incorporates into an RNA-induced silencing complex (RISC), an enzyme complex that serves to target cellular transcripts complementary to the siRNA for specific cleavage and degradation, or translational repression (Hammond <i>et al.</i> , 2000; Nykanen <i>et al.</i> , 2001). MicroRNAs (miRNAs) are endogenous RNAs that trigger gene silencing (Ambros, 2001; Carrington & Ambros, 2003).	
miRNA Pathway	MicroRNAs (miRNAs) are endogenously expressed small ssRNA sequences of ~22 nucleotides in length which naturally direct gene silencing through components shared with the RNAi pathway (Bartel, 2004). Unlike shRNAs, however, the miRNAs are found embedded, sometimes in clusters, in long primary transcripts (pri-miRNAs) of several kilobases in length containing a hairpin structure and driven by RNA Polymerase II (Lee <i>et al.</i> , 2004), the polymerase also responsible for mRNA expression. Drosha, a nuclear RNase III, cleaves the stem-loop structure of the pri-miRNA to generate small hairpin precursor miRNAs (pre-miRNAs) which are ~70 nucleotides in length (Zeng <i>et al.</i> , 2005). The pre-miRNAs are exported from the nucleus to the cytoplasm by exportin-5, a nuclear transport receptor (Lund <i>et al.</i> , 2004; Yi <i>et al.</i> , 2003). Following the nuclear export, the pre-miRNAs are processed by Dicer into a ~22 nucleotides miRNA (mature miRNA) molecule, and incorporated into an miRNA-containing RNA-induced silencing complex (miRISC) (Cullen, 2004).	

# Using miRNA for RNAi Analysis, Continued

Translational Repression versus Target Cleavage	The mature miRNAs regulate gene expression by mRNA cleavage (mRNA is nearly complementary to the miRNA) or translational repression (mRNA is not sufficiently complementary to the miRNA). Target cleavage can be induced artificially by altering the target or the miRNA sequence to obtain complete hybridization (Zeng <i>et al.</i> , 2002). In animals, most miRNAs imperfectly complement their targets and interfere w protein production without directly inducing mRNA degradation (Ambros, 200 Nonetheless, these miRNAs are found associated with the RNAi nuclease AGC (Liu <i>et al.</i> , 2004; Meister <i>et al.</i> , 2004), and at least two miRNAs with close matched to their target sequences, particularly in their 5′ regions, have been shown to cleave cognate mRNAs produced by the BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kits (see below) fully complement their target site and cleave the target mRNA. Sequence analysis showed that the primary cleavage site at the phosphodiester bond in the mRNA found opposite the tenth and eleventh base of the engineered miRNA as predicted for RNAi-mediated cleavage (Elbashir <i>et al.</i> , 2001) similar to siRNA mediated cleavage.		
Using a Vector- Based System to Express Engineered miRNA	A major limitation of using siRNA (diced siRNA or synthetic siRNA) for RNAi analysis in mammalian cells is the transient nature of siRNA. To address this limitation, a number of groups have developed vector-based systems to facilitate expression of engineered short hairpin RNA (shRNA) sequences in mammalian cells using Pol III promoters (Brummelkamp <i>et al.</i> , 2002; Paddison <i>et al.</i> , 2002; Paul <i>et al.</i> , 2002; Sui <i>et al.</i> , 2002; Yu <i>et al.</i> , 2002). However, the use of shRNA vectors for RNAi analysis requires the screening of large number of sequences to identify active sequences and the use of Pol III promoters limits applications such as tissue-specific expression.		
	To overcome the limitations with siRNA and shRNA, we have developed Gateway <sup>®</sup> -adapted expression vectors that enable the expression of engineered miRNA sequences from Pol II promoters. The pcDNA6.2 <sup>™</sup> -GW/ <u>+</u> EmGFP-miR expression vectors (supplied in the BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kits) facilitate the generation of an expression clone containing a ds oligo encoding a pre-miRNA sequence (see page 8). The resulting expression construct may be introduced into dividing mammalian cells for transient expression of the miRNA sequence and initial RNAi screening, if desired. Once initial screening is complete, the pre-miRNA sequence may then be easily and efficiently transferred into the pLenti6/V5-DEST vector (or other suitable destination vector) by Gateway <sup>®</sup> recombination reactions (see page 11).		
	For more information about the BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kits, included components, and generating the expression construct, refer to the BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kit manual.		

# Using miRNA for RNAi Analysis, Continued

Types of miRNA Vectors	The BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kits are supplied with one of the following expression vectors that allow the expression of your engineered pre- miRNA:			
	• pcDNA <sup>™</sup> 6.2-GW/miR			
	Allows expression of the engineered pre-miRNA under the control of the strong, Pol II human CMV (cytomegalovirus) promoter and Herpes Simplex virus (HSV) thymidine kinase (TK) polyadenylation signal			
	• pcDNA <sup>™</sup> 6.2-GW/EmGFP-miR			
	This vector is similar to pcDNA6.2 <sup>™</sup> -GW/miR, except the coding sequence of EmGFP (Emerald Green Fluorescent Protein) is incorporated into the vector such that the pre-miRNA insertion site is in the 3' untranslated (3'UTR) region of the fluorescent protein mRNA. Addition of EmGFP allows tracking of the miRNA expression and provides strong correlation of EmGFP expression with the knockdown of the target gene by your miRNA.			
Advantages of Using Pol II	miRNA vector-based systems that use Pol II promoters for RNAi cassette expression offer the following advantages over traditional siRNA or shRNA expression:			
miRNA Vector- Based Systems	<ul> <li>Enables co-cistronic expression of reporter genes such as GFP (see above) allowing reliable tracking of miRNA expression in mammalian cells</li> </ul>			
	<ul> <li>Allows expression of miRNA from a variety of promoters, including tissue- specific and regulated promoters for <i>in vivo</i> experiments</li> </ul>			
	• Enables expression of multiple miRNAs from a single transcript allowing the knockdown of more than one gene simultaneously (see the BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kit manual for details)			
	• Permits design of predictable RNAi constructs with a high rate of success			
Human CMV Promoter	The BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vectors contain the human cytomegalovirus (CMV) immediate early promoter to allow high-level, constitutive miRNA expression in mammalian cells (Andersson <i>et al.</i> , 1989; Boshart <i>et al.</i> , 1985; Nelson <i>et al.</i> , 1987).			
	We have chosen the human CMV promoter to control vector-based expression of miRNA molecules in mammalian cells for the following reasons:			
	• The promoter is recognized by RNA Polymerase II and controls high-level, constitutive expression of miRNA and co-cistronic reporter genes			
	• The promoter is active in most mammalian cell types			
	<b>Note:</b> Although highly active in most mammalian cell lines, activity of the viral CMV promoter can be down-regulated in some cell lines due to methylation (Curradi <i>et al.</i> , 2002), histone deacetylation (Rietveld <i>et al.</i> , 2002), or both.			

# Using miRNA for RNAi Analysis, Continued

Structure of the Engineered pre- miRNA	The BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vectors are designed to accept engineered pre-miRNA sequences targeting your gene of interest. The engineered pre-miRNA sequence structure is based on the murine miR-155 sequence and the stem-loop structure was optimized to obtain a high knockdown rate. For details on miR-155 and stem-loop optimization, refer to the BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector manual.			
	For optimized knockdown results, we recommend that the ds oligo encoding the engineered pre-miRNA have the following structural features:			
	• Two 4 nucleotides, 5' overhangs complementary to the vector (required for directional cloning)			
	• A 5'G + short 21 nucleotide antisense sequence (mature miRNA) derived from the target gene, followed by			
	• A short spacer of 19 nucleotides to form the terminal loop and			
	<ul> <li>A short sense target sequence with 2 nucleotides removed (Δ2) to create an internal loop</li> </ul>			
	The structural features are depicted in the figure below.			
	TGCT overhang 5'G + antisense Loop sequence Sense ∆2 nt target sequence CAGG overhang			
	For more details on the structure and guidelines to design the oligonucleotide refer to the BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kit manual.			
Pre-miRNA Expression Cassette	The engineered pre-miRNA sequence is cloned into the cloning site of BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vectors that is flanked on either side with sequences from murine miR-155 to allow proper processing of the engineered pre-miRNA sequence (see page 23 for the flanking region sequences).			
The pre-miRNA sequence and adjacent miR-155 flanking regions are demonstrated on the pre-miRNA expression cassette and are shown below. During the Gateware combination reactions, the pre-miRNA expression cassette is transferred vectors.				
	5' miR flanking s'G + antisense target sequence Loop sequence Sense ∆2 nt target sequence region 3' miR flanking region			
	EmGFP 5' miR flanking 5'G + antisense target sequence Loop sequence Sense ∆2 nt 3' miR flanking target sequence region			
	Once the engineered pre-miRNA expression cassette is introduced into the mammalian cells for expression, the pre-miRNA forms an intramolecular stem-loop structure similar to the structure of endogenous pre-miRNA that is then processed by the endogenous Dicer enzyme into a 22 nucleotide mature miRNA.			
	<b>Note:</b> The 21 nucleotides are derived from the target sequence while the 3'-most nucleotide is derived from the native miR-155 sequence (see figure on page 23).			

# The BLOCK-iT<sup>™</sup> Lentiviral Pol II miR RNAi Expression System

Components of	The system includes the following major components
the System	1. The BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kits containing the pcDNA <sup>™</sup> 6.2-GW/miR or pcDNA <sup>™</sup> 6.2-GW/EmGFP-miR vector for producing an expression clone that contains elements required for expression of a double-stranded oligonucleotide encoding an miRNA sequence of interest in mammalian cells using a Pol II promoter. The expression vector containing the pre-miRNA expression cassette may be transfected into mammalian cells for transient RNAi analysis or used to transfer the pre-miRNA expression cassette into the pLenti6/V5-DEST vector (see below) using Gateway <sup>®</sup> Technology. For detailed information about the BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kits and instructions for generating an expression clone, refer to the BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kit manual supplied with the kit.
	2. The pLenti6/V5-DEST vector allows easy transfer of the pre-miRNA expression cassette from the expression clone into a lentiviral destination vector for use with the Lentiviral system components. The destination vector contains the elements required to allow packaging of the expression construct into virions ( <i>e.g.</i> , 5' and 3' LTRs, $\psi$ packaging signal) and a selectable marker to allow generation of stable cell lines. For more information about the pLenti6/V5-DEST vector, see page 65.
	3. The pDONR <sup>™</sup> 221 vector is used as an intermediate to transfer the pre-miRNA expression cassette into the lentiviral expression plasmid (see above) using Gateway <sup>®</sup> Technology. See page 13 for details on pDONR <sup>™</sup> 221 vector.
	4. Gateway <sup>®</sup> BP and LR Clonase <sup>®</sup> II Enzyme Mixes that allow the transfer of the pre-miRNA expression cassette from the expression vectors into the pLenti6/V5-DEST vector using the Rapid BP/LR recombination reaction.
	5. One Shot <sup>®</sup> Stbl3 <sup>™</sup> Competent <i>E. coli</i> to obtain optimal results with lentiviral DNA after transformation.
	6. The ViraPower <sup>™</sup> Packaging Mix that contains an optimized mixture of the three packaging plasmids, pLP1, pLP2, and pLP/VSVG. These plasmids supply the helper functions as well as structural and replication proteins <i>in trans</i> required to produce the lentivirus. For more information about the packaging plasmids, see the <b>Appendix</b> , pages 70–75.
	7. An optimized 293FT producer cell line that stably expresses the SV40 large T-antigen under the control of the human CMV promoter and facilitates optimal production of virus. For more information about the 293FT Cell Line, refer to the 293FT Cell Line manual.

# The BLOCK-iT<sup>™</sup> Lentiviral Pol II miR RNAi Expression System, Continued

tra tai nu hc co all	You will co-transfect the ViraPower <sup>™</sup> Packaging Mix and the pLenti6/V5-DEST expression construct containing the pre-miRNA expression cassette into 293FT cells to produce a replication-incompetent lentivirus, which can then be transduced into the mammalian cell line of interest. Once the lentivirus enters the target cell, the viral RNA is reverse-transcribed, actively imported into the nucleus (Lewis & Emerman, 1994; Naldini, 1999), and stably integrated into the host genome (Buchschacher & Wong-Staal, 2000; Luciw, 1996). Once the lentiviral construct has integrated into the genome, the miRNA is constitutively expressed, allowing you to perform transient RNAi analysis or use Blasticidin selection to generate a stable cell line for long-term knockdown studies.		
Glycoprotein res RN G en	Most retroviral vectors are limited in their usefulness as delivery vehicles by their restricted tropism and generally low titers. In the BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Expression System, this limitation has been overcome by use of the G glycoprotein gene from Vesicular Stomatitis Virus (VSV-G) as a pseudotyping envelope, thus allowing production of a high titer lentivirus with a significantly broadened host cell range (Burns <i>et al.</i> , 1993; Emi <i>et al.</i> , 1991; Yee <i>et al.</i> , 1994).		
Features of the pLenti6/V5-DEST	he pLenti6/V5-DEST vector contains the following elements: Rous Sarcoma Virus (RSV) enhancer/promoter for Tat-independent production of viral mRNA in the producer cell line (Dull <i>et al.</i> , 1998) Modified HIV-1 5' and 3' Long Terminal Repeats (LTR) for viral packaging and reverse transcription of the viral mRNA (Dull <i>et al.</i> , 1998; Luciw, 1996) Note: The U3 region of the 3' LTR is deleted (ΔU3) and facilitates self-inactivation of the 5' LTR after transduction to enhance the biosafety of the vector (Dull <i>et al.</i> , 1998) HIV-1 psi (Ψ) packaging sequence for viral packaging (Luciw, 1996) HIV Rev response element (RRE) for Rev-dependent nuclear export of unspliced viral mRNA (Kjems <i>et al.</i> , 1991; Malim <i>et al.</i> , 1989) Human CMV promoter for high-level, constitutive expression of the miRNA from an RNA Polymerase II-dependent promoter Two recombination sites, <i>att</i> R1 and <i>att</i> R2, for recombinational cloning of the miRNA of interest from the pcDNA <sup>™</sup> 6.2-GW/±EmGFP-miR expression clone using Gateway <sup>®</sup> Technology Chloramphenicol resistance gene (Cm <sup>R</sup> ) located between the two <i>att</i> R sites for counterselection The <i>ccd</i> B gene located between the <i>att</i> R sites for negative selection Blasticidin resistance gene (Izumi <i>et al.</i> , 1991; Kimura <i>et al.</i> , 1994; Takeuchi <i>et al.</i> , 1958; Yamaguchi <i>et al.</i> , 1965) for selection in <i>E. coli</i> and mammalian cells Ampicillin resistance gene for selection in <i>E. coli</i>		

# Gateway<sup>®</sup> Recombination Reactions

Introduction	The Gateway <sup>®</sup> 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 a DNA sequence of interest into multiple vector systems.			
	Review the information in this section to briefly familiarize yourself with the Gateway <sup>®</sup> recombination reactions. For details, refer to the Gateway <sup>®</sup> Technology with Clonase <sup>®</sup> II manual available from at <u>www.invitrogen.com</u> or by contacting Technical Support (see page 79).			
Gateway <sup>®</sup> Vectors	Each of the vectors supplied in the BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Expression System is Gateway <sup>®</sup> -adapted, <i>i.e.</i> contains the appropriate <i>att</i> sites that allow site specific recombination to facilitate the transfer of heterologous DNA sequences between vectors.			
Recombination	Two recombination reactions constitute the basis of the Gateway <sup>®</sup> Technology:			
Reactions	BP Reaction			
	Facilitates recombination of an <i>attB</i> substrate ( <i>attB</i> -PCR product or a linearized <i>attB</i> expression clone) with an <i>attP</i> substrate (donor vector) to create an <i>attL</i> -containing entry clone. This reaction is catalyzed by Gateway <sup>®</sup> BP Clonase <sup>®</sup> II enzyme mix.			
	LR Reaction			
	Facilitates recombination of an <i>att</i> L substrate (entry clone) with an <i>att</i> R substrate (destination vector) to create an <i>att</i> B-containing expression clone. This reaction is catalyzed by Gateway <sup>®</sup> LR Clonase <sup>®</sup> II enzyme mix.			
Note	If you are an experienced user of Gateway <sup>®</sup> Technology and wish to perform the Rapid BP/LR recombination reaction, we recommend that you review the information on page 20 before proceeding with the experiment.			
	<b>Do not</b> use the standard recombination reaction conditions to perform the Rapid BP/LR recombination reaction.			

# Gateway<sup>®</sup> Recombination Reaction, Continued

Pre-miRNA Expression	Since the pcDNA <sup>™</sup> 6.2-GW/miR and pcDNA <sup>™</sup> 6.2-GW/EmGFP-miR expression vectors contain <i>att</i> B sites, the expressions vectors containing the pre-miRNA sequence <b>cannot</b> be transferred directly into the pLenti6/V5-DEST destination vector using a single recombination reaction. To transfer your pre-miRNA expression cassette from pcDNA <sup>™</sup> 6.2-GW/ <u>+</u> EmGFP-miR expression clone into the pLenti6/V5-DEST vector, you need to perform the two Gateway <sup>®</sup> recombination reactions as follows:
	<ol> <li>Generate an entry clone by performing a BP recombination reaction between the <i>att</i>B substrate (pcDNA<sup>™</sup>6.2-GW/miR or pcDNA<sup>™</sup>6.2- GW/EmGFP-miR expression clone) and <i>att</i>P substrate (pDONR<sup>™</sup>221 vector) using Gateway<sup>®</sup> BP Clonase<sup>®</sup> II Enzyme Mix.</li> </ol>
	<ol> <li>Perform an LR recombination reaction between the resulting entry clone (<i>att</i>L substrate) and pLenti6/V5-DEST vector (<i>att</i>R substrate) using Gateway<sup>®</sup> LR Clonase<sup>®</sup> II Enzyme Mix.</li> </ol>
	The standard BP and LR recombination reaction requires more than 2 days for completion. See below for details on expressing the miRNA from pLenti6/V5-DEST destination vector using the Rapid BP/LR Recombination Reaction.
Rapid BP/LR Recombination Reaction	We have developed a Rapid BP/LR recombination reaction to transfer the pre- miRNA expression cassettes into the destination vector that allows the completion of the entire BP and LR reaction in a day. In the Rapid BP/LR Recombination Reaction, instead of isolating the entry clone after BP reaction, the completed BP reaction is transferred directly into the LR reaction to generate expression clones.
	For <b>Rapid BP/LR Recombination Reactions</b> , perform a BP recombination reaction between the pcDNA <sup>™</sup> 6.2-GW/miR or pcDNA <sup>™</sup> 6.2-GW/EmGFP-miR expression clone and pDONR <sup>™</sup> 221 donor vector using Gateway <sup>®</sup> BP Clonase <sup>®</sup> II Enzyme Mix, then perform an LR recombination reaction between the resulting entry clone (pENTR <sup>™</sup> 221/miR) and pLenti6/V5-DEST vector using Gateway <sup>®</sup> LR Clonase <sup>®</sup> II Enzyme Mix (see below) to produce a lentiviral expression clone.
	attB attB attB pcDNA <sup>TM</sup> 6.2- GW/EmGFP-miR linearized expression clone attP attP attP pDONR <sup>TM</sup> 221 vector Gateway <sup>*</sup> BP Clonase <sup>*</sup> II vector Gateway <sup>*</sup> BP Clonase <sup>*</sup> II vector Gateway <sup>*</sup> BP Clonase <sup>*</sup> II vector
	attB attB
	Lentiviral expression clone

# Gateway<sup>®</sup> Recombination Reaction, Continued

Features of pDONR <sup>™</sup> 221	The pDONR <sup>™</sup> 221 vector contains the following elements:		
	• <i>rrn</i> B T1 and T2 transcription terminators for protection of the cloned gene or miRNA from expression by vector-encoded promoters		
	• Two recombination sites, <i>att</i> P1 and <i>att</i> P2, for recombinational cloning of the gene of interest from a Gateway <sup>®</sup> expression clone or <i>att</i> B PCR product		
	• <i>ccdB</i> gene located between the two <i>attP</i> sites for negative selection		
	Chloramphenicol resistance gene located between the two <i>att</i> P sites for counterselection		
	• Kanamycin resistance gene for selection in <i>E. coli</i>		
	• pUC origin for replication and maintenance of the plasmid in <i>E. coli</i>		

For a map of pDONR<sup>™</sup>221, see page 68.

### **Green Fluorescent Protein**

Description	The BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector with EmGFP contains the Emerald Green Fluorescent Protein (EmGFP) derived from <i>Aequorea victoria</i> GFP within the pre-miRNA expression cassette. After transferring the pre-miRNA expression cassette into pLenti6/V5-DEST, you may produce lentiviruses that simultaneously express the EmGFP protein and miRNA, allowing you to visually track the cells in which knockdown is occurring or sort the cells using a flow cytometer.	
Green Fluorescent Protein (GFP)	Green Fluorescent Protein (GFP) is a naturally occurring bioluminescent protein derived from the jellyfish <i>Aequorea victoria</i> (Shimomura <i>et al.</i> , 1962). GFP emits fluorescence upon excitation, and the gene encoding GFP contains all of the necessary information for posttranslational synthesis of the luminescent protein. GFP is often used as a molecular beacon because it requires no species-specific cofactors for function, and the fluorescence is easily detected using fluorescence microscopy and standard filter sets. GFP can function as a reporter gene downstream of a promoter of interest and upstream of one or more pre-miRNAs.	
GFP and Spectral Variants	Modifications have been made to the wild-type GFP to enhance its expression in mammalian systems. These modifications include amino acid substitutions that correspond to the codon preference for mammalian use, and mutations that increase the brightness of the fluorescence signal, resulting in "enhanced" GFP (Zhang <i>et al.</i> , 1996). Mutations have also arisen or have been introduced into GFP that further enhance and shift the spectral properties of GFP such that these proteins will emit fluorescent color variations (reviewed in Tsien, 1998). The Emerald GFP (EmGFP) is a variant of enhanced GFP.	
Note	We have observed reduced EmGFP expression from miRNA-containing vectors due to processing of the transcripts. In most cases, EmGFP expression should remain detectable.	

