

BioModule[™] BLOCK-iT[™] Units with Pol II miR RNAi Expression System

Gateway[®]-adapted expression vector and lentiviral destination vector for high-level expression of microRNA in mammalian cells

Catalog nos. WFGE07, WFGE08

Version B 28 December 2010 25-0883

User Manual

Table of Contents

Table of Contents	iii
Experienced Users Procedure to Generate Expression Clone	v
Kit Contents and Storage	vi
Introduction	1
Overview	1
Using miRNA for RNAi Analysis	7
BioModule™ BLOCK-iT™ Unit with Pol II miR RNAi Expression Vector	12
BioModule [™] BLOCK-iT [™] Unit with Lentiviral Pol II miR RNAi Expression System	18
293FT Cell Line	22
Gateway [®] Recombination Reactions	23
Green Fluorescent Protein	26
Biosafety Features of the Lentiviral System	28
Experimental Outline	30
Methods	
General Information	35
Cloning miRNA	
Designing the Single-Stranded DNA Oligos	37
Generating the Double-Stranded Oligo	43
Checking the Integrity of the ds Oligo	47
Performing the Ligation Reaction	51
Transforming One Shot [®] TOP10 Competent <i>E. coli</i>	53
Analyzing Transformants	54
Chaining pre-miRNAs	57
Removing EmGFP Coding Sequence	59
Transfection and RNAi Analysis	60
Plasmid DNA Purification	61
Transfecting Cells	64
Detecting Fluorescence	69
Generating a Stable Cell Line	70
Expected Results for miRNA Expression	72
Transferring the Pre-miRNA Expression Cassette to Destination Vectors	
Creating Entry Clones for Use with Destination Vectors	75
Performing the Rapid BP/LR Recombination Reaction	
Transforming One Shot [®] Stbl3 [™] Competent <i>E. coli</i>	
Analyzing Transformants	89

Table of Contents, Continued

Lentivirus Production, Transduction, and RNAi Analysis	92
Lentiviral Plasmid DNA Purification	93
Growth and Maintenance of the 293 FT Cell Line	96
Producing Lentivirus in 293FT Cells	
Titering Your Lentiviral Stock	
Transduction and Analysis	114
Expected Results for Lentiviral Expression	119
Troubleshooting	
Appendix I	
Recipes	
Blasticidin	
Performing the Standard BP Recombination Reaction	
Performing the Standard LR Recombination Reaction	
Map of pcDNA [™] 6.2-GW/EmGFP-miR	143
Map of pcDNA [™] 6.2-GW/EmGFP-miR-neg control plasmid	144
Features of pcDNA™6.2-GW/EmGFP-miR	
Map of pcDNA [™] 1.2/V5-GW/ <i>lacZ</i>	
Map and Features of pLenti6/V5-DEST	147
Map of pLenti6/V5-GW/lacZ	
Map and Features of pDONR [™] 221	
Map and Features of pLP1	
Map and Features of pLP2	
Map and Features of pLP/VSVG	
Map of pCMVSPORT6TAg.neo	
Map of pENTR [™] -gus	
Appendix II	
Accessory Products	
Technical Service	
Product Qualification	
Purchaser Notification	
Gateway [®] Clone Distribution Policy	
References	

Experienced Users Procedure to Generate Expression Clone

Introduction A quick procedure t

A quick procedure to generate an expression clone is provided for experienced users of the BioModule[™] BLOCK-iT[™] RNAi Kits. If you are performing this procedure for the first time or for additional protocols, refer to this manual.

Step	Action			
Design single-stranded DNA oligos	Follow the guidelines on page 37 to design single-stranded DNA oligos encoding the pre-miRNA of interest.			
Anneal the single-	1. Set up the following annealing reaction	· ·		
stranded oligos to	200 µM top strand oligo	5 µl		
generate a ds oligo	200 μM bottom strand oligo	5 µl		
	10X Oligo Annealing Buffer	2 µl		
	DNase/RNase-free water	<u>8 µl</u>		
	Total volume	20 µl		
	2. Heat the reaction mixture at 95°C for 4	minutes.		
	3. Remove the sample and set on the labo cool to room temperature for 5-10 min			
	4. Spin down the sample in a microcentri	fuge for 5 seconds. Mix gently.		
	5. Dilute the ds oligo mixture 5,000-fold b 100-fold and 50-fold dilutions: the first the second into 1X Oligo Annealing Bu	into DNase/RNase-free water and		
Clone the ds oligo into	1. Set up the following ligation reaction.			
pcDNA [™] 6.2-GW/EmGFP- miR Vector	5X Ligation Buffer	4 µl		
IIIIK VECIOI	pcDNA [™] 6.2-GW/EmGFP-miR (5 ng/µ	l), linearized 2 μl		
	ds oligo (10 nM; 1:5,000 dilution)	4 µl		
	DNase/RNase-Free water	9 µl		
	<u>T4 DNA Ligase (1 U/μl)</u>	<u> </u>		
	Total volume	20 µl		
	2. Mix reaction well and incubate for 5 m	inutes at room temperature.		
	3. Place reaction on ice and proceed to tra	insform <i>E. coli,</i> below.		
Transform One Shot [®] TOP10 Chemically	 Add 2 μl of the ligation reaction into a Competent <i>E. coli</i> and mix gently. 	Add 2 μ l of the ligation reaction into a vial of One Shot [®] TOP10 Chemically Competent <i>E. coli</i> and mix gently.		
Competent E. coli	2. Incubate on ice for 5-30 minutes.			
	3. Heat-shock the cells for 30 seconds at 42°C without shaking. Immediat transfer the tube to ice.			
	. Add 250 μl of room temperature S.O.C. Medium.			
	5. Incubate at 37°C for 1 hour with shakir	ng.		
	 Spread 20-100 μl of bacterial culture on containing 50 μg/ml spectinomycin an 	· · ·		

vectors, and lentiviral expression, see detailed protocols in this manual.

Kit Contents and Storage

Types of Kits

This manual is supplied with the following products.

Product	Catalog no.
BioModule [™] BLOCK-iT [™] Unit with Pol II miR RNAi	WFGE07
Expression Vector	
BioModule [™] BLOCK-iT [™] Unit with Lentiviral Pol II	WFGE08
miR RNAi Expression System	

Kit ComponentsThe BioModule™ BLOCK-iT™ Unit with Pol II miR RNAi Expression Vector and
BioModule™ BLOCK-iT™ Unit with Lentiviral Pol II miR RNAi Expression
System (referred to as BioModule™ BLOCK-iT™ RNAi Units) include the
following components. For detailed contents, see the following pages.

The BioModule[™] BLOCK-iT[™] RNAi Units are shipped as described below. Upon receipt, store each item as detailed below.

Box	Component	Catalog no.		Shipping	Storage
		WFGE07	WFGE08		
1	BLOCK-iT [™] Pol II miR RNAi Expression Vector Kit	\checkmark	\checkmark	Dry ice	-20°C
2	One Shot [®] TOP10 Chemically Competent <i>E. coli</i>	\checkmark	\checkmark	Dry ice	-80°C
3	PureLink™ HQ Mini Plasmid DNA Purification Kit	\checkmark	\checkmark	Room temperature	Room temperature
4	Lipofectamine [™] 2000 Reagent	\checkmark		Blue ice	4°C (do not freeze)
5	OptiMEM [®] I Reduced Serum Medium	\checkmark	\checkmark	Room temperature	2 to 8°C (in the dark)
6	Spectinomycin	\checkmark	\checkmark	Dry ice	-20°C
7	Blasticidin	\checkmark		-20°C	-20°C
8	Kanamycin	\checkmark	\checkmark	Dry ice	-20°C
9	Ampicillin		\checkmark	Room temperature	4°C
10	LB Broth (1X), liquid	\checkmark	\checkmark	Room temperature	Room temperature
11	4% E-Gel [®] Starter Pak	\checkmark	\checkmark	Room temperature	Room temperature
12	1.2% E-Gel [®] Agarose Gels		\checkmark	Room temperature	Room temperature
13	TrackIt [™] 10 bp DNA Ladder	\checkmark	\checkmark	Room temperature	Room temperature

Kit Components, continued

Box	Component	Catalo	og no.	<u>Shipping</u>	<u>Storage</u>
		<u>WFGE07</u>	<u>WFGE08</u>		
14	TrackIt [™] 10 bp DNA 1 Kb Plus DNA Ladder		\checkmark	Room temperature	Room temperature
15	pLenti6/V5-DEST Gateway® Vector Kit		\checkmark	Blue ice	-20°C
16	pDONR™221		\checkmark	Room temperature	-20°C
17	Gateway [®] BP Clonase™ II Enzyme Mix		\checkmark	Dry ice	-20°C
18	Gateway [®] LR Clonase [™] II Enzyme Mix		\checkmark	Dry ice	-20°C
19	One Shot [®] Stbl3 [™] Chemically Competent <i>E. coli</i>		\checkmark	Dry ice	-80°C
20-21	ViraPower [™] Bsd Lentiviral Support Kit:		\checkmark		
	ViraPower [™] Packaging Mix			Blue ice	-20°C
	Lipofectamine [™] 2000 Reagent			Blue ice	+4°C (do not freeze)
	Blasticidin			Room temperature	-20°C
22	293FT Cell Line		\checkmark	Dry ice	Liquid nitrogen
23	S.N.A.P.™ MidiPrep Kit		\checkmark	Room temperature	Room temperature

BLOCK-iT[™] Pol II miR RNAi Expression Vector Kit

The following reagents are included with the BLOCK-iT[™] Pol II miR RNAi Expression Vector Kit (Box 1). **Store the kit at -20°C.**

Reagent	Composition	Amount
pcDNA [™] 6.2-GW/EmGFP-	5 ng/μl in:	4 x 10 μl
miR, linearized	10 mM Tris-HCl, pH 8.0	
	1 mM EDTA, pH 8.0	
10X Oligo Annealing Buffer	100 mM Tris-HCl, pH 8.0	250 µl
	10 mM EDTA, pH 8.0	
	1 M NaCl	
DNase/RNase-Free Water		3 x 1.5 ml
5X Ligation Buffer	250 mM Tris-HCl, pH 7.6	80 µl
	50 mM MgCl ₂	
	5 mM ATP	
	5 mM DTT	
	25% (w/v) polyethylene glycol-8000	
T4 DNA Ligase	1 (Weiss) U/μl in	20 µl
	10 mM Tris-HCl, pH 7.5	
	50 mM KCl	
	1 mM DTT	
	50% (v/v) glycerol	
EmGFP forward sequencing primer	100 ng/μl in TE Buffer, pH 8.0	20 µl
miRNA reverse sequencing primer	100 ng/μl in TE Buffer, pH 8.0	20 µl
miR- <i>lacZ</i> positive double- stranded (ds) control oligo	10 nM in 1X Oligo Annealing Buffer	50 µl
pcDNA [™] 1.2/V5-GW/lacZ control plasmid	500 ng/μl in TE Buffer, pH 8.0	20 µl
pcDNA [™] 6.2-GW/EmGFP- miR-neg control plasmid	500 ng/μl in TE Buffer, pH 8.0	20 µl

Unit Definition of T4 DNA Ligase Primer Sequences	One (Weiss) unit of T4 DNA Ligase catalyzes the exchange of 1 nmol ³² P-labeled pyrophosphate into $[\gamma/\beta^{-32}P]$ ATP in 20 minutes at 37°C (Weiss <i>et al.</i> , 1968). One unit is equal to approximately 300 cohesive-end ligation units. The table below provides the sequence and the amount of primers included in the kit.		
	Primer	Sequence	Amount
	EmGFP forward sequencing primer	5'-ggcatggacgagctgtacaa-3'	2 μg (323 pmol)
	miRNA reverse sequencing primer	5'-CTCTAGATCAACCACTTTGT-3'	2 μg (332 pmol)
<i>LacZ</i> Control Oligo Sequences	<i>lacZ</i> top and bottom str	niR- <i>lacZ</i> positive ds control oligo are lister rand oligos are annealed and are supplie d oligo that is ready-to-use in the ligation	ed in the kit as a

51).

 LacZ DNA Oligo
 Sequence

 Top strand
 5' - TGCTGAAATCGCTGATTTGTGTAGTCGTTTTGGCCACTGACGACTACACAATCAGCGATTT-3'

 Bottom strand
 5' - CCTGAAATCGCTGATGTGTGTGTGTGTCGTCAGTCGGCCAAAACGACTACACAAATCAGCGATTTC-3'

One Shot[®] TOP10 Reagents

The following reagents are included in the One Shot[®] TOP10 Chemically Competent *E. coli* kit (Box 2). Transformation efficiency is $\ge 1 \times 10^9$ cfu/µg plasmid DNA. **Store Box 2 at -80°C.**

Reagent	Composition	Amount
S.O.C. Medium	2% Tryptone	6 ml
(may be stored at +4°C or	0.5% Yeast Extract	
room temperature)	10 mM NaCl	
	2.5 mM KCl	
	10 mM MgCl ₂	
	10 mM MgSO ₄	
	20 mM glucose	
TOP10 cells		21 x 50 µl
pUC19 Control DNA	10 pg/µl in 5 mM Tris-HCl, 0.5 mM EDTA, pH 8	50 µl

F⁻ mcrA ∆(mrr-hsdRMS-mcrBC) ϕ 80lacZ∆M15 ∆lacX74 recA1 araD139 ∆(ara-leu)7697 galU galK rpsL (Str^R) endA1 nupG

PureLink[™] HQ Mini Plasmid Purification Kit

The components included in the PureLink[™] HQ Mini Plasmid Purification Kit (Box 3) are listed below.

Upon receipt, store all components at room temperature, except store the Resuspension Buffer after addition of RNase at +4°C.

Component	Amount
Resuspension Buffer	120 ml
Lysis Buffer	120 ml
Neutralization/Binding Buffer	2×85 ml
Wash Buffer	16 ml
Elution Buffer (10 mM Tris-HCl, pH 8.5)	15 ml
RNase A	12 mg
Spin Columns	100
Collection Tubes (2 ml)	100
Elution Tubes (1.7 ml)	100

Lipofectamine[™] 2000 Reagent and Opti-MEM[®] I Reduced Serum Medium **Lipofectamine**[™] **2000 Reagent** (Box 4) is a transfection reagent with proprietary formulation. **Store at 4°C. Do not freeze.**

Note: 0.75 ml Lipofectamine[™] 2000 Reagent is included in the ViraPower[™] Bsd Lentiviral Support Kit for Cat. no. WFGE08.

Opti-MEM® I Reduced Serum Medium (Box 5) is used for dilution of the lipid: DNA complexes during transfection. 100 ml Opti-MEM® I Reduced Serum Medium is included with each BioModule[™] BLOCK-iT[™] RNAi Unit. The formulation of the medium is described below.

Opti-MEM[®] I Reduced Serum Medium is a modification of Eagle's Minimal Essential Medium, buffered with HEPES and sodium bicarbonate, and supplemented with hypoxanthine, thymidine, sodium pyruvate, L-glutamine or GLUTAMAX, trace elements and growth factors. The protein level is minimal (15 µg/ml); insulin and transferrin are the only protein supplements. Phenol red is included at a reduced concentration as a pH indicator.

LB Broth and Antibiotics

The following antibiotics and medium are included with the BioModuleTM BLOCK-iTTM RNAi Units (Boxes 6-10).

Antibiotic	Composition	Amount
Spectinomycin	100 mg/ml in deionized water	1 ml
Kanamycin (only with WFGE08)	25 mg/ml in deionized water	1 ml
Ampicillin (only with WFGE08)	Powder	200 mg
Blasticidin	Powder	50 mg
LB Broth (1X), liquid	Formulation per one liter: 10 g SELECT Peptone 140 5 g SELECT Yeast Extract 5 g sodium chloride	500 ml

E-Gel[®] Agarose Gels and DNA Ladders

The following E-Gel[®] agarose gels and DNA Ladders are included with the BioModule[™] BLOCK-iT[™] RNAi Units (Boxes 11-14). **Store at room temperature.**

Item	Composition	Amount
E-Gel [®] 4% Starter Pak	See page 15 for details	6 gels and one E-Gel [®] PowerBase™
E-Gel [®] 1.2% 18 Pak (only with WFGE08)	See page 15 for details	18 gels
TrackIt [™] 10 bp DNA Ladder	0.5 μg/μl in:	20 applications
	10 mM Tris-HCl, pH 7.5	
	10 mM EDTA, pH 8.0	
	0.06% XCFF	
	0.4% Orange G	
	5% glycerol	
TrackIt [™] 1 Kb Plus DNA Ladder (only with WFGE08)	0.1 μg/μl in: 10 mM Tris-HCl, pH 7.5 10 mM EDTA, pH 8.0 0.06% XCFF 0.6% tartrazine 5% glycerol 5 mM NaCl	100 applications

Vectors

The following vectors are included with the BioModule[™] BLOCK-iT[™] RNAi Units (Boxes 15 and 16). For details on pcDNA[™]6.2-GW/EmGFP-miR vector, see page viii. **Store the vectors at -20°C.**

Reagent	Composition	Amount
pLenti6/V5-DEST Gateway® Vector	Lyophilized in TE Buffer, pH 8.0	6 µg
pLenti6/V5-GW/lacZ control	Lyophilized in TE Buffer, pH 8.0	10 µg
pDONR [™] 221 Vector	Lyophilized in TE Buffer, pH 8.0	6 µg

Gateway[®] Clonase[™] The following reagents are included with the Gateway[®] Clonase[™] II Enzyme Mixes (Boxes 17-18).

Store Box 17 and 18 at -20°C for up to 6 months. For long-term storage, store at -80°C.

BP Clonase[™] II

Reagent	Composition	Amount
Gateway [®] BP Clonase [™] II Enzyme Mix	Proprietary	40 µl
Proteinase K Solution	$2 \mu g/\mu l$ in:	40 µl
	10 mM Tris-HCl, pH 7.5	
	20 mM CaCl ₂	
	50% glycerol	
PEG Solution	30% PEG 8000	1 ml
	30 mM MgCl ₂	
pEXP7-tet Positive Control	50 ng/μl in TE Buffer, pH 8.0	20 µl

LR Clonase[™] II

Reagent	Composition	Amount
Gateway [®] LR Clonase [™] II Enzyme Mix	Proprietary	40 µl
Proteinase K Solution	$2 \mu g/\mu l$ in:	40 µl
	10 mM Tris-HCl, pH 7.5	
	20 mM CaCl ₂	
	50% glycerol	
pENTR [™] -gus Positive Control	50 ng/μl in TE Buffer, pH 8.0	20 µl

Note: The pENTR[™]-gus control included with the LR Clonase[™] II Enzyme Mix may be used as a positive control for the LR recombination reaction **only** (page 139).

One Shot[®] Stbl3[™] Chemically Competent *E. coli*

The following reagents are included with the One Shot[®] Stbl3[™] Chemically Competent *E. coli* kit (Box 19). Transformation efficiency is $\geq 1 \times 10^8$ cfu/µg plasmid DNA. **Store Box 19 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 ₂	
	10 mM MgSO ₄	
	20 mM glucose	
Stbl3 [™] Cells		21 x 50 μl
pUC19 Control DNA	10 pg/μl in 5 mM Tris-HCl, 0.5 mM EDTA, pH 8	50 µl

Genotype of Stbl3[™] Cells

F⁻ mcrB mrr hsdS20(r_B⁻, m_B⁻) recA13 supE44 ara-14 galK2 lacY1 proA2 rpsL20(Str^R) xyl-5 λ ⁻ leu mtl-1

ViraPower[™] Bsd Lentiviral Support Kit

The following reagents are included with the ViraPower[™] Bsd Lentiviral Support Kit (Boxes 20-21). Store the ViraPower[™] Packaging Mix and Blasticidin at -20°C. Store Lipofectamine[™] 2000 Reagent at +4°C.

Important: Do not freeze Lipofectamine[™] 2000 Reagent.

Reagent	Composition	Amount
ViraPower [™] Packaging Mix	Contains a mixture of the pLP1, pLP2, and pLP/VSVG plasmids, lyophilized in TE Buffer, pH 8.0	195 µg
Lipofectamine [™] 2000 Reagent	Proprietary	0.75 ml
Blasticidin	Powder	50 mg

293FT Cell Line The 293FT Cell Line (Box 22) is used for the production of lentiviral stocks. The 293FT Cell Line is supplied as one vial containing 3 x 10⁶ frozen cells in 1 ml of Freezing Medium. Upon receipt, store in liquid nitrogen until use. For instructions to thaw, culture, and maintain the 293FT Cell Line, see page 96. S.N.A.P.[™] The following items are included in the S.N.A.P.[™] MidiPrep Kit (Box 23). Store the kit at room temperature, except store the following components at +4°C after opening the kit: Resuspension Buffer

- Precipitation Salt
- RNase A

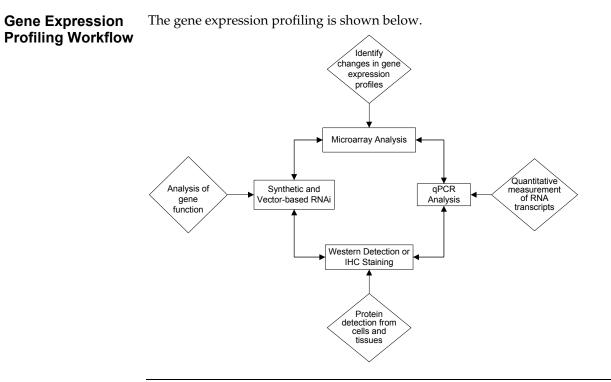
Item	Composition	Quantity
S.N.A.P. [™] MidiPrep		
Column A (Filtering)		20
Column B (Binding)		20
Resuspension Buffer	50 mM Tris-HCl, pH 8.0, 10 mM EDTA	80 ml
RNase A, lyophilized		6 mg
Lysis Buffer	0.2 M NaOH, 1% SDS	80 ml
Precipitation Salt	3 M Potassium acetate, pH 5.2	80 ml
Binding Buffer	7.5 M Guanidine-HCl	2 x 120 ml
Wash Buffer	5 M Guanidine-HCl, 50 mM MOPS, pH 7.0	100 ml
4X Final Wash	400 mM NaCl	80 ml



- Some reagents in the units may be provided in excess of the amount needed.
- Individual documentations detailing general use are included with some of the products supplied in the BioModule[™] BLOCK-iT[™] RNAi Units. To use the products specifically with the BioModule[™] BLOCK-iT[™] RNAi Units, follow the recommended protocols in this manual.

Introduction

Overview	
Introduction	The BioModule [™] BLOCK-iT [™] Unit with Pol II miR RNAi Expression Vector (Cat. no. WFGE07) facilitates the expression of microRNA (miRNA) for use in RNA interference (RNAi) analysis of a target gene in mammalian cells. The kit includes a Gateway [®] -adapted expression vector designed for transient or stable expression of your miRNA and allows for efficient recombination with suitable destination vectors allowing tissue-specific, regulated, or lentiviral expression of the miRNA in mammalian cells using Gateway [®] Technology.
	The BioModule [™] BLOCK-iT [™] Unit with Lentiviral Pol II miR RNAi Expression System (Cat. no WFGE08) combines Invitrogen's BLOCK-iT [™] Pol II miR RNAi and ViraPower [™] Lentiviral technologies to facilitate creation of a replication- incompetent lentivirus that delivers a miRNA sequence of interest to dividing or non-dividing mammalian cells for RNAi analysis.
	The BioModule [™] BLOCK-iT [™] Unit with Pol II miR RNAi Expression Vector and BioModule [™] BLOCK-iT [™] Unit with Lentiviral Pol II miR RNAi Expression System (referred to as BioModule [™] BLOCK-iT [™] RNAi Units) provide qualified reagents and validated protocols to express your miRNA in mammalian cells for RNAi analysis.
	For details on the BLOCK-iT [™] Pol II miR RNAi, Gateway [®] , and ViraPower [™] Lentiviral technologies, see page 5.
BioModule [™] Units for Gene Expression Profiling	The BioModule [™] Transfection and Control Unit with BLOCK-iT [™] Technology is one of the several BioModule [™] Units available from Invitrogen (page 160) for gene expression profiling. Each of the BioModule [™] Units for gene expression profiling includes high-quality reagents and validated protocols with relevant controls for each step of the workflow (see next page). Each unit is designed to provide an integrated workflow that allows you to perform various steps seamlessly during expression analysis.
	Gene expression profiling comprises multiple steps employing various technologies such as microarray analysis or quantitative PCR (qPCR) for analysis at the nucleic acid level; western immunodetection and immunohistochemistry for analysis at the protein level; and RNAi for functional analysis.
	Continued on next page



System Components	The BioModule [™] BLOCK-iT [™] Unit with Pol II miR RNAi Expression Vector includes:
	• pcDNA [™] 6.2-GW/EmGFP-miR expression vector for production of an expression clone containing a double-stranded oligonucleotide (ds oligo) encoding a pre-miRNA sequence for expression in mammalian cells using an RNA Polymerase II (Pol II) promoter, the human cytomegalovirus (CMV) immediate early promoter.
	• Reagents for production of an expression clone containing a double- stranded oligonucleotide (ds oligo) encoding a pre-miRNA (need to order oligos separately).
	• Positive and negative controls for the generation and use of the expression clone.
	• Reagents for plasmid DNA purification and Lipofectamine [™] 2000 Reagent for efficient transfection into mammalian cells.
	The BioModule ^{m} BLOCK-iT ^{m} Unit with Lentiviral Pol II miR RNAi Expression System includes:
	• A BioModule [™] BLOCK-iT [™] Unit with Pol II miR RNAi Expression Vector for production of an expression clone containing a double-stranded oligonucleotide (ds oligo) encoding a pre-miRNA sequence (see above).
	 The pDONR[™]221 vector is used as an intermediate to transfer the pre- miRNA expression cassette (see page 11 for details) into the lentiviral expression plasmid (see below) using Gateway[®] Technology.
	• A pLenti6/V5-DEST destination vector into which the pre-miRNA cassette from the expression clone is transferred using Gateway [®] 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.
	 Gateway[®] BP and LR Clonase[™] II Enzyme Mixes that facilitate the transfer of the pre-miRNA expression cassette from the expression vector into the pLenti6/V5-DEST destination vector.
	 Components of the ViraPower[™] Lentiviral System for production of a replication-incompetent lentivirus that stably expresses the miRNA of interest in both dividing and non-dividing mammalian cells.
	 Reagents for lentiviral plasmid DNA purification and Lipofectamine[™] 2000 Reagent for efficient transfection into mammalian cells.
	For more information about the BLOCK-iT [™] Pol II miR RNAi Technology, ViraPower [™] Lentiviral Technology, and Gateway [®] Technology, see page 5.
	For details on system components, see page 12.

Advantages of the BioModule [™]	Use of the BioModule [™] BLOCK-iT [™] RNAi Units for vector-based expression of miRNA provides the following advantages:
BLOCK-iT [™] RNAi Units	• The BLOCK-iT [™] Pol II miR RNAi Expression Vector provides a rapid and efficient way to clone a ds oligo duplex encoding a desired miRNA target sequence into a vector containing a Pol II promoter for use in RNAi analysis.
	• Gateway [®] -adapted vector for easy transfer of the miRNA of interest from one expression vector (pcDNA [™] 6.2-GW/EmGFP-miR) into other destination vectors.
	• Enables co-cistronic expression of the pre-miRNA of interest with EmGFP to allow visualization of cells expressing the miRNA.
	• Efficiently delivers the miRNA of interest to mammalian cells in culture or <i>in vivo</i> .
	Lentiviral System
	• Generates a replication-incompetent lentivirus that effectively transduces both dividing and non-dividing mammalian cells, thus broadening the potential of RNAi applications beyond those of other traditional retroviral systems (Naldini, 1998).
	• 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 [™] Pol II miR RNAi Technology	The BLOCK-iT [™] 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 [™] Pol II miR RNAi Expression Vector is specially 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 10 for details).
	A variety of BLOCK-iT [™] RNAi products are available from Invitrogen to facilitate RNAi analysis in mammalian and invertebrate systems. For more information about any of the BLOCK-iT [™] RNAi products, see the RNAi Central application portal at www.invitrogen.com/rnai or contact Technical Service (page 163).

ViraPower [™] Lentiviral Technology	The ViraPower [™] 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 [™] system developed by Cell Genesys (Dull <i>et al.</i> , 1998), the ViraPower [™] Lentiviral Technology possesses features which enhance its biosafety while allowing highlevel expression in a wider range of cell types than traditional retroviral systems. The main components of the ViraPower [™] Lentiviral Expression System include:
	• A pLenti-based expression vector (<i>e.g.</i> pLenti6/V5-DEST) into which the sequence of interest is 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.
	• The ViraPower [™] Packaging Mix, an optimized mixture of the three packaging plasmids required for production of the lentivirus.
	• An optimized 293FT producer cell line to facilitate optimal production of virus.
	For more information about these components, see page 12. For more information about the biosafety features of the system, see page 28.
Gateway [®] Technology	The Gateway [®] Technology is a universal cloning method that takes advantage of the site-specific recombination properties of bacteriophage lambda (Landy, 1989) to provide a rapid and highly efficient way to move your DNA sequence of interest into multiple vector systems.
	To express your miRNA of interest in mammalian cells using the BioModule™ BLOCK-iT™ RNAi Units and Gateway® Technology, simply:
	1. Clone a double-stranded oligonucleotide encoding your miRNA sequence of interest into the pcDNA [™] 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.
	 To transfer your pre-miRNA expression cassette into a destination vector including pLenti6/V5-DEST vector, generate an entry clone by performing a BP recombination reaction between the pcDNA[™]6.2-GW/EmGFP-miR expression clone and pDONR[™]221 donor vector.
	 Then perform an LR recombination reaction between the resulting entry clone (pENTR[™]221/miR) and any destination vector including pLenti6/V5-DEST. See page 80 for more details.
	For detailed information about the Gateway [®] Technology, refer to the Gateway [®] Technology with Clonase [™] II manual which is available from our web site (www.invitrogen.com) or by contacting Technical Service (page 163).
	Continued on next nage

Purpose of this Manual	This manual provides an overview of the pathway by which miRNA facilitates gene knockdown in mammalian cells and includes instructions and guidelines to:
	• Design the appropriate single-stranded oligonucleotides representing the target gene.
	• Clone the ds oligo into pcDNA [™] 6.2-GW/EmGFP-miR expression vector.
	• Purify high-quality plasmid DNA for mammalian transfection.
	• Transfect your pcDNA [™] 6.2-GW/EmGFP-miR construct into mammalian cells for transient RNAi analysis or to generate stable cell lines.
	• Perform a Rapid BP/LR recombination reaction with a destination vector of choice to generate an expression clone containing the pre-miRNA sequence of interest.
	Lentiviral expression
	• Co-transfect the pLenti6/V5-GW/EmGFP-miR expression construct and the ViraPower [™] Packaging Mix into the 293FT Cell Line to produce a lentiviral stock.
	• Titer the lentiviral stock.
	 Transduce the lentiviral construct into mammalian cells and perform "transient" RNAi analysis
	• Generate a stably transduced cell line, if desired.
	Sections on troubleshooting and examples of expected results are also included.