# Green Fluorescent Protein, Continued

EmGFP	The EmGFP variant has been described in a published review (Tsien, 1998) and is summarized below. The amino acid mutations are represented by the single letter abbreviation for the amino acid in the consensus GFP sequence, followed by the codon number and the single letter amino acid abbreviation for the substituted amino acid.		
	Fluorescent Protein	GFP Mutations*	
	EmGFP	S65T, S72A, N149K, M153T, I167T	
	*Mutations listed are as described in the literature. When examining the actual sequence the vector codon numbering starts at the first amino acid <b>after</b> the initiation methionine the fluorescent protein, so that mutations appear to be increased by one position. For example, the S65T mutation actually occurs in codon 66 of EmGFP.		
EmGFP Fluorescence	The EmGFP from the pcDNA <sup>™</sup> 6.2-GW/EmGFP-miR expression vector has the following excitation and emission wavelengths, as published in the literature (Tsien, 1998):		
	Excitation (nm) E	mission (nm)	
	487	509	
Filter Sets for Detecting EmGFP Fluorescence	The EmGFP can be detected with standard FITC filter sets. However, for optimized detection of the fluorescence signal, you may use a filter set which is optimized for detection within the excitation and emission ranges for the fluorescent proceeding the filter set for fluorescence microscopy and the manufacturer are listed below.		
	Filter Set	Manufacturer	
	-	Omega (www.omegafilters.com)	

# **Biosafety Features of the System**

Introduction	The lentiviral and packaging vectors supplied in the BLOCK-iT <sup>™</sup> Lentiviral RNAi Expression System are third-generation vectors based on lentiviral vectors developed by Dull <i>et al.</i> , 1998. This third-generation lentiviral system includes a significant number of safety features designed to enhance biosafety and to minimize relation to the wild-type, human HIV-1 virus. These safety features are discussed below.
Biosafety Features of the BLOCK-iT <sup>™</sup> Inducible H1 Lentiviral RNAi Expression System	<ul> <li>The BLOCK-iT<sup>™</sup> Lentiviral Pol II miR RNAi System includes the following key safety features:</li> <li>The pLenti6/V5-DEST expression vector contains a deletion in the 3' LTR (ΔU3) that does not affect generation of the viral genome in the producer cell line, but results in "self-inactivation" of the lentivirus after transduction of the target cell (Yee <i>et al.</i>, 1987; Yu <i>et al.</i>, 1986; Zufferey <i>et al.</i>, 1998). Once integrated into the transduced target cell, the lentiviral genome is no longer capable of producing packageable viral genome.</li> </ul>
	<ul> <li>The number of genes from HIV-1 that are used in the system has been reduced to three (<i>i.e. gag, pol,</i> and <i>rev</i>).</li> <li>The VSV-G gene from Vesicular Stomatitis Virus is used in place of the HIV-1 envelope (Burns <i>et al.</i>, 1993; Emi <i>et al.</i>, 1991; Yee <i>et al.</i>, 1994).</li> <li>Genes encoding the structural and other components required for packaging the viral genome are separated onto four plasmids. All four plasmids have been engineered not to contain any regions of homology with each other to prevent undesirable recombination events that could lead to the generation of a replication-competent virus (Dull <i>et al.</i>, 1998).</li> <li>Although the three packaging plasmids allow expression <i>in trans</i> of proteins required to produce viral progeny (<i>e.g.</i>, gal, pol, rev, env) in the 293FT producer cell line, none of them contain LTRs or the Ψ packaging sequence. This means that none of the HIV-1 structural genes are actually present in the packaged viral genome, and thus, are never expressed in the transduced target cell. No new replication-competent virus can be produced.</li> <li>The lentiviral particles produced in this system are replication-incompetent and only carry the gene of interest. No other viral species are produced.</li> <li>Expression of the <i>gag</i> and <i>pol</i> genes from pLP1 has been rendered Rev-dependent by virtue of the HIV-1 RRE in the gag/pol mRNA transcript. Addition of the RRE prevents gag and pol expression in the absence of Rev (Dull <i>et al.</i>, 1998).</li> <li>A constitutive promoter (RSV promoter) has been placed upstream of the 5' LTR in the pLenti6/V5-DEST expression vector to offset the requirement for Tat in the efficient production of viral RNA (Dull <i>et al.</i>, 1998).</li> </ul>

### Biosafety Features of the System, Continued

### **Biosafety Level 2**



Despite the inclusion of the safety features discussed on the previous page, the lentivirus produced with this system can still pose some biohazardous risk since it can transduce primary human cells. For this reason, we highly recommend that you treat lentiviral stocks generated using this System as Biosafety Level 2 (BL-2) organisms and strictly follow all published BL-2 guidelines with proper waste decontamination. Furthermore, exercise extra caution when creating lentivirus that express shRNA targeting human genes involved in controlling cell division (*e.g.*, tumor suppressor genes).

For more information about the BL-2 guidelines and lentivirus handling, refer to the document, "Biosafety in Microbiological and Biomedical Laboratories," 5<sup>th</sup> Edition, published by the Centers for Disease Control (CDC). This document may be downloaded at the following address:

http://www.cdc.gov/od/ohs/biosfty/bmbl5/bmbl5toc.htm

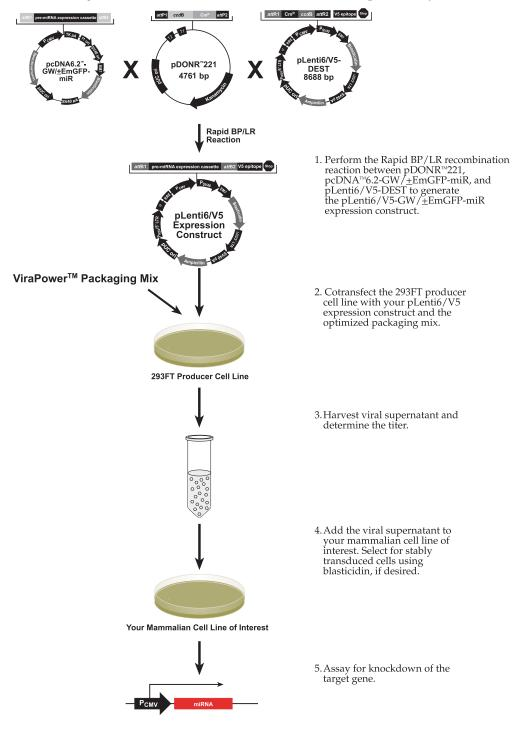


Handle all lentiviruses in compliance with established institutional guidelines. Since safety requirements for use and handling of lentiviruses may vary at individual institutions, we recommend consulting the health and safety guidelines and/or officers at your institution prior to use of the BLOCK-iT<sup>™</sup> Lentiviral Pol II miR RNAi Expression System.

### **Experimental Outline**

**Flow Chart** 

The diagram below describes the general steps required to express an shRNA of interest using the BLOCK-iT<sup>™</sup> Lentiviral Pol II miR RNAi Expression System.



### Methods

### Cloning miRNA

Introduction	You will need to clone your miRNA sequence of interest contained within the engineered pre-miRNA into the pcDNA <sup>™</sup> 6.2-GW/miR or pcDNA <sup>™</sup> 6.2-GW/EmGFP-miR expression vector using the BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kits to generate an expression clone prior to expressing the miRNA sequence of interest from pLenti6/V5-DEST. After generating the expression clone, you will transfer the pre-miRNA expression cassette from the expression clone into the destination vector, pLenti6/V5-DEST using a Rapid BP/LR recombination reaction (see next page).			
	General guidelines for cloning are provided below.			
Using pcDNA6.2- GW/ <u>+</u> EmGFP-miR	To generate an expression clone in pcDNA <sup>™</sup> 6.2-GW/miR or pcDNA <sup>™</sup> 6.2- GW/EmGFP-miR:			
	<ul> <li>Design and synthesize two complementary oligonucleotides containing the miRNA target sequence according to specified guidelines</li> </ul>			
	Anneal the oligonucleotides to create a double-stranded oligonucleotide			
	<ul> <li>Clone the double-stranded oligonucleotide into pcDNA<sup>™</sup>6.2-GW/miR or pcDNA<sup>™</sup>6.2-GW/EmGFP-miR using an optimized 5-minute ligation procedure</li> </ul>			
	• Transform competent <i>E. coli</i> and select for expression clones			
	For detailed instructions and guidelines for generating an expression clone, refer to the BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kit manual. This manual is supplied with the kits and is also available at <u>www.invitrogen.com</u> or by contacting Technical Support (see page 79).			

Introduction	Since the resulting pcDNA <sup>™</sup> 6.2-GW/miR and pcDNA <sup>™</sup> 6.2-GW/EmGFP-miR expression vectors contain <i>att</i> B sites, the expression vectors containing the pre- miRNA expression cassette <b>cannot</b> be used directly with the pLenti6/V5-DEST destination vector to perform the LR recombination reaction.
	To express your miRNA sequence in mammalian cells using the BLOCK-iT <sup>TM</sup> Lentiviral Pol II miR RNAi Expression System, you need to first generate an entry clone containing <i>att</i> L sites by performing a BP recombination reaction, then use the entry clone in an LR recombination reaction with pLenti6/V5-DEST vector to generate a lentiviral expression clone.
	The transfer of the miRNA sequence into the pLenti6/V5-DEST vector can be performed using the standard BP and LR recombination reactions or Rapid BP/LR recombination reactions as described below. See page 11 for an overview of the Gateway <sup>®</sup> recombination reactions.
	To ensure that you obtain the best possible results, we recommend that you read this section, the sections entitled <b>Performing the Rapid BP/LR Recombination Reaction</b> (page 25), and <b>Transforming One Shot</b> <sup>®</sup> <b>Stbl3<sup>™</sup> Competent</b> <i>E. coli</i> (page 29) before beginning.
Choosing a Suitable Protocol	Based on your experimental needs, you may choose between the standard or Rapid BP/LR recombination reactions as described in the table below:

If You Wish to	Then Choose	Described
Generate the expression clones using a fast protocol but obtain at least 10% fewer expression clones than the standard protocol	Rapid BP/LR Recombination Protocol	In this section.
Maximize the number of expression clones generated and isolate entry clones for future use	Standard BP and LR Protocols	On pages 57–62.

Rapid BP/LR	To express your miRNA sequence in mammalian cells using the BLOCK-iT <sup>™</sup>
Recombination	Lentiviral Pol II miR RNAi Expression System, perform the Rapid BP/LR
Reaction	recombination reactions as follows:
	Perform a BP recombination reaction between the pcDNA <sup>™</sup> 6.2-GW/miR or pcDNA <sup>™</sup> 6.2-GW/EmGFP-miR expression clone and pDONR <sup>™</sup> 221 donor vector using Gateway <sup>®</sup> BP Clonase <sup>®</sup> II Enzyme Mix. Then perform a LR recombination reaction between the resulting entry clone (pENTR <sup>™</sup> 221/miR) and pLenti6/V5-DEST vector using Gateway <sup>®</sup> LR Clonase <sup>®</sup> II Enzyme Mix to produce a lentiviral expression clone. See page 11 for an overview on Gateway <sup>®</sup> recombination reactions.

Experimental	To generate an expression clone, you will:
Outline	1. Perform the BP recombination reaction using the <i>att</i> B-expression clone with miR of interest and <i>att</i> P-containing pDONR <sup>™</sup> 221 vector to produce a pENTR <sup>™</sup> 221/miR entry clone.
	2. Mix an aliquot of the BP reaction (containing the entry clone) with the <i>att</i> R-containing pLenti6/V5-DEST vector and perform the LR recombination reaction to produce a lentiviral expression clone.
	3. Transform the reaction mixture into a suitable <i>E. coli</i> host (see page 29).
	4. Select for lentiviral expression clones (see page 23 for a diagram of the recombination region of expression clones in pLenti6/V5-DEST).
Substrates for the	To perform a BP recombination reaction, you need the following substrates:
Recombination Reactions	• Linearized <i>att</i> B-containing expression clones (see the next page for guidelines to linearize <i>att</i> B expression clones)
	• <i>att</i> P-containing donor (pDONR <sup>™</sup> 221) vector (see below)
	To perform an LR recombination reaction, you need the following substrates:
	• Supercoiled <i>att</i> L entry vector (pENTR <sup>™</sup> 221/miR)
	• Supercoiled <i>att</i> R destination vector (pLenti6/V5-DEST)
Donor Vectors	The BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Kits includes the pDONR <sup>™</sup> 221 vector. For a map and a description of the features of pDONR <sup>™</sup> 221, see the <b>Appendix</b> , page 69. You may use other donor vectors, if desired.
Important	The pLenti6/V5-DEST vector is supplied as a supercoiled plasmid. Although the Gateway <sup>®</sup> Technology manual previously recommended using a linearized destination vector for more efficient LR recombination, further testing has found that linearization of pLenti6/V5-DEST is <b>not</b> required to obtain optimal results for any downstream application.

Propagating the Donor and Destination Vectors	To propagate and maintain the pDONR <sup>™</sup> 221 and pLenti6/V5-DEST vectors, use One Shot <sup>®</sup> ccdB 2 T1 <sup>R</sup> Chemically Competent <i>E. coli</i> (see page 77) for transformation. The One Shot <sup>®</sup> ccdB Survival 2 T1 <sup>R</sup> <i>E. coli</i> strain is resistant to CcdB effects and can support the propagation of plasmids containing the ccdB gene.
	Donor Vector
	To maintain the integrity of pDONR <sup>™</sup> 221, select for transformants in media containing 50 µg/mL kanamycin and 15 µg/mL chloramphenicol.
	Destination Vector
	To maintain integrity of pLenti6/V5-DEST, select for transformants in media containing $50-100 \ \mu$ g/mL ampicillin and $15-30 \ \mu$ g/mL chloramphenicol.
	<b>Note: Do not</b> use general <i>E. coli</i> cloning strains including Stbl3 <sup>TM</sup> , TOP10, or DH5 $\alpha$ for propagation and maintenance, as these strains are sensitive to CcdB effects.
Linearizing Expression	For best results, linearize the expression clone using <i>Eag</i> I or <i>Bsr</i> D I (see the guidelines below).
Clones	<ol> <li>Linearize 1–2 μg of the expression clone with a restriction enzyme (<i>Eag</i> I or <i>Bsr</i>D I) that does not digest within the region of interest and is located outside the <i>att</i>B region.</li> </ol>
	<ol> <li>Ethanol precipitate the DNA after digestion by adding 0.1 volume of 3 M sodium acetate followed by 2.5 volumes of 100% ethanol.</li> </ol>
	3. Pellet the DNA by centrifugation. Wash the pellet twice with 70% ethanol.
	<ol> <li>Dissolve the DNA in TE Buffer, pH 8.0 to a final concentration of 50–150 ng/μL.</li> </ol>
	Continued on next nace

Recombination Region of pLenti6/V5-DEST		The recombination region of the lentiviral expression clone resulting from pLenti6/V5-DEST × pENTR <sup>™</sup> 221/miR is shown below. The pENTR <sup>™</sup> 221/miR entry clone is obtained by transferring the pre-miRNA expression cassette from pcDNA <sup>™</sup> 6.2-GW/miR into the pDONR <sup>™</sup> 221 vector.			
		Features of the Recombination Region:			
		Shaded regions correspond to DNA sequences transferred from the pENTR <sup>™</sup> 221/miR entry clone into the pLenti6/V5-DEST vector by recombination. Non-shaded regions are derived from the pLenti6/V5-DEST vector.			
	<b>Note:</b> The DNA sequences transferred from the pENTR <sup>™</sup> 221/miR entry clone contain the pre-miRNA expression cassette.				
2251	ТССТААСААС ТС	CAAT CMV forward priming site TATA CCGCCCCAT TGACGCAAAT GGGCGGTAGG CGTGTACGGT GGGAGGTCTA TATAAGCAGA GCTCGTTTAG			
2331		► Dional start ATCGCCTGG AGACGCCATC CACGCTGTTT TGACCTCCAT AGAAGACACC GACTCTAGAG GATCCACTAG			
		attB1			
2411	TCCAGTGTGG TG	GGAATTCTG CAGATATCAA CAAGTTTGTA CAAAAAAGCA GGCTTTAAAG GGAGGTAGTG AGTCGACCAG GTCTATAGTT GTTCAAACAT GTTTTTCGT CCGAAATTTC CCTCCATCAC TCAGCTGGTC			
	5	5' miR flanking region 3' miR flanking region			
2491		ggettgetg aaggetgta <mark>t getg pie-miRNA</mark> caggaea caaggeetgt tactageaet eacatggaae cegaaegae tteegaeata ega <mark>e dsoligo gtee</mark> tgt gtteeggaea atgategtga gtgtaeettg			
2561		ATCTGGCCG CACTCGAGAT ATCTAGAC CCA GCT TTC TTG TAC AAA GTG GTT GAT ATC CAG TAGACCGGC GTGAGCTCTA TAGATCTG GGT CGA AAG AAC ATG TTT CAC CAA CTA TAG GTC			
		attB2			
2632		GGC CGC TCG AGT CTA GAG GGC CCG CGG TTC GAA GGT AAG CCT ATC CCT AAC CCT CTC CCG GCG GCG AGC TCA GAA TTG GGA GAG			
		V5 epitope V5 (C-term) reverse priming site			
2698		GAT TCT ACG CGT ACC GGT TAG TAA TGA GTTT CTA AGA TGC GCA TGG CCA ATC ATT ACT 			



Since the pLenti6-V5-GW/miR expression construct expresses a pre-miRNA sequence that is processed to form a mature miRNA and not a protein, the V5 epitope is not expressed.

Recombination Region of pLenti6/V5-DEST, continued		The recombination region of the lentiviral expression clone resulting from pLenti6/V5-DEST × pENTR <sup>™</sup> 221/EmGFP-miR is shown below. The pENTR <sup>™</sup> 221/EmGFP-miR entry clone is obtained by transferring the pre-miRNA expression cassette from pcDNA <sup>™</sup> 6.2-GW/EmGFP-miR into the pDONR <sup>™</sup> 221 vector.			
	Features of the Recombination Region:				
		Shaded regions correspond to DNA sequences transferred from the pENTR <sup>™</sup> 221/EmGFP-miR entry clone into the pLenti6/V5-DEST vector by recombination. Non-shaded regions are derived from the pLenti6/V5-DEST vector. Bases 2,473 and 3,192 indicate the coding sequence of EmGFP.			
<b>Note:</b> The DNA sequences transferred from the pENTR <sup>™</sup> 221/miR entry clone contain the pre-miRNA expression cassette including EmGFP coding sequence.					
		CAAT CMV forward priming site TATA			
2251	TCGTAACAAC T	CCGCCCCAT TGACGCAAAT GGGCGGTAGG CGTGTACGGT GGGAGGTCTA TATAAGCAGA GCTCGTTTAG			
2331	1	► ptional start ATCGCCTGG AGACGCCATC CACGCTGTTT TGACCTCCAT AGAAGACACC GACTCTAGAG GATCCACTAG			
		attB1 2473			
2411	TCCAGTGTGG T	GGAATTCTG CAGATATCAA CAAGTTTGTA CAAAAAAGCA GGCTTTAAAA CC <b>ATG</b> GTG AGC AAG			
		GTCTATAGTT GTTCAAACAT GTTTTTTCGT CCGAAATTTT GG TAC CAC TCG TTC			
	EmGFP coding seque	nce			
2485	GGC CCG EmGFP	GGC ATG GAC GAG CTG TAC AAG TAA GCT AAGCACTTCG TGGCCGTCGA TCGTTTAAAG CCG TAC CTG CTC GAC ATG TTC ATT CGA TTCGTGAAGC ACCGGCAGCT AGCAAATTTC			
	Gly	Gly Met Asp Glu Leu Tyr Lys ***			
		5' miR flanking region 3' miR flanking region			
3226		GTCGACCAG TGGATCCTGG AGGCTTGCTG AAGGCTGTA <mark>T GCTG pre-miRNA</mark> CAGGACA CAAGGCCTGT CAGCTGGTC ACCTAGGACC TCCGAACGAC TTCCGACATA CGA <mark>C dsoligo GTCC</mark> TGT GTTCCGGACA			
3296	TACTAGCACT C	ACATGGAAC AAATGGCCCA GATCTGGCCG CACTCGAGAT ATCTAGAC CCA GCT TTC TTG TAC			
	ATGATCGTGA G	TGTACCTTG TTTACCGGGT CTAGACCGGC GTGAGCTCTA TAGATCTG GGT CGA AAG AAC ATG			
3369		attB2 GAT ATC CAG CAC AGT GGC GGC CGC TCG AGT CTA GAG GGC CCG CGG TTC GAA GGT CTA TAG GTC GTG TCA CCG CCG GCG AGC TCA GAT CTC CCG GGC GCC AAG CTT CCA			
	NAC COM NUC				
3432		CCT AAC CCT CTC CTC GGT CTC GAT TCT ACG CGT ACC GGT TAG TAA TGA GTTT GGA TTG GGA GAG GAG CCA GAG CTA AGA TGC CGA TGG CCA ATC ATT ACT			
		/5 (C-term) reverse priming site V5 epitope			



Since the pLenti6-V5-GW/EmGFP-miR expression construct expresses a premiRNA sequence that is processed to form a mature miRNA and not a protein, the V5 epitope is not expressed.

# Performing the Rapid BP/LR Recombination Reaction

Introduction	Follow the guidelines and instructions in this section to perform the Rapid BP/LR recombination reaction using the expression clone containing your pre-miRNA expression cassette, pDONR <sup>™</sup> 221, and pLenti6/V5-DEST vector.
	If you wish to perform the standard BP recombination reaction followed by the standard LR recombination reaction, see page 57 for details.
Rapid BP/LR Protocol	The Rapid BP/LR protocol is used to transfer a gene from one expression clone into another destination vector in 2 consecutive steps: a BP reaction using a donor vector followed by an LR recombination reaction using a destination vector without purification of the intermediate entry clone.
	<b>Note:</b> The Rapid BP/LR protocol allows you to generate expression clones more rapidly than the standard BP and LR protocols provided on page 57. Fewer expression clones are obtained (~10% of the total number of expression clones) using the Rapid BP/LR protocol. If you wish to maximize the number of expression clones generated, <b>do not</b> use this protocol. Use the standard BP and LR recombination protocols on page 57.
Experimental	To perform the Rapid BP/LR protocol, you will:
Outline	<ol> <li>Perform a BP recombination reaction using the linearized expression clone containing your pre-miRNA sequence and pDONR<sup>™</sup>221 to generate the entry clone, pENTR<sup>™</sup>221/miR.</li> </ol>
	<ol> <li>Use a small aliquot of the BP reaction mix to perform the LR recombination reaction using the pLenti6/V5-DEST destination vector to generate the lentiviral expression clone, pLenti6/V5-GW/miR.</li> </ol>
	3. Perform Proteinase K treatment.
Recommended <i>E. coli</i> Host	For optimal results, use Stbl3 <sup>™</sup> E. coli for transformation as this strain is particularly well-suited for use in cloning unstable DNA such as lentiviral DNA, which contains direct repeats. One Shot <sup>®</sup> Stbl3 <sup>™</sup> Chemically Competent <i>E. coli</i> are included in the kit for transformation. For instructions, see Transforming One Shot <sup>®</sup> Stbl3 <sup>™</sup> Competent <i>E. coli</i> , page 29.
Important	<b>Do not</b> transform the BP or LR recombination reaction into <i>E. 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.