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). RNAi 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</i> <i>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).

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 with protein production without directly inducing mRNA degradation (Ambros, 2004). Nonetheless, these miRNAs are found associated with the RNAi nuclease AGO2 (Liu <i>et al.</i> , 2004; Meister <i>et al.</i> , 2004), and at least two miRNAs with close matches to their target sequences, particularly in their 5' regions, have been shown to cleave cognate mRNAs (Yekta <i>et al.</i> , 2004; Yu <i>et al.</i> , 2005).
	The engineered miRNAs produced by the BLOCK-iT TM Pol II miR RNAi Expression Vector Kit (see below) fully complements their target site and cleave the target mRNA. Sequence analysis showed that the primary cleavage site at the phosphodiester bond in the mRNA is found opposite the tenth and eleventh bases 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	Use of siRNA (diced siRNA or synthetic siRNA) for RNAi analysis in mammalian cells is limited by their transient nature. 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 [®] -adapted expression vectors that enable the expression of engineered miRNA sequences from Pol II promoters. The pcDNA6.2 [™] -GW/EmGFP-miR expression vector (supplied in the BioModule [™] Unit with BLOCK-iT [™] Pol II miR RNAi Expression Vector) facilitates the generation of an expression clone containing a ds oligo encoding a pre-miRNA sequence (page 10). 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 [®] recombination reactions (page 23).

miRNA Vector	The pcDNA [™] 6.2-GW/EmGFP-miR Expression Vector included with the unit 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. 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	Using miRNA vector-based systems that use Pol II promoters for RNAi cassette expression offer the following advantages over traditional siRNA or shRNA expression:
	 Enables co-cistronic expression of reporter genes such as GFP (see above) allowing reliable tracking of miRNA expression in mammalian cells
	• Allows expression of miRNA from a variety of promoters, including tissue- specific and regulated promoters for <i>in vivo</i> experiments
	• Enables expression of multiple miRNAs from a single transcript allowing the knockdown of more than one gene simultaneously
	• Permits design of predictable RNAi constructs with a high rate of success
Human CMV Promoter	The pcDNA [™] 6.2-GW/EmGFP-miR expression vector contains 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
	Note: 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.

Design of the Engineered PremiRNA

The engineered pre-miRNA sequence structure is based on the murine miR-155 sequence (Lagos-Quintana *et al.*, 2002). The 5' and 3' flanking regions derived from the miR-155 transcript were inserted in the vector to preserve as much as possible of the miR-155 structure. We optimized the stem-loop structure and a 2 nucleotide internal loop results in higher knockdown rate than the 5 nucleotide/3 nucleotide internal loop found in native miR-155 molecule. An *Msc* I site was incorporated in the terminal loop to aid in sequence analysis. Below the changes are shown made to the native miR-155 to form an engineered pre-miRNA directed against *lacZ* (targeting sequence in bold).

native miR-155		optim	<u>Msc</u> l		
5′-UG	UGUGA	UUGGCC	UG	UU	UUGGCC
CUG UUAAU GACAAUUA 3 '- G^		: : U —	→ !!!!!!	JCGCUGAU GUGUAGU	: A
	internal loop	terminal loop		internal loop	terminal loop

Structure of the Engineered premiRNA The pcDNA[™]6.2-GW/EmGFP-miR expression vector is 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, see above.

For optimized knockdown results, we recommend that the ds oligo encoding the engineered pre-miRNA have the following structural features:

- Two 4 nucleotide, 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
- A short sense target sequence with 2 nucleotides removed ($\Delta 2$) to create an internal loop

The structural features are depicted in the figure below.



For more details on the structure and guidelines to design the oligonucleotides, refer to page 37.

noted as
iteway® ed
anking
he r stem- then miRNA.
nucleotide
s driven s to stronic

BioModule[™] BLOCK-iT[™] Unit with Pol II miR RNAi Expression Vector

Description of the System	 The BioModule[™] BLOCK-iT[™] Unit with Pol II miR RNAi Expression Vector facilitates the generation of an expression construct that permits high-level expression of a pre-miRNA in mammalian cells for RNAi analysis of a target gene. The kit contains the following major components: The pcDNA[™] 6.2-GW/EmGFP-miR linearized plasmid into which a ds oligo encoding the pre-miRNA is cloned to generate an expression clone that contains the elements required for expression of the miRNA in mammalian cells. The pcDNA[™] 6.2-GW/EmGFP-miR vector is supplied linearized with 4-nucleotide 5′ overhangs on each strand to facilitate directional cloning of the ds oligo insert. The resulting expression clone containing the pre-miRNA expression cassette may be transfected into mammalian cells for transient or stable RNAi analysis, or used to transfer the pre-miRNA expression cassette into a suitable destination vector using Gateway[®] Technology. 			
	• T4 DNA Ligase and an optimized ligation buffer to allow 5-minute room temperature ligation of the ds oligo insert into the pcDNA [™] 6.2-GW/EmGFP-miR vector.			
	 One Shot[®] TOP10 Chemically Competent <i>E. coli</i> for high efficiency transformation of the ligation reaction. 			
	• PureLink [™] HQ Mini Plasmid Purification Kit for high-quality plasmid DNA purification suitable for transfection into mammalian cells.			
	• Lipofectamine [™] 2000 Reagent and Opti-MEM [®] Reduced Serum Medium for efficient transfection of your construct into mammalian cells.			
	• E-Gel [®] agarose gels and DNA ladders for checking integrity of the ds oligo by gel electrophoresis.			
Controls	The BioModule [™] BLOCK-iT [™] Unit with Pol II miR RNAi Expression Vector also includes a negative control plasmid and a ds positive control oligo.			
	 The pcDNA[™]6.2-GW/EmGFP-miR-neg control plasmid contains an insert that can form a hairpin structure that is processed into mature miRNA, but is predicted not to target any known vertebrate gene. Thus, this plasmid serves as a suitable negative control for pre-miRNA experiments with pcDNA[™]6.2-GW/EmGFP-miR expression vector. The negative control sequence without 5' overhangs is shown below (page 144 for a map): 5' -GAAATGTACTGCGCGTGGAGACGTTTTGGCCACTGACTGA			
	 The miR-lacZ positive double-stranded (ds) control oligo serves as a positive control during the miRNA expression vector generation. Use this oligo to generate a pcDNA[™]6.2-GW/EmGFP-miR-lacZ expression clone. 			
	 Co-transfecting the resulting pcDNA[™]6.2-GW/EmGFP-miR-<i>lacZ</i> expression clone targeting the <i>lacZ</i> gene and the pcDNA[™]1.2/V5-GW/<i>lacZ</i> reporter plasmid supplied with the kit into mammalian cells provide a means to assess the RNAi response in your cell line by assaying for knockdown of β-galactosidase. 			

BioModule[™] BLOCK-iT[™] Unit with Pol II miR RNAi Expression

Vector, Continued

Generating an miRNA Expression Vector Using the Kit		ing the reagents supplied in the unit, you will perform the following steps to nerate an expression clone in pcDNA™6.2-GW/EmGFP-miR:		
	1.	 Design and synthesize two complementary single-stranded DNA oligonucleotides, with one encoding the miRNA of interest. 		
	2.	Anneal the single-stranded oligonucleotides to generate a double-stranded oligo (ds oligo).		
	3.	Clone the ds oligo into the linearized pcDNA [™] 6.2-GW/EmGFP-miR vector.		
	4.	Transform the ligation reaction into One Shot [®] TOP10 Chemically Competent <i>E. coli</i> and select for spectinomycin-resistant transformants.		
	5.	Use the pcDNA [™] 6.2-GW/EmGFP-miR expression construct for transient RNAi analysis in mammalian cells, isolate stable cell lines expressing the miRNA, or perform a Gateway [®] recombination reaction with a suitable Gateway [®] destination vector to generate a different expression clone.		
Features of the pcDNA [™] 6.2-	The •	e pcDNA [™] 6.2-GW/EmGFP-miR Vector contains the following features: Human CMV promoter for high-level, constitutive expression of the miRNA		
GW/EmGFP-miR Vector		from a RNA Polymerase II-dependent promoter		
Vector	•	5' and 3' miR flanking regions for formation of an engineered pre-miRNA		
	•	Cloning site containing 4-nucleotide 5' overhangs on each DNA strand for directional cloning of the ds oligo encoding the pre-miRNA of interest		
	•	Two recombination sites, <i>att</i> B1 and <i>att</i> B2 sites, flanking the pre-miRNA expression cassette for recombinational cloning of the pre-miRNA expression cassette into a Gateway [®] destination vector		
	•	Herpes Simplex virus (HSV) thymidine kinase (TK) polyadenylation signal for termination and polyadenylation of the transcript		
	•	An EmGFP coding sequence for co-cistronic expression with the pre-miRNA.		
	•	Spectinomycin resistance gene for selection in E. coli		
	•	pUC origin for high-copy maintenance of the plasmid in E. coli		
	•	Blasticidin resistance gene for selection in <i>E. coli</i> and mammalian cells to generate cell lines stably expressing the miRNA		

BioModule[™] BLOCK-iT[™] Unit with Pol II miR RNAi Expression Vector, Continued

Lipofectamine [™] 2000 Reagent	 Lipofectamine[™] 2000 is a proprietary formulation for the delivery of plasmid DNA and RNAi oligomers to mammalian cells for RNAi analysis (Gitlin <i>et al.</i>, 2002; Yu <i>et al.</i>, 2002). Using Lipofectamine[™] 2000 to transfect eukaryotic cells offers the following advantages: Highest transfection efficiency in many cell types and formats (<i>e.g.</i> 96-well).
	Refer to the Cell Lines database at <u>www.invitrogen.com</u> for a list of cell types successfully transfected.
	 Nucleic acid-Lipofectamine[™] 2000 complexes can be added directly to cells in culture medium, in the presence or absence of serum.
	• It is not necessary to remove complexes or change/add medium after transfection, but complexes may be removed after 4-6 hours.
Opti-MEM [®] I Reduced Serum	Opti-MEM [®] I Reduced Serum Medium is a versatile, chemically-defined medium used for diluting the lipid and nucleic acid during transfection.
Medium	The Opti-MEM [®] I Reduced Serum Medium is a multi-purpose medium proven to be useful in reducing serum requirements for a wide variety of cell lines and applications and has been effective in the growth and maintenance of adherent and non-adherent cell lines. When supplemented with 2-4% fetal bovine serum or alternative sera, Opti-MEM [®] I Reduced Serum Medium supports proliferative rates and maximal cell densities comparable to, and in some cases superior to, conventional media supplemented with 10% fetal bovine serum. Relatively non- fastidious cell lines may be maintained in long-term culture with even more substantial serum reduction. If using adherent cell lines and less than 2% serum supplementation or in an agitated system such as in roller bottles, the medium should be further supplemented with 100 mg/l CaCl ₂ .
	The versatility of Opti-MEM [®] I Reduced Serum Medium in the propagation of various cell types makes this medium the optimal choice for many cell culture requirements. For details on using the medium for cell culture, download the Opti-MEM [®] I Reduced Serum Medium manual from www.invitrogen.com.
PureLink [™] HQ Mini Plasmid Purification Kit	The PureLink [™] HQ Mini Plasmid Purification Kit is designed for the isolation of high-quality plasmid DNA that is suitable for restriction enzyme digestion, PCR, sequencing, bacterial cell transformation, and mammalian cell transfection. Use the kit to isolate high-quality plasmid DNA for mammalian transfection. Using the kit, plasmid DNA can be isolated from varying amounts of bacterial cells. The high quality of the isolated plasmid DNA is demonstrated by its low genomic DNA contamination, high supercoiled to nicked forms ratio, and reliable performance in demanding downstream applications such as mammalian cell transfection.
	Bacterial cells are lysed, the lysate is then neutralized and conditions are adjusted for subsequent binding. After clarification by centrifugation, the lysate is processed through the PureLink [™] spin column. The DNA binds to the silica-based membrane in the column, and impurities are removed by a single wash step. The DNA is then eluted in Elution Buffer or water.

BioModule[™] BLOCK-iT[™] Unit with Pol II miR RNAi Expression Vector, Continued

E-Gel [®] System	The E-Gel [®] agarose gel electrophoresis system is a complete bufferless system for agarose gel electrophoresis of DNA samples. Use the E-Gel [®] agarose system to verify the integrity of your ds oligos after annealing and prior to cloning the ds oligo into the pcDNA [™] 6.2-GW/EmGFP-miR Expression Vector.				
	The major components	of the E-Gel® system are	::		
	• E-Gel [®] pre-cast aga	rose gels			
	Electrophoresis bas	se			
	packaged inside a dry, an ion generating syste ethidium bromide for I	disposable, UV-transpar em (TAE buffer system),	gels that include electrodes rent cassette. Each gel contains a pH balancing system, and [®] agarose gels run in a specially se [™] v.4.		
		nnects directly to an elect	er supply in one device. The trical outlet using the adaptor		
E-Gel [®] System	E-Gel [®]				
Specifications	Cassette Dimensions:	8 cm × 10 cm x 0.6	o cm thick		
	Gel Thickness:	3 mm			
	Gel Volume:	20 ml			
	No. of Wells:	12			
	Run length:	5.8 cm			
	Sample volume:	20 µl			
	Separation Range:	100 bp-5 kb (1.2%	gel)		
		20-500 bp (4% gel)		
	Each well is 4.1 mm wi	de and the space betwee	en wells is 1 mm.		
	E-Gel [®] PowerBase [™] v.4	4			
	Dimensions:	12.5 cm x 13 cm x	13.5 cm		
	Weight:	1.19 lbs (540 g) wi			
	Safety:	UL listed and CE	-		
	Temperature:	Ambient 15°C to	40°C		
	Built-in Features:	Alarm, light LED			
	the PowerBase [™] . Use o	nly UL Listed Class 2 Di	with an adaptor included with rect Plug-in Adaptor included ed by the adaptor are shown in		
	Country	Input	Output		
	U.S. and Canada	110-120 V AC, 60 Hz	12 V DC, 880 mA		
	Europe	220-240 V AC, 50 Hz	12 V DC, 880 mA		
	·		Continued on next page		

BioModule[™] BLOCK-iT[™] Unit with Pol II miR RNAi Expression

Vector, Continued

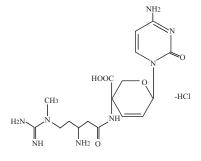
The TrackIt [™] 10 bp and 1 Kb Plus DNA Ladders are suitable for sizing double- stranded DNA fragments on agarose gels. The TrackIt [™] DNA Ladders are formulated with unique tracking dyes, Xylene Cyanol FF (XCFF), tartrazine, or Orange G, that allow you to visually track DNA migration during electrophoresis and also indicate when maximum resolution is achieved.			
The important features of the ladder are listed below:			
• 10 bp Ladder consists of 33 fragments ranging in the size of 10-200 bp and an additional fragment at 1668 bp			
• 1 Kb Plus Ladder consists of 12 bands in 1000 bp increments ranging in size from 1-12 kb, a 1650 bp fragment, and vector DNA fragments from 100-850 bp			
• 10 bp Ladder includes a 100 bp reference band that is ~2-fold brighter for easy band size determination			
• 1 Kb Plus Ladder includes an orientation doublet at 1650 bp and 2000 bp for easy reference			
Formulated with unique tracking dyes			
• Designed for use with E-Gel [®] agarose gels and TBE or TAE agarose gels			
Supplied in a ready-to-load format			
• Visualized with ethidium bromide or SYBR [®] Green staining			
Spectinomycin is an antibiotic isolated from <i>Streptomyces spectabilis</i> which inhibits protein synthesis (elongation) by interfering with peptidyl tRNA translocation in prokaryotic cells. Resistance is conferred by expression of the adenylyltransferase <i>aad</i> A gene that prevents binding of Spectinomycin. The formula for Spectinomycin (dihydrochloride pentahydrate) is $C_{14}H_{24}N_2O_7.2HCl.5H_2O$ and the molecular weight is 495.35. Recommended concentration for selection in bacteria is 50 µg/ml.			

BioModule[™] BLOCK-iT[™] Unit with Pol II miR RNAi Expression Vector, Continued

Blasticidin

Blasticidin S HCl is a nucleoside antibiotic isolated from *Streptomyces griseochromogenes* which inhibits protein synthesis in both prokaryotic and eukaryotic cells (Takeuchi *et al.*, 1958; Yamaguchi *et al.*, 1965). Resistance is conferred by expression of either one of two Blasticidin S deaminase genes: *bsd from Aspergillus terreus* (Kimura *et al.*, 1994) or *bsr* from *Bacillus cereus* (Izumi *et al.*, 1991). These deaminases convert Blasticidin S to a non-toxic deaminohydroxy derivative (Izumi *et al.*, 1991). See page 134 for handling and preparing Blasticidin.

The formula for Blasticidin S is $C_{17}H_{26}N_8O_5$ -HCl, and the molecular weight is 458.9. Recommended concentration is 100 µg/ml for bacterial cells and 2-10 µg/ml for mammalian cells. The structure of Blasticidin is shown below.



LB Broth (1X)

The LB Broth is supplied ready-to-use as a liquid at 1X concentration. The formulation per one liter at 1X: 10 g SELECT Peptone 140, 5 g SELECT Yeast Extract, 5 g sodium chloride. Add the desired antibiotic at the required concentration to the LB Broth and use for growth of bacterial cultures.

Note: The resulting LB Broth is low salt LB medium suitable for use with Blasticidin as well as other antibiotics.

BioModule[™] BLOCK-iT[™] Unit with Lentiviral Pol II miR RNAi Expression System

Introduction	The BioModule [™] BLOCK-iT [™] Unit with Lentiviral Pol II miR RNAi Expression System facilitates highly efficient, <i>in vitro</i> or <i>in vivo</i> delivery of a miRNA sequence to dividing and non-dividing mammalian cells using a replication- incompetent lentivirus.
Components of the System	 The system includes the following major components The BioModule[™] BLOCK-iT[™] Unit with Pol II miR RNAi Expression Vector containing the pcDNA[™]6.2-GW/EmGFP-miR expression vector for production of an expression clone that allows expression of your miRNA sequence of interest in mammalian cells using a Pol II promoter.
	 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, ψ packaging signal) and a selectable marker to allow generation of stable cell lines. For more information about the pLenti6/V5-DEST vector, see page 147.
	• The pDONR [™] 221 vector is used as an intermediate to transfer the pre-miRNA expression cassette into the lentiviral expression plasmid (see above) using Gateway [®] Technology. See page 25 for details on pDONR [™] 221 vector.
	• Gateway [®] BP and LR Clonase [™] 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.
	• One Shot [®] Stbl3 [™] Competent <i>E. coli</i> to obtain optimal results with lentiviral DNA after transformation.
	• The ViraPower [™] 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 pages 152-156.
	• 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, see page 22.
	 PureLink[™] HQ Mini Plasmid Purification Kit for high-quality plasmid DNA purification suitable for transfection into mammalian cells and S.N.A.P.[™] MidiPrep Kit for plasmid DNA purification from lentiviral vectors.
	• Lipofectamine [™] 2000 Reagent and Opti-MEM [®] Reduced Serum Medium for efficient transfection of your construct into mammalian cells.
	• E-Gel [®] agarose gels and DNA ladders for checking integrity of the ds oligo by gel electrophoresis and for analysis of restriction digests to validate lentiviral expression clones.

BioModule[™] BLOCK-iT[™] Unit with Lentiviral Pol II miR RNAi Expression System, Continued

System Overview	You will co-transfect the ViraPower [™] 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.		
VSV Envelope Glycoprotein	Most retroviral vectors are limited in their usefulness as delivery vehicles by their restricted tropism and generally low titers. In the BioModule TM BLOCK-iT TM Unit with 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 Vector	 The 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 (U3) 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[™]6.2-GW/EmGFP-miR expression clone using Gateway[®] Technology Chloramphenicol resistance gene (Cm^R) located between the two <i>att</i>R sites for counterselection 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> 		

BioModule[™] BLOCK-iT[™] Unit with Lentiviral Pol II miR RNAi Expression System, Continued

Gateway [®] Clonase [™] II Enzyme Mixes	The BP and LR Clonase [™] II enzyme mixes combine the proprietary enzyme formulation and 5X Clonase Reaction Buffer previously supplied as separate components in Clonase [™] enzyme mixes into an optimized single-tube format for easier set-up of the BP or LR recombination reaction. The LR Clonase [™] II Enzyme catalyzes the <i>att</i> L x <i>att</i> R Gateway [®] recombination reaction while the BP Clonase [™] II Enzyme catalyzes the <i>att</i> B x <i>att</i> P Gateway [®] recombination reaction. Use the protocol provided on page 80 to perform the recombination reactions using the Rapid protocol or page 133 using the standard protocol. BP and LR Clonase [™] II Enzyme Mixes are supplied with the kit or available
	separately from Invitrogen.
One Shot [®] Stbl3 [™] Chemically Competent <i>E. coli</i>	The Stbl3 TM <i>E. coli</i> strain is derived from the HB101 <i>E. coli</i> strain and is recommended for use when cloning unstable inserts such as lentiviral DNA containing direct repeats. The transformation efficiency of One Shot [®] Stbl3 TM Chemically Competent cells is greater than 1×10^8 cfu/µg DNA.
	Note: One Shot [®] Stbl3 TM cells require IPTG to induce expression from the <i>lac</i> promoter. If blue/white screening is required to select for transformants, make sure that selective plates contain 50 μ g/ml X-gal and 1 mM IPTG.
ViraPower [™] Packaging Mix	The ViraPower [™] Packaging Mix 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 pages 152-156.
S.N.A.P. [™] MidiPrep Kit	The S.N.A.P. [™] MidiPrep Kit allows isolation of highly pure plasmid DNA that is suitable for transfection, manual or automated sequencing, PCR, restriction mapping, ligation, and transformation. The S.N.A.P. [™] resin is supplied in a column and preferentially binds supercoiled plasmid DNA. Use the S.N.A.P. [™] MidiPrep Kit to prepare plasmid DNA from lentiviral vectors.

BioModule[™] BLOCK-iT[™] Unit with Lentiviral Pol II miR RNAi Expression System, Continued

Antibiotics	Kanamycin Sulfate
	Kanamycin is effective as a bacteriocidal agent by inhibiting ribosomal translocation and eliciting miscoding. Resistance is conferred by the KanR- <i>Tn</i> 5 gene product (aminoglycoside phosphotransferase) that modifies the antibiotic and prevents interaction with ribosomes.
	The formula for Kanamycin Sulfate is $C_{18}H_{36}N_4O_{11}$. H_2SO_4 and the molecular weight is 583. Recommended concentration for use is 50 µg/ml.
	Ampicillin
	Ampicillin is a semi-synthetic penicillin derived from the penicillin nucleus, 6-amino-penicillanic acid and causes cell death by inhibiting cell wall biosynthesis. Resistance is conferred by β -lactamase cleavage of the β -lactam ring (<i>bla</i> gene).
	The formula for Ampicillin is $C_{16}H_{18}N_3O_4S_2Na$ and molecular weight is 371.4. Recommended concentration for use is 100 μ g/ml.
	For details on Spectinomycin and Blasticidin, see page 16.
Note	For details on PureLink™ HQ Mini Plasmid Purification Kit, Lipofectamine™ 2000 Reagent, Opti-MEM® Reduced Serum Medium, LB Broth, E-Gel® agarose gels, and DNA ladders, see page 12.

293FT Cell Line

Introduction	The 293FT Cell Line is derived from the 293F Cell Line (see below) and stably expresses the SV40 large T antigen from the pCMVSPORT6TAg.neo plasmid. Expression of the SV40 large T antigen is controlled by the human cytomegalovirus (CMV) promoter and is high-level and constitutive. For more information about pCMVSPORT6TAg.neo, see page 158.
Use of the Cell Line	Studies have demonstrated maximal virus production in human 293 cells expressing SV40 large T antigen (Naldini <i>et al.</i> , 1996), making the 293FT Cell Line a particularly suitable host for generating lentiviral constructs using the ViraPower [™] Lentiviral Expression System including BioModule [™] BLOCK-iT [™] Unit with Lentiviral Pol II miR RNAi Expression System.
Parental Cell Lines	The 293 Cell Line is a permanent line established from primary embryonal human kidney transformed with sheared human adenovirus type 5 DNA (Graham <i>et al.</i> , 1977; Harrison <i>et al.</i> , 1977). The E1A adenovirus gene is expressed in these cells and participates in transactivation of some viral promoters, allowing these cells to produce very high levels of protein.
	The 293-F Cell Line available from Invitrogen (Catalog no. 11625) is a fast- growing variant of the 293 cell line, and was originally obtained from Robert Horlick at Pharmacopeia.
Antibiotic Resistance	293FT cells stably express the neomycin resistance gene from pCMVSPORT6TAg.neo and should be maintained in medium containing 500 μg/ml Geneticin [®] . Expression of the neomycin resistance gene in 293FT cells is controlled by the SV40 enhancer/promoter. Geneticin [®] is available separately from Invitrogen (page 160).

Gateway[®] Recombination Reactions

Introduction	The Gateway [®] Technology is a universal cloning method that takes advantage of the site-specific recombination properties of bacteriophage lambda (Landy, 1989) to provide a rapid and highly efficient way to move your DNA sequence of interest into multiple vector systems.
	Review the information in this section to briefly familiarize yourself with the Gateway [®] recombination reactions. For details, refer to the Gateway [®] Technology with Clonase [™] II manual available from our web site at <u>www.invitrogen.com</u> or by contacting Technical Service (page 163).
Gateway [®] Vectors	Each of the vectors supplied in the BioModule [™] BLOCK-iT [™] RNAi Units is Gateway [®] -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 Reactions	Two recombination reactions constitute the basis of the Gateway [®] Technology: BP Reaction
	Facilitates recombination of an <i>att</i> B substrate (<i>att</i> B-PCR product or a linearized <i>att</i> B expression clone) with an <i>att</i> P substrate (donor vector) to create an <i>att</i> L-containing entry clone. This reaction is catalyzed by BP Clonase TM II enzyme mix.
	LR Reaction
	Facilitates recombination of an $attL$ substrate (entry clone) with an $attR$ substrate (destination vector) to create an $attB$ -containing expression clone. This reaction is catalyzed by LR Clonase TM II enzyme mix.
Note	If you are an experienced user of Gateway [®] Technology and wish to perform the Rapid BP/LR recombination reaction, we recommend that you review the information on this page before proceeding with the experiment.
	Do not use the standard recombination reaction conditions to perform the Rapid BP/LR recombination reaction.

Gateway[®] Recombination Reaction, Continued

Pre-miRNA Expression	Since the pcDNA [™] 6.2-GW/EmGFP-miR expression vector contains <i>att</i> B sites, the expression vector containing the pre-miRNA sequence cannot be transferred directly into any destination vector including pLenti6/V5-DEST using a single recombination reaction.
	To transfer your pre-miRNA expression cassette from pcDNA [™] 6.2- GW/EmGFP-miR expression clone into a destination vector, you need to perform the two Gateway [®] recombination reactions as follows:
	 Generate an entry clone by performing a BP recombination reaction between the <i>att</i>B substrate (pcDNA[™]6.2-GW/EmGFP-miR expression clone) and <i>att</i>P substrate (pDONR[™]221 vector) using BP Clonase[™] II Enzyme Mix.
	 Perform an LR recombination reaction between the resulting entry clone (<i>att</i>L substrate) and a destination vector including pLenti6/V5-DEST vector (<i>att</i>R substrate) using LR Clonase[™] II Enzyme Mix.
	The standard BP and LR recombination reaction requires more than 2 days for completion. See below for details on expressing the miRNA from destination vector using the Rapid BP/LR recombination reaction.
Rapid BP/LR Recombination Reaction	To develop a faster Gateway [®] recombination reaction protocol that transfers the pre-miRNA expression cassette into the destination vector, we have developed a Rapid BP/LR recombination reaction 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 within a day.
	For Rapid BP/LR Recombination Reactions , perform a BP recombination reaction between the pcDNA [™] 6.2-GW/EmGFP-miR expression clone and pDONR [™] 221 donor vector using BP Clonase [™] II Enzyme Mix, then perform an LR recombination reaction between the resulting entry clone (pENTR [™] 221/miR) and a destination vector including pLenti6/V5-DEST vector using LR Clonase [™] II Enzyme Mix (see below) to produce an expression clone.
	Linearized miR expression clone + pDONR TM 221 vector + destination vector destination vector + destination + destination vector + destination + destination vector + destination + destination vector + destination

Gateway[®] Recombination Reaction, Continued

Features of pDONR [™] 221	The pDONR [™] 221 vector (included with Cat. no. WFGE08) 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 [®] expression clone or <i>att</i> B PCR product
	• <i>ccd</i> B gene located between the two <i>att</i> P 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 [™] 221, see page 150.

Green Fluorescent Protein

Description	The pcDNA [™] 6.2-GW/EmGFP-miR expression vector 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 a destination vector, 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 such variant of enhanced GFP.
Note	We have observed reduced EmGFP expression from miRNA-containing vectors when compared to non-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 after the initiation methionine of 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 [™] 6.2-GW/EmGFP-miR expression vector has the following excitation and emission wavelengths, as published in the literature (Tsien, 1998):			
	Excitation (nm) <u>E</u>	Emission (nm)		
	487	509		
Filter Sets for Detecting EmGFP Fluorescence	detection of the fluore for detection within the	etected with standard FITC filter sets. However, for optimal escence signal, you may use a filter set which is optimized he excitation and emission ranges for the fluorescent protein. escence microscopy and the manufacturer are listed below:		
	Filter Set	<u>Manufacturer</u>		
	Omega XF100	Omega (www.omegafilters.com)		

Biosafety Features of the Lentiviral System

Introduction	The lentiviral and packaging vectors supplied in the BioModule [™] BLOCK-iT [™] Unit with Lentiviral Pol II miR 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 its biosafety and to minimize its relation to the wild-type, human HIV-1 virus. These safety features are discussed below.
Biosafety Features of the BioModule [™] BLOCK-iT [™] Unit with Lentiviral Pol II miR RNAi Expression System	 The BioModule[™] BLOCK-iT[™] Unit with Lentiviral Pol II miR RNAi Expression System includes the following key safety features: 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.
	 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>). 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). 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).
	• 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.
	 The lentiviral particles produced in this system are replication-incompetent and only carry the gene of interest. No other viral species are produced. 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).
	• 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).
	Continued on next page

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 miRNA targeting human genes involved in controlling cell division (*e.g.* tumor suppressor genes).

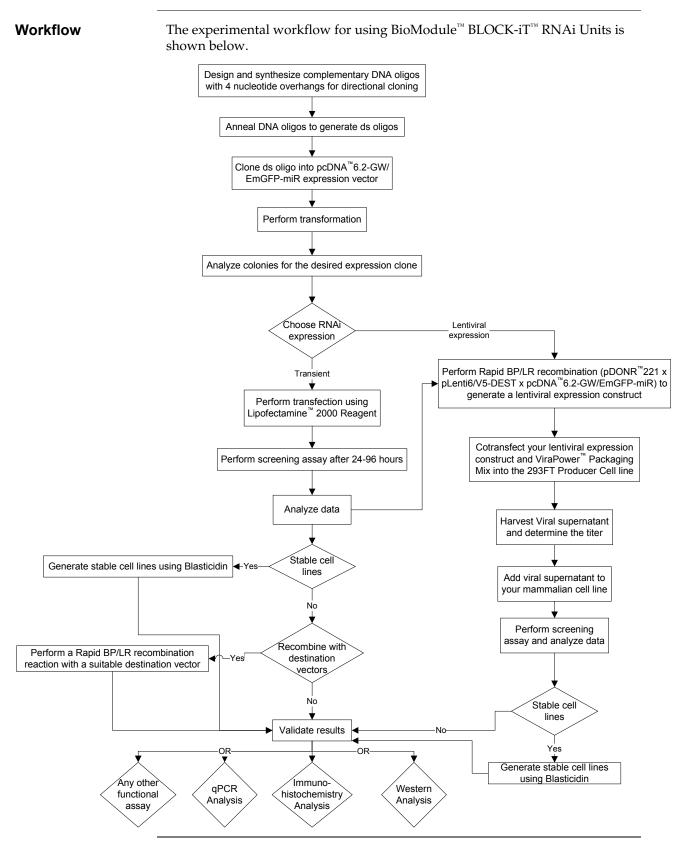
For more information about the BL-2 guidelines and lentivirus handling, refer to the document, "Biosafety in Microbiological and Biomedical Laboratories", 4th Edition, published by the Centers for Disease Control (CDC). This document may be downloaded from the Web at the following address:

http://www.cdc.gov/od/ohs/biosfty/bmbl4/bmbl4toc.htm

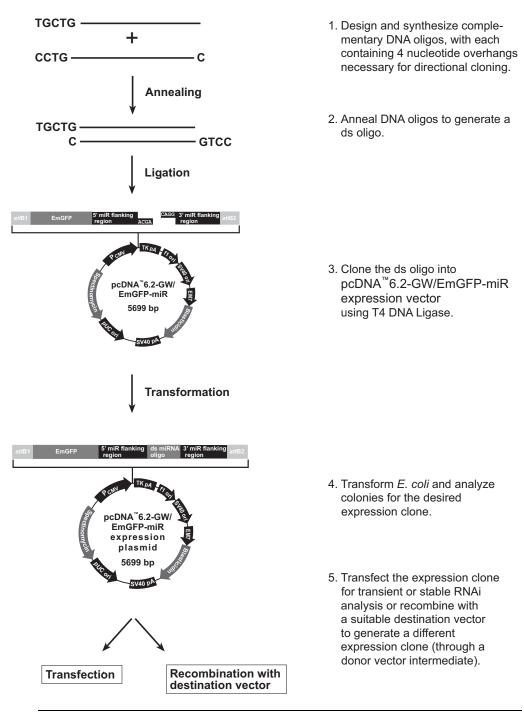


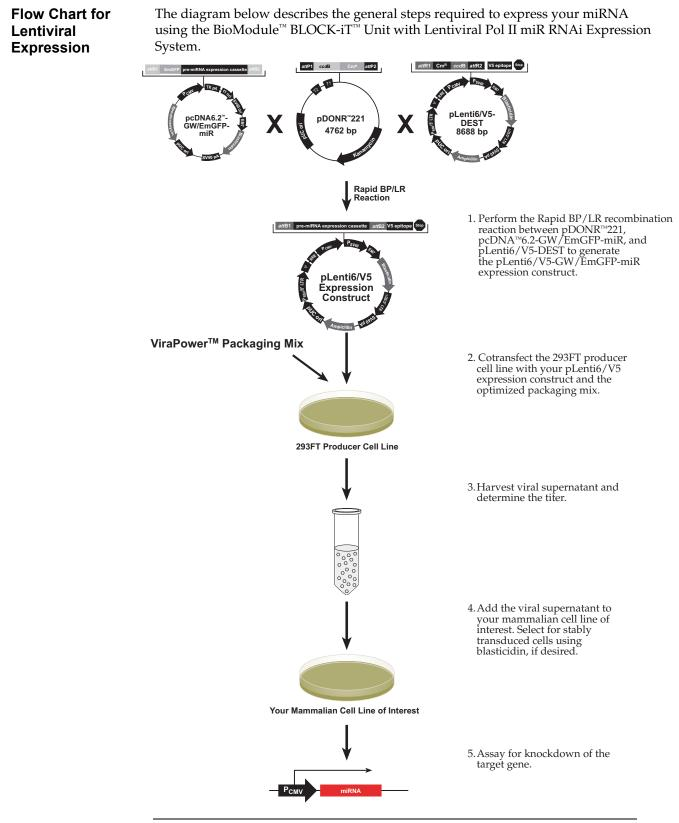
Handle all lentiviruses in compliance with established institutional guidelines. Since safety requirements for use and handling of lentiviruses may vary at individual institutions, we recommend consulting the health and safety guidelines and/or officers at your institution prior to use of the BioModule[™]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

Experimental Outline



Flow Chart for Cloning and Expression The figure below illustrates the major steps necessary to produce a pcDNA[™]6.2-GW/EmGFP-miR expression clone using the BioModule[™] BLOCK-iT[™] Unit with Pol II miR RNAi Expression Vector.





Materials Needed

Materials supplied with the BioModule[™] BLOCK-iT[™] RNAi Units and User Supplied materials are listed below. Ordering information is on page 160.

Step	Supplied in the kit	User Supplied	
Generating ds Oligo	 10X Oligo Annealing Buffer DNase/RNase-Free Water	• Synthesize your "top strand" and "bottom strand" single-stranded oligo (200 µM in water or TE Buffer)	
		• 95°C water bath or heat block	
Checking the Integrity of ds Oligo	 4% E-Gel[®] E-Gel[®] PowerBase[™] TrackIt[™] 10 bp DNA Ladder 	<i>Optional:</i> Sample Buffer	
Ligation Reaction	 pcDNA[™]6.2-GW/EmGFP-miR, linearized 5X Ligation Buffer DNase/RNase-Free Water T4 DNA Ligase (1 U/µl) 	Double-stranded oligo of interest	
Transformation	 One Shot[®] TOP10 Chemically Competent <i>E. coli</i> S.O.C. Medium and LB Broth pUC19 positive control Spectinomycin Ampicillin 	 42°C water bath 37°C shaking and non-shaking incubator 	
Plasmid DNA Purification	PureLink [™] HQ Mini Plasmid Purification Kit	64 ml 100% ethanol	
Transfection	 Lipofectamine[™] 2000 Reagent Opti-MEM[®] I Reduced Serum Medium 	 Mammalian cell line of interest Appropriate tissue culture plates and supplies 	
Rapid BP/LR Recombination Reaction	 pDONR[™]221 vector pLenti6/V5-DEST BP Clonase[™] II enzyme mix LR Clonase[™] II enzyme mix Proteinase K solution One Shot[®] Stbl3[™] Competent <i>E. coli</i> 	 TE Buffer, pH 8.0 (10 mM Tris-HCl, pH 8.0, 1 mM EDTA) If you are not performing lentiviral expression, then you will need: Destination vector of choice Appropriate competent cells 	

Step	Supplied in the kit	User Supplied
Lentiviral Plasmid DNA Purification	S.N.A.P.™ MidiPrep Kit	 240 ml 95% ethanol Sterile water or TE
Lentivirus Production	 ViraPower[™] Packaging Mix 293FT cells Lipofectamine[™] 2000 Reagent Opti-MEM[®] I Reduced Serum Medium 	 Fetal bovine serum (FBS) Complete growth medium for 239FT Cell Appropriate tissue culture plates and supplies
Transduction	• Blasticidin	 Mammalian cell line of choice Complete culture medium for your cell line 6 mg/ml Polybrene[®], if desired Appropriate tissue culture plates and supplies

Methods

General Information

Using Appropriate Protocols	This manual supports protocols to express your miRNA sequence in pcDNA [™] 6.2-GW/EmGFP-miR expression vector for RNAi analysis in mammalian cells and express your miRNA sequence in mammalian cells with lentivirus-based delivery using the two BioModule [™] BLOCK-iT [™] RNAi Units.			
	Be sure to use protocols that are appropriate for the experiment that you wish to perform as shown in the table below:			
	Experiment	Choose Protocols on		
	To express your miRNA sequence in pcDNA™6.2-GW/EmGFP-miR	page 36 (cloning miRNA)		
	Perform transient or stable RNAi analysis in mammalian cells	page 60 (transfection and RNAi analysis)		
	Transfer the pre-miRNA expression cassette into a destination vector including pLenti6/V5-DEST	page 74 (transfer pre-miRNA expression cassette to destination vectors)		
	To express your miRNA sequence in dividing or non-dividing mammalian cells using lentivirus-based delivery	page 92 (lentivirus production, transduction, and RNAi analysis)		
Description of the second seco	 The BioModule[™] BLOCK-iT[™] RNAi Units are CMV promoter-based vector to express miRN analysis. Although the kit has been designed representing a particular target sequence in thof the kit for RNAi analysis assumes that user gene silencing, vector-based production of misystems, and cloning. We highly recommend knowledge of the RNAi pathway and lipid-m The BioModule[™] BLOCK-iT[™] Unit with Lentridesigned to help you create a lentivirus to del sequence in mammalian cells for RNAi analysis mediated transfection, Gateway[®] Technology. For more information on the following topics, RNAi pathway and expression of miRNA references (Brummelkamp <i>et al.</i>, 2002; Cu Sharp, 2002; Sui <i>et al.</i>, 2002; Yu <i>et al.</i>, 2002; Retrovirus biology and the retroviral repl Wong-Staal, 2000 and Luciw, 1996. 	IA in mammalian cell lines for RNAi to help you express miRNA he simplest, most direct fashion, use is are familiar with the principles of IRNA, transfection in mammalian that users possess a working rediated transfection. viral Pol II miR RNAi Expression is liver and express an miRNA sis. We highly recommend that users sue culture techniques, lipid- , and the RNAi pathway. , refer to published references: A in mammalian cells: see published llen, 2004; Kim, 2005; McManus & ; Zeng <i>et al.</i> , 2002) ication cycle: see Buchschacher and		
	1999.			

• General Molecular Biology handbooks, such as *Current Protocols in Molecular Biology* (Ausubel *et al.*, 1994).

Cloning miRNA

Experimental Outline

The experimental steps necessary to clone the miRNA of interest into pcDNA[™]6.2-GW/EmGFP-miR expression vector are outlined below. We recommend performing the steps as described below to achieve the best results.

Step	Action	Page no.
1	Designing and synthesizing the single-stranded oligos.	37
2	Anneal the single-stranded oligo to generate the double- stranded oligo. Dilute the oligo to the required concentration.	43
3	Check the integrity of the double-stranded oligo by gel electrophoresis.	47
4	Clone the double-stranded oligo into pcDNA [™] 6.2- GW/EmGFP-miR expression vector.	51
5	Transform the ligation mixture into One Shot [®] TOP10 Chemically Competent <i>E. coli</i> .	53
6	Analyze transformants by sequencing.	54
7	Optional:	
	Perform chaining of miRNA.	57
	Remove EmGFP Coding Sequence.	59

Designing the Single-Stranded DNA Oligos

Introduction

To use the BioModule[™] BLOCK-iT[™] RNAi Units, you need to design two singlestranded DNA oligonucleotides; one encoding the target pre-miRNA ("top strand" oligo) and the other its complement ("bottom strand" oligo). You will then anneal the top and bottom strand oligos to generate a double-stranded oligonucleotide (ds oligo) suitable for cloning into the pcDNA[™]6.2-GW/EmGFPmiR vector.

The design of the single-stranded oligonucleotides (ss oligos) is critical to the success of both the cloning procedure and ultimately, the RNAi analysis. General guidelines are provided in this section to help you choose the target sequence and to design the ss oligos. Note, that simply following these guidelines does not guarantee that the pre-miRNA will be effective in knocking down the target gene. For a given target gene, you may need to generate and screen multiple pre-miRNA sequences to identify one that is active in gene knockdown studies.



We **strongly** recommend using **Invitrogen's RNAi Designer**, an online tool to help you design and order pre-miRNA sequences for any target gene of interest. The RNAi Designer incorporates the guidelines provided in this manual as well as other design rules into a proprietary algorithm to design pre-miRNA sequences that are compatible for use in cloning into pcDNA[™]6.2-GW/EmGFPmiR. Success rates exceeding 70% have been achieved with the RNAi Designer (i.e. more than 70% of designed miRNAs reduce target gene expression by at least 70%). To use the RNAi Designer, go to <u>www.invitrogen.com/rnaidesigner</u>.

Note: Invitrogen's RNAi Designer is the only online tool for miRNA design at the moment this manual is printed. Other online RNAi designers not intended for miRNA will not necessarily design active pre-miRNA sequences.

Features of Pre- miRNA Insert	When designing the oligos encoding the pre-miRNA, consider that a pre-miRNA insert contains the following features (from 5' to 3' end):
	 5 nucleotides (TGCTG) derived from the endogenous miR-155, an endogenous murine miRNA that is the basis of the miRNA vector system developed in the laboratory of David Turner (U.S. Patent Publication No. 2004/0053876). This also provides a four nucleotide 5' overhang, compatible with a 4 nucleotide overhang in the provided linearized pcDNA[™]6.2-GW/EmGFP-miR to clone the double-stranded oligo.
	• Reverse complement of the 21-nucleotide target sequence (mature miRNA sequence). When transcribed, this is the core sequence that will target your gene of interest, and needs to be antisense to the targeted messenger RNA.
	• 19 nucleotides derived from miR-155 to form a terminal loop with an engineered <i>Msc</i> I site to aid in sequence analysis.
	• Nucleotides 1-8 and 11-21 of the sense target sequence. Note that nucleotides 9 and 10 are removed to form a short internal loop in the mature miRNA, which results in more efficient knockdown.
	• 4 nucleotides derived from endogenous miR-155. This also constitutes the four nucleotide 5' overhang, compatible with a 4 nucleotide overhang in the provided linearized pcDNA [™] 6.2-GW/EmGFP-miR to clone the double-stranded oligo.
	Upon transcription, the mature miRNA sequence and its complement form a stem of the pre-miRNA with a short internal loop, separated by a larger terminal loop. The folded pre-miRNA structure of miR- <i>lacZ</i> is shown below (<i>lacZ</i> targeting sequence in bold)
	UG UU UUGGCC CUG AAAUCGCUGAU GUGUAGUC GUU \ A GACUUUAGCGACUA CACAUCAGCAG / AG^ UCAGUC

Choosing the Target Sequence

When performing RNAi analysis on a particular gene, your choice of target sequence can significantly affect the degree of gene knockdown observed. We recommend following the guidelines below when choosing your target sequence. These are general recommendations only; exceptions may occur.

Length: The target sequence should be 21 nucleotides in length.

Complexity:

- Make sure that the target sequence does **not** contain runs of more than three of the same nucleotide.
- Choose a sequence with low to moderate GC content (~30-50% GC content is suggested).
- Do not choose a target sequence that is a known site for RNA-protein interaction.
- Avoid the following restriction sites, which may be used for optional, advanced features later.

Restriction site	Sequence	Advanced Feature	Page
Msc I	TGGCCA	Alternate sequencing protocol	55
BamH I	GGATCC	miRNA chaining	57
Bgl II	AGATCT	miRNA chaining	57
Sal I	GTCGAC	miRNA chaining	57
Xho I	CTCGAG	miRNA chaining	57
Dra I	TTTAAA	Removal of EmGFP	59

Homology: Make sure that the target sequence does **not** contain significant homology to other genes as this can increase off-target RNAi effects.

Orientation: Choose a target sequence encoding the **sense** sequence of the target mRNA.

Generating the Top Oligo		generate the top oligo sequence, combine these elements (from 5' end to end):
Sequence	1.	5' TGCTG
	2.	Reverse complement of the 21-nucleotide sense target sequence. This is the Mature miRNA Sequence.
	3.	GTTTTGGCCACTGACTGAC (terminal loop).
	4.	Nucleotides 1-8 (5'-3') of sense target sequence.
	_	

5. Nucleotides 11-21 (5'-3') of sense target sequence.

Generating the	To generate the bottom oligo sequence, perform the following steps:
Bottom Oligo	1. Remove 5' TGCT from top oligo sequence (new sequence starts with G).
Sequence	2. Take the reverse complement of the sequence from step 1.
	3. Add CCTG to the 5' end of the sequence from step 2.
Note	 We recommend using Invitrogen's RNAi Designer at <u>www.invitrogen.com/rnaidesigner</u>, which automatically applies the design rules, and produces a high rate of knockdown success. It is not necessary to add 5' phosphates to your single stranded oligos during synthesis. The phosphate groups necessary for ligation are present in the linearized pcDNA[™]6.2-GW/EmGFP-miR.
Example of ss Oligo Design	The diagram below illustrates the required features of the top strand and bottom strand single-stranded oligos as discussed in this section. This particular example lists the sequences of top and bottom strand oligos encoding an miRNA targeting the $lacZ$ gene. These ss oligos were annealed to generate the miR- $lacZ$ positive ds control oligo supplied in the kit.
Top strand oligo	derived from antisense target sequence miR-155 (Mature miRNA Sequence) derived from miR-155 (nucleotides 1-8 and 11-21) 5' - TGCTGAAATCGCTGATTTGTGTGTGTGTGTGTGGCCACTGACCGACTGACGACTACACATCAGCGATTT-3' sequence for s' loop
Bottom strand oligo	derived from miR-155 reverse complement of top strand oligo sequence (minus 5' overhang) 5' - CCTGAAATCGCTGATGTGTGTGTGTCGTCAGTCGGCGAAAACGACTACACAAATCAGCGATTTC-3' sequence for 5' overhang Annealing
ds oligo	↓ 5' -TGCTGAAATCGCTGATTTGGTGTGTGTCGTTTTGGCCACTGACGACTACACATCAGCGATTT-3' 3' -CTTTAGCGACTAAACACATCAGCAAAACCGGTGACTGACT
- Solution of the second secon	We generally purchase unpurified, desalted single-stranded oligos using Invitrogen's custom primer synthesis service (see <u>www.invitrogen.com</u> for more information). The ss oligos obtained anneal efficiently and provide optimal cloning results. Note however, that depending on which supplier you use, the purity and quality of the ss oligos may vary. If you obtain variable annealing and

purity and quality of the ss oligos may vary. If you obtain variable annealing and cloning results using unpurified, desalted oligos, purchase oligos that are HPLC or PAGE-purified.

Cloning Site and Recombination Region of pcDNA[™]6.2-GW/EmGFP-miR Use the diagram on the next page to help you design suitable DNA oligonucleotides to clone into pcDNA[™]6.2-GW/EmGFP-miR after annealing. Note the following features in the diagram next page:

- The pcDNA[™]6.2-GW/EmGFP-miR vector is supplied linearized between nucleotides 1518 and 1519. The linearized vector contains 4 nucleotide overhangs derived from miR-155 sequences. Note that the annealed double-stranded (ds) oligo **must** contain specific 4 nucleotide 5' overhangs on each strand as indicated.
- The light shaded region corresponds to those DNA sequences that will be transferred from the initial pre-miRNA expression vector into the Gateway[®] destination vector (*e.g.* pLenti6/V5-DEST Gateway[®] Vector) following recombination. The dark shaded region represents the EmGFP coding sequence.

Note: Following recombination with a Gateway[®] destination vector, the resulting expression clone will contain a pre-miRNA expression cassette consisting of the EmGFP coding sequence, 5' miR flanking region, miRNA sequence, and the 3' miR flanking region.

The complete sequence of pcDNA[™]6.2-GW/EmGFP-miR is available for downloading from <u>www.invitrogen.com</u> or by contacting Technical Service (page 163). For a map of pcDNA[™]6.2-GW/EmGFP-miR, see page 143.

Cloning Site and Recombination Region of pcDNA[™]6.2-GW/EmGFP-miR, continued

531	CCATTGACGC	CAAT AAATGGGCGG TTTACCCGCC		CGGTGGGAGG GCCACCCTCC	TATA TCTATATAAG AGATATATTC	3' end of CMV promoter CAGAGCTCTC GTCTCGAGAG
591	Puta transcripti TGGCTAACTA ACCGATTGAT	onal start GAGAACCCAC	TGCTTACTGG ACGAATGACC	CTTATCGAAA GAATAGCTTT	TTAATACGAC AATTATGCTG	
651		CTGGCTAGTT GACCGATCAA	AAGCTATCAA TTCGATAGTT		attB1 CAAAAAAGCA GTTTTTTCGT	GGCTTTAAAA CCGAAATTTT
		EmGFP co	ding sequence			
	I			EmGFP for	ward sequencing p	primer site
711	CC ATG GTG GG TAC CAC Met Val	AGC AAG GG TCG TTC CC Ser Lys Gl	G EmGFP	CCG TAC (GAC GAG CTG CTG CTC GAC Asp Glu Leu	TAC AAG TAA ATG TTC ATT Tyr Lys ***
1433		CTTCGTGGCC	GTCGATCGTT	- <i>DI</i> - DI - DI - DI - DI - DI - DI - DI - DI	<i>– Jes</i> – Tagtgagtcg	H H B B B B C C A C C A G G A T
	CGATTCGT	GAAGCACCGG	CAGCTAGCAA	ATTTCCCTCC	ATCACTCAGC	TGGTCACCTA
	5	' miR flanking regi	on		3' miR flank	ing region
1491		TGCTGAAGGC ACGACTTCCG	TGTA <mark>TGCTG P</mark> ACATACGAC	re-miRNA CA ds oligo GT	GGACACAAGG	
					CCTGTGTTCC	GGACAATGAT
			- Bg/ II			GGACAAIGAI
1541	GCACTCACAT CGTGAGTGTA	GGAACAAATG CCTTGTTTAC	Bg/ II	_	GAGATATCTA CTCTATAGAT	GACCCAGCTT CTGGGTCGAA
1541	CGTGAGTGTA	CCTTGTTTAC miRNA reverse	= /6g GCCCAGATCT CGGGTCTAGA	GGCCGCACTC	GAGATATCTA CTCTATAGAT	GACCCAGCTT CTGGGTCGAA
1541	CGTGAGTGTA	CCTTGTTTAC	= /6g GCCCAGATCT CGGGTCTAGA	GGCCGCACTC	GAGATATCTA CTCTATAGAT start TK polyar	GACCCAGCTT

Generating the Double-Stranded Oligo

Introduction	Once you have synthesized the appropriate complementary single-stranded DNA oligos, you will anneal equal amounts of each single-stranded oligo to generate a double-stranded oligo (ds oligo). Guidelines and instructions are provided in this section.
Experimental Outline	 To generate the double-stranded oligo, you will: Calculate the amount of single-stranded oligos needed. Anneal the oligonucleotides to create a double-stranded oligonucleotide. Dilute the double-stranded oligo to a final concentration of 10 nM.
Single-Stranded Oligos	 Before beginning, make sure that you have synthesized the single-stranded oligos with the appropriate sequences required for cloning into the pcDNA[™]6.2-GW/EmGFP-miR vector and for annealing. See the figure below for an illustration. Top strand oligo: Make sure that this oligo contains the sequence TGCTG at the 5' end. Bottom strand oligo: Make sure that this oligo contains the sequence CCTG at the 5' end, has a C at the 3' end, and is complementary to the top strand. Top strand oligo 5'-TGCTG → 3' + Bottom strand oligo 5'-CCTG → C-3' Annealing
	ds oligo 5'-TGCTG → 3' 3'-C → GTCC-5'
Amount of DNA Oligo to Anneal	You will anneal equal amounts of the top and bottom strand oligos to generate the ds oligos. We perform the annealing reaction at a final single-stranded oligo concentration of 50 μM . Annealing at concentrations below 5 μ M significantly reduce the efficiency. Note that the annealing step is not 100% efficient.
Resuspending the Oligos	If your single-stranded oligos are supplied lyophilized, resuspend them in water or TE Buffer to a final concentration of 200 μ M before use.
	Continued on next page

Generating the Double-Stranded Oligo, continued

Materials Needed	You will need the following materials:				
	•	Your "top strand" single-stranded oligo	(200 µM in wate	r or TE Buffer)	
	•	Your "bottom strand" single-stranded ol	igo (200 μM in w	vater or TE Buffer)	
	•	10X Oligo Annealing Buffer (supplied w	ith the kit, Box 1)	
	•	DNase/RNase-Free Water (supplied wit	h the kit, Box 1)		
	•	0.5 ml sterile microcentrifuge tubes			
	•	95°C water bath or heat block			
Annealing Procedure		ow this procedure to anneal your single- o. Note that the final concentration of the	0	6	
	1.	In a 0.5 ml sterile microcentrifuge tube, s reaction at room temperature.	set up the followi	ing annealing	
		Reagent	Amount		
		Top strand DNA oligo (200 μM)	5 µl		
		Bottom strand DNA oligo (200 μ M)	5 µl		
		10X Oligo Annealing Buffer	2 µl		
		DNase/RNase-Free Water	8 µl		
			1		

2. Incubate the reaction at 95°C for 4 minutes.

Total volume

3. Remove the tube containing the annealing reaction from the water bath or the heat block and set on your laboratory bench.

20 µl

- 4. Allow the reaction mixture to cool to room temperature for 5-10 minutes. The single-stranded oligos will anneal during this time.
- 5. Place the sample in a microcentrifuge and centrifuge briefly (~5 seconds). Mix gently.
- 6. Remove 1 μl of the annealing mixture and dilute the ds oligo as directed in **Diluting the ds Oligo**, next page.
- 7. Store the remainder of the 50 μM ds oligo mixture at -20°C (stable for at least a year).

Generating the Double-Stranded Oligo, continued

Diluting the ds Oligo	To clone your ds oligo into pcDNA [™] 6.2-GW/EmGFP-miR, you must dilute the 50 µM stock to a final concentration of 10 nM (<i>i.e.</i> 5,000-fold dilution). We generally perform 100-fold and 50-fold serial dilutions, the first into DNase/RNase-free water and the second into the Oligo Annealing Buffer supplied with the kit. Follow the procedure below to dilute the ds oligo.				
	 Dilute the 50 μM ds oligo mixture (from Annealing Procedure, Step 5, previous page) 100-fold into DNase/RNase-free water (<i>i.e.</i> 1 μl of 50 μM ds oligo into 99 μl of DNase/RNase-free water) to obtain a final concentration of 500 nM. Vortex to mix thoroughly. 				
	2. Dilute the 500 nM ds oligo mixture (from Step 1) 50-fold into Oligo Annealing Buffer as follows to obtain a final concentration of 10 nM.				
	500 nM ds oligo 1 μl				
	10X Oligo Annealing Buffer 5 μl				
	DNase/RNase-free water 44 µl				
	Total volume 50 µl				
	3. Vortex to mix thoroughly. Store the remaining 500 nM ds oligo stock at -20°C.				
	4. Aliquot the 10 nM ds oligo stock and store at -20°C.				
Q Important	The undiluted ds oligos are 5,000-fold more concentrated than the working concentration. When performing the dilutions, be careful not to cross-contaminate the different ds oligo stocks. Remember to wear gloves and change pipette tips after every manipulation.				
Storing the ds Oligo	Once you have diluted your ds oligo, you should have three stocks of annealed ds oligo. Use each stock as follows:				
	• 50 μM ds oligo (undiluted): Use this stock to prepare new diluted ds oligo stocks if existing stocks become denatured or cross-contaminated.				
	• 500 nM ds oligo (100-fold dilution): Use this stock for gel analysis (see Checking the Integrity of the ds Oligo , page 47).				
	• 10 nM ds oligo (5,000-fold dilution): Use this stock for cloning (see Ligation Procedure , page 52).				
	Store the three ds oligo stocks at -20°C.				

Generating the Double-Stranded Oligo, continued



When using the diluted ds oligo stock solutions (*i.e.* 100-fold or 5,000-fold diluted stocks), thaw the solutions on ice. **Do not** heat or allow the ds oligo solutions to reach greater than room temperature as this causes the ds oligos to melt. The concentration of the oligos in the diluted solutions is not high enough to permit re-annealing and instead favors the formation of intramolecular hairpin structures. These intramolecular hairpin structures **will not clone** into pcDNA[™]6.2-GW/EmGFP-miR.

If your diluted ds oligo stock solution is heated, discard the ds oligo solution and prepare new diluted stocks using the procedure on the previous page.

Note: If the 50 μ M ds oligo solution (undiluted stock) becomes heated, the oligos are sufficiently concentrated and may be re-annealed following the annealing procedure on page 44.