# Performing Performing the Rapid BP/LR Recombination

Reaction, Continued

Positive Control	We recommend using the pcDNA <sup>™</sup> 6.2-GW/miR-neg Control Plasmid supplied with the BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Kits as a positive control for the Rapid BP/LR protocol. Dilute the supplied control plasmid 1:10 in sterile water to obtain a final concentration of 50 ng/µL. <b>Do not</b> use the pEXP7-tet supplied with the Gateway <sup>®</sup> BP Clonase <sup>®</sup> II Enzyme Mix or pENTR <sup>™</sup> -gus supplied with the Gateway <sup>®</sup> LR Clonase <sup>®</sup> II Enzyme Mix as positive controls for the Rapid protocol due to the presence of incompatible selection markers.
Gateway <sup>®</sup> Clonase <sup>®</sup> II Enzyme Mixes	The Gateway <sup>®</sup> BP and Gateway <sup>®</sup> LR Clonase <sup>®</sup> II enzyme mixes combine the proprietary enzyme formulation and 5X Clonase Reaction Buffer previously supplied as separate components in Gateway <sup>®</sup> Clonase <sup>®</sup> enzyme mixes into an optimized single-tube format for easier set-up of the BP or LR recombination reaction. The Gateway <sup>®</sup> LR Clonase <sup>®</sup> II Enzyme catalyzes the <i>attL</i> x <i>attR</i> Gateway <sup>®</sup> recombination reaction. Use the protocol provided on page 27 to perform the recombination reactions using the Rapid protocol or page 57 using the standard protocol. Gateway <sup>®</sup> BP and LR Clonase <sup>®</sup> II Enzyme Mixes are supplied with the kit and are also available separately (see page 77). Note: You may perform the BP or LR recombination reaction with the Gateway <sup>®</sup> BP or LR Clonase <sup>®</sup> enzyme mix, if desired. To use the Gateway <sup>®</sup> BP or LR Clonase <sup>®</sup> enzyme mix, follow the protocol provided in this manual.
Converting Femtomoles (fmol) to Nanograms (ng)	Use the following formula to convert femtomoles (fmol) of DNA to nanograms (ng) of DNA required for BP reaction: $ng = (fmol)(N)(\frac{660 \text{ fg}}{fmol})(\frac{1 \text{ ng}}{10^6 \text{ fg}})$ where N is the size of the DNA in bp. For an example, see below. In this example, you need to use 50 fmol of an <i>att</i> B-PCR product in the BP reaction. The <i>att</i> B-PCR product is 2.5 kb in size. Calculate the amount of <i>att</i> B-PCR product required for the reaction (in ng) by using the above equation: $(50 \text{ fmol})(2500 \text{ bp})(\frac{660 \text{ fg}}{fmol})(\frac{1 \text{ ng}}{10^6 \text{ fg}}) = 82.5 \text{ ng of PCR product required}$

# Performing Performing the Rapid BP/LR Recombination

Reaction, Continued

Materials Needed	<ul> <li>Linearized expression clone (50–150 ng/µL in TE Buffer, pl</li> <li>TE Buffer, pH 8.0 (10 mM Tris-HCl, pH 8.0, 1 mM EDTA), s</li> <li>Sterile 0.5 mL microcentrifuge tubes</li> <li><i>Components supplied with the kits</i></li> <li>pDONR<sup>™</sup>221 vector</li> <li>pLenti6/V5-DEST vector</li> <li>pcDNA<sup>™</sup>6.2-GW/±EmGFP-miR-neg control (if desired, sup BLOCK-iT<sup>™</sup> Pol II miR RNAi Expression Vector Kits, Box 1</li> <li>Gateway<sup>®</sup> BP Clonase<sup>®</sup> II enzyme mix (Box 9; store at -20°C immediately before use)</li> <li>Gateway<sup>®</sup> LR Clonase<sup>®</sup> II enzyme mix (Box 10; store at -20°C immediately before use)</li> <li>2 µg/µL Proteinase K solution (supplied with Gateway<sup>®</sup> BI</li> </ul>	see page 7 pplied with ) C until °C until	n	
Setting Up the	<ul><li>enzymes; thaw and keep on ice until use)</li><li>1. Add the following components to sterile 0.5 mL microcentr</li></ul>			
Rapid BP/LR Recombination	room temperature and mix. Component	Sample	Positive	
Reaction	<b>r</b>		Control	
	Linearized <i>att</i> B expression clone from page 22, (20–50 fmol)	1–7 µL	-	
	pcDNA <sup>™</sup> 6.2-GW/ <u>+</u> EmGFPmiR-neg control (diluted to 50 ng/μL)	_	2 μL	
	pDONR <sup>™</sup> 221 vector (150 ng/µL)	1 µL	1 µL	
	TE Buffer, pH 8.0	to 8 µL	5 µL	
	<ol> <li>Remove the Gateway<sup>®</sup> BP Clonase<sup>®</sup> II enzyme mix from −20°C and thaw on ice (~ 2 minutes).</li> </ol>			
	3. Vortex the Gateway <sup>®</sup> BP Clonase <sup>®</sup> II enzyme mix briefly tw time).	ice (2 seco	nds each	
	<ul> <li>time).</li> <li>4. To the samples above, add 2 μL of Gateway<sup>®</sup> BP Clonase<sup>®</sup> I</li> </ul>	I enzyme 1	nix.	
	<ul> <li>time).</li> <li>To the samples above, add 2 µL of Gateway<sup>®</sup> BP Clonase<sup>®</sup> I Mix well by pipetting up and down.</li> <li>Reminder: Return Gateway<sup>®</sup> BP Clonase<sup>®</sup> II enzyme mix to -20</li> </ul>	I enzyme 1	nix.	

### Performing Performing the Rapid BP/LR Recombination

Reaction, Continued

Setting Up the Rapid BP/LR Recombination Reaction, Continued 6. Transfer 3 µL from each BP reaction from Step 5 to clean, sterile 0.5 mL microcentrifuge tubes. This reaction mix contains the resulting **entry clone**, pENTR<sup>™</sup>221/miR with or without EmGFP.

**Note:** Save the remaining BP reaction mix at  $-20^{\circ}$ C for up to 1 week. You can treat the samples with Proteinase K and transform the reaction mix into One Shot<sup>®</sup> TOP10 Chemically Competent *E. coli* as described on page 60 to check the efficiency of the BP reaction and to isolate entry clones for future use. For transformation of the BP reaction only, you can use any *E. coli* including TOP10.

 Add the following components to the tubes containing 3 µL BP reaction from Step 6 at room temperature and mix.

Component	Sample	Positive Control
pLenti6/V5-DEST vector (150 ng/µL)	1 µL	1 μL
TE Buffer, pH 8.0	4 μL	$4 \ \mu L$

- 8. Remove the Gateway<sup>®</sup> LR Clonase<sup>®</sup> II enzyme mix from −20°C and thaw on ice (~ 2 minutes).
- 9. Vortex the Gateway<sup>®</sup> LR Clonase<sup>®</sup> II enzyme mix briefly twice (2 seconds each time).
- 10. To the samples above, add 2 µL of Gateway<sup>®</sup> LR Clonase<sup>®</sup> II enzyme mix. Mix well by pipetting up and down.

Reminder: Return Gateway  $^{\! \mathrm{\$}}$  LR Clonase  $^{\! \mathrm{\$}}$  II enzyme mix to –20°C immediately after use.

11. Incubate the reaction at 25°C for 2–4 hours.

**Note:** The incubation time may be extended from 4 hours to overnight, if more colonies are required.

- 12. Add 1  $\mu$ L of the Proteinase K solution to each reaction. Incubate for 10 minutes at 37°C.
- 13. Proceed to **Transforming One Shot**<sup>®</sup> **Stbl3<sup>™</sup> Competent** *E. coli*, next page. **Note:** You may store the reaction at -20°C for up to 1 week before transformation, if desired.

# Transforming One Shot<sup>®</sup> Stbl3<sup>™</sup> Competent *E. coli*

Introduction	Follow the instructions in this section to transform the LR recombination reaction nto One Shot <sup>®</sup> Stbl3 <sup>TM</sup> Chemically Competent <i>E. coli</i> (Box 8) included with the kit. The transformation efficiency of One Shot <sup>®</sup> Stbl3 <sup>TM</sup> Chemically Competent <i>E. coli</i> is $\frac{1}{2}$ × 10 <sup>8</sup> cfu/µg plasmid DNA.
Materials Needed	<ul> <li>LR recombination reaction (from Step 13, previous page or Step 7, page 62)</li> <li>LB Medium (if performing the pUC19 control transformation)</li> <li>42°C water bath</li> <li>LB plates containing 100 µg/mL ampicillin (two for each transformation; warm at 37°C for 30 minutes before use)</li> <li>37°C shaking and non-shaking incubator</li> <li>Components supplied with the kit</li> <li>One Shot<sup>®</sup> Stbl3<sup>™</sup> Chemically Competent <i>E. coli</i> (Box 8; one vial per transformation; thaw on ice immediately before use)</li> <li>S.O.C. Medium (Box 8; warm to room temperature)</li> <li>pUC19 positive control (if desired to verify the transformation efficiency; Box 8)</li> </ul>
One Shot <sup>®</sup> Stbl3 <sup>™</sup> Transformation	Jse this procedure to transform the LR recombination reaction into One Shot <sup>®</sup> Stbl3 <sup><math>T</math></sup> Chemically Competent <i>E. coli</i> .
Procedure	. Thaw, on ice, one vial of One Shot <sup>®</sup> Stbl3 <sup>™</sup> chemically competent cells for each transformation.
	Add 2 to 3 µL of the LR recombination reaction (from Step 13, previous page or Step 7, page 62) into a vial of One Shot <sup>®</sup> Stbl3 <sup>™</sup> cells and mix gently. <b>Do not mix</b> <b>by pipetting up and down.</b> For the pUC19 control, add 10 pg (1 µL) of DNA into a separate vial of One Shot <sup>®</sup> cells and mix gently.
	8. Incubate the vial(s) on ice for 30 minutes.
	. Heat-shock the cells for 45 seconds at 42°C without shaking.
	<ol> <li>Remove the vial(s) from the 42°C water bath and place them on ice for 2 minutes.</li> </ol>
	. Add 250 μL of pre-warmed S.O.C. Medium to each vial.
	7. Cap the vial(s) tightly and shake horizontally at 37°C for 1 hour at 225 rpm in a shaking incubator.
	8. Spread 25–100 $\mu$ L of the transformation mix on a pre-warmed selective plate and incubate overnight at 37°C. We recommend plating two different volumes to ensure that at least one plate will have well-spaced colonies. For the pUC19 control, dilute the transformation mix 1:10 into LB Medium ( <i>e.g.</i> , add 100 $\mu$ L of the transformation mix to 900 $\mu$ L of LB Medium) and plate 25–100 $\mu$ L.
	<ol> <li>Store the remaining transformation mix at 4°C. Plate out additional cells the next day, if desired.</li> </ol>

# Transforming One Shot<sup>®</sup> Stbl3<sup>™</sup> Competent *E. coli,* Continued

Expected Results	When using One Shot <sup>®</sup> Stbl3 <sup>™</sup> Chemically Competent cells for transformation, the LR recombination reaction should result in greater than 4,000 colonies if the entire LR 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 chloramphenicol-sensitive <b>and</b> ampicillin- and Blasticidin-resistant. Transformants containing a plasmid with a mutated <i>ccdB</i> gene will be chloramphenicol-, ampicillin-, and Blasticidin-resistant. To check your putative expression clone, test for growth on LB plates containing 30 µg/mL chloramphenicol. A true expression clone should not grow in the presence of chloramphenicol.			
Sequencing	Sequencing the expression construct is not required as transfer of the miRNA cassette from pcDNA <sup>™</sup> 6.2-GW/±EmGFP-miR into the pLenti6/V5-DEST vector preserves the orientation of the cassette. However, if you wish to sequence the pLenti6/V5 expression construct, we recommend using the following primers. Refer to the diagram on page 23 for the location of the primer binding sites in the expression vector.			
	Primer	Sequence		
	CMV Forward	5'-CGCAAATGGGCGGTAGGCGTG-3'		
	V5(C-term) Reverse	5'-ACCGAGGAGAGGGTTAGGGAT-3'		
	<b>Note:</b> For information about a convenient custom primer synthesis service go to <u>www.invitrogen.com</u> or call Technical Support (see page 79).			
Maintaining the Expression Clone	Once you have generated your expression clone, maintain and propagate the expression clone in LB medium containing $100 \ \mu g/mL$ ampicillin.			

## **Producing Lentivirus in 293FT Cells**

Introduction	Before creating a stably transduced cell line expressing your shRNA of interest, first produce a lentiviral stock (containing the packaged pLenti6/V5 expression construct) by co-transfecting the optimized ViraPower <sup>™</sup> Packaging Mix and your pLenti6/V5-GW/miR expression construct into the 293FT Producer Cell Line.		
Experimental	To produce lentivirus in 293FT Cells, you will:		
Outline	1. Grow the 293FT Cells to obtain $6 \times 10^6$ 293FT cells for each sample.		
	2. Prepare plasmid DNA of your expression clone.		
	3. Cotransfect the ViraPower <sup>™</sup> Packaging Mix and pLenti6/V5-GW/miR expression plasmid DNA into 293FT Cells using Lipofectamine <sup>®</sup> 2000.		
	4. Harvest virus-containing supernatants 48–72 hours post-transfection.		
293FT Cell Line	The human 293FT Cell Line is supplied with the BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Kits to facilitate optimal lentivirus production (Naldini <i>et al.</i> , 1996). The 293FT Cell Line, a derivative of the 293F Cell Line, stably and constitutively expresses the SV40 large T-antigen from pCMVSPORT6TAg.neo and must be maintained in medium containing Geneticin <sup>®</sup> . For more information about pCMVSPORT6TAg.neo and how to culture and maintain 293FT cells, refer to the 293FT Cell Line manual. This manual is supplied with the BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Kits and is also available at <u>www.invitrogen.com</u> or by contacting Technical Support (see page 79). <b>Note:</b> The 293FT Cell Line is also available separately (see page 77).		
	The health of your 293FT cells at the time of transfection has a critical effect on the success of lentivirus production. Use of "unhealthy" cells can negatively affect the transfection efficiency, resulting in production of a low titer lentiviral stock. For optimal lentivirus production ( <i>i.e.</i> producing lentiviral stocks with the expected titers), follow the guidelines below to culture 293FT cells before use in transfection:		
	• Make sure that cells are greater than 90% viable.		
	<ul> <li>Subculture and maintain cells as recommended in the 293FT Cell Line manual. Do not allow cells to overgrow before passaging. You will need 6 × 10<sup>6</sup> 293FT cells for each sample.</li> </ul>		

• Use cells that have been subcultured for less than 20 passages.

ViraPower™ Packaging Mix	The pLP1, pLP2, pLP/VSVG plasmids are provided in an optimized mixture to facilitate viral packaging of the pLenti6/V5-GW/miR expression vector following cotransfection into 293FT producer cells.			
	The amount of the packaging mix (195 µg) and Lipofectamine <sup>®</sup> 2000 Reagent (0.75 mL) supplied in the BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Kits is sufficient to perform 20 cotransfections in 10 cm plates using the recommended protocol on page 36.			
	Note: The ViraPower <sup>™</sup> Packaging Mix is available separately (page 77) or as part of the ViraPower <sup>™</sup> Bsd Lentiviral Support Kit (see page 77).			
Plasmid Preparation	Once you have generated the expression clone, you <b>must</b> isolate plasmid DNA for transfection. Plasmid DNA for transfection into eukaryotic cells must be very 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. Isolate plasmid DNA using the PureLink <sup>™</sup> Plasmid Purification Kits (page 77) or CsCl gradient centrifugation.			
	Resuspend the purified pLenti6/V5-GW/miR expression plasmid in sterile water or TE Buffer, pH 8.0, to a final concentration ranging from 0.1–3.0 $\mu$ g/ $\mu$ L. You will need 3 $\mu$ g of the expression plasmid for each transfection.			
	Important: Do not use mini-prep plasmid DNA for transfection.			
Lipofectamine <sup>®</sup> 2000	The Lipofectamine <sup>®</sup> 2000 reagent supplied with the BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Kits is a proprietary, cationic lipid-based formulation suitable for the transfection of nucleic acids into eukaryotic cells (Ciccarone <i>et al.</i> , 1999). Using Lipofectamine <sup>®</sup> 2000 to transfect 293FT cells offers the following advantages:			
	Provides the highest transfection efficiency in 293FT cells			
	<ul> <li>DNA-Lipofectamine<sup>®</sup> 2000 complexes can be added directly to cells in culture medium in the presence of serum</li> </ul>			
	• Removal of complexes or medium change or addition following transfection is not required, although complexes can be removed after 4–6 hours without loss of activity			
	Note: Lipofectamine <sup>®</sup> 2000 is available separately or as part of the ViraPower <sup>™</sup> Bsd Lentiviral Support Kit (see page 77).			
Opti-MEM <sup>®</sup> I	To facilitate optimal formation of DNA-Lipofectamine <sup>®</sup> 2000 complexes, use Opti-MEM <sup>®</sup> I Reduced Serum Medium (see page 77). For more information about Opti-MEM <sup>®</sup> I, see <u>www.invitrogen.com</u> or contact Technical Support (see page 79).			

miR Positive Control	Generate the miR Positive Control using the reagents included in the kit as follows:		
	<ul> <li>Generate the pcDNA<sup>™</sup>6.2-GW/<u>+</u>EmGFP-miR-<i>lacZ</i> expression control using the <i>lacZ</i> double-stranded oligo and pcDNA<sup>™</sup>6.2-GW/<u>+</u>EmGFP-miR expression vector included with the BLOCK-iT<sup>™</sup> Pol II miR RNAi Expression Vector Kit and as described in the expression vector manual.</li> </ul>		
	• Use the pcDNA <sup>™</sup> 6.2-GW/ <u>+</u> EmGFP-miR- <i>lacZ</i> expression control to generate the lentiviral construct with pLenti6-V5-DEST vector using the Rapid BP/LR recombination reaction as described in this manual.		
	• Use the resulting lentiviral expression construct, pLenti6/V5-GW/ <u>+</u> EmGFP- miR- <i>lacZ</i> , to generate a miR control lentiviral stock ( <i>lacZ</i> targeting miRNA).		
	Once generated, the miR control lentivirus may be transduced into mammalian cell lines (see page 47) to express an miRNA targeted to the human <i>lacZ</i> gene, and may be used as a control for the RNAi response in these cell lines.		
pLenti6/V5- GW/ <i>lacZ</i> Positive Control	A pLenti6/V5-GW/ <i>lacZ</i> positive control vector is included with the pLenti6/V5- DEST vector for use as an expression control in the ViraPower <sup>™</sup> Lentiviral Expression System. The β-galactosidase is expressed as a C-terminally tagged fusion protein that may be easily detected by western blot or functional assay. For details on the vector, see page 67.		
	To propagate and maintain the control plasmid:		
	<ol> <li>Use the vector stock solution to transform a <i>recA</i>, <i>endA E</i>. <i>coli</i> strain like Stbl3<sup>™</sup>, TOP10, DH5α<sup>™</sup>-T1<sup>R</sup>, or equivalent. Use 10 ng of plasmid for transformation.</li> </ol>		
	<ol> <li>Select transformants on LB agar plates containing 100 µg/mL ampicillin (for Stbl3<sup>™</sup> cells) or LB agar plates containing 100 µg/mL ampicillin and 50 µg/mL Blasticidin (for TOP10 or DH5α).</li> </ol>		
	3. Prepare a glycerol stock of a transformant containing plasmid for long-term storage. Propagate the plasmid in LB containing 100 $\mu$ g/mL ampicillin.		
	4. Use the pLenti6/V5-GW/ <i>lacZ</i> positive control to generate a control lentiviral stock (expressing the LacZ protein).		
	<ol> <li>Use the pLenti6/V5-GW/<i>lacZ</i> lentiviral control and the pLenti6/V5-GW/<u>±</u>EmGFP-miR-<i>lacZ</i> lentiviral control in cotransduction experiments as a positive control for lentiviral induced RNAi analysis in your system (see page 46 for details).</li> </ol>		

Positive controls (see previous page to generate positive controls; resuspend in sterile water to a concentration of 1 μg/μL) 293FT cells cultured in the appropriate medium ( <i>i.e.</i> D-MEM supplemented with 10% FBS, 2 mM L-glutamine, 0.1 mM MEM Non-Essential Amino Acids, and 1% penicillin/streptomycin). You will need 6 × 10 <sup>6</sup> 293FT cells for each sample. Note: D-MEM already contains 4 mM L-glutamine, which is enough to support cell growth of the 293FT Cell Line. However, since L-glutamine slowly decays over time, supplement the medium with 2 mM L-glutamine to ensure that the concentration of L-glutamine will not get too low over time due to its slow degradation. 293FT cells grow vell in 6 mM L-glutamine, but higher concentrations of L-glutamine may reduce growth. Dpti-MEM <sup>®</sup> I Reduced Serum Medium (pre-warmed; see page 32) Fetal bovine serum (FBS; see page 77) Complete growth medium containing sodium pyruvate ( <i>i.e.</i> D-MEM supplemented with 10% FBS, 2 mM L-glutamine, 0.1 mM MEM Non-Essential Amino Acids, 1% penicillin/streptomycin, and 1 mM MEM Sodium Pyruvate)
<ul> <li>with 10% FBS, 2 mM L-glutamine, 0.1 mM MEM Non-Essential Amino Acids, and 1% penicillin/streptomycin). You will need 6 × 10<sup>6</sup> 293FT cells for each sample.</li> <li>Note: D-MEM already contains 4 mM L-glutamine, which is enough to support cell growth of the 293FT Cell Line. However, since L-glutamine slowly decays over time, supplement the medium with 2 mM L-glutamine to ensure that the concentration of L-glutamine will not get too low over time due to its slow degradation. 293FT cells grow vell in 6 mM L-glutamine, but higher concentrations of L-glutamine may reduce growth.</li> <li>Dpti-MEM<sup>®</sup> I Reduced Serum Medium (pre-warmed; see page 32)</li> <li>Fetal bovine serum (FBS; see page 77)</li> <li>Complete growth medium containing sodium pyruvate (<i>i.e.</i> D-MEM supplemented with 10% FBS, 2 mM L-glutamine, 0.1 mM MEM Non-Essential</li> </ul>
growth of the 293FT Cell Line. However, since L-glutamine slowly decays over time, supplement the medium with 2 mM L-glutamine to ensure that the concentration of L-glutamine will not get too low over time due to its slow degradation. 293FT cells grow well in 6 mM L-glutamine, but higher concentrations of L-glutamine may reduce growth. Dpti-MEM <sup>®</sup> I Reduced Serum Medium (pre-warmed; see page 32) Fetal bovine serum (FBS; see page 77) Complete growth medium containing sodium pyruvate ( <i>i.e.</i> D-MEM supplemented with 10% FBS, 2 mM L-glutamine, 0.1 mM MEM Non-Essential
Fetal bovine serum (FBS; see page 77) Complete growth medium containing sodium pyruvate ( <i>i.e.</i> D-MEM supplemented with 10% FBS, 2 mM L-glutamine, 0.1 mM MEM Non-Essential
Complete growth medium containing sodium pyruvate ( <i>i.e.</i> D-MEM supplemented with 10% FBS, 2 mM L-glutamine, 0.1 mM MEM Non-Essential
supplemented with 10% FBS, 2 mM L-glutamine, 0.1 mM MEM Non-Essential
<b>Note:</b> MEM Sodium Pyruvate provides an extra energy source for the cells and is available separately, see page 77. See note above for L-glutamine concentration.
Sterile, 10 cm tissue culture plates (one each for the lentiviral construct, positive control, and negative control)
Sterile, tissue culture supplies
and 15 mL sterile, capped, conical tubes
Cryovials
ponents supplied with the kits
/iraPower <sup>™</sup> Packaging Mix
Lipofectamine <sup>®</sup> 2000 transfection reagent (store at 4°C and mix gently before use)
Continued on next page

#### Recommended Transfection Conditions

We produce lentiviral stocks in 293FT cells using the **optimized** transfection conditions shown below. The amount of lentivirus produced using these recommended conditions (at a titer of  $1 \times 10^5$  to  $1 \times 10^7$  transducing units (TU)/mL) is generally sufficient to transduce  $1 \times 10^6$  to  $1 \times 10^8$  cells at a multiplicity of infection (MOI) = 1.