Checking the Integrity of the ds Oligo

Introduction	You may verify the integrity of your annealed ds oligo using 4% E-Gel [®] Starter Pak included in the kit (Box 11) for agarose gel electrophoresis. We suggest running an aliquot of the annealed ds oligo and comparing it to an aliquot of each starting single-stranded oligo. Note: 4% E-Gel [®] resolves these fragments much better than regular 4% agarose gels.
Experimental	To perform agarose gel electrophoresis, you will:
Outline	1. Prepare your samples for electrophoresis.
	2. Analyze the samples on 4% E-Gel [®] agarose gels.
	3. Visualize the results.
Materials Needed	You will need the following materials:
	• 4% E-Gel [®] (supplied with the kit)
	 E-Gel[®] PowerBase[™] (supplied with the kit)
	• TrackIt [™] 10 bp DNA Ladder (supplied with the kit)
	• Sample Buffer (page 160)
<u>,</u> Ι,	Follow these recommendations to obtain the best results with E-Gel [®] agarose gels:
	• All wells in the gel must contain sample or water. Avoid introducing bubbles while loading, as bubbles will cause bands to distort.
	• The One-Step Loading method described on the next page is routinely used to load E-Gel [®] gels. Use the Two-Step Loading method, if the One-Step Loading method produces fuzzy or indistinct bands, or the gel was removed from its plastic pouch for more than 30 minutes. Refer to the E-Gel [®] Technical Guide (available from <u>www.invitrogen.com</u>) for details on the Two-Step Loading method.
Preparing Samples	You may prepare DNA samples for E-Gel [®] agarose gels in deionized water or loading buffer (recommended final loading buffer concentration is 10 mM Tris- HCl; 1 mM EDTA, pH 7.5; 0.005% bromophenol blue; and 0.005% xylene cyanol FF). If you wish to use 10X BlueJuice [™] Gel Loading Buffer or TrackIt [™] Loading Buffer (page 160), dilute this buffer 50- to 200-fold to obtain the optimal dye concentration.
	Prepare each sample for analysis in total sample volume of 20 µl as follows:
	 Annealed ds oligo: To 5 μl of the 500 nM stock, add 15 μl deionized water or loading buffer
	 Each starting single-stranded oligo: Dilute the 200 μM stock 400-fold to 500 nM. To 5 μl of the 500 nM stock, add 15 μl deionized water or loading buffer
	 Mix 2 µl TrackIt[™] 10 bp DNA Ladder with 18 µl deionized water. Loading undiluted TrackIt[™] DNA Ladder on an E-Gel[®] will result in loss of resolution.

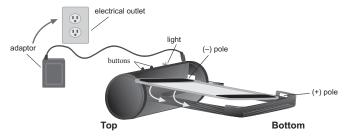
Checking the Integrity of the ds Oligo, Continued

Running E-Gel[®]

Pre-run the E-Gel[®] agarose gel for 2 minutes **with the comb in place** before loading your samples to ensure proper resolution of your DNA fragments.

Each E-Gel[®] cassette is supplied individually wrapped and ready for use. To set up and use an E-Gel[®], follow the instructions below:

1. Plug the E-Gel[®] PowerBase[™] v.4 into an electrical outlet using the adaptor plug.



- 2. Open the package containing the gel and insert the gel (with the comb in **place**) into the apparatus right edge first. Press firmly at the top and bottom to seat the gel in the base. You should hear a snap when it is in place. The Invitrogen logo should be located at the bottom of the base, close to the positive pole. See the above diagram. A **steady, red light** illuminates when the E-Gel[®] is correctly inserted (Ready Mode).
- 3. **Press and hold** either button until the **red** light turns to a **flashing green light**. This indicates the start of the 2-minute pre-run.
- 4. At the end of the pre-run, current automatically shuts off. The **flashing green** light changes to a **flashing red** light and the PowerBase[™] **beeps rapidly**.
- 5. **Press and release** either button to stop the beeping. The light changes from a **flashing red** light to a **steady red** light.
- 6. Remove the comb from the E-Gel[®] using both hands to lift the comb gently by rolling the comb slowly towards you. Remove any excess fluid using a pipette.
- 7. Load 20 μl of sample into each well (see previous page for sample preparation).
- 8. Load 20 µl of the diluted TrackIt[™] 10 bp DNA Ladder (see previous page) on the well. Load 20 µl of water into any remaining empty wells.
- 9. Proceed immediately to electrophoresis, next page.

Checking the Integrity of the ds Oligo, Continued

Electrophoresis of E-Gels [®]	1.	Choose the 30-minute run for single-comb gels on the E-Gel [®] PowerBase [™] v.4. For the 30-minute run, press and release the 30-min button to start the 30-minute electrophoresis run. The light changes to a steady green light.
		Note: The actual running time of the E-Gel [®] gel may vary between 30-33 minutes for single-comb gels.
	2.	Current through the E-Gel [®] gel automatically shuts off at the end of each run. The E-Gel [®] PowerBase [™] v.4 signals the end of the run with a flashing red light and rapid beeping .
	3.	Press and release either button to stop the beeping. The light turns to a steady red light.
	4.	At the end of the run, remove the gel cassette from the power unit and analyze your results using a UV transilluminator.
Note	E-0	Gel® agarose gels can only be used once. Do not re-use them.
What You Should See		hen analyzing an aliquot of the annealed ds oligo reaction by agarose gel ctrophoresis, we generally see the following:
	•	A detectable higher molecular weight band representing annealed ds oligo running around 60 - 70 bp.
	•	A faintly detectable lower molecular weight band representing unannealed oligos that form hairpins, running at around 30-35 bp.
	nex	r an example of expected results obtained from agarose gel analysis, see the xt page. If the band representing ds oligo is weak or if you do not see a band, Troubleshooting , page 121 for tips to troubleshoot your annealing reaction.

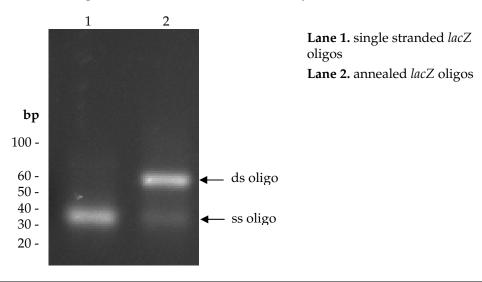
Checking the Integrity of the ds Oligo, continued

Example of Expected Results

In this experiment, *lacZ* control oligos (see page ix for the sequence of each DNA oligo) were annealed (50 μ M final concentration) using the reagents supplied in the kit and following the procedure in this manual to generate the *lacZ* ds control oligo. The annealing reaction was diluted 100-fold in water to a concentration of 500 nM. Aliquots of the diluted ds oligo (5 μ l; 2.5 pmol) and the corresponding single-stranded oligos (mixed but not annealed; 2.5 pmol per oligo) were analyzed on a 4% E-Gel[®].

Results: The *lacZ* oligo annealing reaction shows a clearly detectable, higher molecular weight band that differs in size from each component single-stranded oligo. Remaining unannealed ss oligos are also weakly detectable.

Note: The agarose gel is non-denaturing; therefore, the single-stranded oligos do not resolve at the expected size due to formation of secondary structure.



Performing the Ligation Reaction

Introduction	After generating your ds oligo and diluting to the appropriate concentration, clone the ds oligo into the pcDNA [™] 6.2-GW/EmGFP-miR vector and transform your ligation reaction into competent TOP10 <i>E. coli</i> . It is important to have everything you need set up and ready to use to ensure that you obtain the best results. We suggest that you read the sections entitled Performing the Ligation Reaction (pages 51-52) and Transforming One Shot [®] TOP10 Competent <i>E. coli</i> (page 53) before beginning.
	Note: If you want to perform miRNA chaining, refer to page 57.
Important	You will use T4 DNA Ligase and a 5X Ligation Buffer supplied with the kit to facilitate ligation of your ds oligo with the linearized pcDNA [™] 6.2-GW/EmGFP-miR vector. When performing the ligation reaction, note the following:
	• The T4 DNA Ligase and the 5X Ligation Buffer supplied with the kit have been optimized to permit ligation of the ds oligo into the pcDNA [™] 6.2-GW/EmGFP-miR vector in 5 minutes at room temperature . T4 DNA Ligase preparations and reaction buffers available from other manufacturers may not be appropriate for use in this application.
	 Traditional ligation reactions are performed at 16°C overnight. This is not recommended for this application. Follow the ligation procedure on the next page.
Amount of ds Oligo to Use	For optimal results, use approximately a 15:1 molar ratio of ds oligo insert: vector for ligation.
Ligation Controls	We recommend a ligation with DNase/RNase-Free Water instead of oligo as negative control for the ligation reaction.
	We recommend including the miR- <i>lacZ</i> positive double-stranded (ds) control oligo supplied with the kit as a positive control in your ligation experiment. The miR- <i>lacZ</i> positive ds control oligo is supplied ready-to-use as a 10 nM stock in 1X Oligo Annealing Buffer. See page ix for the sequence of each strand of the <i>lacZ</i> ds control oligo.
	Note: Once you have cloned the <i>lacZ</i> ds control oligo into pcDNA ^{III} 6.2-GW/EmGFP-miR, you may use the resulting expression clone as a positive control for the RNAi response in your mammalian cell line. Simply co-transfect the expression clone and the pcDNA ^{<math>III1.2/V5-GW/lacZ reporter plasmid supplied with the kit into your mammalian cell line and assay for knockdown of β-galactosidase expression.</math>}
Important	Reminder: When using the 10 nM ds oligo stock solution for cloning, thaw the solution on ice. Do not thaw the ds oligo by heating or the ds oligo duplexes may melt and form intramolecular hairpin structures. After use, return the tube to -20°C storage.

Performing the Ligation Reaction, continued

Materials Needed	You will need the following materials:			
	• Double-stranded oligo of interest (10 nM in 1X C on ice before use)	ligo Annea	aling Buffe	r; thaw
	 pcDNA[™]6.2-GW/EmGFP-miR, linearized (5 ng/μl, supplied with the kit, Box 1; thaw on ice before use) 			kit,
	• 5X Ligation Buffer (supplied with the kit, Box 1)			
	• DNase/RNase-Free Water (supplied with the kit	t, Box 1)		
	• T4 DNA Ligase (1 U/ μ l, supplied with the kit, Be	ox 1)		
Ligation Procedure).	
	Reagent	Sample	Positive control	Negative control
	5X Ligation Buffer	4 µl	4 µl	4 µl
	pcDNA [™] 6.2-GW/EmGFP-miR, linearized (5 ng/µl)	2 µl	2 µl	2 µl

	-	-		
miR-ds oligo (10 nM; <i>i.e.</i> 1:5,000 dilution; page 45)	4 µl			
miR- <i>lacZ</i> positive ds control oligo (10 nM)		4 µl		
DNase/RNase-Free Water	9 µl	9 µl	13 µl	
T4 DNA Ligase (1 U/μl)	1 µl	1 µl	1 µl	
Total volume	20 µl	20 µl	20 µl	1

2. Mix reaction well by pipetting up and down.

Note: The presence of PEG and glycerol (supplied by the Ligation Buffer and the T4 DNA Ligase) will make the reaction mixture viscous. Be sure to mix the reaction thoroughly by pipetting up and down. **Do not vortex.**

3. Incubate for 5 minutes at room temperature.

Note: Extending the incubation time may result in a higher yield of colonies. Do not exceed 2 hours.

4. Place the reaction on ice and proceed to **Transforming One Shot**[®] **TOP10 Competent** *E. coli*, next page.

Note: You may store the ligation reaction at -20°C overnight.

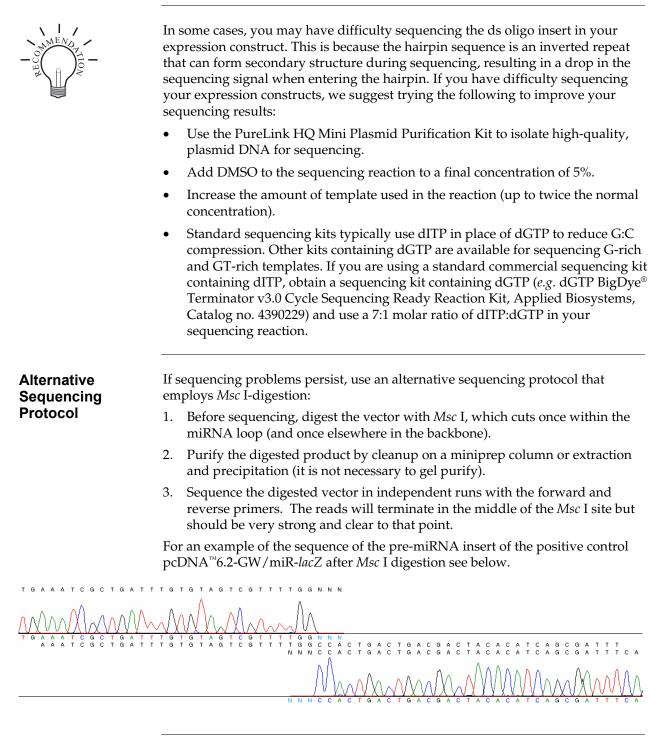
Transforming One Shot[®] TOP10 Competent *E. coli*

Introduction	int inc	nce you have performed the ligation reaction, transform your ligation mixture to competent <i>E. coli</i> . One Shot [®] TOP10 Chemically Competent <i>E. coli</i> (Box 2) are cluded with the kit to facilitate transformation. Follow the guidelines and structions provided in this section.
	No	te: One Shot [®] TOP10 <i>E. coli</i> possess a transformation efficiency of 1 x 10 ⁹ cfu/μg DNA.
Materials Needed	Yo	ou will need the following materials:
	٠	Ligation reaction (from Step 3, previous page)
	•	One Shot [®] TOP10 Chemically Competent <i>E. coli</i> (supplied with the kit, Box 2; one vial per transformation; thaw on ice immediately before use)
	•	S.O.C. Medium (supplied with the kit, Box 2; warm to room temperature)
	•	pUC19 positive control (supplied with the kit, Box 2; if desired)
	•	42°C water bath
	•	LB plates containing 50 μ g/ml spectinomycin (two for each transformation; warm at 37°C for 30 minutes before use). See page 133 for recipe.
	•	LB plates containing 100 μ g/ml ampicillin (if transforming pUC19 control)
	•	37°C shaking and non-shaking incubator
	for	te: Low salt LB agar plates containing $100 \ \mu\text{g/ml}$ Blasticidin can also be used to select transformants. Be sure to use low salt agar plates and check pH carefully for Blasticidin work efficiently. For more information on Blasticidin and recipes, see page 134.
One Shot [®] TOP10 Transformation Procedure	Ch	e this procedure to transform your ligation reaction into One Shot [®] TOP10 nemically Competent <i>E. coli</i> . For a positive control, transform 10 pg (1 μl) of JC19 plasmid into a vial of One Shot [®] TOP10 chemically competent <i>E. coli</i> .
	1.	Add 2 μl of the ligation reaction (from Step 3, previous page) into a vial of One Shot [®] TOP10 chemically competent <i>E. coli</i> and mix gently. Do not mix by pipetting up and down.
	2.	Incubate on ice for 5 to 30 minutes.
		Note: Longer incubations seem to have a minimal effect on transformation efficiency. The length of the incubation is at the user's discretion.
	3.	Heat-shock the cells for 30 seconds at 42°C without shaking.
	4.	Immediately transfer the tubes to ice.
	5.	Add 250 µl of room temperature S.O.C. Medium.
	6.	Cap the tube tightly and shake horizontally (200 rpm) at 37°C for 1 hour.
	7.	Spread 50-200 μ l from each transformation on a pre-warmed LB agar plate containing 50 μ g/ml spectinomycin 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, plate 20-100 μ l of the transformation reaction on LB plates containing 100 μ g/ml ampicillin.
	8.	An efficient ligation reaction may produce several hundred colonies.
	9.	Proceed to analyzing transformants, next page.

Analyzing Transformants

Analyzing	To analyze positive clones, we recommend that you:
Transformants	1. Pick 5-10 spectinomycin-resistant colonies and culture them overnight in LB or SOB medium containing 50 μ g/ml spectinomycin.
	Note: Low salt LB containing 100 μ g/ml Blasticidin can also be used to grow transformants. See page 134.
	 Isolate plasmid DNA using the PureLink[™] HQ Mini Plasmid Purification Kit included with the BioModule[™] units. See page 61 for protocol details.
	3. Sequence each pcDNA [™] 6.2-GW/EmGFP-miR expression construct (see below) to confirm the following:
	a. The presence and correct orientation of the ds oligo insert.
	b. The sequence of the ds oligo insert.
	Note: Because of the small size of the ds oligo insert, we do not recommend using restriction enzyme analysis to screen transformants.
	 sequence of the ds oligo insert. When screening transformants, we find that up to 20% of the clones contain mutated inserts (generally 1 or 2 bp deletions within the ds oligo). The reason for this is not known, but may be due to triggering of repair mechanisms within <i>E. coli</i> as a result of the inverted repeat sequence within the ds oligo insert. Note: Expression clones containing mutated ds oligo inserts generally elicit a poor RNAi response in mammalian cells. Identify expression clones with the correct ds oligo sequence and use these clones for your RNAi analysis.
Sequencing	To facilitate sequencing of your pcDNA [™] 6.2-GW/EmGFP-miR expression clones, use the EmGFP forward sequencing primer and miRNA reverse sequencing primer supplied with the kit (Box 1). See the diagram on page 41 for the location of the priming sites.
Note	If you download the sequence for pcDNA [™] 6.2-GW/EmGFP-miR from our web site, note that the overhang sequences will be shown already hybridized to their complementary sequences (<i>e.g.</i> TGCT will be shown hybridized to ACGA and CAGG will be shown hybridized to GTCC).

Analyzing Transformants, continued



Analyzing Transformants, continued

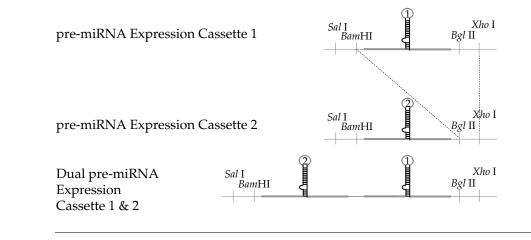
Long-Term Storage	Once you have identified the correct expression clone, be sure to purify the colony and make a glycerol stock for long-term storage. We recommend that you store a stock of plasmid DNA at -20°C.
	 Streak the original colony out for a single colony on an LB plate containing 50 μg/ml spectinomycin.
	2. Isolate a single colony and inoculate into 1-2 ml of LB containing 50 μ g/ml spectinomycin.
	3. Grow until the culture reaches stationary phase.
	4. Mix 0.85 ml of culture with 0.15 ml of sterile glycerol and transfer to a cryovial.
	5. Store the glycerol stock at -80°C.
	Note: Low salt LB containing 100 μ g/ml Blasticidin can also be used to grow transformants.
What to Do Next	 Once you have obtained your pcDNA[™]6.2-GW/EmGFP-miR expression clone, you have the following options: Transfect the expression clone directly into the mammalian cell line of interest to perform transient RNAi analysis (see Transfecting Cells, page 64).
	• Transfect the expression clone directly into the mammalian cell line of interest and isolate stable transfectants that knock down the gene of interest constitutively (see Generating a Stable Cell Line , page 70).
	• Perform miRNA chaining to express multiple pre-miRNAs from one single construct (see Chaining multiple pre-miRNAs , next page).
	• Remove the EmGFP coding sequence from your pcDNA [™] 6.2-GW/EmGFP- miR expression clone (see Removing EmGFP Coding Sequence , page 59).
	• Perform an LR recombination reaction with your expression construct and a suitable Gateway [®] destination vector to generate an expression clone in an alternative backbone (see Performing the Rapid BP/LR Recombination Reaction , page 80).
	• Perform an LR recombination reaction with your expression construct and a pLenti5/V5-DEST destination vector to generate an expression clone to perform lentiviral expression (see Performing the Rapid BP/LR Recombination Reaction , page 80).

Chaining pre-miRNAs

Introduction

miRNAs are sometimes expressed in clusters in long primary transcripts driven by RNA Pol II (Lee *et al.*, 2004). Our vectors support chaining of miRNAs to express them in one primary transcript, thus ensuring co-cistronic expression of multiple miRNAs. In the final construct, the original pattern of restriction sites is regenerated, making the construct amenable to multiples rounds of chaining. The figure below shows the principle of chaining two miRNAs, derived from two different miRNA vectors, into one miRNA expression vector.

Note: Chaining together miRNAs targeting different genes usually results in slightly reduced knockdown of each gene. Chaining different miRNAs targeting the same gene or repeating one miRNA can enhance knockdown. Due to increased processing, EmGFP expression is attenuated by miRNA chaining.



Restriction Strategy Two strategies of restriction digestions are possible:

- a. A combination of *Bam*H I and *Xho* I to excise the pre-miRNA insert, and *Bgl* II and *Xho* I to digest the pre-miRNA expression vector used as backbone
- b. A combination of *Sal* I and *Bgl* II to excise the pre-miRNA insert, and *Sal* I and *Bam*H I to digest the pre-miRNA expression vector used as backbone

The procedure for the first strategy is described on the next page. For the second strategy, change the restriction enzymes used in the procedure.

Chaining pre-miRNAs, continued

Materials Needed	will need the following materials: Appropriate restriction enzymes 2% E-Gel® agarose gel PureLink™ Quick Gel Extraction Kit (page 160) One Shot® TOP10 Competent Cells (page 160)	
Procedure for Chaining	ow is a protocol for chaining miRNAs.	
	Insert: Digest 2 μ g pcDNA ^{M} 6.2-GW/EmGFP-miR-1 with 10 units <i>Bam</i> H 10 units <i>Xho</i> I for 2 hours at 37° C.	I and
	Backbone: Digest 1 μ g pcDNA TM 6.2-GW/EmGFP-miR-2 with 10 units <i>Bg</i> and 10 units <i>Xho</i> I for 2 hours at 37° C.	<i>l</i> II
	Run fragments on 2% E-Gel® or other high percentage agarose gels.	
	Excise the backbone and insert fragments from the gel. Purify the fragments from the PureLink [™] Quick Gel Extraction Kit from Invitrogen or equiva	
	Ligate the purified backbone and insert fragment at a 1:4 molar ratio, usi DNA ligase from Invitrogen or equivalent.	ing T4
	Transform competent cells, such as <i>E. coli</i> . One Shot [®] TOP10 as described page 53.	l on
	Analyze resulting clones as described on pages 54-56	
	Test construct for both miRNAs by transfecting cells as described on pag	șe 60.

Removing EmGFP Coding Sequence

Introduction	Depending on your experiment, you may not wish to express EmGFP from pre-miRNA expression construct. If you have previously established a pcDNA ^{TM} 6.2-GW/EmGFP-miR clone that works well, you can remove the EmGFP coding sequence by <i>Dra</i> I digestion and self-ligation of the vector, forming a pcDNA ^{TM} 6.2-GW/ miR clone expressing the same pre-miRNA. The section describes a procedure for removing the EmGFP coding sequence from pcDNA ^{TM} 6.2-GW/EmGFP-miR.					
Materials Needed	Yo	u will need the following materials:				
	٠	Appropriate restriction enzymes				
	•	 PureLink[™] Quick Gel Extraction Kit (page 160) 				
	•	One Shot [®] TOP10 Competent Cells (page 160)				
Procedure for Removing EmGFP		slow is a protocol for removing the EmGFP Coding Sequence from pcDNA [™] 6.2- W / EmGFP-miR.				
	1.	Digest 1 μ g pcDNA TM 6.2-GW/EmGFP-miR with 10 units <i>Dra</i> I for 2 hours at 37°C.				
	2.	Run fragments on 0.8% E-Gel [®] or other low percentage agarose gels.				
	3.	Excise the vector fragment from the gel and purify the fragment using the PureLink [™] Quick Gel Extraction Kit from Invitrogen or equivalent.				
	4.	Ligate the purified backbone and insert fragment at a 1:4 molar ratio, using T4 DNA ligase.				
	5.	Transform competent cells, such as <i>E. coli</i> . One Shot [®] TOP10 as described on page 53.				
	6.	 Analyze resulting clones by restriction analysis with <i>Dra</i> I; no 750 bp fragment should be visible. 				

Transfection and RNAi Analysis

Experimental Outline

The experimental steps necessary to transfect the pcDNA[™]6.2-GW/EmGFPmiR expression clone into a mammalian cell line of choice and perform RNAi analysis are outlined below. We recommend performing the steps as described below to achieve the best results.

Step	Action	Page no.
1	Purify plasmid DNA from the verified expression clone using PureLink™ HQ Mini Plasmid DNA Purification Kit.	61
2	Culture the mammalian cell of choice to the required cell density.	64
3	Perform transient transfection using Lipofectamine [™] 2000 Reagent.	67
4	Perform analysis of the EmGFP fluorescent protein from the expression clone using fluorescent microscopy to determine the expression of your miRNA.	69
5	Assay for target gene knockdown using qPCR, Western analysis, immunohistochemistry, or any other functional assay.	
6	Optional:	
	Generate stable cell lines that constitutively express your miRNA.	70
	Transfer the pre-miRNA expression cassette into appropriate Gateway [®] destination vectors to allow for expression of the miRNA in viral systems or using tissue-specific promoters.	74

Plasmid DNA Purification

Introduction	 Once you have obtained your expression clone, isolate plasmid DNA for transfection. Plasmid DNA for transfection into eukaryotic cells must be very clean and free from contamination with phenol or sodium chloride. Contaminants will kill the cells, and salt will interfere with lipid complexing, decreasing transfection efficiency. The PureLink[™] HQ Mini Plasmid Purification Kit (Box 3) is included with each BioModule[™] BLOCK-iT[™] RNAi Units for isolating high quality plasmid DNA. Note: If you are using the BioModule[™] BLOCK-iT[™] Unit with Lentiviral Pol II miR RNAi Expression System, do not use the S.N.A.P.[™] MidiPrep Kit for plasmid DNA purification at this step. Use the PureLink[™] HQ Mini Plasmid Purification Kit. 			
Experimental				
Outline	1. Grow <i>E. coli</i> cells.			
	2. Prepare bacterial lysates.			
	3. Bind the plasmid DNA from the lysate to the PureLink ^{TM} Spin Column.			
	4. Perform washing to remove impurities.			
	5. Elute the plasmid DNA in Elution Buffer.			
Materials Needed	You will need the following materials:			
	 PureLink[™] HQ Mini Plasmid Purification Kit (Box 3, supplied with the kit) 64 ml 100% ethanol 			
	Sterile microcentrifuge tubes			
	Microcentrifuge			
Before Starting	 Prepare the Resuspension Solution with the RNase A supplied in the kit. Resuspend the lyophilized RNase A (12 mg) in 200 µl of Resuspension Solution, and then add the resuspended mixture to the remaining Resuspension Solution for a final concentration of 0.1 mg/ml RNase A. After mixing, store the Resuspension Solution with RNase A at +4°C. Stable for up to 6 months. 			
	• Prepare the Wash Buffer with ethanol. Add 64 ml of 96–100% ethanol to the entire volume of Wash Buffer (16 ml) to obtain a total volume of 80 ml.			
	• Before each use, check the Neutralization/Binding Buffer and Lysis Buffer for a white salt precipitate. If present, place each buffer in a 37°C water bath for 5 minutes or until the salts redissolve and the solution clears.			
	Do not shake the Lysis Buffer, as this can lead to foaming.			

Plasmid DNA Purification, Continued

Column Capacity	Each column has a DNA binding capacity of up to 60 µg plasmid DNA.				
	For cell volumes $>2 \times 10^{9}$ cells, prepare cells as separate lysates of $\le 1-2 \times 10^{9}$ cells each as described below, and load lysates consecutively on the same column as described in Binding DNA , Step 3, below.				
Preparing	To prepare the bacterial cell lysate:				
Bacterial Cell Lysates	1.	In a microcentrifuge tube, pellet 1–3 ml $(1–2 \times 10^9)$ of <i>E. coli</i> cells from overnight cultures by centrifugation in a tabletop centrifuge at $1,500 \times g$ for 15 minutes. Remove the culture media completely.			
	2.	Completely resuspend the pellet in 240 μ l Resuspension Solution, prepared with RNase A as described on the previous page.			
	3.	Add 240 μ l Lysis Buffer to the above solution. Mix gently by inverting the tube 4–8 times.			
	4.	Incubate for 3-5 minutes at room temperature. Do not exceed 5 minutes.			
	5.	Add 340 μ l Neutralization/Binding Buffer, and immediately mix gently by inverting the tube 4-8 times.			
	6.	Centrifuge for 10 minutes at maximum speed in a tabletop centrifuge to clarify the cell lysate.			
	7.	Proceed to Binding DNA , below.			
Binding DNA	1.	Place a PureLink [™] spin column inside a 2-ml collection tube.			
	2.	Pipette or decant the supernatant from Step 6, above, into the spin column.			
	3.	Centrifuge the column at room temperature at $10,000-14,000 \times g$ for 1 minute. Discard the flowthrough, and place the column back in the tube.			
		If you are loading multiple samples on the same column, repeat Steps 2–3 for each lysate preparation.			
	4.	Add 650 μl Wash Buffer, prepared with ethanol as described on the previous page, to the column.			
	5.	Centrifuge the column at room temperature at $10,000-14,000 \times g$ for 1 minute. Discard the flowthrough from the collection tube, and place the column back in the tube.			
	6.	Centrifuge the column at maximum speed for 1–3 minutes to remove the residual wash buffer.			
	7.	Proceed to Eluting DNA, next page.			
		Continued on next page			

Plasmid DNA Purification, Continued

Eluting DNA	 Place the spin column in a clean 1.7-ml elution tube supplied with the kit. Add the following volume of Elution Buffer (10 mM Tris-HCl, pH 8.5) or sterile, distilled water (pH >7.0): Add 50 μl Elution Buffer or water to the center of the column if the expected DNA yield is <30 μg. Add 100 μl Elution Buffer or water to the center of the column if the expected DNA yield is >30 μg. Incubate the column at room temperature for 1 minute, then centrifuge at maximum speed for 1 minute. The elution tube contains your purified DNA. Remove and discard the column. Store the plasmid DNA at -20°C.
	Determine the quantity and quality of the DNA as described below.
Estimating DNA Yield and Quality	DNA Yield You can estimate the quantity of the purified plasmid DNA using UV absorbance at 260 nm or Quant-iT [™] DNA Assay Kits. <i>UV Absorbance</i>
	 Dilute an aliquot of the plasmid DNA sample in 10 mM Tris-HCl, pH 7.5. Mix well. Transfer to a cuvette (1-cm optical path length).
	 Measure the A₂₆₀ of the solution using a spectrophotometer blanked against 10 mM Tris-HCl, pH 7.5.
	 Calculate the amount of DNA using the following formula:
	DNA (μ g) = A ₂₆₀ × 50 μ g/(1 A ₂₆₀ x 1 ml) x dilution factor x total sample volume (ml) Assumption: For dsDNA, A ₂₆₀ = 1 for a 50 μ g/ml solution measured in a cuvette
	with an optical path length of 1 cm.
	Quant-iT™ DNA Assay Kits
	The Quant-iT DNA Assay Kits (page 160) provide a rapid, sensitive, and specific method for DNA quantitation with minimal interference from RNA, protein, or other common contaminants that affect UV absorbance readings.
	The kit contains a state-of-the-art quantitation reagent, pre-diluted standards for standard curve, and a ready-to-use buffer. The assay is performed in a microtiter plate format and is designed for reading in standard fluorescent microplate readers. Follow manufacturer's recommendations to perform the assay.
	DNA Purity
	Typically, DNA isolated using the PureLink TM HQ Mini Plasmid Purification Kit has an $A_{260}/A_{280} > 1.80$ when samples are diluted in Tris-HCl (pH 7.5). An A_{260}/A_{280} of >1.80 indicates that the DNA is reasonably clean of proteins that could interfere with downstream applications. Absence of contaminating genomic DNA and RNA may be confirmed by agarose gel electrophoresis.

Transfecting Cells

Introduction	This section provides general guidelines to transfect your pcDNA [™] 6.2- GW/EmGFP-miR expression construct into the mammalian cell line of interest to perform transient RNAi analysis. Performing transient RNAi analysis is useful to:
	Quickly test multiple miRNA sequences to a particular target gene
	Quickly screen for an RNAi response in your mammalian cell line
	• Test the effect of gene knockdown on your particular transient assay
	Once you have tested various miRNA target sequences using transient transfection, you may use the most efficient miRNA expression clone for further transient assays, generate stable transfectants, or transfer the optimal miRNA expression cassettes into suitable destination vectors for use in other RNAi applications (<i>e.g.</i> use of alternative promoters and/or viral transduction).
Experimental	To perform transfection, you will:
Outline	1. Grow the mammalian cell line of choice.
	2. Transfect the plasmid DNA into mammalian cells using Lipofectamine [™] 2000.
	3. Harvest cells after 18-48 hours to perform transient knockdown experiments.
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	Transfection efficiency
	Transcription rate of the target gene of interest
	Stability of the target protein
	Growth characteristics of your mammalian cell line
	 Activity of the promoter driving the miRNA expression cassette
	Take these factors into account when designing your RNAi experiments.
Lipofectamine [™] 2000 Reagent	Lipofectamine [™] 2000 Reagent (Ciccarone <i>et al.,</i> 1999) is included with BioModule [™] Units to transfect plasmid DNA into eukaryotic cells and offers the following advantages:
	• Provides the highest transfection efficiency in many mammalian cell types.
	• DNA-Lipofectamine [™] 2000 complexes can be added directly to cells in culture medium in the presence of serum.
	• Removal of complexes, medium change, or medium addition following transfection are not required, although complexes can be removed after 4-6 hours without loss of activity.
	An optimized transfection protocol for mammalian cells is described on page 67. However, depending on your cell type, you may need to optimize the transfection conditions such as cell number, amount of the transfection reagent and plasmid DNA, and the time period to assay for target gene knockdown to obtain the best results.
	Cell-type specific transfection protocols are available at <u>www.invitrogen.com/rnai</u> .

Opti-MEM [®] I	To facilitate optimal formation of DNA-Lipofectamine [™] 2000 complexes, we recommend using Opti-MEM [®] I Reduced Serum Medium supplied with the kit.
Positive and Negative Controls	If you have performed the positive control reaction and have cloned the miR-lacZ positive ds control oligo supplied with the kit into pcDNA TM 6.2-GW/EmGFP-miR, we recommend using the resulting pcDNA TM 6.2-GW/EmGFP-miR-lacZ expression construct as a positive control to assess the RNAi response in your cell line. Simply co-transfect the pcDNA TM 6.2-GW/EmGFP-miR-lacZ expression construct and the pcDNA TM 1.2/V5-GW/lacZ reporter plasmid supplied with the kit into your mammalian cells and assay for knockdown of β -galactosidase expression 24-48 hours post-transfection using Western blot analysis or activity assay. For more information about the pcDNA TM 1.2/V5-GW/lacZ reporter plasmid, recommendations for transfection, and methods to assay for β -galactosidase activity, see below.
	As negative control , perform parallel transfections with the pcDNA [™] 6.2-GW/EmGFP-miR-neg control plasmid.
pcDNA [™] 1.2/V5- GW/ <i>lacZ</i> Reporter Plasmid	The pcDNA [™] 1.2/V5-GW/ <i>lacZ</i> reporter plasmid is supplied with the kit for use as a positive control to assay for the RNAi response in your mammalian cell line. In this vector, β-galactosidase is expressed as a C-terminally tagged fusion protein under the control of the human cytomegalovirus (CMV) promoter (Andersson <i>et al.</i> , 1989; Boshart <i>et al.</i> , 1985; Nelson <i>et al.</i> , 1987). See page 146 for more information. The pcDNA [™] 1.2/V5-GW/ <i>lacZ</i> vector is supplied as 500 ng/µl of plasmid DNA in
	TE Buffer, pH 8.0. Dilute the stock as necessary for use in transfection (see below). If you wish to propagate the plasmid, transform a <i>recA</i> , <i>end</i> A <i>E</i> . <i>coli</i> strain such as TOP10. Use 10 ng of plasmid for transformation and select on LB agar plates containing 100 μ g/ml ampicillin.
Transfecting the <i>LacZ</i> -Containing Reagents	To perform RNAi analysis using the <i>lacZ</i> control reagents, you will co-transfect the pcDNA [™] 1.2/V5-GW/ <i>lacZ</i> reporter plasmid and pcDNA [™] 6.2-GW/EmGFP-miR- <i>lacZ</i> expression construct that you have generated into your mammalian cell line. For optimal results, we recommend using 6-fold more expression construct DNA than reporter plasmid DNA in the co-transfection.
	For example, use 600 ng of pcDNA [™] 6.2-GW/miR- <i>lacZ</i> DNA and 100 ng of pcDNA [™] 1.2/V5-GW/ <i>lacZ</i> DNA when transfecting cells plated in a 24-well format.
Assaying for β-galactosidase Expression	If you perform RNAi analysis using the control expression clone containing the <i>lacZ</i> ds oligo (pcDNA TM 6.2-GW/EmGFP-miR- <i>lacZ</i>), you may assay for β -galactosidase expression by western blot analysis using β -gal Antiserum (Cat. no. R901-25), by activity assay using cell-free lysates (Miller, 1972) and FluoReporter [®] <i>lacZ</i> /Galactosidase Quantitation Kit (Cat. no. F-2905), or by staining the cells for activity using the β -Gal Staining Kit (Cat. no. K1465-01) for fast and easy detection of β -galactosidase expression. For an example of results, see page 72.



The β -galactosidase protein expressed from the pcDNA^{\square}1.2/V5-GW/*lacZ* control plasmid is fused to a V5 epitope and is approximately 119 kDa in size. If you are performing Western blot analysis, you may also use the Anti V5 Antibodies available from Invitrogen (*e.g.* Anti-V5-HRP Antibody; Cat. no. R961-25 or Anti-V5-AP Antibody, Cat. no. R962-25) for detection. For more information, refer to our Web site (<u>www.invitrogen.com</u>) or call Technical Service (page 163).

- We recommend using Opti-MEM[®] I Reduced Serum Medium to dilute Lipofectamine[™] 2000 and plasmid DNA before complexing.
- **Do not** add antibiotics to media during transfection as this causes cell death.
- Use low-passage cells, and make sure that cells are healthy and greater than 90% viable before transfection.
- Maintain the same seeding conditions between experiments.
- **Transfect cells at 90-95% confluence.** Transfecting cells at a lower density allows a longer time interval to elapse between transfection and assay time, and minimizes the loss of cell viability due to cell overgrowth. Depending on the nature of the target gene, transfecting cells at higher densities may be suitable with optimization of conditions.
- Test serum-free media for compatibility with Lipofectamine[™] 2000 Reagent since some serum-free formulations (*e.g.* CD293, SFM II, VP-SFM) may inhibit cationic lipid-mediated transfection.

Materials NeededYou will need the following materials:•Mammalian cell line of interest cultured in the appropriate growth medium•Lipofectamine™ 2000 Reagent (supplied with the kit, store at 4°C until use)•Opti-MEM® I Reduced Serum Medium (supplied with the kit, pre-warm to
37°C before use)•pcDNA™6.2-GW/EmGFP-miR plasmid DNA•Appropriate tissue culture plates and supplies•Appropriate controls



Transfection Protocol	Use the following procedure to transfect plasmid DNA into mammalian cells in a 24-well format . All amounts and volumes are given on a per well basis.				
	For other formats, see Scaling Up or Down Transfections (next page). Prepare complexes using a DNA (µg) to Lipofectamine [™] 2000 (µl) ratio of 1:2 to 1:3 for most cell lines. Transfect cells at high cell density for high efficiency, high expression levels, and to minimize cytotoxicity.				
	Use this procedure as a starting point; optimize transfections as described in Optimizing Transfection (see below), especially if you are transfecting a mammalian cell line for the first time.				
	1. Adherent cells: One day before transfection, plate 0.5-2 x 10 ⁵ cells in 500 μl growth medium without antibiotics such that cells will be 90-95% confluent at the time of transfection.				
	Suspension cells: Just prior to preparing complexes, plate 4-8 x 10^5 cells in 500 µl of growth medium without antibiotics.				
	2. For each transfection sample, prepare complexes as follows:				
	a. Dilute DNA (0.8 µg) in 50 µl Opti-MEM [®] I Reduced Serum Medium without serum. Mix gently.				
	b. Mix Lipofectamine [™] 2000 gently before use, then dilute the appropriate amount (2 µl) in 50 µl Opti-MEM [®] I Medium. Incubate for 5 minutes at room temperature.				
	 Note: Combine diluted Lipofectamine[™] 2000 with diluted DNA within 30 minutes. c. After the 5 minute incubation, combine the diluted DNA with diluted Lipofectamine[™] 2000 (total volume = 100 µl). Mix gently and incubate for 20 minutes at room temperature (solution may appear cloudy). 				
	Note: Complexes are stable for 6 hours at room temperature.3. Add 100 µl of complexes to each well containing cells and medium. Mix gently by rocking the plate back and forth.				
	 Incubate cells at 37°C in a CO₂ incubator for 18-48 hours prior to testing for expression. Medium may be changed after 4-6 hours. 				
Optimizing Transfection	To obtain the highest transfection efficiency and low cytotoxicity, optimize transfection conditions by varying cell density as well as DNA and Lipofectamine [™] 2000 concentrations.				
	 Make sure that cells are greater than 90% confluent 				
	 Vary DNA (μg): Lipofectamine[™] 2000 (μl) ratios from 1:0.5 to 1:5 				
	 Depending on the nature of the target gene, transfecting cells at higher densities may also be considered when optimizing conditions. 				

Scaling Up or
DownTo transfect cells in different tissue culture formats, vary the amounts of
Lipofectamine[™] 2000, DNA, cells, and medium used in proportion to the relative
surface area, as shown in the table.
With automated, high-throughput systems, a complexing volume of 50 µl is

recommended for transfections in 96-well plates. **Note:** You may perform rapid 96-well plate transfections by plating cells directly into the transfection mix. Prepare complexes in the plate and directly add cells at twice the cell density as in the basic protocol in a 100 µl volume. Cells will adhere as usual in the presence of complexes.

Culture vessel	Surf. Area per well ¹	Vol. of plating medium	DNA (µg) in media volume (µl)	Lipofectamine [™] 2000 (µl) in media volume (µl)
96-well	0.3 cm ²	100 µl	0.2 μg in 25 μl	0.5 µl in 25 µl
24-well	2 cm^2	500 µl	0.8 μg in 50 μl	2.0 µl in 50 µl
12-well	4 cm^2	1 ml	1.6 µg in 100 µl	4.0 μl in 100 μl
6-well	10 cm ²	2 ml	4.0 μg in 250 μl	10 µl in 250 µl
60-mm	20 cm ²	5 ml	8.0 µg in 0.5 ml	20 µl in 0.5 ml
10-cm	60 cm ²	15 ml	24 µg in 1.5 ml	60 µl in 1.5 ml

¹Surface areas may vary depending on the manufacturer.

Detecting Fluorescence

Introduction	You can perform analysis of the EmGFP fluorescent protein from the express clone in either transiently transfected cells or stable cell lines (see next page). Once you have transfected your expression clone into mammalian cells, you detect EmGFP protein expression directly in cells by fluorescence microscop other methods that use light excitation and detection of emission. See below recommended fluorescence microscopy filter sets. The EmGFP expression is essentially 100% correlated with the expression of			ext page). cells, you may nicroscopy or ee below for
	miRNA.	Tis essentially 100%		ession of your
Filters for Use with EmGFP	The EmGFP can be detected with standard FITC filter sets. However, for optimal 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 protein such as the Omega XF100 filter set for fluorescence microscopy.			
	The spectral characteris	stics of EmGFP are li	sted in the table below:	
	Fluorescent Protein	Excitation (nm)	<u>Emission (nm)</u>	
	EmGFP	487	509	
	For information on obta (www.omegafilters.com (<u>www.chroma.com</u>).		s, contact Omega Optica ology Corporation	al, Inc.
Fluorescence Microscope	fluorescence microscop	e with FITC filter or	nGFP in cells using an ir Omega XF100 filter (ava n culture or a flow cytor	ailable from
Color Camera		Ve recommend using	is compatible with the n a digital camera or high	
Detecting Transfected Cells	fluorescence. Medium	can be removed and due to the medium. continue growing th		ng viewing to
What You Should See	signal that should be ea fluorescence signal of Em	asy to detect above the GFP from miRNA-cont	beled and emit a green f ne background fluoresce taining vectors is reduced non-miRNA containing vec	nce. Note: The due to
	functional miRNA. How the miRNA and demon	wever, cells with red astrate knockdown si wn are generally low	the highest knockdown luced fluorescence may since the expression level er than that required to ts, page 72.	still express s required to

Generating a Stable Cell Line

Introduction	Once you have determined that the miRNA in your pcDNA [™] 6.2-GW/EmGFP-miR expression clone is functional, you may wish to establish stable cell lines that constitutively express your miRNA. As negative control, establish cell lines expressing pcDNA [™] 6.2-GW/EmGFP-miR-neg control plasmid.
Blasticidin Selection	The pcDNA [™] 6.2-GW/EmGFP-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 are stably transfected with the pcDNA [™] 6.2-GW/EmGFP-miR construct.
	For more information about how to prepare and handle Blasticidin, see page 134.
Experimental	To generate stable cell lines, you will:
Outline	1. Grow the mammalian cell line of choice.
	2. Determine antibiotic sensitivity for your mammalian cell line.
	3. Perform transfection using Lipofectamine [™] 2000 Reagent.
	4. On the next day, replace the medium with fresh, complete medium containing Blasticidin.
	5. Continue to replace the medium with fresh, complete medium containing Blasticidin every 3-4 days until Blasticidin-resistant colonies are visible.
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 untransfected mammalian cell line (<i>i.e.</i> perform a kill curve experiment). Typically, concentrations ranging from 2-10 μ g/ml Blasticidin are sufficient to kill most untransfected mammalian cell lines. We recommend that you test a range of concentrations (see protocol below) to ensure that you determine the minimum concentration necessary for your cell line.
	 Plate cells at approximately 25% confluence. Prepare a set of 6 plates. Allow cells to adhere overnight.
	2. The next day, substitute culture medium with medium containing varying concentrations of Blasticidin (<i>e.g.</i> 0, 2, 4, 6, 8, 10 μ g/ml Blasticidin).
	 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.
	Continued on next page

Generating a Stable Cell Line, continued

Materials Needed	You will need the following materials:				
	 Mammalian cell line of interest (make sure that cells are healthy and > 90% viable before beginning) 				
	• pcDNA [™] 6.2-GW/EmGFP-miR expression clone				
	• pcDNA [™] 6.2-GW/EmGFP-miR-neg control plasmid				
	• Lipofectamine [™] 2000 Reagent (supplied with the kit, store at 4°C until use)				
	 Opti-MEM[®] I Reduced Serum Medium (supplied with the kit, pre-warm to 37°C before use) 				
	• Blasticidin (5 to 10 mg/ml)				
	Appropriate tissue culture dishes and supplies				
Guidelines for Transfection and Selection	Guidelines are provided below to transfect your pcDNA [™] 6.2-GW/EmGFP-miR expression clone and pcDNA [™] 6.2-GW/EmGFP-miR-neg control plasmid into the mammalian cell line of choice and to select for stable cell lines using Blasticidin.				
	 Perform transfection of your pcDNA[™]6.2-GW/EmGFP-miR expression construct and pcDNA[™]6.2-GW/EmGFP-miR-neg control plasmid into cells using the transfection protocol on page 67. Use separate wells for separate constructs. 				
	2. Four to six hours after transfection, remove the medium and replace with fresh growth medium. Incubate the cells overnight at 37°C.				
	3. The following day (Day 2), trypsinize and replate cells into a larger-sized tissue culture format in fresh complete medium containing the appropriate concentrations of Blasticidin.				
	Example: If transfecting cells in a 6-well format, trypsinize and replate cells into 10 cm tissue culture plates in medium containing Blasticidin.				
	4. Replace medium with fresh medium containing Blasticidin every 3-4 days until Blasticidin- resistant colonies can be identified (generally 10-14 days after selection).				
	5. Pick at least 10 Blasticidin-resistant colonies per construct and expand each clone.				
	6. Assay for target gene knockdown, compare to uninduced cells and cells stably transfected with pcDNA [™] 6.2-GW/EmGFP-miR-neg control plasmid.				
The Next Step	Once you have assessed the expression of your miRNA of interest using EmGFP fluorescence, we recommend that you assay for gene knockdown using qPCR, Western analysis, immunohistochemistry, or any other functional assay.				
	A variety of BioModule [™] Units that include qualified reagents and validated protocols are available from Invitrogen to perform validation experiments (page 160).				

Expected Results for miRNA Expression

20% -10% -0% -

Ctrl

lacZ

luc

Neg

Introduction	Examples of results obtained with pcDNA [™] 6.2-GW/ EmGFP-miR expression vector for miRNA expression are shown below.		
Knockdown of Reporter Gene	In this experiment, pcDNA [™] 6.2-GW/ EmGFP-miR expression vector containing ds oligo encoding miRNA targeting the <i>lacZ</i> , luciferase reporter gene or a negative control (neg) were generated following the recommended protocols and using the reagents supplied in the BioModule [™] BLOCK-iT [™] RNAi Units. Note that the miR- <i>lacZ</i> positive ds control oligo and negative control vectors used in this experiment are supplied with the kit.		
	GripTite [™] 293 MSR cells (Cat. no. R795-07) were grown to 90% confluence. Individual wells in a 24-well plate were transfected using Lipofectamine [™] 2000 Reagent with 100 ng of the pcDNA [™] 1.2/V5-GW/ <i>lacZ</i> reporter plasmid and co- transfected with 300 ng of the <i>lacZ</i> , luc or neg pre-miRNA expression vectors as indicated. Non-specific plasmid DNA was added to a total of 500 ng DNA. Cell lysates were prepared 48 hours after transfection and assayed for β-galactosidase activity using the FluoReporter [®] <i>lacZ</i> /Galactosidase Quantitation Kit (Cat. no. F-2905).		
	Results: Potent and specific inhibition of β -galactosidase activity is evident from the <i>lacZ</i> -derived miRNA and not from the luciferase-derived or negative control miRNA for the pcDNA ^{\mathbb{M}} 6.2-GW/EmGFP-miR vector.		
	110% 100% 90% 100% 1		

Continued on next page

Expected Results for miRNA Expression, Continued

Knockdown of HeLa cells transfected with pcDNA[™]6.2-GW/EmGFP-miR vector containing lacZ- (top panel) or lamin A/C-directed (bottom panel) miRNA inserts were Endogenous Lamin A/C fixed and stained four days after transfection. A single field of cells is shown in each row to reveal nuclei (Hoechst dye), EmGFP (auto-fluorescence), lamin A/C (immunofluorescence using mouse monoclonal antibody, sc-7292, from Santa Cruz Biotechnology), and the overlay of EmGFP and lamin A/C signal. Results: The lamin A/C-miRNA transfected cells in which EmGFP fluorescence is detectable show markedly reduced lamin A/C staining, indicating tight correlation between EmGFP expression and lamin knockdown. In the control *lacZ*-miRNA transfected wells, lamin signals in non-EmGFP and EmGFP expressing cells are similar, indicating no effect of the *lacZ*-miRNA on lamin A/C expression. EmGFP + Nuclei EmGFP Lamin A/C Lamin A/C **Detection:** Hoechst dye auto-fluorescence immunofluorescence overlay lacZ miRNA Lamin A/C miRNA

73

Transferring the Pre-miRNA Expression Cassette to Destination Vectors

Experimental Outline

The experimental steps necessary to transfer the pre-miRNA expression cassette from pcDNA[™]6.2-GW/EmGFP-miR expression clone into other Gateway[®] destination vectors are outlined below. We recommend performing the steps as described below to achieve the best results.

Step	Action	Page no.
1	Generate an expression clone.	36
2	Perform the Rapid BP/LR Gateway [®] recombination reactions to transfer the pre-miRNA expression cassette from an expression clone into a destination vector.	80
3	Transform the recombination reaction mix into One Shot [®] TOP10 Chemically Competent <i>E. coli</i> or One Shot [®] Stbl3 ^{M} Chemically Competent <i>E. coli</i> (for lentiviral destination vectors).	87
4	Analyze transformants and purify plasmid DNA for transfection.	54
5	Perform transient transfection using Lipofectamine [™] 2000 Reagent into a mammalian cell line of choice.	67
6	Perform analysis of the EmGFP fluorescent protein from the expression clone using fluorescent microscopy to determine the expression of your miRNA.	69
7	Assay for target gene knockdown using qPCR, Western analysis, immunohistochemistry, or any other functional assay.	
8	Optional:	
	Perform lentiviral expression of your miRNA, if you used a lentiviral destination vector.	92

Creating Entry Clones for Use with Destination Vectors

Introduction	The pcDNA [™] 6.2-GW/EmGFP-miR expression vector contains <i>att</i> sites to facilitate the transfer of the pre-miRNA expression cassette into appropriate Gateway [®] destination vectors to allow for expression of the miRNA in viral systems or using tissue-specific promoters. The pre-miRNA is transcribed by RNA Polymerase II (Pol II); the pre-miRNA expression cassette can be transferred to other Gateway [®] adapted destination vectors utilizing Pol II promoters to control expression of the pre-miRNA.
Important	Since the pcDNA [™] 6.2-GW/EmGFP-miR expression vector contains <i>att</i> B sites, the expression vector containing the pre-miRNA expression cassette cannot be used directly with a destination vector to perform the LR recombination reaction.
Transferring the Cassette	To transfer your pre-miRNA expression cassette into other destination vectors, you need to first generate an entry clone containing <i>att</i> L sites by performing a BP recombination reaction, then use the resulting entry clone in an LR recombination reaction with a destination vector to generate a new miRNA expression clone.
	The transfer of the miRNA sequence into the destination vector can be performed using the standard BP and LR recombination reactions or Rapid BP/LR recombination reactions as described on the next page.
	See below for an overview of the Gateway [®] recombination reactions and page 79 for the recombination region.
Gateway [®] Recombination Reactions	Two recombination reactions constitute the basis of the Gateway [®] Technology as described below. You will perform both Gateway [®] recombination reactions to transfer the pre-miRNA expression cassette from pcDNA [™] 6.2-GW/EmGFP-miR expression vector to a new destination vector as outlined below.
	BP Reaction
	Facilitates recombination of an <i>att</i> B substrate (linearized <i>att</i> B expression clone) with an <i>att</i> P substrate (donor vector) to create an <i>att</i> L-containing entry clone. This reaction is catalyzed by BP Clonase TM II enzyme mix.
	You will recombine pcDNA [™] 6.2-GW/EmGFP-miR expression clone (<i>att</i> B substrate) with an <i>att</i> P substrate (pDONR [™] 221 donor vector) first to form an entry clone.
	LR Reaction
	Facilitates recombination of an $attL$ substrate (entry clone) with an $attR$ substrate (destination vector) to create an $attB$ -containing expression clone. This reaction is catalyzed by LR Clonase TM II enzyme mix.
	The resulting entry clone (<i>att</i> L substrate) from the BP reaction is then recombined with the destination vector (<i>att</i> R substrate) to form a new miRNA expression clone.
	Continued on next page

Creating Entry Clones for Use with Destination Vectors,

Continued

Choosing aBased on your experimental needs, you may choose between the standard orSuitable ProtocolRapid BP/LR recombination reactions as described in the table below:

If You Wish to	Then Choose	Described
Generate the expression clone using a fast protocol but obtain fewer (~10% of the total number of clones) expression clones than the standard protocol	Rapid BP/LR Recombination Protocol	On page 80.
Maximize the number of expression clones generated and isolate entry clones for future use	Standard BP and LR Protocols	On page 135.

Substrates for the To perform a BP recombination reaction, you need the following substrates: Recombination • Linearized attB-containing expression clones (see page 78 for guidelines to Reactions linearize *att*B expression clones) *att*P-containing donor (pDONR[™]221) vector (see below) • To perform an LR recombination reaction, you need the following substrates: Supercoiled *att*L entry vector (pENTR[™]221/miR) Supercoiled *att*R destination vector (*e. g.*, pLenti6/V5-DEST) **Donor Vector** A large variety of donor vectors are available from Invitrogen. We recommend using the pDONR[™]221 vector. The pDONR[™]221 vector is supplied with the BioModule[™] BLOCK-iT[™] Unit with Lentiviral Pol II miR RNAi Expression System (Cat. no. WFGE08).

Creating Entry Clones for Use with Destination Vectors, Continued

Appropriate Destination Vectors	tion The various Gateway [®] vectors have widely different transcriptional and to		technical tom entiviral	
	A list of Gateway [®] destination vectors that are compatible with the pcDNA [™] 6.2- GW/EmGFP-miR expression vector is shown below. For more information on these destination vectors, visit www.invitrogen.com or contact Technical Service (page 163).			
	Destination Vector Catalog No.			
	Standard Destination Vectors			
	pLenti6/V5-DEST (included with WFGE08) V496-10			

Standard Destination Vectors	
pLenti6/V5-DEST (included with WFGE08)	V496-10
pLenti6/UbC/V5-DEST	V499-10
pEF-DEST51	12285-011
pT-REx [™] -DEST30	12301-016
pEF5/FRT/V5-DEST [™] (Flp-In [™])	V6020-20
N-terminal reporter tag vectors, e.g.:	
pcDNA [™] 6.2/nGeneBLAzer [™] -DEST	12578-068, 12578-050
pcDNA [™] 6.2/N-YFP-DEST	V358-20
MultiSite Gateway [®] Destination Vectors	
pDEST [™] /R4-R3	12567-023
pLenti6/R4R2/V5-DEST	K591-10

Transferring the pre-miRNA expression cassette from the pcDNA[™]6.2-GW/EmGFP-miR to the pLenti6/BLOCK-iT[™]-DEST destination vector will not yield a functional miRNA expression vector because this vector does not carry a Pol II promoter upstream of the *att*R1 site. **Expression of the pre-miRNA** requires the destination vector to supply a Pol II promoter.

For lentiviral expression, transfer to pLenti6/V5-DEST as described on page 80.

Resuspending the Donor and Destination Vector

Important

The donor and destination vectors are supplied as 6 μ g of supercoiled plasmid, lyophilized in TE Buffer, pH 8.0. To use, simply resuspend each plasmid DNA in 40 μ l sterile water to a final concentration of 150 ng/ μ l.