Condition	Amount
Tissue culture plate size	10 cm (one per lentiviral construct)
Number of 293FT cells to transfect	$6 \times 10^{6}$ cells (see <b>Recommendation</b> on page 31 to prepare cells for transfection)
Amount of ViraPower <sup>™</sup> Packaging Mix	9 μg (9 μL of 1 μg/μL stock)
Amount of pLenti6/V5-GW/miR expression plasmid	3 µg
Amount of Lipofectamine <sup>®</sup> 2000 Reagent to use	36 µL

**Note:** You may produce lentiviral stocks using other tissue culture formats, but keep in mind that optimization will be necessary to obtain the expected titers.

The recommended procedure to co-transfect 293FT cells differs from the traditional Lipofectamine<sup>®</sup> 2000 transfection procedure in that you will:

- First prepare DNA:Lipofectamine<sup>®</sup> 2000 complexes and add them to plates containing growth media, then
- Add the 293FT cells to the media containing DNA:Lipofectamine<sup>®</sup> 2000 complexes and allow the cells to attach and transfect overnight (see details on the next page).

Using this procedure, we consistently obtain lentiviral stocks with titers that are **3 to 4-fold higher** than lentiviral stocks generated using the traditional Lipofectamine<sup>®</sup> 2000 transfection procedure (*i.e.* plating cells first followed by transfection with DNA:Lipofectamine<sup>®</sup> 2000 complexes). You may use the traditional Lipofectamine<sup>®</sup> 2000 transfection procedure, if desired, but keep in mind that the viral titer obtained may be lower (see **Alternative Transfection Procedure**, page 37).



Transfection Procedure	Follow the procedure below to cotransfect 293FT cells. Include a negative control (no DNA, no Lipofectamine <sup>®</sup> 2000) in your experiment to help evaluate results. You will need $6 \times 10^6$ 293FT cells for each sample.					
	1. <b>For each transfection sample,</b> prepare DNA-Lipofectamine <sup>®</sup> 2000 confollows:					
		a. In a sterile 5 mL tube, combine 9 μg of the ViraPower <sup>™</sup> Packaging Mix and 3 μg of pLenti6/V5-GW/miR expression plasmid DNA (12 μg total) in 1.5 mL of Opti-MEM <sup>®</sup> I Medium without serum. Mix gently.				
		b. In a separate sterile 5 mL tube, mix Lipofectamine <sup>®</sup> 2000 gently before use, then dilute 36 µL in 1.5 mL of Opti-MEM <sup>®</sup> I Medium without serum. Mix gently and incubate for 5 minutes at room temperature.				
		c. After the 5 minute incubation, combine the diluted DNA with the diluted Lipofectamine <sup>®</sup> 2000. Mix gently.				
		d. Incubate for 20 minutes at room temperature to allow the DNA- Lipofectamine <sup>®</sup> 2000 complexes to form. The solution may appear cloudy, but this will not impede the transfection.				
	2.	While DNA-lipid complexes are forming, trypsinize and count the 293FT cells. Resuspend the cells at a density of $1.2 \times 10^6$ cells/mL in growth medium containing serum (or Opti-MEM <sup>®</sup> I Medium containing serum).				
	3.	Add the DNA-Lipofectamine <sup>®</sup> 2000 complexes to a 10 cm tissue culture plate containing 5 mL of growth medium containing serum (or Opti-MEM <sup>®</sup> I Medium containing serum). <b>Do not add antibiotics to the medium.</b>				
	4.	Add 5 mL of the 293FT cell suspension ( $6 \times 10^6$ total cells) to the plate containing media and DNA-Lipofectamine <sup>®</sup> 2000 complexes and mix gently by rocking the plate back and forth. Incubate the cells overnight at 37°C in a CO <sub>2</sub> incubator.				
	5.	The next day, remove the media containing the DNA-Lipofectamine <sup>®</sup> 2000 complexes and replace with complete culture medium containing sodium pyruvate (see page 34).				
		<b>Note:</b> Expression of the VSV G glycoprotein causes 293FT cells to fuse, resulting in the appearance of multinucleated syncitia. This morphological change is normal and does not affect lentivirus production.				
	6.	Harvest virus-containing supernatants 48–72 hours posttransfection by removing medium to a 15 mL sterile, capped, conical tube.				
		<b>Note:</b> Minimal differences in viral yield are observed whether supernatants are collected 48 or 72 hours posttransfection.				
		<b>Caution:</b> Remember that you are working with infectious virus at this stage. Follow the recommended guidelines for working with BL-2 organisms (see pages 17 and 40 for more information).				
	7.	Centrifuge at 3000 rpm for 5 minutes at 4°C to pellet cell debris. Perform filtration step, if desired (see <b>Note</b> on the next page).				
	8.	Pipet viral supernatants into cryovials in 1 mL aliquots. Store viral stocks at –80°C. Proceed to <b>Titering Your Lentiviral Stock</b> , page 38.				
		Continued on next page				

Alternative Transfection Procedure	An alternative transfection procedure is provided below to cotransfect 293FT cells. Note that use of this procedure generally results in production of lentiviral stocks with a slightly lower titer that those produced when using the recommended <b>Transfection Procedure</b> , previous page.			
	1. The day before transfection, plate 293FT cells in a 10 cm tissue culture plate such that they will be 90–95% confluent on the day of transfection ( <i>i.e.</i> $6 \times 10^6$ cells in 10 mL of growth medium containing serum).			
	2. On the day of transfection, remove the culture medium from the 293FT cells and replace with 5 mL of growth medium containing serum (or Opti-MEM <sup>®</sup> I Medium containing serum). <b>Do not include antibiotics in the medium.</b>			
	3. Prepare DNA-Lipofectamine <sup>®</sup> 2000 complexes as instructed in the recommended <b>Transfection Procedure</b> , Step 1, previous page.			
	<ol> <li>Add the DNA-Lipofectamine<sup>®</sup> 2000 complexes dropwise to each plate of cells. Mix gently by rocking the plate back and forth. Incubate the cells overnight at 37°C in a CO₂ incubator.</li> </ol>			
	5. Follow Steps 5–8 as instructed in the recommended <b>Transfection Procedure</b> , previous page.			
Note	If you plan to use your lentiviral construct for <i>in vivo</i> applications, filter your viral supernatant through a sterile, 0.45 $\mu$ m low protein binding filter after the low-speed centrifugation step (see Step 7, previous page) to remove any remaining cellular debris. We recommend using Millex <sup>®</sup> -HV 0.45 $\mu$ m PVDF filters (Millipore, Cat. no. SLHV033RB) for filtration.			
	If you wish to concentrate your viral stock to obtain a higher titer, perform the filtration step first before concentrating your viral stock.			
Long-Term Storage	Place lentiviral stocks at –80°C for long-term storage. Repeated freezing and thawing is not recommended as it may result in loss of viral titer. When stored properly, viral stocks of an appropriate titer should be suitable for use for up to one year. After long-term storage, re-titer your viral stocks before transducing the mammalian cell line of interest.			
Scaling Up Virus Production	It is possible to scale up the cotransfection experiment to produce a larger volume of lentivirus, if desired. For example, we have scaled up the cotransfection experiment from a 10 cm plate to a T-175 cm <sup>2</sup> flask and harvested up to 30 mL of viral supernatant. To scale up your cotransfection, increase the number of cells plated and the amounts of DNA, Lipofectamine <sup>®</sup> 2000, and medium used in proportion to the difference in surface area of the culture vessel.			

## **Titering Your Lentiviral Stock**

Introduction	Before transducing a mammalian cell line of interest and expressing the miRNA for RNAi analysis, we highly recommend determining the titer of your lentiviral stock. While this procedure is not required for some applications, it is necessary if:			
	• You wish to control the number of integrated copies of the lentivirus			
	• You wish to generate reproducible gene knockdown results			
Titering Methods	You can determine the titer of your lentiviral stock using any of the following methods:			
	• Blasticidin selection (usually takes 2 weeks to determine the titer)			
	• EmGFP detection (usually takes 4 days post-transduction to determine the titer), if the miRNA sequence was cloned into pcDNA <sup>™</sup> 6.2-GW/EmGFP-miR vector			
Experimental	To determine the titer of a lentiviral stock, you will:			
Outline	1. Prepare 10-fold serial dilutions of your lentiviral stock			
	2. Transduce the different dilutions of lentivirus into a mammalian cell line of choice in the presence of Polybrene <sup>®</sup>			
	3. Select for stably transduced cells using the appropriate selection agent			
	4. Stain and count the number of antibiotic-resistant colonies in each dilution			
Factors Affecting	A number of factors can influence lentiviral titers including:			
Viral Titer	• The characteristics of the cell line used for titering (see the next page).			
	• The age of your lentiviral stock. Viral titers may decrease with long-term storage at -80°C. If your lentiviral stock has been stored for longer than 6 months, titer or re-titer your lentiviral stock prior to use in an RNAi experiment.			
	• Number of freeze/thaw cycles. Viral titers can decrease as much as 10% with each freeze/thaw cycle.			
	• Improper storage of the lentiviral stock. Lentiviral stocks should be aliquotted and stored at -80°C (see page 37 for recommended storage conditions).			

Continued on next page

Polybrene® is a registered trademark of Abbott Laboratories

Selecting a Cell Line	<ul> <li>You may titer your lentiviral stock using any mammalian cell line of choice. Generally, we recommend using the same mammalian cell line to titer your lentiviral stock as you will use to perform your expression studies. However, in some instances, you may wish to use a different cell line to titer your lentivirus (<i>e.g.</i>, if you are performing RNAi studies in a non-dividing cell line or a primary cell line). In these cases, we recommend that you choose a cell line with the following characteristics to titer your lentivirus:</li> <li>Grows as an adherent cell line</li> <li>Easy to handle</li> <li>Exhibits a doubling time in the range of 18–25 hours</li> <li>Non-migratory</li> <li>We generally use the HT1080 human fibrosarcoma cell line (ATCC, Cat. no. CCL-121) for titering purposes.</li> <li>Important: You may use other cell lines, including HeLa and NIH/3T3, to titer your lentivirus. However, note that the titer obtained when using HeLa cells or NIH/3T3 cells is approximately 10-fold lower than the titer obtained when using HT1080 cells.</li> </ul>
Note	The titer of a lentiviral construct may vary depending on which cell line is chosen (see <b>Selecting a Cell Line</b> , above). If you have more than one lentiviral construct, we recommend that you titer all of the lentiviral constructs using the same mammalian cell line.
Blasticidin Selection	The pLenti6/V5-GW/miR expression construct contains the Blasticidin resistance gene ( <i>bsd</i> ) (Kimura <i>et al.</i> , 1994) to allow for Blasticidin selection (Takeuchi <i>et al.</i> , 1958; Yamaguchi <i>et al.</i> , 1965) of mammalian cells that have stably transduced the lentiviral construct. If you are using the BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Kits, Blasticidin is supplied with the kit. Blasticidin is also available separately or as part of the ViraPower <sup>™</sup> Bsd Lentiviral Support Kit (see page 77 for ordering information). For more information about how to prepare and handle Blasticidin, and determine the Blasticidin sensitivity, refer to the <b>Appendix</b> , page 63.
Using Polybrene <sup>®</sup> During Transduction	Transduction of lentivirus into mammalian cells may be enhanced if cells are transduced in the presence of hexadimethrine bromide (Polybrene <sup>®</sup> ). For best results, we recommend performing transduction in the presence of Polybrene <sup>®</sup> . Note however, that some cells are sensitive to Polybrene <sup>®</sup> ( <i>e.g.</i> , primary neurons). Before performing any transduction experiments, you may want to test your cell line for sensitivity to Polybrene <sup>®</sup> . If your cells are sensitive to Polybrene <sup>®</sup> ( <i>e.g.</i> , exhibit toxicity or phenotypic changes), do not add Polybrene <sup>®</sup> during transduction. In this case, cells should still be successfully transduced.

Preparing and Storing	Follow the instructions below to prepare Polybrene <sup>®</sup> (Sigma-Aldrich, Cat. no. H9268):			
Polybrene <sup>®</sup>	1. Prepare a 6 mg/mL stock solution in deionized, sterile water.			
	2. Filter-sterilize and dispense 1 mL aliquots into sterile microcentrifuge tubes.			
	3. Store at -20°C for long-term storage. Stock solutions may be stored at -20°C for up to 1 year. Do not freeze/thaw the stock solution more than 3 times as this may result in loss of activity.			
	<b>Note:</b> The working stock may be stored at 4°C for up to 2 weeks.			
Materials Needed	• Your Lenti6/V5-DEST lentiviral stock (store at -80°C until use)			
	Adherent mammalian cell line of choice			
	Complete culture medium for your cell line			
	• 6 mg/mL Polybrene <sup>®</sup> , if desired			
	6-well tissue culture plates			
	<ul> <li>Crystal violet (Sigma-Aldrich<sup>®</sup>, Cat. no. C3886; prepare a 1% crystal violet solution in 10% ethanol), if you are using Blasticidin selection for titering</li> </ul>			
	• Inverted fluorescence microscope and appropriate filters for EmGFP visualization (see page 15 for filters), if you are using EmGFP titering method			
	Phosphate-Buffered Saline (PBS; page 77)			
	Components supplied with the kits			
	• Blasticidin (10 mg/mL stock), if you are using Blasticidin selection for titering			
CAUTION	Remember that you will be working with media containing infectious virus. Follow the recommended Federal and institutional guidelines for working with BL-2 organisms.			
	• Perform all manipulations within a certified biosafety cabinet.			
	Treat media containing virus with bleach.			
	• Treat used pipets pipette tipe, and other tissue culture supplies with block			

- Treat used pipets, pipette tips, and other tissue culture supplies with bleach and dispose of as biohazardous waste.
- Wear gloves, a laboratory coat, and safety glasses or goggles when handling viral stocks and media containing virus.

### Transduction and Titering Procedure

Follow the procedure below to determine the titer of your lentiviral stock using the mammalian cell line of your choice.

**Note:** You will use **at least** one 6-well plate for every lentiviral stock to be titered (one mock well plus five dilutions).

- If you have generated a lentiviral stock of the pLenti6-V5-GW/miR-*lacZ* control construct with or without EmGFP, perform titering using the Blasticidin or EmGFP method.
- If you generated a lentiviral stock of the pLenti6-V5-GW/*lacZ* control construct, use the Blasticidin titering method.
- 1. The day before transduction (Day 1), trypsinize and count the cells, plating them in a 6-well plate such that they will be 30–50% confluent at the time of transduction. Incubate cells at 37°C overnight.

**Example:** When using HT1080 cells, we plate  $2 \times 10^5$  cells per well in a 6-well plate.

2. On the day of transduction (Day 2), thaw your lentiviral stock and prepare 10-fold serial dilutions ranging from 10<sup>-2</sup> to 10<sup>-6</sup>. For each dilution, dilute the lentiviral construct into complete culture medium to a final volume of 1 mL. **DO NOT** vortex.

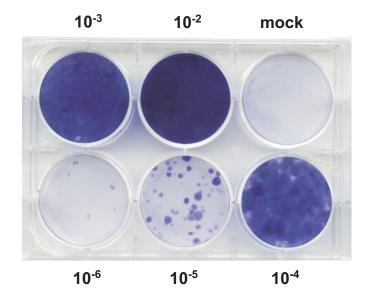
**Note:** You may prepare a wider range of serial dilutions (10<sup>-2</sup> to 10<sup>-8</sup>), if desired.

- 3. Remove the culture medium from the cells. Mix each dilution gently by inversion and add to one well of cells (total volume = 1 mL).
- 4. Add Polybrene<sup>®</sup> (if desired) to each well to a final concentration of 6 μg/mL. Swirl the plate gently to mix. Incubate at 37°C overnight.
- 5. The following day (Day 3), remove the media containing virus and replace with 2 mL of complete culture medium.
- 6. The following day (Day 4), proceed to Steps 7–8 for the **EmGFP titering method** or proceed to Steps 9–14 for the **Blasticidin titering method**.
- Determine the titer by flow cytometry on Day 4 for titering EmGFP. For each viral dilution well of the 6 well plate, trypsinize and resuspend the cells in complete media at a concentration of 10–500 cells/µL.
- 8. Using a flow cytometry system, determine the percentage of GFP-positive cells for each dilution (see next page). Determine the titer using the formula described on the next page.
- 9. For Blasticidin selection, remove the medium on Day 4 and replace with complete culture medium containing the appropriate amount of Blasticidin to select for stably transduced cells.
- 10. Replace medium with fresh medium containing Blasticidin every 3-4 days.
- 11. After 10–12 days of selection (day 14–16), you should see no live cells in the mock well and discrete Blasticidin-resistant colonies in one or more of the dilution wells. Remove the medium and wash the cells twice with PBS.
- 12. Add crystal violet solution (1 mL for 6-well dish; 5 mL for 10 cm plate) and incubate for 10 minutes at room temperature.
- 13. Remove the crystal violet stain and wash the cells with PBS. Repeat wash.
- 14. Count the blue-stained colonies and determine your lentiviral stock titer.

Preparing Cells for Flow Cytometry	If you have used the EmGFP titering method, prepare cells for flow cytometry according to the established protocols in use at your flow cytometry facility. Refer to page 15 for spectral characteristics of EmGFP. The steps below provide general guidelines, but other methods may be suitable.				
		4 post-transduction, dissoc ssociation buffer.	iate the cells from the plate b	y using trypsin or	
	the cell 1% FBS	l pellet in flow cytometry bu	ve residual media componer offer such as calcium/magne analysis on your flow cytom out may be done, if desired.	sium-free PBS with	
	in calciv of cells;	um/magnesium-free PBS. How	pmetry, use 2% formaldehyde or rever, these fixatives may increa ed, mock-transduced cells as a r	se autofluorescence	
	negativ				
Calculating Lentiviral Titer	Calculate the EmGFP lentivirus titers from the dilutions at which the percentage of EmGFP-positive cells fall within the range of 1–30% (Sastry <i>et al.</i> , 2002; White <i>et al.</i> , 1999) to avoid analyzing dilution samples containing multiple integrated lentiviral genomes, which may result in underestimating the viral titer, or dilution samples containing too few transduced cells, which will give inaccurate results. The titer is expressed as transducing units (TU)/mL.				
	Use the following formula to calculate the titer: $[F \times C/V] \times D$				
	C = total n V = volum	<ul> <li>F = frequency of GFP-positive cells (percentage obtained divided by 100)</li> <li>C = total number of cells in the well at the time of transduction</li> <li>V = volume of inoculum in mL</li> <li>D = lentivirus dilution</li> </ul>			
	An example for calculating the lentiviral titer is provided below. An EmGFP lentiviral stock was generated using the protocol on the previous page. The following data were generated after performing flow cytometry analysis:				
		Lentivirus Dilution	% EmGFP Positive Cells		
		10-2	91.5%		
		10-3	34.6%		
		10-4	4.4%		
	of EmGFP- positive ce	positive cells falls into the c lls is 4.4/100 = 0.044, multip 1 (the volume of inoculum)	is used to calculate the titer s lesired range of 1–30%. The f lied by 2 × 10 <sup>5</sup> (the number c . Thus the calculation is as fo × 200,000)/1] × 10 <sup>4</sup>	requency of EmGFP- of cells in the well)	

The lentiviral titer for this example is  $8.8 \times 10^7 \text{ TU/mL}$ .