Creating Entry Clones for Use with Destination Vectors, Continued

Propagating the Donor and Destination Vectors	Donor Vector If you wish to propagate and maintain the pDONR [™] 221 vector supplied with cat. no WFGE08, we recommend using One Shot [®] <i>ccdB</i> Survival T1 ^R Chemically Competent <i>E. coli</i> or Library Efficiency [®] DB3.1 [™] Competent <i>E. coli</i> from Invitrogen (page 160) for transformation. The <i>ccdB</i> Survival T1 ^R and DB3.1 [™] <i>E. coli</i> strains are resistant to CcdB effects and can support the propagation of plasmids containing the <i>ccdB</i> gene. To maintain the integrity of the vector, select for transformants in media containing 50 µg/ml kanamycin and 15 µg/ml chloramphenicol.
	Destination Vector
	If you wish to propagate and maintain the pLenti6/V5-DEST vector supplied with Cat. no WFGE08, we recommend using Library Efficiency [®] DB3.1 [™] Competent <i>E. coli</i> from Invitrogen (page 160) for transformation. The DB3.1 [™] <i>E. coli</i> strain is resistant to CcdB effects and can support the propagation of plasmids containing the <i>ccdB</i> gene. To maintain integrity of the vector, select for transformants in media containing 50-100 µg/ml ampicillin and 15-30 µg/ml chloramphenicol. Do not use One Shot [®] <i>ccdB</i> Survival T1 ^R Chemically Competent <i>E. coli</i> to propagate lentiviral vectors as it is more susceptible to recombination and results in lower yields.
	Note: Do not use general <i>E. coli</i> cloning strains including Stbl3 [™] , TOP10, or DH5α for propagation and maintenance, as these strains are sensitive to CcdB effects.
Linearizing Expression Clones	We recommend that you linearize the pcDNA ^{m} 6.2-GW-EmGFP-miR expression clone using <i>Eag</i> I or <i>Bsr</i> D I.
	1. 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 pre-miRNA region of interest and is located outside the <i>att</i> B region.
	2. Ethanol precipitate the DNA after digestion by adding 0.1 volume of 3 M sodium acetate followed by 2.5 volumes of 100% ethanol.
	3. Pellet the DNA by centrifugation. Wash the pellet twice with 70% ethanol.
	 Dissolve the DNA in TE Buffer, pH 8.0 to a final concentration of 50-150 ng/µl.
Recombination Region	The recombination region of the lentiviral expression clone resulting from pLenti6/V5-DEST x pENTR [™] 221/EmGFP-miR entry clone is shown on the next page.
	For details on the recombination region of other destination vectors, refer to the manual supplied with each destination vector.
	Continued on next page

Creating Entry Clones for Use with Destination Vectors, Continued

Regior	 Combination egion of the recombination region of the lentiviral expression clone resulting from pLenti6/V5-DEST x pENTR[™]221/EmGFP-miR entry clone is shown below. The pENTR[™]221/EmGFP-miR entry clone is obtained by transferring the premiRNA expression cassette from pcDNA[™]6.2-GW/EmGFP-miR into pDONR[™]221. Features of the Recombination Region: Shaded regions correspond to those DNA sequences transferred from the pENTR[™]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 2473 and 3192 indicate the coding sequence of EmGFP. Note: The DNA sequences transferred from the pENTR[™]221/miR entry clone contain pre-miRNA expression cassette including EmGFP coding sequence. 		miR entry clone is shown below. is obtained by transferring the pre- 5.2-GW/EmGFP-miR into equences transferred from the the pLenti6/V5-DEST vector by rived from the pLenti6/V5-DEST ding sequence of EmGFP. pENTR [™] 221/miR entry clone contain the	
2251	TCGTAACAAC TCC		IV forward priming site	GGGAGGTCTA TATAAGCAGA GCTCGTTTAG
2331	TGAACCGTCA GA		CGCTGTTT TGACCTCCAT	AGAAGACACC GACTCTAGAG GATCCACTAG
			<i>att</i> B1	2473
2411	TCCAGTGTGG TGC		AGTTT <mark>GTA CAAAAAAGCA</mark> TCAAACAT GTTT <mark>TTTCGT</mark>	
	EmGFP coding sequence	EmGFP forward sequencing pr	0102	Met Val Ser Lys
2485	GGC CCG EmGFP Gly	GGC ATG GAC GAG CTG CCG TAC CTG CTC GAC Gly Met Asp Glu Leu	ATG TTC ATT CGA	AAGCACTTCG TGGCCGTCGA TCGTTTAAAG TTCGTGAAGC ACCGGCAGCT AGCAAATTTC
		5	5' miR flanking region	3' miR flanking region
3226		ICGACCAG TGGATCCTGG AG AGCTGGTC ACCTAGGACC TC		
3296		CATGGAAC AAATGGCCCA GA GTACCTTG TTTACCGGGT CT.		TAGATCTG GGT CGA AAG AAC ATG
3369				attB2 A GAG GGC CCG CGG TTC GAA GGT T CTC CCG GGC GCC AAG CTT <u>CCA</u>
3432	TTC GGA TAG G	GA TTG GGA GAG GAG CCA	GAG CTA AGA TGC GC.	T ACC GGT TAG TAA TGA GTTT A TGG CCA ATC ATT ACT
	V5	(C-term) reverse priming site	┘ V5 epitope	



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

Introduction	Follow the guidelines and instructions in this section to perform the Rapid BP/LR recombination reaction using the pcDNA [™] 6.2-GW/EmGFP-miR vector containing the pre-miRNA expression cassette, pDONR [™] 221, and a destination vector. See below for details on the Rapid protocol.			
	If you wish to perform the standard BP recombination reaction followed by the standard LR recombination reaction, see page 135.			
Rapid BP/LR Protocol	The Rapid BP/LR protocol is used to transfer a pre-miRNA expression cassette from an expression clone into a destination vector in 2 steps - a BP reaction using a donor vector followed by an LR reaction using a destination vector without purification of the intermediate entry clone.			
	Note: Using this protocol allows you to generate expression clones more rapidly than the standard BP and LR protocols provided in the Gateway [®] Technology with Clonase [™] II manual. 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, do not use this protocol. Use the standard BP and LR protocols on page 135.			
Note	This protocol is for experienced users of the Gateway [®] Technology. If you are unfamiliar with the Gateway [®] system, refer to the Gateway [®] Technology with Clonase [™] II manual available at www.invitrogen.com.			
Experimental	To perform the Rapid BP/LR protocol, you will:			
Outline	 Perform a BP recombination reaction using the linearized expression clone containing your pre-miRNA sequence and pDONR[™]221 to generate the entry clone. 			
	2. Use a small aliquot of the BP reaction mix to perform the LR recombination reaction using the destination vector to generate the new miRNA expression clone.			
	3. Perform Proteinase K treatment.			
	4. Transform the reaction mixture into a suitable <i>E. coli</i> host.			
	5. Select for expression clones.			
	Based on the destination vector that you are using, you need to perform the appropriate LR recombination reaction as described below.			
•	If you are using standard destination vectors (see page 77), perform the LR recombination reaction with LR Clonase™ II enzyme mix as described on page 84,			
	If you are using MultiSite Gateway [®] destination vectors (see page 77), perform the LR recombination reaction with LR Clonase[™] Plus enzyme mix as described on page 85.			

<i>E. coli</i> Host	Once you have performed the Rapid BP/LR recombination reaction, you will transform the recombination reaction into competent <i>E. coli</i> and select for the appropriate transformants. You may use any <i>recA</i> , <i>endA E. coli</i> strain including TOP10, DH5 α^{TM} , or equivalent for transformation.		
	DO NOT transform the 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. See page 160 for ordering information on competent cells.		
	Important: When performing the LR recombination reaction with the lentiviral destination vectors, transformation into the Stbl3 TM <i>E. coli</i> strain is recommended for optimal results (see below).		
Recommended <i>E. coli</i> Host for pLenti6/V5-DEST	For optimal results with pLenti6/V5-DEST, we recommend using Stbl3 [™] E. <i>coli</i> for transformation as this strain is particularly well-suited for use in cloning unstable DNA such as lentiviral DNA containing direct repeats. One Shot [®] Stbl3 [™] Chemically Competent <i>E. coli</i> are included with Cat. no WFGE08 for transformation. For instructions, see Transforming One Shot[®] Stbl3[™] Competent <i>E. coli</i> , page 87.		
Positive Control	We recommend using the pcDNA [™] 6.2-GW/miR-neg Control Plasmid supplied with the kit 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.		
	Do not use the pEXP7-tet supplied with the BP Clonase [™] II Enzyme Mix or pENTR [™] -gus supplied with the LR Clonase [™] II Enzyme Mix as positive controls for the Rapid protocol due to the presence of incompatible selection markers.		
Converting Femtomoles (fmol)	Use the following formula to convert femtomoles (fmol) of DNA to nanograms (ng) of DNA required for BP reaction:		
to Nanograms (ng)	$ng = (fmol)(N)(\frac{660 fg}{fmol})(\frac{1 ng}{10^6 fg})$		
	where N is the size of the DNA in bp. An example using <i>att</i> B-PCR product is described below.		
	In this example, you need to use 50 fmol of an <i>att</i> B-PCR in the BP reaction. The <i>att</i> B-PCR is 2.5 kb in size. Calculate the amount of <i>att</i> B-PCR required for the reaction (in ng) by using the above equation:		
	$(50 \text{ fmol})(2500 \text{ bp})(\frac{660 \text{ fg}}{\text{fmol}})(\frac{1 \text{ ng}}{10^6 \text{ fg}}) = 82.5 \text{ ng of product required}$		
	Continued on next page		

Materials Needed	You will need the following materials:
	• Linearized expression clone (50-150 ng/µl in TE Buffer, pH 8.0, see page 78)
	 pDONR[™]221 vector (supplied with Cat. no WFGE08, resuspend to 150 ng/µl in sterile water)
	 Destination vector including pLenti6/V5-DEST (150 ng/µl in TE Buffer, pH 8.0)
	 BP Clonase[™] II enzyme mix (supplied with the kit, store at -20°C until immediately before use)
	• LR Clonase [™] II enzyme mix (supplied with the kit for standard destination vectors) or LR Clonase [™] Plus enzyme mix (for MultiSite Gateway [®] destination vectors); store at -20°C until immediately before use
	 2 μg/μl Proteinase K solution (supplied with Clonase[™] enzymes; thaw and keep on ice until use)
	• TE Buffer, pH 8.0 (10 mM Tris-HCl, pH 8.0, 1 mM EDTA)
	• Appropriate competent cells (One Shot [®] Stbl3 [™] Competent <i>E. coli</i> are supplied with Cat. no. WFGE08 for transforming the pLenti6/V5-DEST constructs)
	Sterile 0.5 ml microcentrifuge tubes

Continued

Setting Up the Rapid BP/LR Recombination Reaction

Follow this procedure to perform the Rapid BP/LR reaction between your linearized expression clone, pDONR[™]221 vector, and the destination vector.

1. Add the following components to sterile 0.5 ml microcentrifuge tubes at room temperature and mix.

Component	Sample	Positive Control
Linearized <i>att</i> B expression clone from page 78, (75-180 ng)	1-7 µl	
pcDNA [™] 6.2-GW/EmGFP-miR-neg Control Plasmid (50 ng/µl)		2 µl
pDONR [™] 221 vector (150 ng/µl)	1 µl	1 µl
TE Buffer, pH 8.0	to 8 µl	5 μl

- 2. Remove the BP Clonase[™] II enzyme mix from -20°C and thaw on ice (~ 2 minutes).
- 3. Vortex the BP Clonase[™] II enzyme mix briefly twice (2 seconds each time).
- To the sample above, add 2 µl BP Clonase[™] II enzyme mix. Mix well by pipetting up and down.

Reminder: Return BP Clonase[™] II enzyme mix to -20°C immediately after use.

5. Incubate the reaction at 25°C for 1 hour.

Important: Unlike the standard BP reaction, **do not** add Proteinase K to the samples. Instead, proceed immediately to the next step.

6. Depending on the destination vector that you will use, perform the appropriate LR reaction as described on the following pages using an aliquot of the BP reaction mix that contains the resulting **entry clone**.

Note: Save the remaining BP reaction mix at -20°C for up to 1 week. You can treat the samples with Proteinase K and transform the reaction mix into One Shot[®] TOP10 Chemically Competent *E. coli* to check the efficiency of the BP reaction. This also allows you to isolate entry clones for future use.

LR Reaction for
Standard
Destination
Vectors

Use this LR recombination reaction for standard destination vectors. For LR recombination reaction with MultiSite Gateway[®] destination vectors, see next page.

- 1. Transfer 3 µl from each of the BP reaction from Step 5, previous page to clean, sterile 0.5 ml microcentrifuge tubes.
- 2. Add the following components to the microcentrifuge tubes containing 3 μ l BP reaction at room temperature and mix.

Component	Sample	Positive Control
Standard destination vector (150 ng/µl)	1 µl	1 µl
TE Buffer, pH 8.0	4 μl	4 μl

- 3. Remove the LR Clonase[™] II enzyme mix from -20°C and thaw on ice (~ 2 minutes).
- 4. Vortex the LR Clonase[™] II enzyme mix briefly twice (2 seconds each time).
- 5. To the samples above, add 2 µl LR Clonase[™] II enzyme mix. Mix well by pipetting up and down.

Reminder: Return LR Clonase[™] II enzyme mix to -20°C immediately after use.

6. 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.

7. Add 1 µl of the Proteinase K solution to each reaction. Incubate for 10 minutes at 37°C.

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

8. Transform an appropriate *E. coli* strain as recommended for your destination vector.

For pLenti6/V5-DEST, proceed to Transforming One Shot[®] Stbl3[™] Competent *E. coli*, next page.

9. Pick transformants and isolate plasmid DNA using PureLink[™] HQ Mini Plasmid Purification Kit. Perform restriction analysis to select a clone that has incorporated the pre-miRNA expression cassette (sequence analysis is not necessary).

LR Reaction for
MultiSite
Gateway®
Destination
Vectors

Use this LR recombination reaction for MultiSite Gateway[®] destination vectors. For LR recombination reaction with standard destination vectors, see previous page.

- 1. Transfer 6 µl from each of the BP reaction from Step 5, page 83, to clean, sterile 0.5 ml microcentrifuge tubes.
- 2. Add the following components to the microcentrifuge tubes containing 6 μ l BP reaction at room temperature and mix.

Component	Sample	Positive Control
MultiSite Gateway [®] Destination vector (60 ng/µl)	1 µl	1 µl
5' pENTR [™] vector (60 ng/μl)	1 µl	1 µl
3' pENTR [™] vector (60 ng/µl)	1 µl	1 µl
5X LR Clonase [™] Plus Buffer	3 µl	3 µl
TE Buffer, pH 8.0	4 µl	4 μl

3. Remove the LR Clonase[™] Plus enzyme mix from -20°C and thaw on ice (~ 2 minutes).

- 4. Vortex the LR Clonase[™] Plus enzyme mix briefly twice (2 seconds each time).
- 5. To the samples above, add 4 µl of LR Clonase[™] Plus enzyme mix. Mix well by pipetting up and down.

Reminder: Return LR Clonase[™] Plus enzyme mix to -20°C immediately after use.

- 6. Incubate the reaction at 25°C overnight.
- 7. Add 1 µl of the Proteinase K solution to each reaction. Incubate for 10 minutes at 37°C.
- 8. Transform an appropriate *E. coli* strain as recommended for your destination vector.

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

 Pick transformants and isolate plasmid DNA using PureLink[™] HQ Mini Plasmid Purification Kit. Perform restriction analysis to select a clone that has incorporated the pre-miRNA expression cassette (sequence analysis is not necessary).

What You ShouldWhen using One Shot® Stbl3™ or TOP10 Chemically Competent *E. coli* for
transformation, the LR recombination reaction should result in greater than 4,000
colonies if the entire LR reaction is transformed and plated.

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 and 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 $30 \ \mu\text{g/ml}$ chloramphenicol. A true expression clone should not grow in the presence of chloramphenicol.
Sequencing the expression construct is not required as transfer of the miRNA cassette from pcDNA [™] 6.2-GW/EmGFP-miR miRNA vector into the destination vector preserves the orientation of the cassette. However, you may sequence your expression construct using appropriate sequencing primers, if desired.
Once you have generated your expression clone, maintain and propagate the expression clone in LB medium containing the appropriate antibiotic.
Once you have obtained your new expression clone, we recommend that you test the clone by transfecting an appropriate mammalian cell line to perform transient RNAi analysis. Refer to the manual supplied with the destination vector to take advantage of the features of the new expression clone.
Once you have assessed the gene knockdown, we recommend that you validate the results using additional methods such as qPCR, Western analysis, immunohistochemistry, or any other functional assay.
A variety of BioModule [™] Units that include qualified reagents and validated protocols are available from Invitrogen to perform validation experiments (page 160).

Transforming One Shot[®] Stbl3[™] Competent *E. coli*

Introduction	fro E. c	low the instructions in this section to transform the LR recombination reaction m a lentiviral destination vector into One Shot [®] Stbl3 [™] Chemically Competent <i>coli</i> (Box 19) included with the kit. The transformation efficiency of One Shot [®] 13 [™] Chemically Competent <i>E. coli</i> is 1 x 10 ⁸ cfu/µg plasmid DNA.
Materials Needed	Yo	u will need the following materials:
	•	LR recombination reaction (page 84)
	•	One Shot [®] Stbl3 [™] Chemically Competent <i>E. coli</i> (supplied with the kit, Box 19; one vial per transformation; thaw on ice immediately before use)
	•	S.O.C. Medium (supplied with the kit, Box 19; warm to room temperature)
	•	pUC19 positive control (if desired to verify the transformation efficiency; supplied with the kit, Box 19)
	•	LB Medium (if performing the pUC19 control transformation)
	•	42°C water bath
	•	LB plates containing 100 μ g/ml ampicillin (two for each transformation; warm at 37°C for 30 minutes before use)
	•	37°C shaking and non-shaking incubator
One Shot [®] Stbl3 [™] Transformation		e this procedure to transform the LR recombination reaction into One Shot [®] l3™ Chemically Competent <i>E. coli</i> .
Procedure	1.	Thaw, on ice, one vial of One ${\rm Shot}^{\rm I\!S}$ ${\rm Stbl3^{\rm TM}}$ chemically competent cells for each transformation.
	2.	Add 2-3 µl of the LR recombination reaction (from Step 7, page 84 or Step 7, page 141) into a vial of One Shot [®] Stbl3 [™] cells and mix gently. Do not mix by pipetting up and down. For the pUC19 control, add 10 pg (1 µl) of DNA into a separate vial of One Shot [®] cells and mix gently.
	3.	Incubate the vial(s) on ice for 30 minutes.
	4.	Heat-shock the cells for 45 seconds at 42°C without shaking.
	5.	Remove the vial(s) from the 42°C water bath and place them on ice for 2 minutes.
	6.	Add 250 µl 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 μ 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 and plate 25-100 μ l.
	9.	Store the remaining transformation mix at +4°C. Plate out additional cells the next day, if desired.

Transforming One Shot[®] Stbl3[™] Competent *E. coli*, Continued

What You Should See	When using One Shot [®] Stbl3 [™] 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 and 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 $30 \ \mu\text{g/ml}$ chloramphenicol. A true expression clone should not grow in the presence of chloramphenicol.

Analyzing Transformants

Introduction	We recommend analyzing the transformants using restriction digestion as described below, even if you have observed a successful LR recombination. This allows you to confirm the presence of the insert as well as ensure the absence of any aberrant lentiviral vector recombination (between the LTRs). Depending on the sequences, lentiviral vectors have a tendency to undergo aberrant recombination (recombine or delete within regions such as between the LTRs), which can occur from time to time, even during vector propagation.
Materials Needed	You will need the following materials:
	Appropriate restriction enzymes
	• LB medium containing 100 µg/ml ampicillin (page 133 for recipe)
	 S.N.A.P.[™] MidiPrep Kit or PureLink[™] HQ Mini Plasmid Kit to isolate plasmid DNA (supplied with the kit)
	• 1.2% E-Gel [®] agarose gels (supplied with the kit)
	• TrackIt [™] 1 Kb Plus DNA Ladder (supplied with the kit)
	• TE Buffer
Restriction Digest Analysis	This protocol allows you to quickly analyze plasmid DNA isolated using S.N.A.P. [™] MidiPrep Kit from 2 transformants. If you wish to use a mini-prep kit for plasmid DNA isolation or screen more transformants, an alternate protocol is included on the next page.
	1. Grow 2 ampicillin- and Blasticidin-resistant colonies overnight in LB medium containing 100 μ g/ml ampicillin.
	2. Isolate plasmid DNA using S.N.A.P. [™] MidiPrep Kit (see page 93 for protocol).
	3. Perform restriction analysis to confirm the presence of the insert. Use a restriction enzyme or a combination of enzymes that cut once in the vector and once in the insert.
	4. Analyze the digests on 1.2% E-Gel [®] agarose gels (supplied with the kit) as follows:
	• To 10 µl of the digest, add 10 µl TE buffer and mix well.
	 Load 20 µl sample on a pre-run 1.2% E-Gel[®] agarose gels (see page 48 for pre-running protocol).
	 Mix 2 μl TrackIt[™] 1 Kb Plus DNA Ladder with 18 μl deionized water and load 20 μl of the diluted ladder. Loading undiluted TrackIt[™] DNA Ladder on an E-Gel[®] will result in loss of resolution.
	• Perform electrophoresis for 30 minutes using a E-Gel [®] PowerBase [™] v.4 as described on page 48.
	• At the end of the run, remove the gel cassette from the power unit and analyze your results using a UV transilluminator. Example of results are shown on page 91.
	5. Use the plasmid DNA from the positive clone for producing the lentivirus (page 105).

Analyzing Transformants, Continued

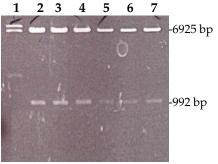
Alternate Protocol	This alternate protocol allow kit for plasmid DNA isolation	vs you to screen more transformants using a mini-prep on.	
	1. Isolate plasmid DNA us	sing PureLink™ HQ Mini Plasmid Kit (page 61).	
	2. Perform restriction dige	st analysis as described on the previous page.	
	this diluted DNA to retr Cells as described on pa	id DNA from the positive clone to 1:500 in TE. Use 1 µl of cansform into One Shot [®] Stbl3 [™] Chemically Competent ge 87.Plate approximately one-tenth of the ates containing 100 µg/ml ampicillin.	
		ter culture, and isolate plasmid DNA using S.N.A.P.™ ed on page 93 for use in lentivirus production (page 105).	
Sequencing	Sequencing the expression construct is not required as transfer of the miRNA cassette from pcDNA [™] 6.2-GW/EmGFP-miR into the pLenti6/V5-DEST vector preserves the orientation of the cassette. If you wish to sequence your pLenti6/V5 expression construct, we recommend using the following primers. Refer to the diagram on page 79 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'	
		vitrogen has a custom primer synthesis service. For more www.invitrogen.com) or call Technical Service (page 160).	
Maintaining the Expression Clone		pur expression clone, maintain and propagate the fum containing 100 μ g/ml ampicillin.	

Analyzing Transformants, Continued

Example of Expected Results with pLenti6/V5-DEST

In this experiment, MAP2 (microtubule-associated protein 2) microRNA was cloned into pcDNA[™]6.2-GW/EmGFP-miR as described in this manual. The resulting expression clone and the pcDNA[™]6.2-GW/EmGFP-miR-neg control plasmid (supplied with the kit) were each recombined with pLenti6/V5-DEST using the Rapid BP/LR Gateway[®] recombination reaction and transformed into One Shot[®] Stbl3[™] Chemically Competent cells as described in this manual to produce lentiviral expression clones.

The transformants were subjected to restriction digestion using *Xho* I and *Spe* I, followed by analysis on a 1.2% E-Gel[®] agarose gel as described in this section. **Results:** The gel analysis shows the correct digestion pattern indicating proper LR recombination with the lentiviral vector. No aberrant lentiviral vector recombination is observed.



Lane 1

Digest from pLenti6/V5-DEST Lanes 2-4 Digest from pcDNA[™]6.2-GW/EmGFP-miRneg control lentiviral expression clone Lanes 5-7 Digest from pcDNA[™]6.2-GW/EmGFP-miR-MAP2 lentiviral expression clone

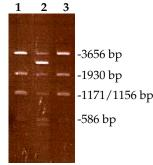
Example of Expected Results with pLenti6/R4R2/V5-DEST

In this experiment, MAP2 (microtubule-associated protein 2) microRNA was cloned into pcDNA[™]6.2-GW/EmGFP-miR as described in this manual. The resulting expression clone and the pcDNA[™]6.2-GW/EmGFP-miR-neg control plasmid (supplied with the kit) were each recombined with MultiSite pLenti6/R4R2/V5-DEST using the Rapid BP/LR Gateway[®] recombination reaction and transformed into One Shot[®] Stbl3[™] Chemically Competent cells as described in this manual to produce lentiviral expression clones.

The transformants were analyzed by restriction digestion using *Xho* I and *Afl* II. The digests were analyzed on a 1.2% E-Gel[®] agarose gel as described in this section.

Results: The gel analysis shows the correct digestion pattern indicating proper LR recombination with the lentiviral vector for lanes 1 and 3. Lane 2 shows additional bands indicating aberrant recombination of the lentiviral vector.

Lanes 1-3: Digest from pcDNA[™]6.2-GW/EmGFP-miR-MAP2 lentiviral expression clone



Lentivirus Production, Transduction, and RNAi Analysis

Experimental Outline

The experimental steps necessary to generate replication-incompetent lentivirus that delivers the miRNA sequence of interest into dividing or nondividing mammalian cells for RNAi analysis are outlined below. We recommend performing the steps as described below to achieve the best results.

Step	Action	Page no.
1	Purify plasmid DNA from the lentiviral expression clone using the S.N.A.P. [™] MidiPrep Kit.	93
2	Thaw and culture 239FT Cells to the required density.	96
3	Produce a lentiviral stock (containing the packaged pLenti6/V5 expression construct) by co-transfecting the optimized ViraPower [™] Packaging Mix and your pLenti6/V5-GW/miR expression construct into the 293FT Producer Cell Line.	101
4	Determine the titer of the lentiviral stock using Blasticidin selection or EmGFP detection.	108
5	Transduce the lentiviral construct into your mammalian cell to express the miRNA of interest.	114
6	Assay for target gene knockdown using qPCR, Western analysis, immunohistochemistry, or any other functional assay.	

Lentiviral Plasmid DNA Purification

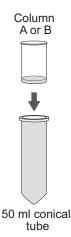
Introduction	Once you have generated your expression clone using pLenti6/V5-DEST, you must isolate plasmid DNA for transfection. Plasmid DNA for transfection into eukaryotic cells must be very clean and free from contamination with phenol and sodium chloride. Contaminants will kill the cells, and salt will interfere with lipid complexing, decreasing transfection efficiency. We recommend isolating lentiviral plasmid DNA using the S.N.A.P. [™] MidiPrep Kit supplied with Cat. no. WFGE08. Important: Do not use mini-prep plasmid DNA for transfection or PureLink [™] HQ Mini			
	Plasmid DNA Purification Kit for lentiviral plasmid DNA isolation.			
Experimental	To purify plasmid DNA from pLenti6/V5-DEST construct, you will:			
Outline	1. Lyse cells using the Lysis Buffer.			
	2. Precipitate protein and genomic DNA. Filter the lysate/precipitate through Column A.			
	3. Add Binding Buffer to the flowthrough and apply to Column B.			
	4. Wash the bound plasmid and dry the resin by centrifugation.			
	5. Elute the plasmid DNA with TE buffer.			
CAUTION	The Binding Buffer and Wash Buffer contain a chaotropic salt. Use gloves and protective eye wear when handling these solutions.			
Materials Needed	You will need the following materials:			
	• S.N.A.P. [™] MidiPrep Kit (Box 23 supplied with Cat. no WFGE08)			
	100 ml overnight culture			
	• 240 ml 95% ethanol for diluting 4X Final Wash			
	• 500 ml sterile bottle for diluting 4X Final Wash			
	Sterile 50 ml conical tubes			
	Centrifuge with rotor and rotor adapters for 50 ml conical tubesTE Buffer			
Before Starting	 Resuspend the entire contents of the RNase A tube (6 mg) in 1 ml of Resuspension Buffer. Add the resulting solution to the remaining Resuspension Buffer. Resuspension Buffer containing RNase A should now be stored at +4°C. 			
	• Add the 80 ml 4X Final Wash to 240 ml 95% ethanol to make 320 ml 1X Final Wash. Store in a 500 ml sterilized bottle.			
	• Check the Lysis Buffer for a white precipitate. If present, place the buffer in a 37°C water bath for 5 minutes or until the solution clears.			

Lentiviral Plasmid DNA Purification, Continued



Use the plasmid DNA isolation protocol described below to maximize the yield from lentiviral plasmids. We recommend using 50-100 ml overnight culture and process the 100 ml bacterial culture as two samples until DNA binding and then process the column as a single sample to obtain the best results.

Lysis and Removal of Precipitate



All centrifugations are at room temperature except as noted.

- 1. Pick 2 colonies and culture in LB medium containing $100 \mu g/ml$ ampicillin for 6-8 hours to obtain a starter culture. Transfer the starter culture to 100 ml LB medium containing $100 \mu g/ml$ ampicillin and culture overnight.
- 2. Transfer 100 ml overnight bacterial culture into two sterile 50 ml conical tubes.
- 3. Centrifuge the culture at 4000 x g for 5-10 minutes at 4°C to pellet the cells. Discard the medium. Process each tubes as two samples.
- 4. Resuspend the cell pellet in each tube in 4 ml Resuspension Buffer with RNase by vortexing or gently pipetting up and down.
- 5. Add 4 ml Lysis Solution to each tube and mix by inverting gently 5-6 times. Incubate for 3 minutes at room temperature.
- 6. Add 4 ml Precipitation Salt to each tube and invert gently 6-8 times. Incubate 5 minutes in ice. Invert tubes twice during incubation to ensure even formation of precipitate within the solution. Do not centrifuge.
- 7. During the incubation, label two, sterile 50 ml conical tubes "A" and one sterile 50 ml tube "B". Place one S.N.A.P.[™] MidiPrep Column A (Filtering) into each 50 ml conical tube labeled "A", and one S.N.A.P.[™] MidiPrep Column B (Binding) into the 50 ml conical tube labeled "B" (see adjacent figure).
- 8. After incubation, transfer each solution from Step 6 onto each Column A and centrifuge for 5 minutes at 3,000 x g. Discard Column A.

Save the filtrate containing the plasmid DNA.

9. Proceed to Plasmid Binding, next page.

Lentiviral Plasmid DNA Purification, Continued

Plasmid Binding	1.	To each filtrate from Step 8, previous page, add 12 ml Binding Buffer and mix by gently inverting twice.
	2.	Transfer the solution from one filtrate to Column B in a 50 ml conical tube.
	3.	Centrifuge for 2 minutes at 1,000 x g. Discard the flowthrough.
		The plasmid DNA from one sample is now bound to Column B.
	4.	Transfer the solution from the second filtrate to Column B in a 50 ml conical tube from Step 3.
	5.	Centrifuge for 2 minutes at 1,000 x g. Discard the flowthrough.
		The plasmid DNA from both samples is now bound to Column B.
	6.	Add 5 ml Wash Buffer to the column and centrifuge for 1 minute at 2,000 x g. Discard the flowthrough.
	7.	Add 5 ml 1X Final Wash Buffer to the column and centrifuge for 2 minutes at $2,000 \ge g$.
	8.	Add 10 ml 1X Final Wash Buffer to the column and centrifuge for 2 minutes at 2,000 x g. Discard the flowthrough.
	9.	Centrifuge the Column B at >4,000 x g for 5 minutes to dry the resin.
Plasmid Elution	1.	To elute the plasmid DNA, transfer Column B to a new, sterile 50 ml conical tube and add 750 μ l TE buffer directly to column resin.
	2.	Incubate for 3 minutes at room temperature.
		Note: Do not elute in less than 750 μl. Plasmid recovery will decrease.
	3.	Centrifuge for 5 minutes at >4,000 x g. The plasmid DNA is now eluted from Column B.
	4.	Remove and discard the column.
	5.	Determine plasmid DNA yield and purity as described on page 63. Ensure that the A_{260}/A_{280} of the purified plasmid DNA is >1.8 to obtain good transfection efficiency.