Expected Results When titering pLenti/V5 lentiviral stocks using HT1080 cells, we generally obtain titers ranging from 5 × 10<sup>5</sup> to 2 × 10<sup>7</sup> transducing units (TU)/mL. For an example of expected results obtained from a typical titering experiment using Blasticidin, see below.
 Note: If the titer of your lentiviral stock is less than 1 × 10<sup>5</sup> TU/mL, we recommend producing a new lentiviral stock. See page 31 and the Troubleshooting section, page 52 for more tips and guidelines to optimize the viral yield.
 Example of Expected Results In this experiment, a pLenti6 lentiviral stock was generated using the protocol on page 36. HT1080 cells were transduced with 10-fold serial dilutions of the lentiviral supernatant (10<sup>-2</sup> to 10<sup>-6</sup> dilutions) or untransduced (mock) following the protocol on page 41. Forty-eight hours post-transduction, the cells were placed under Blasticidin selection (10 µg/mL). After 10 days of selection, the cells were



stained with crystal violet (see plate below), and colonies were counted.

In the plate above, the colony counts were:

- Mock: no colonies
- 10<sup>-2</sup> dilution: confluent; undeterminable
- 10<sup>-3</sup> dilution: confluent; undeterminable
- 10<sup>-4</sup> dilution: confluent; undeterminable
- 10<sup>-5</sup> dilution: 46
- 10<sup>-6</sup> dilution: 5

Thus, the titer of this lentiviral stock is  $4.8 \times 10^6$  TU/mL (*i.e.* average of  $46 \times 10^5$  and  $5 \times 10^6$ ).

# **Transduction and Analysis**

Introduction	Once you have generated lentiviral stocks with suitable titers, you are ready to transduce the lentiviral constructs into a mammalian cell line of choice to inducibly express the miRNA of interest and perform RNAi analysis.			
	<b>Reminder:</b> Remember that each lentiviral construct contains a deletion in the 3' LTR that leads to self-inactivation of the lentivirus after transduction into mammalian cells. Once integrated into the genome, the lentivirus can no longer produce packageable virus.			
Experimental	To perform transduction, you will:			
Outline	1. Determine the Multiplicity of Infection (MOI) and antibiotic sensitivity for your cell line.			
	2. Grow the mammalian cell line of choice.			
	3. Transduce the mammalian cell line of choice with your lentiviral construct in the presence of Polybrene <sup>®</sup> .			
	4. Harvest cells after 48–96 hours to perform transient knockdown experiments or select for stably transduced cells using Blasticidin.			
	5. Expand at least 5 Blasticidin-resistant colonies and analyze each clone to assay for knockdown of the target gene.			
Factors Affecting Gene Knockdown	A number of factors can influence the degree to which expression of your gene of interest is reduced ( <i>i.e.</i> gene knockdown) in an RNAi experiment including:			
Levels	Transduction efficiency			
	MOI used to transduce cells			
	Transcription rate of the target gene of interest			
	Stability of the target protein			
	Growth characteristics of your mammalian cell line			
	Activity of your miRNA in transient transfections			
	Take these factors into account when designing your transduction and RNAi experiments.			
Transient vs. Stable Expression	After transducing your lentiviral construct into the mammalian cell line of choice, you may assay for target gene knockdown in the following ways:			
	• Pool a heterogeneous population of cells and test for gene knockdown directly after transduction ( <i>i.e.</i> "transient" RNAi analysis). Note that you must wait for a minimum of 48–72 hours after transduction before harvesting your cells to allow expressed miRNA molecules to accumulate in transduced cells.			
	• Select for stably transduced cells using Blasticidin. This requires a minimum of 10–12 days after transduction, but allows generation of clonal cell lines that stably express the miRNA sequence.			

Determining Antibiotic Sensitivity for Your Cell Line	Before selecting for stably transduced cells, you must first determine the minimum concentration of Blasticidin required to kill your untransduced mammalian cell line ( <i>i.e.</i> perform a kill curve experiment). For guidelines to perform a kill curve experiment, see page 64. If you titered your lentiviral construct in the same mammalian cell line that you are using to generate a stable cell line, then you may use the same concentration of Blasticidin for selection that you used for titering.	
Multiplicity of Infection (MOI)	To obtain optimal expression of your miRNA and therefore, the highest degree of target gene knockdown, you will need to transduce the lentiviral construct into your mammalian cell line of choice using a suitable MOI. MOI is defined as the number of virus particles per cell and generally correlates with the number of integration events and as a result, expression. Typically, miRNA expression levels increase as the MOI increases.	
Determining the Optimal MOI	<ul> <li>A number of factors can influence determination of an optimal MOI including:</li> <li>The nature of your mammalian cell line (<i>e.g.</i>, non-dividing vs. dividing cell type; see Note, below)</li> <li>The transduction efficiency of your mammalian cell line</li> <li>The nature of your target gene of interest</li> <li>If you are transducing the lentiviral construct into a mammalian cell line of choice for the first time, use a range of MOIs (<i>e.g.</i>, 0, 1, 5, 10, 50) to determine the MOI required to obtain the optimal level of target gene knockdown.</li> </ul>	
Note	In general, lentiviral constructs transduce non-dividing cell types less efficiently than actively dividing cell lines. If you are transducing your lentiviral construct into a non-dividing cell type, you may need to increase the MOI to achieve optimal target gene knockdown.	
Preparing Mammalian Cells	Initiate your mammalian cell line of choice that will be used for titering. Grow the cells in the appropriate medium. Cells should be >95% viable.	

Positive Control	If you have generated two positive control lentiviral constructs (pLenti6/V5-GW/ <u>+</u> EmGFP-miR- <i>lacZ</i> control and pLenti6/V5-GW/ <i>lacZ</i> control constructs) as described on page 33, you may use the controls in cotransduction experiments to verify the lentiviral induced RNAi response in mammalian cells.
	For cotransductions, use a 3:1 MOI ratio of pLenti6/V5-GW/miR- <i>lacZ</i> to pLenti6/V5-GW/ <i>lacZ</i> expression clone. For expected results, see page 50.
	The $\beta$ -galactosidase protein expressed from the pLenti6/V5-GW/ <i>lacZ</i> control lentiviral construct is approximately 121 kDa in size. You may assay for $\beta$ -galactosidase expression by activity assay FluoReporter <sup>®</sup> <i>lacZ</i> /Galactosidase Quantitation Kit or by staining the cells for activity using the $\beta$ -Gal Staining Kit for fast and easy detection of $\beta$ -galactosidase expression, see page 77 for ordering information.
<b>Q</b> Important	Remember that viral supernatants are generated by harvesting spent media containing virus from the 293FT producer cells. Spent media lacks nutrients and may contain some toxic waste products. If you are using a large volume of viral supernatant to transduce your mammalian cell line ( <i>e.g.</i> , 1 mL of viral supernatant per well in a 6-well plate), note that growth characteristics or morphology of the cells may be affected during transduction. These effects are generally alleviated after transduction when the media is replaced with fresh, complete media.
Concentrating Virus	It is possible to concentrate VSV-G pseudotyped lentiviruses using a variety of methods without significantly affecting the transducibility of the lentivirus. If the titer of your lentiviral stock is relatively low (less than $5 \times 10^5$ TU/mL) and your experiment requires that you use a large volume of viral supernatant ( <i>e.g.</i> , a relatively high MOI), you may wish to concentrate your virus before proceeding to transduction. For details and guidelines to concentrate your virus, refer to published reference sources (Yee, 1999).
Materials Needed	• Titered lentiviral stock (store at -80°C until use)
	Mammalian cell line of choice
	Complete culture medium for your cell line
	• 6 mg/mL Polybrene <sup>®</sup> , if desired
	Appropriately sized tissue culture plates for your application
	Components supplied with the kits
	• 10 mg/mL Blasticidin stock (if selecting for stably transduced cells)
	Continued on next page

Transduction Procedure	Follow the procedure below to transduce your mammalian cell line of choice with your lentiviral construct.
	<ol> <li>Plate cells in complete media as appropriate for your application. When determining the density at which to plate cells, remember to take into account the length of time cells will be cultured prior to performing RNAi analysis (<i>e.g.</i>, 48 hours vs. 120 hours).</li> </ol>
	2. On the day of transduction (Day 1), thaw the lentiviral stock and dilute (if necessary) the appropriate amount of virus (at a suitable MOI) into fresh complete medium. Keep the total volume of medium containing virus as low as possible to maximize transduction efficiency. <b>DO NOT</b> vortex.
	<ol> <li>Remove the culture medium from the cells. Mix the medium containing virus gently by pipetting and add to the cells.</li> </ol>
	I. Add Polybrene <sup>®</sup> (if desired) to a final concentration of 6 μg/mL. Swirl the plate gently to mix. Incubate at 37°C overnight.
	<b>Note:</b> If you are transducing cells with undiluted viral stock and are concerned about possible toxicity or growth effects caused by overnight incubation, it is possible to incubate cells for as little as 6 hours prior to changing medium.
	5. The following day (Day 2), remove the medium containing virus and replace with fresh, complete culture medium.
	5. The following day (Day 3), perform one of the following:
	• Harvest the cells and assay for inhibition of your target gene if you are performing transient expression experiments. If you wish to assay the cells at a later time, you may continue to culture the cells or replate them into larger-sized tissue culture formats as necessary.
	• Remove the medium and replace with fresh, complete medium containing the appropriate amount of Blasticidin to select for stably transduced cells. Proceed to Step 7.
	<ol> <li>Replace medium with fresh medium containing Blasticidin every 3–4 days until Blasticidin-resistant colonies can be identified (generally 10–12 days after selection).</li> </ol>
	<ol> <li>Pick at least 5 Blasticidin-resistant colonies (see Note, below) and expand each clone to assay for knockdown of the target gene.</li> </ol>
Note	Integration of the lentivirus into the genome is random. Depending upon the influence of the surrounding genomic sequences at the integration site, you may see varying levels of target gene knockdown from different Blasticidin-resistant clones. Test at least 5 Blasticidin-resistant clones and select the clone that provides the optimal degree of gene knockdown for further studies.

Performing RNAi Analysis	Use any method as appropriate to assay for knockdown of your target gene including functional analysis, immunofluorescence, western blot, or real-time quantitative RT-PCR (qRT-PCR) with the appropriate fluorogenic LUX <sup>™</sup> primers. For more information about LUX <sup>™</sup> primers or to design LUX <sup>™</sup> primers, see <u>www.invitrogen.com/lux</u> .
	If you have used the BLOCK-iT <sup>™</sup> Lentiviral Pol II miR RNAi Expression System with EmGFP, you can detect EmGFP fluorescence using fluorescence microscopy or flow cytometry (see the next page).
Expected Results	When performing RNAi studies using pLenti6/V5 lentiviral constructs, we generally observe significant inhibition of gene expression within 48 to 120 hours after transduction. The degree of gene knockdown depends on the time of assay, stability of the protein of interest, and on the other factors listed on page 44. Note that 100% gene knockdown is generally not observed, but > 80% is possible with optimized conditions
	For an example of results obtained from RNAi experiments using the pLenti6-V5- DEST, see page 50.

## **Detecting Fluorescence**

Introduction	You can perform analysis of the EmGFP fluorescent protein from the expression clone in either transiently transfected cells or stable cell lines, if you used pcDNA <sup>™</sup> 6.2- GW/EmGFP-miR for cloning your miRNA sequence. Once you have transfected your expression clone into mammalian cells, you may detect EmGFP protein expression directly in cells by fluorescence microscopy or other methods that use light excitation and detection of emission. The EmGFP expression is essentially 100% correlated with the expression of your miRNA. See below for recommended fluorescence microscopy filter sets.		
Filters for Use with EmGFP	EmGFP can be detected with standard FITC filter sets. However, for optimal detection of the fluorescence signal, use a filter set which is optimized for detection within the excitation and emission ranges for the fluorescent protein, such as the Omega XF100-2 or XF100-3 filter set for fluorescence microscopy.		
	The spectral characteristics of EmGFP are listed in the table below:		
	Excitation (nm) Emission (nm)		
	487 509		
	For information on obtaining these filter sets, contact Omega Optical, Inc. ( <u>www.omegafilters.com</u> ) or Chroma Technology Corporation ( <u>www.chroma.com</u> ).		
Fluorescence Microscope	You may view the fluorescence signal of EmGFP in cells using an inverted fluorescence microscope with an FITC filter or Omega XF100 filter (available from <u>www.omegafilters.com</u> ) for viewing cells in culture or a flow cytometry system.		
Color Camera	If desired, you may use a color camera that is compatible with the microscope to photograph the cells.		
Detecting Transfected Cells	After transduction, allow the cells to recover for 24 to 48 hours before assaying for fluorescence. Medium can be removed and replaced with PBS during viewing to avoid any fluorescence due to the medium. Be sure to replace PBS with fresh medium if you wish to continue growing the cells. <b>Note:</b> Cells can be incubated further to optimize expression of EmGFP.		
Expected Results	Cells expressing EmGFP will appear brightly labeled and will emit a green fluorescence signal that should be easy to detect above the background fluorescence. <b>Note:</b> The fluorescence signal of EmGFP from miRNA-containing vectors is reduced when compared to non-miRNA containing vectors. Cells with bright fluorescence will demonstrate highest knockdown with a functional miRNA. However, cells with reduced fluorescence may still express the miRNA and demonstrate knockdown since the expression levels required to observe gene knockdown are generally lower than that required to detect EmGFP expression.		

### **Examples of Expected Results**

### Example of Expected Results

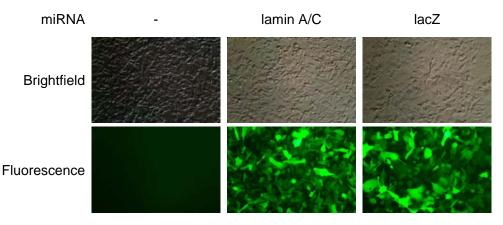
In this experiment, a double-stranded miR-*lacZ* control oligo (directed towards exogenous LacZ) and a double-stranded miR-lamin oligo (directed towards endogenous lamin) were cloned into pcDNA6.2<sup>™</sup>-GW/EmGFP-miR expression vector using the BLOCK-iT<sup>™</sup> Pol II miR RNAi Expression Vector Kit. The miR-*lacZ* and miR-lamin cassettes were transferred into the pLenti6/V5-DEST vector using the Rapid BP/LR recombination reaction to generate the pLenti6-GW/EmGFP-miR-*lacZ* and pLenti6-GW/EmGFP-miR-lamin expression constructs, respectively. Lentiviral stocks were generated and titered in HT1080 cells following the protocols in this manual.

HT1080 cells plated in a 12-well plate were co-transduced with the following lentiviral particles:

- Expressing the *lacZ*-directed miRNA with EmGFP at an MOI of 30 and pLenti6/V5-GW/*lacZ* at an MOI of 10
- Expressing the lamin A/C-directed miRNA with EmGFP at an MOI of 30 and pLenti6/V5-GW/*lacZ* at an MOI of 10.

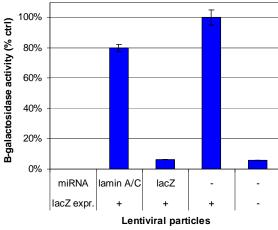
Cells were harvested 48 hours post-transduction and subjected to fluorescence microscopy using the appropriate filters as described in this manual. Fluorescence microscopy results are shown in figure A below and indicate the expression of EmGFP in virtually all cells transduced with the miRNA lentiviruses.

#### **Figure A**



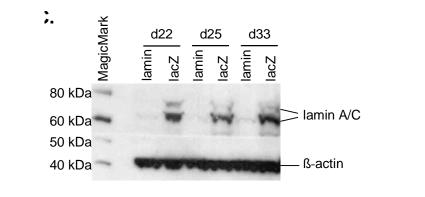
Cell lysates were then prepared from duplicate wells 48 hours (*i.e.* 2 days) after transduction.  $\beta$ -galactosidase activity was determined from equivalent amounts of cell lysate using the FluoReporter<sup>®</sup> *lacZ*/Galactosidase Quantitation Kit (see page 77). Results are shown in figure B, next page.

### Examples of Expected Results, Continued



HeLa cells were transduced with the *lacZ* and lamin miRNA lentiviruses and stable cell lines were generated using Blasticidin selection as described in this manual. Cell lysates were prepared from stably transduced cells after 22–33 days and subjected to western blot analysis. The western blot was cut in half and one half was probed with Anti-Lamin A/C Antibody (1:1,000 dilution, BD Biosciences, Cat. no. 612162) and the other half was probed with Anti- $\beta$ -Actin Antibody (1:5,000 dilution, Abcam, Cat. no. ab6276). The blot was developed using WesternBreeze<sup>®</sup> Chemiluminescent Kit, see page 77. Western blot results are shown in figure C, below. The results show substantial decrease in the lamin A/C protein isoforms (molecular weight of lamin A is 70 kDa and lamin C is 65 kDa) even 33 days post-transduction indicating the reliable and stable transduction of lamin-directed miRNA in lentiviral system.





## Troubleshooting

Introduction	Review the information in this section to troubleshoot your lentiviral expression experiments. To troubleshoot oligo design and cloning experiments, refer to the BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kit manual.

### Rapid BP/LR Reaction and Transformation

The table below lists some potential problems and possible solutions that may help you troubleshoot the Rapid BP/LR recombination and transformation procedures.

Problem	Reason	Solution
Few or no colonies obtained from sample reaction <b>and</b> the transformation control gave colonies	Incorrect antibiotic used to select for transformants	Select for transformants on LB agar plates containing 100 µg/mL ampicillin.
	Rapid BP/LR reaction may not work for your insert	Use the standard BP and LR recombination reactions as described on page 57.
	BP recombination reaction is treated with Proteinase K	<b>Do not</b> treat reaction the BP reaction with Proteinase K before the LR reaction.
	Did not use the suggested Gateway® Clonase® II enzyme mixes or Gateway® Clonase® II	<ul> <li>Make sure to store the Gateway<sup>®</sup> BP and LR Clonase<sup>®</sup> II enzyme mix at -20°C or -80°C.</li> </ul>
	enzyme mixes were inactive	• Do not freeze/thaw the Gateway <sup>®</sup> BP and LR Clonase <sup>®</sup> II enzyme mix more than 10 times.
		• Use the recommended amount of Gateway <sup>®</sup> BP and LR Clonase <sup>®</sup> II enzyme mix (see page 27).
		Test another aliquot of the Gateway <sup>®</sup> Clonase <sup>®</sup> II enzyme mix.
	Not enough LR reaction transformed	Transform 2–3 μL of the LR reaction into One Shot <sup>®</sup> Stbl3 <sup>™</sup> Chemically Competent <i>E. coli</i> .
	Not enough transformation mixture plated	Increase the amount of <i>E. coli</i> plated.
	Did not perform the 1 hour grow-out period before plating the transformation mixture	After the heat-shock step, add S.O.C. Medium and incubate the transformation mixture for 1 hour at 37°C with shaking before plating.
	Too much BP reaction used in the LR reaction	Use 3 $\mu$ L BP reaction for the LR reaction.

## Troubleshooting, Continued

Problem	Reason	Solution
Different sized colonies ( <i>i.e.</i> large and small) appear when using TOP10 <i>E. coli</i> for transformation	Some transformants contain plasmids in which unwanted recombination has occurred between 5' and 3' LTRs	Always use the One Shot <sup>®</sup> Stbl3 <sup>™</sup> Chemically Competent <i>E. coli</i> supplied with the kit for transformation. Stbl3 <sup>™</sup> <i>E. coli</i> are recommended for cloning unstable DNA including lentiviral DNA containing direct repeats and generally give rise to fewer unwanted recombinants.
Few or no colonies obtained from the transformation control	Competent cells stored incorrectly	<ul> <li>Store the One Shot<sup>®</sup> Stbl3<sup>™</sup> Chemically Competent <i>E. coli</i> at -80°C.</li> <li>Thaw a vial of One Shot<sup>®</sup> cells on ice immediately before use.</li> </ul>
	After addition of DNA, competent cells mixed by pipetting up and down	After adding DNA, mix competent cells gently. <b>Do not mix by pipetting up and down.</b>

### Rapid BP/LR Reaction and Transformation, Continued

### Generating the Lentiviral Stock

The table below lists some potential problems and possible solutions that may help you troubleshoot your co-transfection and titering experiments.

Problem	Reason	Solution
Low viral titer	Low transfection efficiency:	
	• Used poor quality expression construct plasmid DNA ( <i>i.e.</i> DNA from a mini-prep)	• <b>Do not</b> use plasmid DNA from a miniprep for transfection. Use PureLink <sup>™</sup> Plasmid Purification Kits or CsCl gradient centrifugation to prepare plasmid DNA.
	• Unhealthy 293FT cells; cells exhibit low viability	• Use healthy 293FT cells under passage 20; do not overgrow.
	<ul> <li>Cells transfected in media containing antibiotics (<i>i.e.</i> Geneticin<sup>®</sup>)</li> </ul>	• <b>Do not</b> add Geneticin <sup>®</sup> in the media during transfection as this reduces transfection efficiency and causes cell death.
	Plasmid DNA:transfection     reagent ratio incorrect	<ul> <li>Use a DNA (in μg):Lipofectamine<sup>®</sup> 2000 (in μL) ratio ranging from 1:2 to 1:3.</li> </ul>
	• 293FT cells plated too sparsely	• Plate cells such that they are 90–95% confluent at the time of transfection OR use the recommended transfection protocol ( <i>i.e.</i> add cells to media containing DNA:lipid complexes; page 36).

Problem	Reason	Solution
Low viral titer	Transfected cells not cultured in media containing sodium pyruvate	One day after transfection, remove media containing DNA:lipid complexes and replace with complete media containing sodium pyruvate. Sodium pyruvate provides an extra energy source for the cells.
	Lipofectamine <sup>®</sup> 2000 Reagent handled incorrectly	<ul> <li>Store at 4°C. Do not freeze.</li> <li>Mix gently by inversion before use. Do not vortex.</li> </ul>
	Viral supernatant harvested too early	Viral supernatants can generally be collected 48–72 hours post-transfection. If many cells are still attached to the plate and look healthy at this point, wait an additional 24 hours before harvesting the viral supernatant.
	Viral supernatant too dilute	Concentrate virus using any method of choice (Yee, 1999).
	Viral supernatant frozen and thawed multiple times	<b>Do not</b> freeze/thaw viral supernatant more than 3 times.
	Poor choice of titering cell line	Use HT1080 cells or another adherent cell line with the characteristics discussed on page 39.
	Target gene is essential for cell viability	Make sure that your target gene is not essential for cell viability or growth by performing a transient transfection with the entry construct containing the miRNA of interest.
	Polybrene <sup>®</sup> not included during titering procedure	Transduce the lentiviral construct into cells in the presence of Polybrene <sup>®</sup> .
No colonies obtained upon titering	Too much Blasticidin used for selection	Determine the Blasticidin sensitivity of your cell line by performing a kill curve experiment. Use the minimum Blasticidin concentration required to kill your untransduced cell line.
	Viral stocks stored incorrectly	Aliquot and store stocks at -80°C. Do not freeze/thaw more than 3 times.
	Polybrene <sup>®</sup> not included during transduction	Transduce the lentiviral construct into cells in the presence of Polybrene <sup>®</sup> .
Titer indeterminable; cells confluent	Too little Blasticidin used for selection	Increase amount of Blasticidin used for selection.
	Viral supernatant not diluted sufficiently	Titer lentivirus using a wider range of 10-fold serial dilutions ( <i>e.g.</i> , $10^{-2}$ to $10^{-8}$ ).

### Generating the Lentiviral Stock, Continued

## Troubleshooting, Continued

# Transduction and RNAi Analysis

The table below lists some potential problems and possible solutions that may help you troubleshoot transduction and knockdown experiment.

Problem	Reason	Solution
Low levels of gene knockdown observed	<ul> <li>Low transduction efficiency:</li> <li>Polybrene<sup>®</sup> not included during transduction</li> <li>Non-dividing cell type used</li> </ul>	<ul> <li>Transduce the lentiviral construct into cells in the presence of Polybrene<sup>®</sup>.</li> <li>Transduce your lentiviral construct into cells using a higher MOI.</li> </ul>
	MOI too low	Transduce your lentiviral construct into cells using a higher MOI.
	Cells harvested and assayed too soon after transduction	Do not harvest cells until at least 48–72 hours after transduction to allow expressed miRNA to accumulate in transduced cells. If low levels of knockdown are observed at 48 hours, culture cells for a longer period of time before assaying for gene knockdown or place cells under Blasticidin selection. <b>Note:</b> Placing cells under Blasticidin selection can improve gene knockdown results by killing untransduced cells.
	Target gene is important for cell viability	Make sure that your target gene is not essential for cell viability or growth.
	Viral stocks not titered	Titer the lentiviral stock using the procedure on page 41 before use.
	Viral stock stored incorrectly	<ul> <li>Aliquot and store stocks at -80°C.</li> <li>Do not freeze/thaw more than 3 times.</li> <li>If stored for longer than 6 months, re-titer stock before use.</li> </ul>
	miRNA with weak activity chosen	Select a different target region. If possible, screen miRNA first by transient transfection of the expression construct to verify its activity, then perform BP/LR recombination with the pLenti6/V5-DEST vector and proceed to generate lentivirus. <b>Note:</b> Generally, transient transfection greatly overexpresses miRNA, so moderately active expression clones may be less active when expressed from a lentiviral construct.

Problem	Reason	Solution
No gene knockdown observed	miRNA with no activity chosen	Select a different target region. If possible, screen miRNA first by transient transfection of the expression construct to verify activity, then perform BP/LR recombination with the pLenti6/V5-DEST vector and proceed to generate lentivirus.
	Viral stocks stored incorrectly	Aliquot and store stocks at -80°C. Do not freeze/thaw more than 3 times.
	MOI too low	Transduce your lentiviral construct into cells using a higher MOI.
Cytotoxic effects observed after transduction	Target gene is essential for cell viability	Make sure that your target gene is not essential for cell viability or growth.
	Large volume of viral supernatant used for transduction	• Remove the "spent" media containing virus and replace with fresh, complete media.
		• Concentrate the virus (Yee, 1999).
	Polybrene <sup>®</sup> used during transduction	Verify the sensitivity of your cells to Polybrene <sup>®</sup> . If cells are sensitive, omit the Polybrene <sup>®</sup> during transduction.
	Too much Blasticidin used for selection	Determine the Blasticidin sensitivity of your cell line by performing a kill curve. Use the minimum Blasticidin concentration required to kill your untransduced cell line.
Non-specific off-target gene knockdown observed	Target sequence contains strong homology to other genes	Select a different target region.
No fluorescence signal detected with expression clone containing EmGFP	Incorrect filters used to detect fluorescence	Use the recommended filter sets for detection of fluorescence (see page 49). Use an inverted fluorescence microscope for analysis. If desired, allow the protein expression to continue for an additional 1–3 days before assaying for fluorescence. <b>Note:</b> The expression levels required to observe gene knockdown are generally lower than that required to detect EmGFP expression. Knockdown may still occur in non-EmGFP positive cells.

## Appendix

## Performing the Standard BP Recombination Reaction

Introduction	General guidelines and instructions are provided below and in the next section to perform a standard BP recombination reaction using the linearized expression clone containing your pre-miRNA expression cassette ( <i>attB</i> substrate) and a donor vector, and to transform the reaction mixture into One Shot® TOP10 Chemically Competent <i>E. coli</i> host to select for entry clones (page 60). Use the standard BP reaction to obtain the maximum number of colonies or generate an entry clone for future use.	
Experimental	To generate an entry clone, you will:	
Outline	<ol> <li>Perform a BP recombination reaction using the linearized <i>att</i>B-containing expression clone and <i>att</i>P-containing pDONR<sup>™</sup>221 vector.</li> </ol>	
	2. Transform the reaction mixture into competent <i>E. coli</i> host.	
	3. Select for entry clones.	
Donor Vector and Expression Clone	See page 22 for details on the donor vector, propagating the donor vector, and for linearizing the expression clone.	
Positive Control	pEXP7-tet is provided as a positive control for the BP reaction. pEXP7-tet is an approximately 1.4 kb linear fragment and contains <i>att</i> B sites flanking the tetracycline resistance gene and its promoter (Tc <sup>r</sup> ). Using the pEXP7-tet fragment in a BP reaction with a donor vector results in entry clones that express the tetracycline resistance gene. The efficiency of the BP recombination reaction can easily be determined by streaking entry clones onto LB plates containing 20 $\mu$ g/mL tetracycline.	
<i>E. coli</i> Host Strain	You may use any <i>recA</i> , <i>endA E</i> . <i>coli</i> strain including TOP10, OmniMAX <sup>TM</sup> 2-T1 <sup>R</sup> or equivalent for transformation. <b>Do not</b> use <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.	

## Performing the Standard BP Recombination Reaction, Continued

Materials Needed	For BP Recombination			
	• Linearized <i>att</i> B expression clone, see page 22			
	• TE Buffer, pH 8.0 (10 mM Tris-HCl, pH 8.0, 1 mM EDTA), see page 77			
	Components supplied with the kits			
	• $pDONR^{TM}221$ vector (150 ng/ $\mu$ L)			
	<ul> <li>Gateway<sup>®</sup> BP Clonase<sup>®</sup> II enzyme mix (keep at -20°C until immediately before use)</li> </ul>			
	• 2 µg/µL Proteinase K solution (keep on ice until use)			
	• pEXP7-tet positive control (50 ng/ $\mu$ L)			
	For transformation			
	• One Shot <sup>®</sup> TOP10 Chemically Competent <i>E. coli</i> (or equivalent; one vial per transformation; thaw on ice before use), see page 77			
	• LB Medium			
	<ul> <li>LB plates containing 50 μg/mL kanamycin (two for each transformation; warm at 37°C for 30 minutes)</li> </ul>			
	• 42°C water bath			
	• 37°C shaking and non-shaking incubator			
	Components supplied with the kits			
	• S.O.C. Medium (warm to room temperature)			
	• Positive control ( <i>e.g.</i> , pUC19, use as a control for transformation if desired)			
	Continued on next page			

## Performing the Standard BP Recombination Reaction, Continued

#### Setting Up the BP Recombination Reaction

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

**Note:** To include a negative control, set up a second sample reaction and omit the Gateway<sup>®</sup> BP Clonase<sup>®</sup> II enzyme mix (see Step 4).

Components	Sample	Positive Control
Linearized <i>att</i> B expression clone from Step 4, page 22 (20–50 fmol)	1–7 µL	_
pDONR <sup>™</sup> 221 vector (150 ng/μL)	1 µL	1 µL
pEXP7-tet positive control (50 ng/µL)	-	2 µL
TE Buffer, pH 8.0	to 8 µL	5 µL

- 2. Remove the Gateway<sup>®</sup> BP Clonase<sup>®</sup> II enzyme mix from -20°C and thaw on ice (~ 2 minutes).
- 3. Vortex the Gateway<sup>®</sup> BP Clonase<sup>®</sup> II enzyme mix briefly twice (2 seconds each time).
- Add 2 µL of Gateway<sup>®</sup> BP Clonase<sup>®</sup> II enzyme mix to the sample and positive control vials. Do not add Gateway<sup>®</sup> BP Clonase<sup>®</sup> II to the negative control vial. Mix well by vortexing briefly twice (2 seconds each time).

**Reminder:** Return Gateway<sup>®</sup> BP Clonase<sup>®</sup> II enzyme mix to  $-20^{\circ}$ C immediately after use.

5. Incubate reactions at 25°C for 1 hour.

**Note:** For most applications, a 1 hour incubation will yield a sufficient number of entry clones. Depending on your needs, the length of the recombination reaction can be extended up to 18 hours. An overnight incubation typically yields 5–10 times more colonies than a 1 hour incubation.

- 6. Add 1 μL of the Proteinase K solution to each reaction. Incubate for 10 minutes at 37°C.
- 7. Proceed to Transformation Protocol, next page.

Note: You may store the BP reaction at  $-20^{\circ}$ C for up to 1 week before transformation, if desired.

## Performing the Standard BP Recombination Reaction, Continued

One Shot <sup>®</sup> TOP10 Transformation Protocol	Use this procedure to transform the BP recombination reaction into One Shot <sup>®</sup> TOP10 Chemically Competent <i>E. coli</i> . If you are using any other competent cell, follow the manufacturer's protocol.	
	1. Thaw, on ice, one vial of One Shot <sup>®</sup> TOP10 chemically competent cells for each transformation.	
	<ol> <li>Add 1 μL of the BP recombination reaction (from, Step 6, previous page) into a vial of One Shot<sup>®</sup> TOP10 cells and mix gently. <b>Do not mix by pipetting up</b> and down. For the pUC19 control, add 10 pg (1 μL) of DNA into a separate vial of One Shot<sup>®</sup> cells and mix gently.</li> </ol>	
	3. Incubate the vial(s) on ice for 30 minutes.	
	4. Heat-shock the cells for 30 seconds at 42°C without shaking.	
	5. Remove the vial(s) from the 42°C bath and place them on ice for 2 minutes.	
	6. Add 250 μL of room temperature S.O.C. medium to each vial.	
	7. Cap the vial(s) tightly and shake horizontally (225 rpm) at 37°C for 1 hour.	
	8. Before plating, dilute the transformation mix 1:10 into LB Medium ( <i>e.g.</i> , remove 20 $\mu$ L of the transformation mix and add to 180 $\mu$ L of LB Medium)	
	9. Spread 20 $\mu$ L and 100 $\mu$ L from each transformation on a prewarmed selective plate and incubate overnight at 37°C.	
	An efficient BP recombination reaction may produce hundreds of colonies (> 1,500 colonies if the entire BP reaction is transformed and plated).	
Verifying Entry Clones	You may verify the entry clones by performing restriction digestion analysis. Sequencing of the entry clone is not required as transfer of the pre-miRNA expression cassette from pcDNA6.2 <sup><math>\text{TM}</math></sup> -GW/ <u>+</u> EmGFP-miR into the pDONR <sup><math>\text{TM}</math></sup> 221 vector preserves the orientation of the cassette.	
	After verifying the entry clone, isolate plasmid DNA using your method of choice. We recommend using PureLink <sup>™</sup> Plasmid Purification Kits (see page 77). For the LR recombination reaction, you will need purified plasmid DNA at a concentration of 50–150 ng/µL in TE, pH 8.0. Proceed to the LR recombination reaction, next page.	
Verifying pEXP7- tet Entry Clones	If you included the pEXP7-tet control in your BP recombination reaction, you may transform One Shot <sup>®</sup> TOP10 competent cells using the protocol on the previous page. The efficiency of the BP reaction may then be assessed by streaking entry clones onto LB agar plates containing 20 µg/mL tetracycline. True entry clones should be tetracycline-resistant.	

## Performing the Standard LR Recombination Reaction

Introduction	After obtaining an entry clone containing the pre-miRNA expression cassette, perform an LR recombination reaction between the entry clone (previous page) and a destination vector (pLenti6/V5-DEST), and transform the reaction mixture into One Shot <sup>®</sup> Stbl3 <sup>™</sup> Competent <i>E. coli</i> to select for expression clones (see page 29). Use the standard LR reaction, if the Rapid BP/LR protocol produces fewer (~20–30) colonies.	
Experimental Outline	<ol> <li>To generate an expression clone, you will:</li> <li>Perform a LR recombination reaction using the <i>att</i>L-containing entry clone (previous page) and <i>att</i>R-containing pLenti6/V5-DEST.</li> <li>Transform the reaction mixture into a One Shot<sup>®</sup> Stbl3<sup>™</sup> Competent <i>E. coli</i>.</li> <li>Select for expression clones.</li> </ol>	
Destination Vector	See page 22 for details on the destination vector and instructions for propagating the destination vector.	
Important	The pLenti6/V5-DEST vector is supplied as a supercoiled plasmid. Although the Gateway <sup>®</sup> Technology manual previously recommended using a linearized destination vector for more efficient LR recombination, further testing has found that linearization of pLenti6/V5-DEST is not required to obtain optimal results for any downstream application.	
Recommended <i>E. coli</i> Host	For optimal results, use Stbl3 <sup>™</sup> E. coli for transformation as this strain is particularly well-suited for use in cloning unstable DNA such as lentiviral DNA containing direct repeats. One Shot <sup>®</sup> Stbl3 <sup>™</sup> Chemically Competent E. coli are included in the kit for transformation. For instructions, see Transforming One Shot <sup>®</sup> Stbl3 <sup>™</sup> Competent E. coli, page 29.	
Positive Control	The pENTR <sup><math>m</math></sup> -gus plasmid is provided with the Gateway <sup>®</sup> LR Clonase <sup>®</sup> II Enzyme Mix for use as a positive control for recombination. Using the pENTR <sup><math>m</math></sup> -gus entry clone in an LR recombination reaction with a destination vector will allow you to generate an expression clone containing the gene encoding $\beta$ -glucuronidase ( <i>gus</i> ) (Kertbundit <i>et al.</i> , 1991).	
Recombination Region of pLenti6/V5-DEST	The recombination region of the expression clone resulting from pLenti6/V5-DEST × entry clone is shown on page 23.	

## Performing the LR Recombination Reaction, Continued

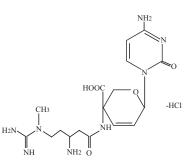
Materials Needed	•	Purified plasmid DNA of your entry clor	ne (50–150 ng	/μL in TE, pH 8.0)		
	•	• TE Buffer, pH 8.0 (10 mM Tris-HCl, pH 8.0, 1 mM EDTA)				
	•	• LB agar plates containing 100 µg/mL ampicillin (two for each transformation; warm at 37°C for 30 minutes before use) to select for expression clones				
	C01	nponents supplied with the kits	150 / 1 :			
	•	• pLenti6/V5-DEST (supplied with the kit, 150 ng/µL in TE, pH 8.0)				
	•	Gateway <sup>®</sup> LR Clonase <sup>®</sup> II enzyme mix (keep at –20°C until immediately before use)				
	•	• 2 $\mu$ g/ $\mu$ L Proteinase K solution (thaw and keep on ice until use)				
	٠	<ul> <li>pENTR<sup>™</sup>-gus positive control (50 ng/µL)</li> </ul>				
	٠	One Shot <sup>®</sup> Stbl3 <sup>™</sup> Chemically Competent	: E. coli			
	•	• S.O.C. Medium (warm to room temperature)				
Setting Up the LR Recombination	1.	Add the following components to 1.5 mI temperature and mix.	. microcentril	fuge tubes at room		
Reaction		<b>Note:</b> To include a negative control, set up a second sample reaction and omit the Gateway LR Clonase <sup>®</sup> II enzyme mix (see Step 4).				
		Component	Sample	<b>Positive Control</b>		
		Entry clone (50–150 ng/reaction)	1–7 µL			
		pLenti6/V5-DEST vector (150 ng/µL)	1 µL	1 µL		
		pENTR <sup>™</sup> -gus (50 ng/µL)		2 µL		
		TE Buffer, pH 8.0	to 8 µL	5 µL		
	2.	Remove the Gateway <sup>®</sup> LR Clonase <sup>®</sup> II en: (~ 2 minutes).	zyme mix fro	m −20°C and thaw on	ı ice	
	3.	Vortex the Gateway <sup>®</sup> LR Clonase <sup>®</sup> II enzyme mix briefly twice (2 seconds each time)				
	4.	Add 2 µL of Gateway <sup>®</sup> LR Clonase <sup>®</sup> II en control vials. Do not add Gateway <sup>®</sup> LR C control vial. Mix well by vortexing briefl	Clonase® II enz	zyme mix to the negat		
		Reminder: Return Gateway® LR Clonase®	II enzyme mix	to -20°C immediately a	fter use.	
	5.	Incubate reactions at 25°C for 1 hour.				
		<b>Note:</b> For most applications, 1 hour will yield a sufficient number of colonies for analysis. For large plasmids ( $\geq$ 10 kb), longer incubation times ( <i>i.e.</i> 18 hour incubation) will yield more colonies and are recommended.				
	6.	Add 1 $\mu$ L of the Proteinase K solution. In	cubate for 10	minutes at 37°C.		
	7.	Proceed to Transforming One Shot® Stb	ol3 <sup>™</sup> Compete	ent <i>E. coli</i> (page 29).		
Expected Results	LŔ	rou use <i>E. coli</i> cells with a transformation e reaction should give > 5,000 colonies if th ted. See page 30 for confirming the expres	e entire LR re	-	and	

## Blasticidin

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

Preparing and Storing Stock Solutions Always wear gloves, mask, goggles, and protective clothing (*e.g.*, a laboratory coat) when handling Blasticidin. Weigh out Blasticidin and prepare solutions in a hood.

Blasticidin may be obtained in 50 mg aliquots (see page 77 for ordering information).

- Blasticidin is soluble in water and acetic acid.
- Prepare a stock solution of 5 to 10 mg/mL Blasticidin in sterile water and filter-sterilize the solution.
- Aliquot in small volumes suitable for one time use 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 week at 4°C and 6–8 weeks at –20°C.
- pH of the aqueous solution should not exceed 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 discard the unused portion.
- Medium containing Blasticidin may be stored at 4°C for up to 2 weeks.

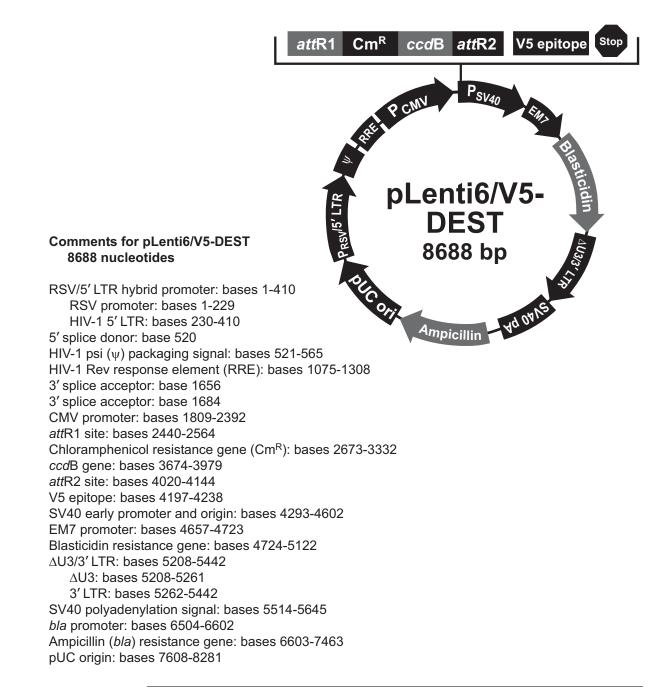
## Blasticidin, Continued

Determining Antibiotic Sensitivity	Since you will be selecting for stably transduced cells using Blasticidin, you must first determine the minimum concentration of Blasticidin required to kill your untransduced mammalian cell line ( <i>i.e.</i> perform a kill curve experiment). Typically, concentrations ranging from 2–10 $\mu$ g/mL Blasticidin are sufficient to kill most untransduced mammalian cell lines. Test a range of concentrations (see protocol below) to ensure that you determine the minimum concentration necessary for your cell line.
	<ol> <li>Plate cells at approximately 25% confluence. Prepare a set of 6 plates. Allow cells to adhere overnight.</li> </ol>
	2. The next day, substitute culture medium with medium containing varying concentrations of Blasticidin ( <i>e.g.</i> , 0, 2, 4, 6, 8, 10 $\mu$ g/mL Blasticidin).
	3. Replenish the selective media every 3–4 days and observe the percentage of surviving cells.
	4. Determine the appropriate concentration of Blasticidin that kills the cells within 10–14 days after addition of antibiotic.

## Map and Features of pLenti6/V5-DEST

### Map of pLenti6/V5-DEST

The map below shows the elements of pLenti6/V5-DEST. DNA from the entry clone replaces the region between bases 2,447 and 4,130. The sequence for pLenti6/V5-DEST is available at <u>www.invitrogen.com</u> or by contacting Technical Support (see page 79).