Growth and Maintenance of the 293FT Cell Line

Introduction	The human 293FT Cell Line is supplied with the BioModule [™] BLOCK-iT [™] Lentiviral Pol II miR RNAi Expression System 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 [®] . General guidelines for handling the cells and protocols for thawing, sub culturing, and freezing the cells are included in this section.
CAUTION	Handle as potentially biohazardous material under at least Biosafety Level 2 containment. This product contains Dimethyl Sulfoxide (DMSO), a hazardous material. Review the Material Safety Data Sheet before handling.
General Cell	Follow the general guidelines below to grow and maintain 293FT cells.
Handling	• Make sure that all solutions and equipment that come in contact with the cells are sterile. Always use proper sterile technique and work in a laminar flow hood.
	• Before starting experiments, be sure to establish the cells and have some frozen stocks. We recommend using early-passage cells for your experiments.
	• For general maintenance of cells, pass 293FT cells when they are > 80% confluent. Avoid overgrowing cells before passaging.
	• Maintain 293FT cells in complete medium containing 500 µg/ml Geneticin [®] .
	• Use trypan blue exclusion to determine cell viability. Log phase cultures should be >90% viable.
	• When thawing or subculturing cells, transfer cells into medium warmed to room temperature.
	• Cells should be at the appropriate confluence and at greater than 90% viability prior to transfection.
Materials Needed	You will need the following materials:
	• 15 ml sterile, conical tubes
	Appropriate sized tissue culture flasks and pipettes
	Complete medium (next page)
	• 50 mg/ml Geneticin [®]
	Phosphate-Buffered Saline (PBS; Invitrogen, Catalog no. 10010-023)
	Reagents for counting cells
	Trypsin/versene (EDTA) solution or other trypsin solution
	Freezing Medium (next page)
	Table-top centrifuge
	Cryovials (if needed)

Media for 293FT Cells		The table below lists the recommended complete medium, freezing medium, and antibiotic concentration required to maintain and culture the 293FT Cell Line. Note: FBS does not need to be heat-inactivated for use with the 293FT Cell Line.				
	The cell culture media are available from Invitrogen individually or you may purchase the BioModule [™] Lentiviral 293 Unit from Invitrogen (page 160) that contains all necessary media, antibiotics, serum, and Trypan Blue for growth, maintenance, and checking cell viability of 293 FT cells.					
		Complete Medium	[Antibiotic]]	Freezing Medium	
	D	-MEM (high glucose)	500 µg/ml	90% co	omplete medium	
	10	% fetal bovine serum (FBS)	Geneticin®	10% DMSO	MSO	
		1 mM MEM Non-Essential mino Acids (NEAA)				
	2 :	mM L-glutamine				
	1:	mM MEM Sodium Pyruvate				
	1%	% Pen-Strep (optional)				
			1			
Preparing Medium	Prepare the complete D-MEM medium containing 10% FBS supplemented with 0.1 mM MEM Non-Essential Amino Acids and 2 mM L-glutamine as described below using the BioModule [™] Lentiviral 293 Unit from Invitrogen (page 160):					
	Perform all steps in a tissue culture hood under sterile conditions.					
	1. Remove 100 ml D-MEM from 1 L D-MEM bottle and replace with 100 ml FBS.					
	2.	To the bottle of medium, add	d the following:		-	
		200 mM L-Glutamine (100X)			10 ml	
		10 mM MEM Non-Essential Amino Acids (100X) 10 ml				
		100 mM MEM Sodium Pyru			10 ml	
	•	Optional: Penicillin-Streptom			10 ml	
	3. Filter sterilize the medium using 0.45μ m filtration device.					
	 Store the complete medium at 4°C until use. The medium is stable for 6 months at 4°C (avoid introducing any contamination into the medium). 					
	5.	To an aliquot of the complete medium with 500 µg/ml Ge		neticin®	• to prepare complete	
					Continued on next page	
					1 6	

Thawing Cells	Freezing Medium. Store	The 293FT Cell Line is supplied in a vial containing 3 x 10 ⁶ cells in 1 ml of Freezing Medium. Store frozen 293FT cells in liquid nitrogen until ready to use. Use the following procedure to thaw 293FT cells to initiate cell culture. Thaw cells			
	÷ .	medium without Geneticin [®] .			
	 Remove the vial of fr 37°C water bath. 	ozen cells from liquid nitrogen and thaw quickly in a			
	vial with 70% ethano	The completely thawed, decontaminate the outside of the ol, and transfer the cells to a sterile 15 ml tube containing ge the cells at 150-200 x g and resuspend the cells in 2 ml ithout Geneticin [®] .			
	3. Transfer the cells to Twithout Geneticin [®] .	0 1			
	 Incubate the flask ov bottom of the flask. 	0 0			
	<i>y</i> 1	 The next day, aspirate off the medium and replace with fresh, complete medium containing 500 µg/ml Geneticin[®]. 			
	6. Incubate the cells and				
	7. Proceed to Subculturing Cells , next page.				
	We recommend subcultu before use in other applic	uring cells for a minimum of 3 passages after thawing cations.			
Subculturing Conditions	Use the following recom procedure to subculture	mended conditions to subculture 293FT cells. For a cells, see below.			
	Parameter	Recommended Condition			
	Cell density	$> 5 \times 10^5$ viable cells/ml (> 80% confluent)			
	Culture vessel	T-75 cm ² to T-162 cm ² disposable sterile T-flasks. Dilute cells in a total working volume of 15-20 ml for T-75 cm ² flasks and 40-50 ml for T-162 cm ² flasks			
	Seeding density	2-5 x 10 ⁴ viable cells/cm ²			
	Incubation conditions	37° C incubator with a humidified atmosphere of $5-10\%$ CO ₂ in air; loosen caps to allow for oxygenation/aeration			

Determining Viability and Cell Density	ollow the procedure below to determine viable and total cell counts using the ypan blue exclusion method. Trypan Blue Stain (0.4%) is included with the ioModule [™] Lentiviral 293 Unit (page 160).	
	rypan blue is a vital dye. The chromophore is negatively charged and does not iteract with the cell unless the membrane is damaged. Therefore, cells that iteract with the cell unless the membrane is damaged. Therefore, cells that iteract with the dye are viable while cells that absorb the dye (blue cells) are non- able.	
	Transfer a small aliquot of the cell suspension to a microcentrifuge tube and dilute the cells such that the total number of cells counted will not be less than 100 or more than 1,000.	
	To 1 ml of the diluted cell suspension, add 100 µl Trypan Blue Stain (0.4%) solution. Gently aspirate to mix.	
	Record the dilution factor. The dilution factor equals the total volume (amount of cell suspension and amount of trypan blue) divided by the amount of cell suspension.	
	Incubate the cells with the trypan blue solution for 1-2 minutes.	
	Count all cells (including the blue cells) using a Coulter Counter or manually using a hemocytometer chamber.	7
	To calculate the total cells per ml in suspension, multiply the total count by the dilution factor.	
	To determine the viability, count only the blue cells. Calculate the % viability	7:
	[1.00 - (Number of blue cells ÷ Number of total cells)] x 100	
	ell viability should be at least 95% for healthy log-phase cultures.	
Subculturing Cells	se this procedure to subculture 293FT cells grown in a T-75 cm ² flask. If you are sing other-sized flasks, scale the reagent volumes accordingly.	ò
	Remove all medium from the flask and wash the cells once with 10 ml PBS to remove excess medium and serum. Serum contains inhibitors of trypsin.)
	Add 2 ml of trypsin/versene (EDTA) solution to the monolayer and incubate 1-5 minutes at room temperature until cells detach. Check the cells under a microscope and confirm that most of the cells have detached. If cells are still attached, incubate a little longer until most of the cells have detached.	j
	Add 8 ml complete medium containing Geneticin [®] and transfer the cell suspension to a 15 ml sterile, conical tube.	
	Determine viable and total cell counts (see procedure above).	
	Seed cells at the recommended density (see table on previous page), diluting in pre-warmed complete medium containing 500 μ g/ml Geneticin [®] . Incubate flasks as recommended (see table on previous page).	
	Maintain cells as adherent monolayer cultures in complete medium containing 500 μg/ml Geneticin [®] .	
	For the transfection protocol, you will need 6 x 10^6 293FT cells for each sample (page 105).	
	<i>a</i>	_

Freezing Cells	Once you have established the cells, we recommend freezing some cells for future use as described below.
	• Freeze cells at a density of at least 3 x 10 ⁶ viable cells/ml.
	• Use a freezing medium composed of 90% complete medium and 10% DMSO. Prepare freezing medium immediately before use . Filter-sterilize the freezing medium and store at +4°C until use. Discard any remaining freezing medium after use.
	Before starting, label cryovials and prepare freezing medium (see above). Keep the freezing medium on ice.
	1. Culture the desired quantity of 293FT cells to 70-90% confluency.
	 Remove the cells from the tissue culture flask(s) following Steps 1-3, Subculturing Cells, page 99.
	3. Determine viable and total cell counts (see procedure on previous page) and calculate the volume of freezing medium required to yield a final cell density of $\ge 3 \times 10^6$ cells/ml.
	4. Prepare the required volume of freezing medium (see above).
	5. Centrifuge the cells suspension (from Step 2) at 250 x g for 5 minutes in a table top centrifuge at room temperature. Carefully aspirate off the medium and resuspend the cell pellet in the pre-determined volume of chilled freezing medium.
	 Dispense aliquots of this suspension (frequently mixing to maintain a homogeneous cell suspension) into cryovials according to manufacturer's specifications.
	 Freeze cells in an automated, controlled-rate freezing apparatus or using a manual method following standard procedures. For ideal cryopreservation, the freezing rate should be decrease at 1°C per minute.
	8. Transfer vials to liquid nitrogen storage.
	Note: You may check the viability and recovery of frozen cells 24 hours after storing cryovials in liquid nitrogen by following the procedure outlined in Thawing Cells , page 98.
Transfection Methods	The 293FT Cell Line is generally amenable to transfection using standard methods including lipid-mediated transfection with Lipofectamine [™] 2000. See page 105 for transfection protocol.
Transient or Stable Transfection	The 293FT Cell Line may be transiently transfected with any plasmid. Make sure that cells are healthy at the time of plating. Overgrowth of cells prior to passaging can compromise their transfection efficiency.
	293FT cells can be used as hosts to generate a stable cell line expressing your gene of interest from most plasmids. The introduced plasmid must contain a selection marker other than neomycin resistance. Stable cell lines can then be generated by transfection and dual selection with Geneticin [®] and the appropriate selection agent.
	Note: Since 293FT cells stably express the SV40 large T antigen, we do not recommend generating stable cell lines with plasmids that contain the SV40 origin of replication.

Producing Lentivirus in 293FT Cells

Introduction	Before you can create a stably transduced cell line expressing your miRNA, you need to produce a lentiviral stock (containing the packaged pLenti6/V5 expression construct) by co-transfecting the optimized ViraPower [™] Packaging Mix and your pLenti6/V5-GW/miR expression construct into the 293FT Producer Cell Line. The following section provides protocols and instructions to generate a lentiviral stock.		
Experimental	To produce lentivirus in 293FT Cells, you will:		
Outline	1. Grow the 293FT Cells to obtain 6×10^6 293FT cells for each sample (page 96).		
	2. Prepare plasmid DNA of your expression clone (page 93).		
	 Cotransfect the ViraPower[™] Packaging Mix and pLenti6/V5-GW/miR expression plasmid DNA into 293FT Cells using Lipofectamine[™] 2000. 		
	4. Harvest virus-containing supernatants 48-72 hours post-transfection.		
	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.		
	• Subculture and maintain cells as recommended on page 96. Do not allow cells to overgrow before passaging. You will need 6 x 10 ⁶ 293FT cells for each sample.		
	• 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 your pLenti6/V5-GW/miR expression vector following cotransfection into 293FT producer cells.		
	The amount of the packaging mix (195 µg) and Lipofectamine [™] 2000 Reagent (0.75 ml) supplied in the BioModule [™] BLOCK-iT [™] Lentiviral Pol II miR RNAi Expression System is sufficient to perform 20 cotransfections in 10 cm plates using the recommended protocol on page 105.		
	To use the ViraPower TM Packaging Mix, resuspend the contents of one tube (195 μ g) in 195 μ l of sterile water to obtain a 1 μ g/ μ l stock.		
	Note: ViraPower [™] Packaging Mix is available separately from Invitrogen (page 160) or as part of the ViraPower [™] Bsd Lentiviral Support Kit (page 160).		

miR Positive Control	You may generate a miR Positive Control using the reagents included in the kit as follows:
	• Generate the pcDNA [™] 6.2-GW/EmGFP-miR- <i>lacZ</i> expression control using the <i>lacZ</i> double-stranded oligo and pcDNA [™] 6.2-GW/EmGFP-miR expression vector included with the BLOCK-iT [™] Pol II miR RNAi Expression Vector Kit and as described on page 36.
	• Use the pcDNA [™] 6.2-GW/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/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 (page 117) 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 ^{M} 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 149.
	To propagate and maintain the control plasmid:
	1. Resuspend the vector in 10 μ l of sterile water to prepare a 1 μ g/ μ l stock solution.
	 Use the stock solution to transform a <i>recA</i>, <i>endA E</i>. <i>coli</i> strain like Stbl3[™], TOP10, DH5α[™]-T1^R, or equivalent. Use 10 ng of plasmid for transformation.
	 Select transformants on LB agar plates containing 100 µg/ml ampicillin (for Stbl3[™] cells) or LB agar plates containing 100 µg/ml ampicillin and 50 µg/ml Blasticidin (for TOP10 or DH5α).
	4. Prepare a glycerol stock of a transformant containing plasmid for long-term storage. Propagate the plasmid in LB containing $100 \ \mu g/ml$ ampicillin.
	5. Use the pLenti6/V5-GW/ <i>lacZ</i> positive control to generate a control lentiviral stock (expressing the LacZ protein).
	6. Use the pLenti6/V5-GW/ <i>lacZ</i> lentiviral control and the pLenti6/V5-GW/EmGFP-miR- <i>lacZ</i> lentiviral control in cotransduction experiments as a positive control for lentiviral induced RNAi analysis in your system (see page 115 for details).

Materials Needed	You will need the following materials:	
	• pLenti6/V5-GW/miR expression cons Resuspend the purified pLenti6/V5-G	struct (0.1-3.0 μ g/ μ l in TE Buffer, pH 8.0). W/miR expression plasmid in TE Buffer, g from 0.1-3.0 μ g/ μ l. You will need 3 μ g of ection.
	• Positive controls (see previous page to sterile water to a concentration of 1 µg	generate positive controls; resuspend in //µl)
	• ViraPower [™] Packaging Mix (supplied sterile water to a concentration of 1 μg	-
	• 293FT cells (page 96), you will need 6	x 10 ⁶ 293FT cells for each sample.
		ent (supplied with the kit; store at +4°C
		n (supplied with the kit, pre-warmed; see
	• Fetal bovine serum (FBS)	
	• Complete growth medium for 239FT C	Cell (see page 97 for a recipe).
	• Sterile, 10 cm tissue culture plates (one controls)	e each for lentiviral construct and
	• Sterile, tissue culture supplies	
	• 5 and 15 ml sterile, capped, conical tub	Des
	Cryovials	
Recommended Transfection Conditions	We produce lentiviral stocks in 293FT cells transfection conditions below. The amoun recommended conditions at a titer of 1×10^{6} is generally sufficient to transduce 1×10^{6} infection (MOI) = 1.	t of lentivirus produced using these 0 ⁵ to 1 x 10 ⁷ transducing units (TU)/ml
	Condition	Amount
	Tissue culture plate size	10 cm (one per lentiviral construct)
	Number of 293FT cells to transfect	6 x 10 ⁶ cells (see Recommendation on page 101 to prepare cells for transfection)
	Amount of ViraPower [™] Packaging Mix	9 μ g (9 μ l of 1 μ g/ μ l stock)
	Amount of pLenti6/V5-GW/miR expression plasmid	3 µg
	Amount of Lipofectamine [™] 2000 Reagent to use	36 μl
	Note: You may produce lentiviral stocks using	other tissue culture formats, but

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



The recommended procedure to co-transfect 293FT cells differs from the traditional Lipofectamine[™] 2000 transfection procedure in that you will:

- First prepare DNA:Lipofectamine[™] 2000 complexes and add them to plates containing growth media, then
- Add the 293FT cells to the media containing DNA:Lipofectamine[™] 2000 complexes, 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[™] 2000 transfection procedure (i.e. plating cells first followed by transfection with DNA:Lipofectamine[™] 2000 complexes). You may use the traditional Lipofectamine[™] 2000 transfection procedure, if desired, but keep in mind that the viral titer obtained may be lower (see **Alternative Transfection Procedure**, page 106).

Transfection Procedure	a n	low the procedure below to cotransfect 293FT cells. We recommend including egative control (no DNA, no Lipofectamine [™] 2000) in your experiment to help luate your results.
	1.	For each transfection sample, prepare DNA-Lipofectamine [™] 2000 complexes as follows:
		a. In a sterile 5 ml tube, dilute 9 μg ViraPower [™] Packaging Mix and 3 μg pLenti6/V5-GW/miR expression plasmid DNA (12 μg total) in 1.5 ml of Opti-MEM [®] I Medium without serum. Mix gently.
		b. In a separate sterile 5 ml tube, mix Lipofectamine [™] 2000 gently before use, then dilute 36 µl in 1.5 ml of Opti-MEM [®] 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 [™] 2000. Mix gently.
		d. Incubate for 20 minutes at room temperature to allow the DNA-Lipid 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 1.2×10^6 cells/ml in growth medium containing serum (or Opti-MEM [®] I Medium containing serum).
	3.	Add the DNA-Lipofectamine [™] 2000 complexes to a 10 cm tissue culture plate containing 5 ml of growth medium containing serum (or Opti-MEM [®] I Medium containing serum). Do not include antibiotics in the medium.
	4.	Add 5 ml of the 293FT cell suspension (6 x 10^6 total cells) to the plate containing media and DNA-Lipofectamine TM 2000 complexes and mix gently by rocking the plate back and forth. Incubate the cells overnight at 37°C in a CO ₂ incubator.
	5.	The next day, remove media containing the DNA-Lipofectamine [™] 2000 complexes and replace with complete culture medium containing sodium pyruvate (<i>i.e.</i> D-MEM containing 10% FBS, 2 mM L-glutamine, 0.1 mM MEM Non-Essential Amino Acids, 1% penicillin/streptomycin, and 1 mM MEM Sodium Pyruvate).
		Note: Expression of the VSV glycoprotein causes 293FT cells to fuse, resulting in the appearance of multinucleated syncitia which is normal and does not affect lentivirus production. See page 107 for an example of expected results after transfection.
	6.	Harvest virus-containing supernatants 48-72 hours post-transfection by removing medium to a 15 ml sterile, capped, conical tube.
		Note: Minimal differences in viral yield are observed whether supernatants are collected 48 or 72 hours post-transfection.
		Caution: Remember that you are working with infectious virus at this stage. See pages 28 and 110.
	7.	Centrifuge at 3000 rpm for 5 minutes at +4°C to pellet cell debris. Perform filtration step, if desired (see Note on the next page).
	8.	Pipet viral supernatants into cryovials in 1 ml aliquots. Store viral stocks at -80°C.
		Continued on next page

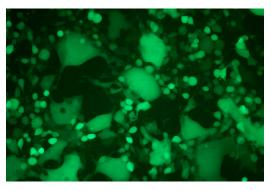
Alternative Transfection Procedure	An alternative transfection procedure is provided below to cotransfect 293FT cells. Note that use of this procedure generally results in production of lentiviral stocks with a slightly lower titer that those produced when using the Transfection Procedure , previous 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> 6×10^6 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 [®] I Medium containing serum). Do not include antibiotics in the medium.			
	3. Prepare DNA-Lipofectamine [™] 2000 complexes as instructed in the Transfection Procedure , Step 1, previous page.			
	 Add the DNA-Lipofectamine[™] 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. 			
	5. Follow Steps 5-8 as instructed in the Transfection Procedure , previous page.			
Note	If you plan to use your lentiviral construct for <i>in vivo</i> applications, we recommend filtering your viral supernatant through a sterile, 0.45 μm 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-HV 0.45 μm PVDF filters (Millipore, Catalog no. SLHVR25LS) 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, we recommend re-titering your viral stocks before transducing your 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 ² flask and harvested up to 30 ml of viral supernatant. If you wish to scale up your cotransfection, remember that you will need to increase the number of cells plated and the amounts of DNA, Lipofectamine [™] 2000, and medium used in proportion to the difference in surface area of the culture vessel.			

Example of Expected Results

In this experiment, 293FT Cells were co-transfected with ViraPower[™] Packaging Mix and a lentiviral expression plasmid expressing the cycle 3 GFP gene (pLenti6/V5-GW/cycle 3 GFP) using the transfection protocol on page 105 with Lipofectamine[™] 2000 Reagent.

48 hours post-transfection, the cells were subjected to fluorescence microscopy to visualize cycle GFP fluorescence using suitable filters.

Result: Cycle 3 GFP expression is observed in most cells indicating an efficient transfection. Multinucleated syncitia (fused producer cells) are seen due to VSV glycoprotein expression.



Titering Your Lentiviral Stock

Introduction	Before proceeding to transduce the mammalian cell line of interest and express 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 Guidelines and protocols are provided in this section.			
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)			
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 the mammalian cell line of choice in the presence of Polybrene ^{®.}			
	3. Based on the titering method used:			
	• Select for stably transduced cells using Blasticidin. Stain and count the number of Blasticidin-resistant colonies in each dilution.			
	• Determine the titer by flow cytometry 4 days post-transduction, if using EmGFP.			
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 for more information).			
	• 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, we recommend titering or re-titering 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 your lentiviral stock. Lentiviral stocks should be aliquotted and stored at -80°C. 			
	Continued on next page			

.

Selecting a Cell Line	 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: Grows as an adherent cell line Easy to handle Exhibits a doubling time in the range of 18-25 hours Non-migratory We generally use the HT1080 human fibrosarcoma cell line (ATCC, Catalog no. CCL-121) for titering purposes. 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.
Note	The titer of a lentiviral construct may vary depending on which cell line is chosen. 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/EmGFP-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 BioModule [™] BLOCK-iT [™] Lentiviral Pol II miR RNAi Expression System, Blasticidin is supplied with the kit. Blasticidin is also available separately from Invitrogen or as part of the ViraPower [™] Bsd Lentiviral Support Kit (see page 161 for ordering information). For more information about how to prepare and handle Blasticidin, and determine the Blasticidin sensitivity, refer to page 134.
Using Polybrene [®] During Transduction	Transduction of lentivirus into mammalian cells may be enhanced if cells are transduced in the presence of hexadimethrine bromide (Polybrene [®]). For best results, we recommend performing transduction in the presence of Polybrene [®] . Note, however, that some cells are sensitive to Polybrene [®] (<i>e.g.</i> primary neurons). Before performing any transduction experiments, you may want to test your cell line for sensitivity to Polybrene [®] . If your cells are sensitive to Polybrene [®] (<i>e.g.</i> exhibit toxicity or phenotypic changes), do not add Polybrene [®] during transduction. In this case, cells should still be successfully transduced.

Preparing and	Follow the instructions below to prepare Polybrene® (Sigma, Catalog no. H9268):			
Storing Polybrene [®]	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.			
	Note: The working stock may be stored at +4°C for up to 2 weeks.			
Materials Needed	You will need the following materials:			
	• pLenit6/V5-GW/EmGFP-miR 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 [®] , if desired			
	6-well tissue culture plates			
	 Blasticidin (10 mg/ml stock) and crystal violet (Sigma, Catalog no. C3886; prepare a 1% crystal violet solution in 10% ethanol), if you are using Blasticidin selection for titering 			
	• Inverted fluorescence microscope and appropriate filters for EmGFP visualization (see page 27 for filters), if you are using EmGFP titering method			
	• PBS (Cat. no. 10010-023)			
Preparing Mammalian Cells	Initiate your mammalian cell line of choice that will be used for titering. Grow cells in the appropriate medium. You will use at least one 6-well plate for every lentiviral stock to be titered (one mock well plus five dilutions). Cells should be >95% viable.			
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 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			

• Wear gloves, a laboratory coat, and safety glasses or goggles when handling viral stocks and media containing virus.

Transduction and	Foll
Titering Procedure	mar

ollow the procedure below to determine the titer of your lentiviral stock using the nammalian cell line of choice.

Note: If you have generated a lentiviral stock of the pLenti6-V5-GW/EmGFP-miR-*lacZ* control construct, perform titering using the Blasticidin or EmGFP method, and if you generated a lentiviral stock of the pLenti6-V5-GW/*lacZ* control construct, use 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 usually plate 2×10^5 cells/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 10⁻² to 10⁻⁶. 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 (10⁻² to 10⁻⁸), 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[®] (if desired) to each well to a final concentration of $6 \mu g/ml$. 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 **EmGFP titering method** or proceed to Steps 9-14 for **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.

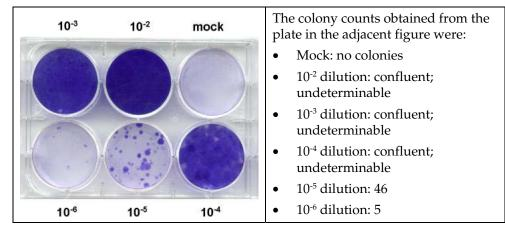
Preparing Cells for Flow Cytometry	If you have used EmGFP titering method, prepare cells for flow cytometry according to the established protocols in use at your flow cytometry facility. Refer to page 27 for spectral characteristics of EmGFP. The steps below provide general guidelines, and other methods may be suitable.			
	1. At day 4 post-transduction, disso trypsin or cell dissociation buffer.	ciate the cells from the plate by usi	ing	
	resuspend the cell pellet in flow c calcium/magnesium free PBS wit analysis on your flow cytometer. analysis, but may be done, if desi cytometry, use 2% formaldehyde or p PBS. However, these fixatives may in	ove residual media components an ytometry buffer such as th 1% FBS at the required density for Fixing the cells is not necessary for red. Note: To fix your cells before flow paraformaldehyde in calcium/magnesis crease autofluorescence of cells, thus i lls as a negative control for flow cytom	for r v ium free t is critical	
		d the lowest dilution of virus (<i>i.e.</i> 7 s, respectively, to set up the param		
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). This is to avoid analyzing dilution samples containing multiple integrated lentiviral genomes, which may result in an underestimate of the viral titer, or dilution samples containing too few transduced cells, which will give inaccurate results. Titer is expressed as transducing units (TU)/ml. Use the following formula to calculate the titer: $[F \times C/V] \times D$ F = frequency of GFP-positive cells (percentage obtained divided by 100) C = total number of cells in the well at the time of transduction V = volume of inoculum in ml D = lentivirus dilution 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%		
	In the above example, the 10^{-4} dilution percentage of EmGFP-positive cells far frequency of EmGFP-positive cells is number of cells in the well) divided b calculation is as follows: [(0.044 × 2)]	alls into the desired range of $1-30\%$ $4.4/100 = 0.044$, multiplied by $2 \times$	₀. The 10⁵ (the	

The lentiviral titer for this example is $8.8\times 10^7\, TU/ml.$

What You Should See	When titering pLenti6/V5 lentiviral stocks using HT1080 cells, we generally obtain titers ranging from 5 x 10^5 to 2 x 10^7 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 1×10^5 TU/ml, we recommend producing a new lentiviral stock. See page 101 and the Troubleshooting section, page 121 for more tips and guidelines to optimize your viral yield.

Example of Expected Results Using Blasticidin Titering Method

In this experiment, a pLenti6 lentiviral stock was generated using the protocol on page 105. HT1080 cells were transduced with 10-fold serial dilutions of the lentiviral supernatant (10^{-2} to 10^{-6} dilutions) or untransduced (mock) following the protocol in this manual. Forty-eight hours post-transduction, the cells were placed under Blasticidin selection ($10 \ \mu g/ml$). After 10 days of selection, the cells were stained with crystal violet (see plate below), and colonies were counted.



Thus, the titer of this lentiviral stock is 4.8 x 10⁶ TU/ml (*i.e.* average of 46 x 10⁵ and 5 x 10⁶).

Transduction and Analysis

Introduction	Once you have generated a lentiviral stock with a suitable titer, you are ready to transduce the lentiviral construct into your mammalian cell line to express the miRNA of interest and perform RNAi analysis. Guidelines are provided below. Reminder: Remember that your 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 [®] .		
	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 Levels	 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: 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.		
	Cautional an and and		

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 134. 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	A number of factors can influence determination of an optimal MOI including the nature of your mammalian cell line (<i>e.g.</i> non-dividing vs. dividing cell type; see Note below), its transduction efficiency, and the nature of your target gene of interest. If you are transducing your lentiviral construct into the mammalian cell line of choice for the first time, we recommend using a range of MOIs (<i>e.g.</i> 0, 1, 5, 10, 50) to determine the MOI required to obtain the optimal degree of target gene knockdown.			
Note	In general, non-dividing cell types transduce lentiviral constructs 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 an optimal degree of target gene knockdown.			
Preparing Mammalian Cells	Initiate your mammalian cell line of choice that will be used for transduction. Grow the cells in the appropriate medium. Cells should be >95% viable.			
Positive Controls	If you have generated two positive control lentiviral constructs (pLenti6/V5-GW/EmGFP-miR- <i>lacZ</i> control and pLenti6/V5-GW/ <i>lacZ</i> control constructs) as described on page 102, 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 119.			
	The β -galactosidase protein expressed from the pLenti6/V5-GW/ <i>lacZ</i> control lentiviral construct is approximately 121 kDa in size. You may assay for β -galactosidase expression by western blot analysis using β -gal Antiserum (Catalog no. R901-25), activity assay FluoReporter [®] <i>lacZ</i> /Galactosidase Quantitation Kit (Catalog no. F-2905), or by staining the cells for activity using the β -Gal Staining Kit (Catalog no. K1465-01) for fast and easy detection of β -galactosidase expression.			