# Map and Features of pLenti6/V5-DEST, Continued

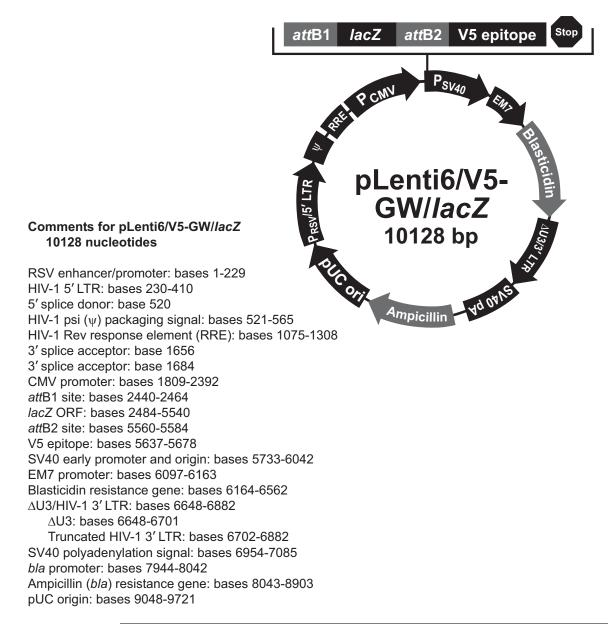
Features of the<br/>VectorThe pLenti6/V5-DEST (8,688 bp) vector contains the following elements. All<br/>features have been functionally tested and the vector is fully sequenced.

Feature	Benefit
Rous Sarcoma Virus (RSV) enhancer/promoter	Allows Tat-independent production of viral mRNA (Dull et al., 1998).
HIV-1 truncated 5' LTR	Permits viral packaging and reverse transcription of the viral mRNA (Luciw, 1996).
5' splice donor and 3' acceptors	Enhances the biosafety of the vector by facilitating removal of the $\Psi$ packaging sequence and RRE such that expression of the gene of interest in the transduced host cell is no longer Rev-dependent (Dull <i>et al.</i> , 1998).
HIV-1 psi (ψ) packaging signal	Allows viral packaging (Luciw, 1996).
HIV-1 Rev response element (RRE)	Permits Rev-dependent nuclear export of unspliced viral mRNA (Kjems <i>et al.</i> , 1991; Malim <i>et al.</i> , 1989).
CMV promoter	Permits high-level, constitutive expression of the gene or miRNA of interest (Andersson <i>et al.</i> , 1989; Boshart <i>et al.</i> , 1985; Nelson <i>et al.</i> , 1987).
attR1 and attR2 sites	Bacteriophage $\lambda$ -derived DNA recombination sequences that permit recombinational cloning of the gene of interest from a Gateway <sup>®</sup> entry clone (Landy, 1989).
Chloramphenicol resistance gene (Cm <sup>R</sup> )	Allows counterselection of the plasmid.
<i>ccd</i> B gene	Permits negative selection of the plasmid.
V5 epitope	Allows detection of the recombinant fusion protein by the Anti-V5 Antibodies (Southern <i>et al.</i> , 1991). <b>This feature is not used when</b> <b>expressing miRNAs.</b>
SV40 early promoter and origin	Allows high-level expression of the selection marker and episomal replication in cells expressing the SV40 large T antigen.
EM7 promoter	Synthetic prokaryotic promoter for expression of the selection marker in <i>E. coli</i> .
Blasticidin (bsd) resistance gene	Permits selection of stably transduced mammalian cell lines (Kimura <i>et al.,</i> 1994).
ΔU3/HIV-1 truncated 3' LTR	Allows viral packaging but self-inactivates the 5' LTR for biosafety purposes (Dull <i>et al.,</i> 1998). The element also contains a polyadenylation signal for transcription termination and polyadenylation of mRNA in transduced cells.
SV40 polyadenylation signal	Allows transcription termination and polyadenylation of mRNA.
<i>bla</i> promoter	Allows expression of the ampicillin resistance gene.
Ampicillin resistance gene (β-lactamase)	Allows selection of the plasmid in <i>E. coli</i> .
pUC origin	Permits high-copy replication and maintenance in <i>E. coli</i> .

## Map of pLenti6/V5-GW/lacZ

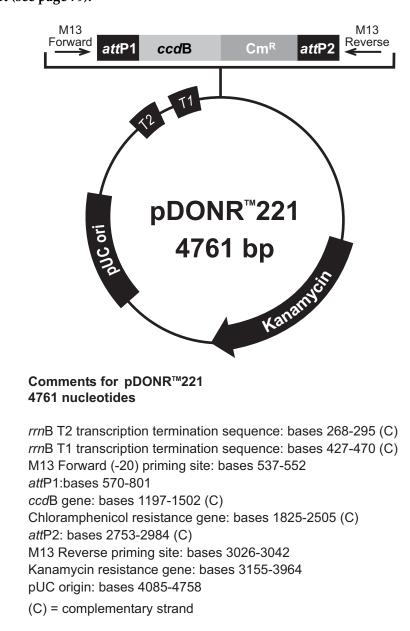
**Description** pLenti6/V5-GW/*lacZ* is a 10,128 bp control vector expressing β-galactosidase, and was generated using the Gateway<sup>®</sup> LR recombination reaction between an entry clone containing the *lacZ* gene and pLenti6/V5-DEST. β-galactosidase is expressed as a C-terminal V5 fusion protein with a molecular weight of approximately 121 kDa.

Map of pLenti6/V5-<br/>GW/lacZThe map below shows the elements of pLenti6/V5-GW/lacZ. The sequence of<br/>the vector is available at <a href="http://www.invitrogen.com">www.invitrogen.com</a> or by contacting Technical<br/>Support (see page 79).



## Map and Features of pDONR<sup>™</sup>221

Map of<br/>pDONR<sup>™</sup>221The map below shows the elements of pDONR<sup>™</sup>221 vector. The sequence of<br/>pDONR<sup>™</sup>221 is available at <u>www.invitrogen.com</u> or by contacting Technical<br/>Support (see page 79).



# Map and Features of pDONR<sup>™</sup>221, Continued

### Features of pDONR<sup>™</sup>221

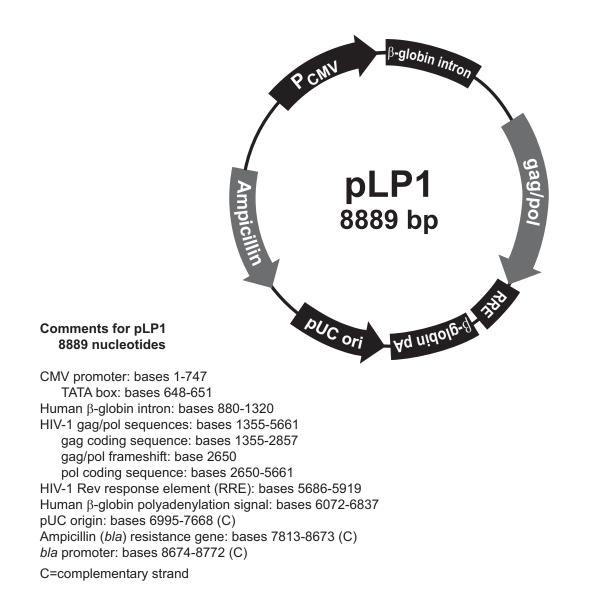
 $pDONR^{T}$  221 (4,761 bp) contains the following elements. All features have been functionally tested and the vectors fully sequenced.

Feature	Benefit
<i>rrn</i> B T1 and T2 transcription terminators	Protects the cloned gene or miRNA from expression by vector-encoded promoters, thereby reducing possible toxicity (Orosz <i>et al.</i> , 1991).
M13 Forward (–20) priming site	Allows sequencing in the sense orientation.
<i>att</i> P1 and <i>att</i> P2 sites	Bacteriophage $\lambda$ -derived DNA recombination sequences that allow recombinational cloning of the gene of interest from a Gateway <sup>®</sup> expression clone or <i>att</i> B PCR product (Landy, 1989).
ccdB gene	Allows negative selection of the plasmid.
Chloramphenicol resistance gene (Cm <sup>R</sup> )	Allows counterselection of the plasmid.
Kanamycin resistance gene	Allows selection of the plasmid in <i>E. coli</i> .
pUC origin	Allows high-copy replication and maintenance in <i>E. coli</i> .

## Map and Features of pLP1

#### pLP1 Map

The figure below shows the features of the pLP1 vector. Note that the *gag* and *pol* genes are initially expressed as a gag/pol fusion protein, which is self-cleaved by the viral protease into individual Gag and Pol polyproteins. The sequence of pLP1 is available at <u>www.invitrogen.com</u> or by contacting Technical Support (see page 79).



## Map and Features of pLP1, Continued

# Features of pLP1

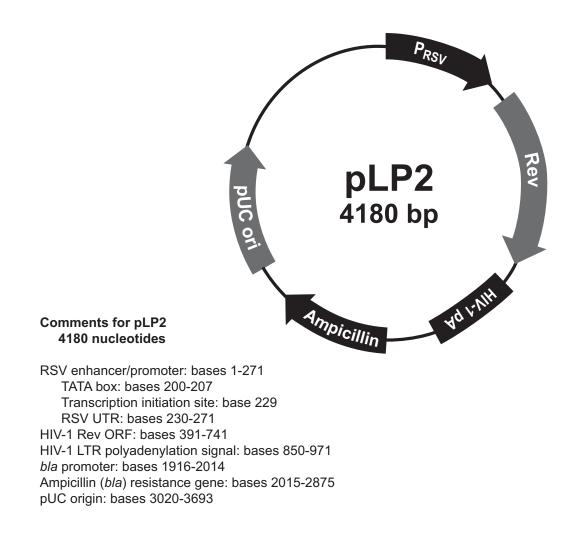
pLP1 (8,889 bp) contains the following elements. All features have been functionally tested.

Feature	Benefit
Human cytomegalovirus (CMV) promoter	Permits high-level expression of the HIV-1 <i>gag</i> and <i>pol</i> genes in mammalian cells (Andersson <i>et al.,</i> 1989; Boshart <i>et al.,</i> 1985; Nelson <i>et al.,</i> 1987).
Human β-globin intron	Enhances expression of the <i>gag</i> and <i>pol</i> genes in mammalian cells.
HIV-1 gag coding sequence	Encodes the viral core proteins required for forming the structure of the lentivirus (Luciw, 1996).
HIV-1 <i>pol</i> coding sequence	Encodes the viral replication enzymes required for replication and integration of the lentivirus (Luciw, 1996).
HIV-1 Rev response element (RRE)	Permits Rev-dependent expression of the <i>gag</i> and <i>pol</i> genes.
Human $\beta$ -globin polyadenylation signal	Allows efficient transcription termination and polyadenylation of mRNA.
pUC origin of replication (ori)	Permits high-copy replication and maintenance in <i>E. coli</i> .
Ampicillin (bla) resistance gene	Allows selection of the plasmid in <i>E. coli</i> .

## Map and Features of pLP2

pLP2 Map

The figure below shows the features of the pLP2 vector. The sequence of pLP2 is available at <u>www.invitrogen.com</u> or by contacting Technical Support (see page 79).



# Map and Features of pLP2, Continued

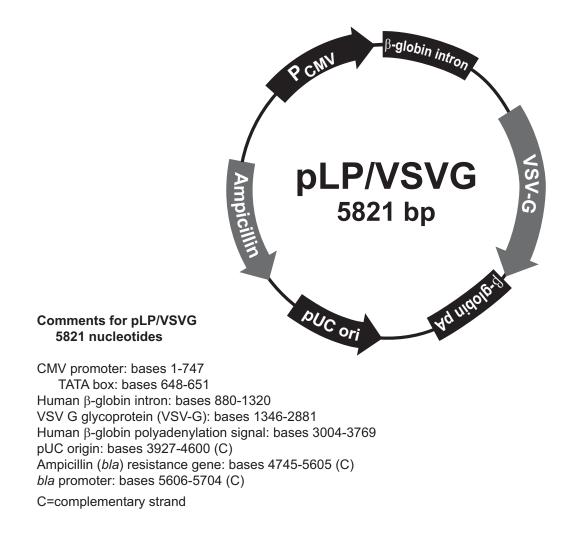
Features of<br/>pLP2pLP2 (4,180 bp) contains the following elements. All features have been<br/>functionally tested.

Feature	Benefit
RSV enhancer/promoter	Permits high-level expression of the <i>rev</i> gene (Gorman <i>et al.,</i> 1982).
HIV-1 Rev ORF	Encodes the Rev protein which interacts with the RRE on pLP1 to induce Gag and Pol expression, and on the pLenti6/V5-DEST expression vector to promote the nuclear export of the unspliced viral RNA for packaging into viral particles.
HIV-1 LTR polyadenylation signal	Allows efficient transcription termination and polyadenylation of mRNA.
Ampicillin (bla) resistance gene	Allows selection of the plasmid in <i>E. coli</i> .
pUC origin of replication ( <i>ori</i> )	Permits high-copy replication and maintenance in <i>E. coli.</i>

## Map and Features of pLP/VSVG

#### pLP/VSVG Map

The figure below shows the features of the pLP/VSVG vector. The sequence of pLP/VSVG is available at <u>www.invitrogen.com</u> or by contacting Technical Support (see page 79).



# Map and Features of pLP/VSVG, Continued

Features of<br/>pLP/VSVGpLP/VSVG (5,821 bp) contains the following elements. All features have been<br/>functionally tested.

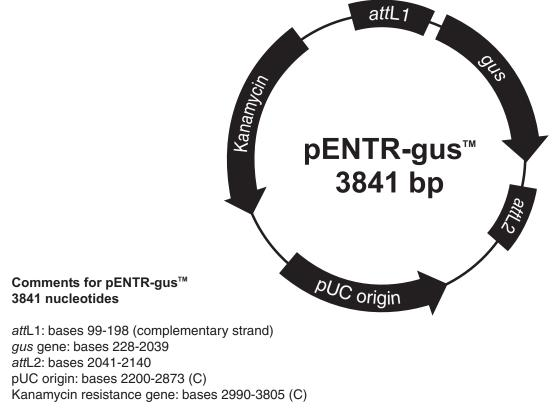
Feature	Benefit
Human CMV promoter	Permits high-level expression of the VSV-G gene in mammalian cells (Andersson <i>et al.</i> , 1989; Boshart <i>et al.</i> , 1985; Nelson <i>et al.</i> , 1987).
Human β-globin intron	Enhances expression of the VSV-G gene in mammalian cells.
VSV G glycoprotein (VSV-G)	Encodes the envelope G glycoprotein from Vesicular Stomatitis Virus to allow production of a pseudotyped retrovirus with a broad host range (Burns <i>et al.</i> , 1993; Emi <i>et al.</i> , 1991; Yee <i>et al.</i> , 1994).
Human $\beta$ -globin polyadenylation signal	Allows efficient transcription termination and polyadenylation of mRNA.
pUC origin of replication (ori)	Permits high-copy replication and maintenance in <i>E. coli</i> .
Ampicillin (bla) resistance gene	Allows selection of the plasmid in <i>E. coli</i> .

## Map of pENTR<sup>™</sup>-gus

Vector

Description pENTR<sup>™</sup>-gus is a 3,841 bp entry clone containing the *Arabidopsis thaliana* gene for β-glucuronidase (*gus*) (Kertbundit *et al.*, 1991). The *gus* gene was amplified using PCR primers containing attB recombination sites. The amplified PCR product was then used in a BP recombination reaction with pDONR201<sup>™</sup> to generate the entry clone. For more information about the BP recombination reaction, refer to the Gateway® Technology with Clonase® II manual which is available at www.invitrogen.com or by contacting Technical Support (see page 79).

Map of Control The figure below summarizes the features of the pENTR<sup>™</sup>-gus vector. **The** sequence for pENTR<sup>™</sup>-gus is available at <u>www.invitrogen.com</u> or by contacting Technical Support (see page 79).



C = complementary strand

## **Additional Products**

#### Accessory Products

Many of the reagents supplied in the BLOCK-iT<sup>™</sup> Lentiviral Pol II miR RNAi Expression System as well as other products suitable for use with the kits are available separately. Ordering information is provided below. For more information, go to <u>www.invitrogen.com</u> or contact Technical Support (see page 79).

Product	Amount	Cat. no.
BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kit	20 constructions	K4935-00
BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vector Kit with EmGFP	20 constructions	K4936-00
Gateway <sup>®</sup> BP Clonase <sup>™</sup> II Enzyme Mix	20 reactions	11789-020
	100 reactions	11789-100
Gateway <sup>®</sup> LR Clonase <sup>™</sup> II Enzyme Mix	20 reactions	11791-020
	100 reactions	11791-100
One Shot <sup>®</sup> Stbl3 <sup>™</sup> Chemically Competent <i>E. coli</i>	$20\times 50~\mu L$	C7373-03
One Shot® TOP10 Chemically Competent E. coli	$20\times 50~\mu L$	C4040-03
OmniMAX <sup>™</sup> 2-T1 <sup>R</sup>	$20\times 50~\mu L$	C854003
Blasticidin	50 mg	R21001
ViraPower <sup>™</sup> Bsd Lentiviral Support Kit	20 reactions	K4970-00
One Shot <sup>®</sup> ccdB Survival <sup>™</sup> 2 T1 <sup>R</sup> Chemically Competent Cells	10 reactions	A10460
ViraPower <sup>™</sup> Lentiviral Packaging Mix	60 reactions	K4975-00
Lipofectamine <sup>®</sup> 2000 Reagent	0.75 mL	11668-027
	1.5 mL	11668-019
pDONR™221 Vector	6 µg	12536-017
Opti-MEM <sup>®</sup> I Reduced Serum Medium	100 mL	31985-062
	500 mL	31985-070
293FT Cell Line	$3 \times 10^{6}$ cells	R700-07
Phosphate-Buffered Saline (PBS), pH 7.4	500 mL	10010-023
	1 L	10010-031
FluoReporter <sup>®</sup> lacZ/Galactosidase Quantitation Kit	1,000 assays	F-2905
β-Gal Staining Kit	1 kit	K1465-01
Ampicillin	200 mg	11593027
TE, pH 8.0	500 mL	AM9849
PureLink <sup>™</sup> HiPure Plasmid DNA Purification MidiPrep Kit	25 reactions	K2100-04
PureLink <sup>™</sup> HQ Mini Plasmid Purification Kit	100 preps	K2100-01
Fetal Bovine Serum (FBS), Certified	500 mL	16000-044
MEM Sodium Pyruvate Solution (100X)	100 mL	11360-070
WesternBreeze® Chemiluminescent Kit Anti-Mouse	1 kit	WB7104

# Additional Products, Continued

BLOCK-iT <sup>™</sup> RNAi Designer	The BLOCK-iT <sup>™</sup> RNAi Designer is an online tool to help you design and order microRNA sequences for any target gene of interest. The RNAi Designer incorporates the guidelines provided in this manual as well as other design rul into a proprietary algorithm to design microRNA sequences that are compatib for use in cloning into the BLOCK-iT <sup>™</sup> Pol II miR RNAi Expression Vectors.	
BLOCK-iT <sup>™</sup> RNAi Products	A large variety of BLOCK-iT <sup>™</sup> RNAi products are available to facilitate RNAi analysis including Stealth <sup>™</sup> RNAi, Validated Stealth <sup>™</sup> RNAi Collection, Validated miRNA Vector Collection, and a large selection of RNAi vectors.	
	For details, visit the RNAi central portal or contact Technical Support (see page 79).	

## **Technical Support**

#### Web Resources



- Visit the Invitrogen website at <u>www.invitrogen.com</u> for:
- Technical resources, including manuals, vector maps and sequences, application notes, SDSs, FAQs, formulations, citations, handbooks, etc.
- Complete technical support contact information
- Access to the Invitrogen Online Catalog
- Additional product information and special offers

### **Contact Us**

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

<b>Corporate Headquarte</b> 5791 Van Allen Way Carlsbad, CA 92008 US Tel: 1 760 603 7200 Tel (Toll Free): 1 800 95 Fax: 1 760 602 6500 E-mail: <u>tech_support@</u>	5A 55 6288	Japanese Headquarters: LOOP-X Bldg. 6F 3-9-15, Kaigan Minato-ku, Tokyo 108-0022 Tel: 81 3 5730 6509 Fax: 81 3 5730 6519 E-mail: jpinfo@invitrogen.com	<b>European Headquarters:</b> Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF, UK Tel: +44 (0) 141 814 6100 Tech Fax: +44 (0) 141 814 6117 E-mail: <u>eurotech@invitrogen.com</u>
SDS	Safety Data Sheet	s (SDSs) are available on our website a	t <u>www.invitrogen.com/sds</u> .
Certificate of Analysis	The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to www.invitrogen.com/support and search for the Certificate of Analysis by product lot number, which is printed on the box.		e available on our website. Go to
Limited Warranty	with high-quality satisfied with our about an Invitrog All Invitrogen pro- certificate of anal- meet those specifi- <u>product.</u> No warr warranty is appli- instructions. The product unless th the order. Invitrogen makes occasional typogr warranty of any k discover an error Representatives. Life Technologie incidental, indire warranty is sole a	to f Life Technologies Corporation) is consistent of Life Technologies Corporation) is consistent of the service. Our goal is to ensure products and our service, contact our Technolucts are warranted to perform according site. The Company will replace, free of ications. This warranty limits the Company is granted for products beyond the cable unless all product components ar Company reserves the right to select the Company agrees to a specified method every effort to ensure the accuracy of if aphical or other error is inevitable. The cind regarding the contents of any publications, please reports a corporation shall have no responsible to consequential loss or damage with a exclusive. No other warranty is material or fitness for the contents of the consequential loss or fitness for the consequent conseque	ure that every customer is 100% d have any questions or concerns nical Support Representatives. ding to specifications stated on the charge, any product that does not <u>bany's liability to only the price of the</u> eir listed expiration date. No e stored in accordance with the method(s) used to analyze a od in writing prior to acceptance of the publications, but realizes that the erefore the Company makes no lications or documentation. If you rt it to our Technical Support <b>bility or liability for any special,</b> <b>hatsoever. The above limited</b> <b>ade, whether expressed or implied,</b>

### **Purchaser Notification**

Limited Use Label License No. 19: Gateway® Cloning Products The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for profit entity). The purchase of this product does not convey a license under any method claims in the foregoing patents or patent applications, or to use this product with any recombination sites other than those purchased from 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 Clonase<sup>™</sup> 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.

Limited Use Label License No. 23: GUS control vector	The GUS positive control vector in these products is claimed in patents and patent applications (See U.S. Patent No. 5,599,670 and Great Britain Patent No. 2,197,653) licensed to Life Technologies by Cambia Biosystems, L.L.C. ("CBL"). Use of the GUS gene is restricted to use as a positive control. Any other use may require a license from CBL.
Limited Use Label License No. 27: RNA Transfection	Use of this product in conjunction with methods for the introduction of RNA molecules into cells may require licenses to one or more patents or patent applications. Users of these products should determine if any licenses are required.
Limited Use Label License No. 51: Blasticidin and the Blasticidin Selection Marker	Blasticidin and the blasticidin resistance gene ( <i>bsd</i> ) 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. 54: ULB ccdB Selection Technology	This product is the subject of one or more of U.S. Patent Numbers 5,910,438, 6,180,407, and 7,176,029 and corresponding foreign patents and is sold under license from the Université Libre de Bruxelles for research purposes only. ccdB selection technology is described in Bernard et al., "Positive Selection Vectors Using the F Plasmid ccdB Killer Gene" Gene 148 (1994) 71-74. 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). For licensing information for use in other than research, please contact: Out Licensing, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008; Phone (760) 603-7200 or e-mail.
Limited Use Label License No: 108 Lentiviral Technology	The Lentiviral Technology (based upon the lentikat <sup>™</sup> system) is licensed from Cell Genesys, Inc., under U.S. Patent Nos. 5,686,279; 5,834,256; 5,858,740; 5,994,136; 6,013,516; 6,051,427; 6,165,782 and 6,218,187 and corresponding patents and applications in other countries for internal research purposes only. Use of this technology for gene therapy applications or bioprocessing other than for non- human research use requires a license from Cell Genesys (Cell Genesys, Inc. 342 Lakeside Drive, Foster City, California 94404). 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), including non-gene therapy research and target validation applications in laboratory animals.

Limited Use Label License No: 109 Retroviral Helper Lines	Retroviral helper cell lines are licensed from Wisconsin Alumni Research Foundation, under U.S. Patents and corresponding patents and applications in other countries for internal research purposes only. Use of these cell lines for Commercial Purposes requires a license from Life Technologies.
Limited Use Label License No. 127: GFP with Heterologous Promoter	This product and its use is the subject of one or more of U.S. Patent Nos. 5,491,084 and 6,146,826, and foreign equivalents. This product is sold under license from Columbia University. Rights to use this product are limited to research use only, and expressly exclude the right to manufacture, use, sell or lease this product for use for measuring the level of toxicity for chemical agents and environmental samples in cells and transgenic animals. No other rights are conveyed. Not for human use or use in diagnostic or therapeutic procedures. Inquiry into the availability of a license to broader rights or the use of this product for commercial purposes should be directed to Columbia Innovation Enterprise, Columbia University, Engineering Terrace-Suite 363, New York, New York 10027.
Limited Use Label License No. 267: Mutant GFP Products	This product and its use is the subject of one or more of U.S. Patent Nos. 6,090,919, 5,804,387, 5,994,077, and foreign equivalents.
Limited Use Label License No. 270: miRNA Vectors	This product is produced and sold under license from the University of Michigan. Use of this product is permitted for research purposes only. Any other use requires a license from the University of Michigan, Office of Technology Transfer, 3003 S. State St., Suite 2071, Ann Arbor, MI 48190-1280.
Limited Use Label License No. 271: miRNA Constructs	This product is produced and sold under license from Galapagos Genomics N.V, for research use only and not for therapeutic or diagnostic use in humans. This product is not sold with license to use this product in conjunction with adenoviral vectors.
Limited Use Label License No. 272: Humanized GFP	This product is the subject of one or more of U.S. Patent Numbers 5,786,464, 5,795,737, 5,874,304, and 6,114,148 and foreign equivalents licensed by Life Technologies Corporation. This product is sold for research use only. Not for therapeutic or diagnostic use in humans.

Limited Use Label License No: 317 LentiVector<sup>®</sup> Technology This product is licensed under U.S. Pat. Nos. 5,817,491; 5,591,624; 5,716,832; 6,312,682; 6,669,936; 6,235,522; 6,924,123 and foreign equivalents from Oxford BioMedica (UK) Ltd., Oxford, UK, and is provided for use in academic and commercial in vitro and in vivo research for elucidating gene function, and for validating potential gene products and pathways for drug discovery and development, but excludes any use of LentiVector® technology for: creating transgenic birds for the purpose of producing useful or valuable proteins in the eggs of such transgenic birds, the delivery of gene therapies, and for commercial production of therapeutic, diagnostic or other commercial products not intended for research use where such products do not consist of or incorporate a lentiviral vector. Information about licenses for commercial uses excluded under this license is available from Oxford BioMedica (UK), Ltd., Medawar Centre, Oxford Science Park, Oxford OX4 4GA UK enquiries@oxfordbiomedica.co.uk or BioMedica Inc., 11622 EI Camino Real #100, San Diego CA 92130-2049 USA. LentiVector is a registered US and European Community trade mark of Oxford BioMedica plc.

Limited Use Label License No. 352: Vectors containing genes for Green Fluorescent Proteins

This product and its use is the subject of one or more of U.S. Patent Nos. 5,777,079, 6,066,476, and 6,319,669 and foreign equivalents. 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. Buyer means an academic entity, or a not for profit entity which is not owned or controlled by a for profit entity (an "NFPE") No rights are conveyed to modify or clone the gene encoding GFP contained in this product. The Buyer cannot sell (a) this product, (b) its components, or (c) materials made by the employment of this product or its components (collectively, "Materials") to any third party, or otherwise transfer Materials to a for profit entity, or any entity owned or controlled by a for profit entity unless such transfer has been approved in writing by Life Technologies Corporation, or otherwise use Materials for Commercial Purposes. The Buyer may transfer information or materials made through the employment of this product to an academic entity or an NFPE for the purposes of scientific collaboration, provided that such transfer is pursuant to the terms of a material transfer agreement ("MTA") that may be obtained as described below. 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; (4) resale of the product or its components, whether or not such product or its components are resold for use in research or (5) any use by a for-profit entity or an entity owned or controlled by a for profit entity. 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 this product, or any of its components was used in the manufacture of such product. If the Buyer 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, or to obtain the MTA discussed above, contact Licensing Department, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008. Phone (760) 603-7200 or outlicensing@lifetech.com.

### Limited Use Label License No. 358: Research Use Only

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

# Gateway<sup>®</sup> 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 <sup>®</sup> Technology.
Gateway <sup>®</sup> Entry Clones	Invitrogen understands that Gateway <sup>®</sup> 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 <sup>®</sup> Expression Clones	Invitrogen also understands that Gateway <sup>®</sup> 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 <sup>®</sup> expression clones for a nominal fee (\$10 per clone) payable to Invitrogen.
Additional Terms and Conditions	We would ask that such distributors of Gateway <sup>®</sup> entry and expression clones indicate that such clones may be used only for research purposes, that such clones incorporate the Gateway <sup>®</sup> Technology, and that the purchase of Gateway <sup>®</sup> Clonase <sup>®</sup> from Invitrogen is required for carrying out the Gateway <sup>®</sup> recombinational cloning reaction. This should allow researchers to readily identify Gateway <sup>®</sup> containing clones and facilitate their use of this powerful technology in their research. Use of Invitrogen's Gateway <sup>®</sup> Technology, including Gateway <sup>®</sup> 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

Ambros, V. (2001) MicroRNAs: Tiny Regulators with Great Potential. Cell 107, 823-826

Ambros, V. (2004) The functions of animal microRNAs. Nature 431, 350-355

- Anandalakshmi, R., Pruss, G. J., Ge, X., Marathe, R., Mallory, A. C., Smith, T. H., and Vance, V. B. (1998) A Viral Suppressor of Gene Silencing in Plants. Proc. Natl. Acad. Sci. USA 95, 13079-13084
- 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
- Bartel, D. P. (2004) MicroRNAs: genomics, biogenesis, mechanism, and function. Cell 116, 281-297
- Bernstein, E., Caudy, A. A., Hammond, S. M., and Hannon, G. J. (2001) Role for a Bidentate Ribonuclease in the Initiation Step of RNA Interference. Nature 409, 363-366
- 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
- Brummelkamp, T. R., Bernards, R., and Agami, R. (2002) A System for Stable Expression of Short Interfering RNAs in Mammalian Cells. Science 296, 550-553
- Buchschacher, G. L., Jr., and Wong-Staal, F. (2000) Development of Lentiviral Vectors for Gene Therapy for Human Diseases. Blood 95, 2499-2504
- Burns, J. C., Friedmann, T., Driever, W., Burrascano, M., and Yee, J.-K. (1993) Vesicular Stomatitis Virus G Glycoprotein Pseudotyped Retroviral Vectors: Concentration to a Very High Titer and Efficient Gene Transfer into Mammalian and Nonmammalian Cells. Proc. Natl. Acad. Sci. USA 90, 8033-8037
- Carrington, J. C., and Ambros, V. (2003) Role of MicroRNAs in Plant and Animal Development. Science 301, 336-338
- Ciccarone, V., Chu, Y., Schifferli, K., Pichet, J.-P., Hawley-Nelson, P., Evans, K., Roy, L., and Bennett, S. (1999) Lipofectamine<sup>™</sup> 2000 Reagent for Rapid, Efficient Transfection of Eukaryotic Cells. Focus 21, 54-55
- Cogoni, C., and Macino, G. (1997) Isolation of Quelling-Defective (qde) Mutants Impaired in Posttranscriptional Transgene-Induced Gene Silencing in *Neurospora crassa*. Proc. Natl. Acad. Sci. USA 94, 10233-10238
- Cogoni, C., and Macino, G. (1999) Gene Silencing in *Neurospora crassa* Requires a Protein Homologous to RNA-Dependent RNA Polymerase. Nature 399, 166-169
- Cogoni, C., Romano, N., and Macino, G. (1994) Suppression of Gene Expression by Homologous Transgenes. Antonie Van Leeuwenhoek 65, 205-209

Cullen, B. R. (2004) Derivation and function of small interfering RNAs and microRNAs. Virus Res 102, 3-9.

Cullen, B. R. (2004) Transcription and processing of human microRNA precursors. Mol Cell 16, 861-865

- Curradi, M., Izzo, A., Badaracco, G., and Landsberger, N. (2002) Molecular Mechanisms of Gene Silencing Mediated by DNA Methylation. Mol. Cell. Biol. 22, 3157-3173
- Dull, T., Zufferey, R., Kelly, M., Mandel, R. J., Nguyen, M., Trono, D., and Naldini, L. (1998) A Third-Generation Lentivirus Vector with a Conditional Packaging System. J. Virol. 72, 8463-8471
- Elbashir, S. M., Harborth, J., Lendeckel, W., Yalcin, A., Weber, K., and Tuschl, T. (2001) Duplexes of 21-Nucleotide RNAs Mediate RNA Interference in Cultured Mammalian Cells. Nature 411, 494-498
- Emi, N., Friedmann, T., and Yee, J.-K. (1991) Pseudotype Formation of Murine Leukemia Virus with the G Protein of Vesicular Stomatitis Virus. J. Virol. 65, 1202-1207
- Gorman, C. M., Merlino, G. T., Willingham, M. C., Pastan, I., and Howard, B. H. (1982) The Rous Sarcoma Virus Long Terminal Repeat is a Strong Promoter When Introduced into a Variety of Eukaryotic Cells by DNA-mediated Transfection. Proc. Natl. Acad. Sci. USA 79, 6777-6781
- Hammond, S. M., Bernstein, E., Beach, D., and Hannon, G. J. (2000) An RNA-Directed Nuclease Mediates Genetic Interference in *Caenorhabditis elegans*. Nature 404, 293-296
- 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
- Jones, A. L., Thomas, C. L., and Maule, A. J. (1998) *De novo* Methylation and Co-Suppression Induced by a Cytoplasmically Replicating Plant RNA Virus. EMBO J. 17, 6385-6393
- Kertbundit, S., Greve, H. d., Deboeck, F., Montagu, M. V., and Hernalsteens, J. P. (1991) *In vivo* Random βglucuronidase Gene Fusions in *Arabidopsis thaliana*. Proc. Natl. Acad. Sci. USA *88*, 5212-5216
- Ketting, R. F., Fischer, S. E., Bernstein, E., Sijen, T., Hannon, G. J., and Plasterk, R. H. (2001) Dicer Functions in RNA Interference and in Synthesis of Small RNA Involved in Developmental Timing in *C. elegans*. Genes Dev. 15, 2654-2659
- Kim, V. N. (2005) MicroRNA biogenesis: coordinated cropping and dicing. Nat Rev Mol Cell Biol 6, 376-385.
- 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
- Kjems, J., Brown, M., Chang, D. D., and Sharp, P. A. (1991) Structural Analysis of the Interaction Between the Human Immunodeficiency Virus Rev Protein and the Rev Response Element. Proc. Natl. Acad. Sci. USA 88, 683-687
- Landy, A. (1989) Dynamic, Structural, and Regulatory Aspects of Lambda Site-specific Recombination. Ann. Rev. Biochem. 58, 913-949
- Lee, Y., Kim, M., Han, J., Yeom, K. H., Lee, S., Baek, S. H., and Kim, V. N. (2004) MicroRNA genes are transcribed by RNA polymerase II. Embo J 23, 4051-4060

- Lewis, P. F., and Emerman, M. (1994) Passage Through Mitosis is Required for Oncoretroviruses but not for the Human Immunodeficiency Virus. J. Virol. *68*, 510-516
- Li, W. X., and Ding, S. W. (2001) Viral Suppressors of RNA Silencing. Curr. Opin. Biotechnol. 12, 150-154
- Liu, J., Carmell, M. A., Rivas, F. V., Marsden, C. G., Thomson, J. M., Song, J. J., Hammond, S. M., Joshua-Tor, L., and Hannon, G. J. (2004) Argonaute2 is the catalytic engine of mammalian RNAi. Science 305, 1437-1441
- Luciw, P. A. (1996) in *Fields Virology* (Fields, B. N., Knipe, D. M., Howley, P. M., Chanock, R. M., Melnick, J. L., Monath, T. P., Roizman, B., and Straus, S. E., eds), 3rd Ed., pp. 1881-1975, Lippincott-Raven Publishers, Philadelphia, PA
- Lund, E., Guttinger, S., Calado, A., Dahlberg, J. E., and Kutay, U. (2004) Nuclear export of microRNA precursors. Science *303*, 95-98.
- Malim, M. H., Hauber, J., Le, S. Y., Maizel, J. V., and Cullen, B. R. (1989) The HIV-1 Rev Trans-activator Acts Through a Structured Target Sequence to Activate Nuclear Export of Unspliced Viral mRNA. Nature 338, 254-257
- McManus, M. T., and Sharp, P. A. (2002) Gene Silencing in Mammals by Small Interfering RNAs. Nature Rev. Genet. 3, 737-747
- Meister, G., Landthaler, M., Patkaniowska, A., Dorsett, Y., Teng, G., and Tuschl, T. (2004) Human Argonaute2 mediates RNA cleavage targeted by miRNAs and siRNAs. Mol Cell 15, 185-197
- Naldini, L. (1998) Lentiviruses as Gene Transfer Agents for Delivery to Non-dividing Cells. Curr. Opin. Biotechnol. 9, 457-463
- Naldini, L. (1999) in *The Development of Human Gene Therapy* (Friedmann, T., ed), pp. 47-60, Cold Spring Harbor Laboratory Press, Cold Spring Harbor, NY
- Naldini, L., Blomer, U., Gage, F. H., Trono, D., and Verma, I. M. (1996) Efficient Transfer, Integration, and Sustained Long-Term Expression of the Transgene in Adult Rat Brains Injected with a Lentiviral Vector. Proc. Natl. Acad. Sci. USA *93*, 11382-11388
- Napoli, C., Lemieux, C., and Jorgensen, R. (1990) Introduction of a Chalcone Synthase Gene into *Petunia* Results in Reversible Co-Suppression of Homologous Genes *in trans*. Plant Cell 2, 279-289
- 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
- Nykanen, A., Haley, B., and Zamore, P. D. (2001) ATP Requirements and Small Interfering RNA Structure in the RNA Interference Pathway. Cell 107, 309-321
- Orosz, A., Boros, I., and Venetianer, P. (1991) Analysis of the Complex Transcription Termination Region of the *Escherichia coli rrnB* Gene. Eur. J. Biochem. 201, 653-659

- Paddison, P. J., Caudy, A. A., Bernstein, E., Hannon, G. J., and Conklin, D. S. (2002) Short Hairpin RNAs (shRNAs) Induce Sequence-Specific Silencing in Mammalian Cells. Genes Dev. 16, 948-958
- Paul, C. P., Good, P. D., Winer, I., and Engelke, D. R. (2002) Effective Expression of Small Interfering RNA in Human Cells. Nat. Biotechnol. 20, 505-508
- Rietveld, L. E., Caldenhoven, E., and Stunnenberg, H. G. (2002) In vivo Repression of an Erythroid-Specific Gene by Distinct Corepressor Complexes. EMBO J. 21, 1389-1397
- Romano, N., and Macino, G. (1992) Quelling: Transient Inactivation of Gene Expression in *Neurospora crassa* by Transformation with Homologous Sequences. Mol. Microbiol. *6*, 3343-3353
- Sastry, L., Johnson, T., Hobson, M. J., Smucker, B., and Cornetta, K. (2002) Titering Lentiviral vectors:comparison of DNA, RNA and marker expression methods. Gene Ther. *9*, 1155-1162
- Shimomura, O., Johnson, F. H., and Saiga, Y. (1962) Extraction, Purification and Properties of Aequorin, a Bioluminescent Protein from the Luminous hHydromedusan, Aequorea. Journal of Cellular and Comparative Physiology 59, 223-239
- Smith, C. J., Watson, C. F., Bird, C. R., Ray, J., Schuch, W., and Grierson, D. (1990) Expression of a Truncated Tomato Polygalacturonase Gene Inhibits Expression of the Endogenous Gene in Transgenic Plants. Mol. Gen. Genet. 224, 477-481
- 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
- Sui, G., Soohoo, C., Affar, E. B., Gay, F., Shi, Y., Forrester, W. C., and Shi, Y. (2002) A DNA Vector-Based RNAi Technology to Suppress Gene Expression in Mammalian Cells. Proc. Natl. Acad. Sci. USA 99, 5515-5520
- 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
- Tsien, R. Y. (1998) The Green Fluorescent Protein. Annu. Rev. Biochem. 67, 509-544
- van der Krol, A. R., Mur, L. A., Beld, M., Mol, J. N., and Stuitje, A. R. (1990) Flavonoid Genes in *Petunia*: Addition of a Limited Number of Gene Copies May Lead to a Suppression of Gene Expression. Plant Cell 2, 291-299
- Voinnet, O., Pinto, Y. M., and Baulcombe, D. C. (1999) Suppression of Gene Silencing: A General Strategy Used by Diverse DNA and RNA Viruses of Plants. Proc. Natl. Acad. Sci. USA *96*, 14147-14152
- White, S. M., Renda, M., Nam, N. Y., Klimatcheva, E., Y.Zhu, Fisk, J., Halterman, M., Rimel, B. J., Federoff, H., Pandya, S., Rosenblatt, J. D., and Planelles, V. (1999) Lentivirus vectors using human and simian imunodeficiency virus elements. J Virology 73, 2832-2840

- 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
- Yee, J.-K., Miyanohara, A., LaPorte, P., Bouic, K., Burns, J. C., and Friedmann, T. (1994) A General Method for the Generation of High-Titer, Pantropic Retroviral Vectors: Highly Efficient Infection of Primary Hepatocytes. Proc. Natl. Acad. Sci. USA *91*, 9564-9568
- Yee, J. K. (1999) in *The Development of Human Gene Therapy* (Friedmann, T., ed), pp. 21-45, Cold Spring Harbor Laboratory Press, Cold Spring Harbor, NY
- Yee, J. K., Moores, J. C., Jolly, D. J., Wolff, J. A., Respess, J. G., and Friedmann, T. (1987) Gene Expression from Transcriptionally Disabled Retroviral Vectors. Proc. Natl. Acad. Sci. USA 84, 5197-5201
- Yekta, S., Shih, I. H., and Bartel, D. P. (2004) MicroRNA-directed cleavage of HOXB8 mRNA. Science 304, 594-596
- Yi, R., Qin, Y., Macara, I. G., and Cullen, B. R. (2003) Exportin-5 mediates the nuclear export of premicroRNAs and short hairpin RNAs. Genes Dev 17, 3011-3016
- Yu, J. Y., DeRuiter, S. L., and Turner, D. L. (2002) RNA Interference by Expression of Short-interfering RNAs and Hairpin RNAs in Mammalian Cells. Proc. Natl. Acad. Sci. USA *99*, 6047-6052
- Yu, S. F., Ruden, T. v., Kantoff, P. W., Garber, C., Seiberg, M., Ruther, U., Anderson, W. F., Wagner, E. F., and Gilboa, E. (1986) Self-Inactivating Retroviral Vectors Designed for Transfer of Whole Genes into Mammalian Cells. Proc. Natl. Acad. Sci. USA *83*, 3194-3198
- Yu, Z., Raabe, T., and Hecht, N. B. (2005) MicroRNA122a Reduces Expression of the Post-Transcriptionally Regulated Germ Cell Transition Protein 2 (Tnp2) Messenger RNA (mRNA) by mRNA Cleavage. Biol Reprod 18
- Zeng, Y., Wagner, E. J., and Cullen, B. R. (2002) Both Natural and Designed MicroRNAs Can Inhibit the Expression of Cognate mRNAs When Expressed in Human Cells. Mol Cell 9, 1327-1333
- Zeng, Y., Yi, R., and Cullen, B. R. (2005) Recognition and cleavage of primary microRNA precursors by the nuclear processing enzyme Drosha. Embo J 24, 138-148
- Zhang, G., Gurtu, V., and Kain, S. (1996) An Enhanced Green Fluorescent Protein Allows Sensitive Detection of Gene Transfer in Mammalian Cells. Biochem. Biophys. Res. Comm. 227, 707-711
- Zufferey, R., Dull, T., Mandel, R. J., Bukovsky, A., Quiroz, D., Naldini, L., and Trono, D. (1998) Selfinactivating Lentivirus Vector for Safe and Efficient *in vivo* Gene Delivery. J. Virol. 72, 9873-9880

©2010 Life Technologies Corporation. All rights reserved.

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.



Corporate Headquarters 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