Q 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 their transducibility. If the titer of your lentiviral stock is relatively low (less than 5×10^5 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	 You will need the following materials before starting: Your 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®, if desired Appropriately sized tissue culture plates for your application 10 mg/ml Blasticidin stock (if selecting for stably transduced cells) 			

Transduction Procedure		llow the procedure below to transduce the mammalian cell line of choice with ur lentiviral construct.
	1.	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).
	2.	On the day of transduction (Day 1), thaw your 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. DO NOT vortex.
	3.	Remove the culture medium from the cells. Mix the medium containing virus gently by pipetting and add to the cells.
	4.	Add Polybrene [®] (if desired) to a final concentration of 6 μ g/ml. Swirl the plate gently to mix. Incubate at 37°C overnight.
		Note: 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.
	6.	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.
	7.	Replace medium with fresh medium containing Blasticidin every 3-4 days until Blasticidin-resistant colonies can be identified (generally 10-12 days after selection).
	8.	Pick at least 5 Blasticidin-resistant colonies (see Note on the next page) and expand each clone to assay for knockdown of the target gene.
		Continued on next page

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. We recommend testing at least 5 Blasticidin-resistant clones and selecting the clone that provides the optimal degree of gene knockdown for further studies.				
Performing RNAi Analysis	You may use any method as appropriate to assay for knockdown of your target gene including functional analysis, immunofluorescence, western blot, or qRT-PCR with the appropriate LUX [™] primers. For more information about LUX [™] primers, see our Web site (<u>www.invitrogen.com/lux</u>). A variety of BioModule [™] Units that include qualified reagents and validated protocols are available from Invitrogen to perform validation experiments (page 160).				
	You can also detect EmGFP fluorescence using fluorescence microscopy or flow cytometry (see page 69).				
What You Should See	When performing RNAi studies using pLenti6/V5 lentiviral constructs, we generally observe inhibition of gene expression within 48-120 hours after transduction. The degree of gene knockdown depends on the time of assay, stability of the protein of interest, and on other factors listed on page 114. 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 next page.				

Expected Results for Lentiviral Expression

Example of Expected Results

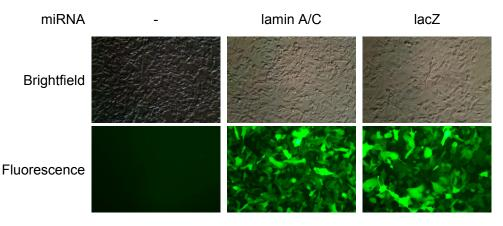
In this experiment, a double-stranded miR-lacZ control oligo (directed towards exogenous LacZ) and a double-stranded miR-lamin oligo (directed towards endogenous lamin) were each cloned into pcDNA6.2[™]-GW/EmGFP-miR expression vector using the BLOCK-iT[™] Pol II miR RNAi Expression Vector Kit. The miR-lacZ 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.

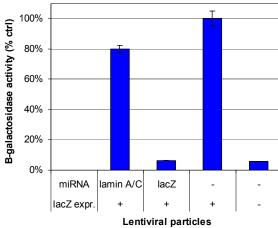
Figure A



Cell lysates were then prepared from duplicate wells 48 hours (*i.e.* 2 days) after transduction. β -galactosidase activity was determined from equivalent amounts of cell lysate using the FluoReporter[®] *lacZ*/Galactosidase Quantitation Kit (Cat. no. F-2905) from Invitrogen. Results are shown in figure B, next page.

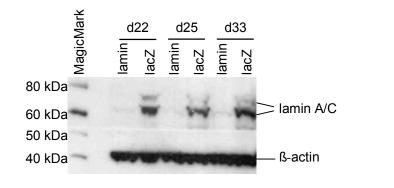
Expected Results for Lentiviral Expression, Continued

Example of Expected Results, continued
Figure B depicts a dramatic decrease in β -galactosidase activity only in cells cotransduced with the *lacZ*-directed miRNA and not lamin-directed miRNA.



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 (1:1000 dilution, BD Biosciences, Cat. no. 612162) and the other half was probed with Anti-β-Actin Antibody (1:5000 dilution, Abcam, Cat. no. ab6276). The blot was developed using WesternBreeze[®] Chemiluminescent Kit available from Invitrogen. 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 expression, lentiviral expression, and RNAi experiments.

Annealing Reaction

The table below lists some potential problems and possible solutions that may help you troubleshoot the annealing reaction.

Problem	Reason	Solution
Weak band representing ds oligo observed on an agarose gel	Single-stranded oligos designed incorrectly	Verify that the sequence of the bottom strand oligo is complementary to the sequence of the top strand oligo. If not, re-synthesize the bottom strand oligo.
	Allowed oligos to cool at +4°C instead of room temperature during annealing procedure	After heating to 95°C, anneal the oligos by setting the microcentrifuge tube at room temperature for 5-10 minutes (see the procedure on page 44).
	Did not anneal equal amounts of top and bottom strand oligo	Anneal equal amounts of the top and bottom strand oligo using the procedure on page 44.
No band representing ds oligo observed on an agarose gel	Single-stranded oligos designed incorrectly	Verify that the sequence of the bottom strand oligo is complementary to the sequence of the top strand oligo. If not, re-synthesize the bottom strand oligo.
	Used the wrong single-stranded oligos	Make sure that you mix single-stranded oligos with complementary sequence.

Ligation and Transformation Reactions

The table below lists some potential problems and possible solutions that may help you troubleshoot the ligation and transformation procedures.

Problem	Reason	Solution
Few spectinomycin- resistant colonies obtained on the selective plate	Single-stranded oligos designed incorrectly	Make sure that each single-stranded oligo contains the 4 nucleotides on the 5' end required for cloning into pcDNA [™] 6.2- GW/EmGFP-miR:
		Top strand oligo: include TGCT on the 5' end.
		Bottom strand oligo: include CCTG on the 5' end.
	ds oligos were degraded	Store the 10 nM ds oligo stock in 1X Oligo Annealing Buffer.
		Avoid repeated freeze/thaw cycles. Aliquot the 10 nM ds oligo stock and store at -20°C.

Problem	Reason	Solution
Few spectinomycin- resistant colonies obtained on the selective plate, continued	ds oligos stored incorrectly	Store the ds oligo stocks at -20°C.
	500 nM ds oligo stock solution diluted into water instead of 1X Oligo Annealing Buffer	To dilute the 50 μ M ds oligo reaction:
		Dilute the 50 µM stock 100-fold into DNase/RNase-free water to generate a 500 nM stock.
		Dilute the 500 nM stock 50-fold into 1X Oligo Annealing Buffer to generate a 10 nM stock. Use the 10 nM stock for cloning.
	10 nM ds oligo stock solution heated above room temperature prior to use	Thaw ds oligo stock solution on ice or at +4°C prior to use.
		Important: Diluted ds oligos will melt and form intramolecular hairpins if heated above room temperature. These hairpins will not clone into pcDNA [™] 6.2-GW/EmGFP-miR.
	Incorrect vector: insert ratio used in ligation reaction	Dilute the 50 µM ds oligo mixture as instructed on page 45 to generate a 10 nM
	Forgot to dilute annealed ds oligo 1:5,000 before use	stock. Use the 10 nM ds oligo stock for cloning.
	Annealed ds oligo diluted incorrectly	
	Ligation reaction not adequately mixed or incorrectly mixed prior to incubation	Mix the ligation reaction well by pipetting up and down.
		Note: Flicking the tube is not adequate to mix the reagents.
		Do not vortex the ligation reaction.
	Did not use the 5X Ligation Buffer supplied with the kit	Use the T4 DNA Ligase and 5X Ligation Buffer supplied with the kit for ligation as these reagents have been optimized to facilitate 5-minute ligation at room temperature.
		Important: Other T4 DNA Ligase and ligation buffers may not support 5-minute, room temperature ligation.
	Ligation reaction not incubated for long enough	Extend the incubation time of the ligation reaction up to 2 hours at room temperature.
	Ligation reaction incubated overnight at 16°C	The ligation conditions used to clone the ds oligo into pcDNA [™] 6.2-GW/EmGFP-miR differ from traditional ligation conditions. Incubate the ligation reaction at room temperature for 5 minutes.

Ligation and Transformation Reactions, continued

Problem	Reason	Solution
Few spectinomycin- resistant colonies obtained on the selective plate, continued	Not enough transformation mixture plated	Increase the amount of the transformation mixture plated.
	Selective plates contained too much spectinomycin	Use LB agar plates containing 50 µg/ml spectinomycin for selection.
	Did not use the competent cells supplied with the kit	Use the One Shot [®] TOP10 Chemically Competent <i>E. coli</i> supplied with the kit; trans- formation efficiency is > 1 x 10^9 cfu/µg DNA.
	Not enough of the ligation reaction transformed	Increase the amount of ligation reaction transformed.
	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 bacterial culture for 1 hour at 37°C with shaking before plating.
Many clones contain inserts with sequence	Poor quality single-stranded oligos used	Use high-quality single-stranded oligos.
mutations	Oligo preparation contains mutated sequences	Use mass spectrometry to check for peaks of the wrong mass.
	Oligo preparation contains contaminants	Order HPLC or polyacrylamide gel (PAGE)- purified oligos.
		Order oligos from Invitrogen's custom primer synthesis service (see <u>www.invitrogen.com</u> for details).
	Did not use the competent cells supplied with the kit	Use One Shot [®] TOP10 Chemically Competent <i>E. coli</i> supplied with the kit; transformation efficiency is $> 1 \times 10^9$ cfu/µg DNA.
Poor sequencing results	Loss of sequencing signal in the hairpin region due to secondary	Use high-quality, purified plasmid DNA for sequencing.
	structure formation	Add DMSO to the sequencing reaction to a final concentration of 5%.
		Increase the amount of template used for sequencing (up to twice the normal amount).
		Use a 7:1 molar ratio of dITP:dGTP in your sequencing reaction.
		Cut your construct with <i>Msc</i> I and purify before sequencing (see page 55).
No colonies obtained on the selective plate	Used the wrong antibiotic for selection	Select for transformants on LB agar plates containing 50 μ g/ml spectinomycin.

Ligation and Transformation Reactions, continued

Transfection and RNAi Analysis

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

Problem	Reason	Solution
Low levels of gene knockdown observed due to low transfection efficiency	Antibiotics added to the media during transfection Cells not healthy or too sparse at the time of transfection Not enough plasmid DNA transfected Not enough Lipofectamine [™] 2000 Reagent used	Do not add antibiotics to the media during transfection. Use healthy cells. Plate cells such that they will be 90-95% confluent at the time of transfection. Increase the amount of plasmid DNA transfected. Optimize the transfection conditions for your cell line by varying the amount of Lipofectamine [™] 2000 Reagent used.
Low levels of gene knockdown observed (other causes)	Didn't wait long enough after transfection before assaying for gene knockdown	Repeat the transfection and wait for a longer period of time after transfection before assaying for gene knockdown. Perform a time course of expression to determine the point at which the highest degree of gene knockdown occurs.
	ds oligo insert in your pcDNA™6.2-GW/EmGFP-miR construct contains mutations	When analyzing spectinomycin-resistant transformants, sequence the ds oligo insert to verify its sequence. Select constructs containing the correct ds oligo insert for use in RNAi analysis.
	miRNA sequence not optimal due to the target region selected	Select a different target region.

Problem	Reason	Solution
Cytotoxic effects observed after transfection	Too much Lipofectamine [™] 2000 Reagent used	Optimize the transfection conditions for your cell line by varying the amount of Lipofectamine [™] 2000 Reagent used.
	Plasmid DNA not pure	Prepare purified plasmid DNA for transfection. We recommend using the PureLink [™] HQ Mini Plasmid Purification Kit as described on page 61.
	Targets an essential gene	Make sure that your target gene is not essential for cell viability or growth.
No gene knockdown	miRNA with no activity chosen	Select a different target region.
observed	pre-miRNA designed incorrectly	Follow the guidelines on pages 37 to select the target sequence and design the single-stranded oligos.
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	Be sure to use the recommended filter sets for detection of fluorescence and use an inverted fluorescence microscope for analysis. If desired, allow the protein expression to continue for additional days before assaying for fluorescence.
		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.

Transfection and RNAi Analysis, continued

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 and 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 135.
control gave colonies	BP recombination reaction is treated with Proteinase K	Do not treat the BP reaction with Proteinase K before the LR reaction.
	Did not use the suggested Clonase [™] II enzyme mixes or Clonase [™] II enzyme mixes were inactive	• Make sure to store the BP and LR Clonase [™] II enzyme mix at -20°C or -80°C.
		 Do not freeze/thaw the BP and LR Clonase[™] II enzyme mix more than 10 times.
		• Use the recommended amount of BP and LR Clonase [™] II enzyme mix (page 80).
		• Test another aliquot of the Clonase [™] II enzyme mix.
	Not enough LR reaction transformed	Transform 2-3 μl of the LR reaction into One Shot [®] Stbl3 [™] 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 the recommended amount of BP reaction for the LR reaction.
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 [®] Stbl3 [™] Chemically Competent <i>E. coli</i> supplied with the kit for transformation of lentiviral constructs. Stbl3 [™] <i>E. coli</i> are recommended for cloning unstable DNA including lentiviral DNA containing direct repeats and generally give rise to fewer unwanted recombinants.

Problem	Reason	Solution
Few or no colonies obtained from the	Competent cells stored incorrectly	• Store the One Shot [®] Chemically Competent <i>E. coli</i> at -80°C.
transformation control		• Thaw a vial of One Shot [®] cells on ice immediately before use.
	After addition of DNA, competent cells mixed by pipetting up and down	After adding DNA, mix competent cells gently. Do not mix by pipetting up and down.

Rapid BP/LR Reaction and Transformation, continued

Generating the 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) Unhealthy 293FT cells; cells exhibit low viability Cells transfected in media containing antibiotics (<i>i.e.</i> Geneticin[®]) Plasmid DNA:transfection reagent ratio incorrect 293FT cells plated too sparsely 	 Do not use plasmid DNA from a miniprep for transfection. Use S.N.A.P.[™] MidiPrep Kit to prepare plasmid DNA. Use healthy 293FT cells under passage 20; do not overgrow. Do not add Geneticin[®] in the media during transfection as this reduces transfection efficiency and causes cell death. Use a DNA (in μg):Lipofectamine[™] 2000 (in μl) ratio ranging from 1:2 to 1:3. 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 105).
	Transfected cells not cultured in media containing sodium pyruvate Lipofectamine [™] 2000 Reagent	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. • Store at +4°C. Do not freeze.
	handled incorrectly	 Store at +4°C. Do not freeze. Mix gently by inversion before use. Do not vortex.

Problem	Reason	Solution
Low viral titer	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	Do not 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 109.
	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 [®] not included during titering procedure	Transduce the lentiviral construct into cells in the presence of Polybrene [®] .
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 [®] not included during transduction	Transduce the lentiviral construct into cells in the presence of Polybrene [®] .
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

Transduction and RNAi Analysis

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

Problem	Reason	Solution
Low levels of gene knockdown observed	 Low transduction efficiency: Polybrene[®] not included during transduction 	• Transduce the lentiviral construct into cells in the presence of Polybrene [®] .
	 Non-dividing cell type used 	 Transduce your lentiviral construct into cells using a higher MOI.
	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.
		Note: 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 111 before use.
	Viral stock stored incorrectly	• Aliquot and store stocks at -80°C.
		• Do not freeze/thaw more than 3 times.
		• If stored for longer than 6 months, re-titer stock before use.
	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.
		Note: 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 its 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	Target gene is essential for cell viability	Make sure that your target gene is not essential for cell viability or growth.
transduction	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 [®] used during transduction	Verify the sensitivity of your cells to Polybrene [®] . If cells are sensitive, omit the Polybrene [®] 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	Be sure to use the recommended filter sets for detection of fluorescence and use an inverted fluorescence microscope for analysis. If desired, allow the protein expression to continue for additional 1-3 days before assaying for fluorescence.
		Note: 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.

Transduction and RNAi Analysis, continued

Troubleshooting, Continued

E-Gel[®] Analysis

The table below provides solutions to some problems that you may encounter with E-Gel[®] agarose gels.

Problem	Cause	Solution	
No current	Copper contacts in the base are damaged due to improper use	Make sure the copper contacts in the base are intact.	
	Expired or defective gel cassette	Use fresh gel cassette. Use properly stored gels before the specified expiration date.	
	E-Gel [®] cassette is not inserted properly into a base	Remove cassette and reinsert; a steady red light illuminates on the base when the cassette is correctly inserted and power is on.	
	Incorrect adaptor used	Use only UL Listed Class 2 Direct Plug-in Adaptor included with the E-Gel [®] PowerBase [™] .	
Poor resolution or smearing of	Sample is overloaded	Load the recommended amount of sample on the gel (page 47). Less DNA is required since E-Gel [®] agarose gels are thinner.	
bands	Very low volume of sample loaded or sample was not	Avoid introducing bubbles while loading the samples. Bubbles will cause band distortion.	
	loaded properly	Load the recommended sample volume based on the gel type and loading method.	
		For proper band separation, we recommend keeping sample volumes uniform. Load deionized water or TE into any empty wells.	
	Gel was not electrophoresed immediately after sample	For best results, run the gel within 15 minutes of sample loading.	
	loading	If you cannot run the gel immediately after sample loading, use the Two-Step Loading method (refer to the E-Gel [®] Technical Guide available at www.invitrogen.com).	
	Expired gel used	Use properly stored gels before the expiration date.	
	Longer electrophoresis run time or high current during the run	Longer run times cause an increase in the current, resulting in poor band migration or a melted gel. Do not run the gel longer than the recommended time for each gel type.	
Sample	Sample is overloaded	Load the recommended sample volume per well.	
leaking from		Use the Two-Step Loading method.	
the wells	Wells damaged during comb removal	Remove the comb gently without damaging the wells.	
Failure Mode indicated by a steady red light and continuous rapid beeping	Defective cassette	Disconnect the base and replace gel cassette with a fresh gel cassette. Press and release the power button to return to Ready Mode.	
	Cold cassette or improper operating conditions	Use a cassette stored at room temperature. Avoid storing gel cassettes at 4°C. Use E-Gel [®] PowerBase [™] and E-Gel [®] Base at room temperature (20-25°C).	

Troubleshooting, Continued

Plasmid Purification

Review the information provided in the table below to troubleshoot your plasmid purification experiments using the PureLinkTM HQ Mini Plasmid Kit or S.N.A.P.TM MidiPrep Kit.

Problem	Cause	Solution	
Low yield	Incomplete lysis or too much cell lysate has clogged the column	Decrease cell culture volume used. Use up to 100 ml of an overnight culture, density 2×10^9 cells/ml (OD ₆₀₀ < 3.0 per ml). For more dense cultures, decrease volume of culture.	
		Decrease cell culture density. Use LB medium instead of TB (Terrific Broth).	
		Ensure complete resuspension of the bacterial cell pellet.	
	Poor quality of starting material.	Check the growth conditions of the cell culture to ensure plasmid propagation. Use a high copy number plasmid if possible.	
	Incorrect elution conditions	Ensure that the DNA elution was performed with the Elution Buffer supplied in the kit or sterile water (pH >7.0)	
	Genomic DNA present	Perform all mixing steps very gently to avoid damaging DNA. Genomic DNA must be intact to be efficiently removed.	
Plasmid resistant to restriction enzyme digestion	Denatured, supercoiled plasmid present (migrates below native, supercoiled plasmid)	Do not incubate longer than 3 minutes in Lysis Buffer.	

Appendix I

Recipes			
Preparing Antibiotic Stock Solutions	 Prepare the following antibiotic stock solutions. Store all stock solutions at 4°C for up to 2 weeks. For long-term storage, store at -20°C in small aliquots to avoid repeated freezing and thawing. Spectinomycin (10 mg/ml) To 100 µl Spectinomycin solution (100 mg/ml) supplied with the kit, add 900 µl sterile deionized water to produce a 10 mg/ml stock solution. Filter-sterilize. Ampicillin (50 mg/ml) To a sterile microcentrifuge tube, add 50 mg Ampicillin powder supplied with the kit. Add 1 ml sterile deionized water to the tube and mix well to produce a 50 mg/ml stock solution. Filter sterilize. Kanamycin (25 mg/ml) Use the 25 mg/ml solution supplied with the kit as the stock solution. Blasticidin (10 mg/ml) See next page for details. 		
Preparing LB Broth with Antibiotics		e appropriate antibiotic as described in the table supplied with the kit and antibiotic stock solution Recipe To 20 ml LB Broth (1X), add 40 µl ampicillin stock solution (50 mg/ml). To 20 ml LB Broth (1X), add 40 µl kanamycin stock solution (25 mg/ml). To 20 ml LB Broth (1X), add 100 µl Spectinomycin Stock solution (10 mg/ml). To 20 ml LB Broth (1X), add 200 µl	
LB Agar Plates	 Adjust the pH of the solution Autoclave on liquid cycle f Allow solution to cool to 55 10 cm plates. 	Blasticidin stock solution (10 mg/ml). LB Agar (page 160) in 900 ml deionized water. on to 7.0 with NaOH and bring volume up to 1 L. or 20 minutes at 15 psi. 5°C. Add appropriate antibiotic and pour into l store at +4°C. Plates containing antibiotic are	

Blasticidin

Handling Blasticidin	Always wear gloves, mask, goggles, and protective clothing (<i>e.g.</i> a laboratory coat) when handling Blasticidin. Weigh out Blasticidin and prepare solutions in a hood.			
Preparing and Storing Stock Solutions	Blasticidin may be obtained separately from Invitrogen (see page 160) in 50 mg aliquots. Blasticidin is soluble in water. Sterile water is generally used to prepare stock solutions of 5 to 10 mg/ml.			
	• Dissolve Blasticidin in sterile water and filter-sterilize the solution.			
	• Aliquot in small volumes suitable for one time use (see next to last point below) and freeze at -20°C for long-term storage or store at +4°C for short-term storage.			
	• Aqueous stock solutions are stable for 1-2 weeks at +4°C and 6-8 weeks at -20°C.			
	• pH of the aqueous solution should be 7.0 to prevent inactivation of Blasticidin.			
	• Do not subject stock solutions to freeze/thaw cycles (do not store in a frost- free freezer).			
	• Upon thawing, use the required amount and store the thawed stock solution at +4°C for up to 2 weeks.			
	• Medium containing Blasticidin may be stored at +4°C for up to 2 weeks.			
Determining Antibiotic Sensitivity	Since you will be selecting for stable cells using Blasticidin, 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 μ g/ml Blasticidin are sufficient to kill most untransduced mammalian cell lines. We recommend that you test a range of concentrations (see protocol below) to ensure that you determine the minimum concentration necessary for your cell line.			
	1. Plate cells at approximately 25% confluence. Prepare a set of 6 plates. Allow cells to adhere overnight.			
	2. The next day, substitute culture medium with medium containing varying concentrations of Blasticidin (<i>e.g.</i> 0, 2, 4, 6, 8, 10 μ 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.			

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 138). Use the standard BP reaction, if you wish to obtain the maximum number of colonies or generate an entry clone for future use.		
Experimental	To generate an entry clone, you will:		
Outline	1. Perform a BP recombination reaction using the linearized <i>att</i> B-containing expression clone and <i>att</i> P-containing pDONR [™] 221 vector.		
	2. Transform the reaction mixture into competent <i>E. coli</i> host.		
	3. Select for entry clones.		
Donor Vector and Expression Clone	See page 78 for details on the donor vector, resuspending the donor vector, and propagating the donor vector. See page 78 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 ^r). 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 μ 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 ^{m} 2-T1 ^{R} or equivalent for transformation. Do not 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	You will need the following materials:			
	• Linearized <i>att</i> B expression clone (page 78)			
	 pDONR[™]221 vector (supplied with Cat. no. WFGE08; resuspend to 150 ng/µl with water) 			
	• BP Clonase [™] II enzyme mix (supplied with the kit; keep at -20°C until immediately before use)			
	• TE Buffer, pH 8.0 (10 mM Tris-HCl, pH 8.0, 1 mM EDTA)			
	 2 µg/µl Proteinase K solution (supplied with the BP Clonase[™] II enzyme mix; thaw and keep on ice until use) 			
	• pEXP7-tet positive control (50 ng/µl; supplied with the kit)			
	For transformation			
	• One Shot [®] TOP10 Chemically Competent <i>E. coli</i> (one vial per transformation; thaw on ice before use)			
	• S.O.C. Medium (supplied with the competent cells, warm to room temperature)			
	• Positive control (<i>e.g.</i> pUC19 supplied with the competent cells, use as a control for transformation if desired)			
	• LB Medium			
	 LB plates containing 50 μg/ml kanamycin (two for each transformation; warm at 37°C for 30 minutes) 			
	• 42°C water bath			
	• 37°C shaking and non-shaking incubator			
	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 BP Clonase[™] II enzyme mix (see Step 4).

Components	Sample	Positive Control
Linearized <i>att</i> B expression clone from Step 4, page 78 (20-50 fmol)	1-7 µl	
pDONR™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 BP Clonase[™] II enzyme mix from -20°C and thaw on ice (~ 2 minutes).

- 3. Vortex the BP Clonase[™] II enzyme mix briefly twice (2 seconds each time).
- Add 2 µl of BP Clonase[™] II enzyme mix to the sample and positive control vials. Do not add BP Clonase[™] II to the negative control vial. Mix well by vortexing briefly twice (2 seconds each time).

Reminder: Return BP Clonase[™] II enzyme mix to -20°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 μ l 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°C for up to 1 week before transformation, if desired.

Performing the Standard BP Recombination Reaction, Continued

One Shot [®] TOP10 Transformation Protocol			
	Thaw, on ice, one vial of One Shot [®] TOP10 Chemically Competent cells for each transformation.		
	2. Add 1 µl of the BP recombination reaction (from, Step 6, previous page) into a vial of One Shot [®] TOP10 cells and mix gently. Do not mix by pipetting up and down. For the pUC19 control, add 10 pg (1 µl) of DNA into a separate vial of One Shot [®] cells and mix gently.		
	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 μl of the transformation mix and add to 180 μl of LB Medium)		
	 Spread 20 µl and 100 µl from each transformation on a prewarmed selective plate and incubate overnight at 37°C. 		
	An efficient BP recombination reaction may produce hundreds of colonies (> 1500 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 [™] -GW/EmGFP-miR into the pDONR [™] 221 vector preserves the orientation of the cassette.		
	Once you have verified the entry clone, isolate plasmid DNA using PureLink [™] HQ Mini Plasmid Purification Kits (page 160). 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 [®] TOP10 Competent cells using the above protocol. Assess the efficiency of the BP reaction by streaking entry clones onto LB agar plates containing 20 μ g/ml tetracycline. True entry clones should be tetracycline-resistant.		

Performing the Standard LR Recombination Reaction

Introduction	Once you have obtained an entry clone containing the pre-miRNA expression cassette, you will perform an LR recombination reaction between the entry clone (previous page) and a destination vector or pLenti6/V5-DEST, and transform the reaction mixture into competent <i>E. coli</i> to select for expression clones (see page 87). Use the standard LR reaction, if the Rapid BP/LR protocol produces fewer (~20-30) colonies.		
Experimental	To generate an expression clone, you will:		
Outline	 Perform a LR recombination reaction using the <i>att</i>L-containing entry clone (previous page) and <i>att</i>R-containing destination vector including pLenti6/V5-DEST. 		
	2. Transform the reaction mixture into competent <i>E. coli</i> .		
	3. Select for expression clones.		
Destination Vector	See page 78 for details on the destination vector, resuspending the destination vector, and propagating the destination vector.		
Recommended <i>E. coli</i> Host	For optimal results, we recommend using Stbl3 [™] <i>E. coli</i> for transformation as this strain is particularly well-suited for use in cloning unstable DNA such as lentiviral DNA containing direct repeats. One Shot [®] Stbl3 [™] Chemically Competent <i>E. coli</i> are included with Cat. no. WFGE08 for transformation. For instructions, see Transforming One Shot [®] Stbl3 [™] Competent <i>E. coli</i> , page 87.		
	You can use One Shot [®] TOP10 Competent Cells, if you are using non-lentiviral destination vectors.		
Positive Control	The pENTR TM -gus plasmid is provided with the LR Clonase TM II Enzyme Mix for use as a positive control for recombination. Using the pENTR TM -gus entry clone in an LR recombination reaction with a destination vector will allow you to generate an expression clone containing the gene encoding β -glucuronidase (<i>gus</i>) (Kertbundit <i>et al.</i> , 1991).		

Performing the Standard LR Recombination Reaction, Continued

Materials Needed	You will need the following materials:
	• Purified plasmid DNA of your entry clone (50-150 ng/µl in TE, pH 8.0)
	 Destination vector (150 ng/µl in TE, pH 8.0); pLenti6/V5-DEST is supplied with Cat. no. WFGE08
	 LR Clonase[™] II enzyme mix (supplied with the kit, keep at -20°C until immediately before use)
	• TE Buffer, pH 8.0 (10 mM Tris-HCl, pH 8.0, 1 mM EDTA)
	• 2 µg/µl Proteinase K solution (supplied with the LR Clonase [™] II enzyme mix; thaw and keep on ice until use)
	 pENTR[™]-gus positive control (50 ng/µl; supplied with the LR Clonase[™] II enzyme mix)
	• Competent <i>E. coli</i> or One Shot [®] Stbl3 [™] Chemically Competent <i>E. coli</i> (supplied with WFGE08, for use with pLenti6/V5-DEST)
	• S.O.C. Medium (supplied with the competent cells, warm to room temperature)
	• 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
~	
Note	Depending on the destination vector that you will use, perform the appropriate LR reaction as described on the following pages using plasmid DNA from the resulting entry clone .

Performing the Standard LR Recombination Reaction, Continued

LR Recombination
Reaction for
Standard
Destination
Vectors

Use this LR recombination reaction for standard destination vectors. For LR recombination reaction with MultiSite Gateway[®] destination vectors, see next page.

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 LR Clonase[™] II enzyme mix (see Step 4).

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

- Remove the LR Clonase[™] II enzyme mix from -20°C and thaw on ice (~ 2 minutes).
- 3. Vortex the LR Clonase[™] II enzyme mix briefly twice (2 seconds each time).
- Add 2 µl of LR Clonase[™] II enzyme mix to the sample and positive control vials. Do not add LR Clonase[™] II enzyme mix to the negative control vial. Mix well by vortexing briefly twice (2 seconds each time).

Reminder: Return LR Clonase[™] II enzyme mix to -20°C immediately after use.

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

Note: For most applications, 1 hour will yield a sufficient number of colonies for analysis. Depending on your needs, the length of the recombination reaction can be extended up to 18 hours. For large plasmids (\geq 10 kb), longer incubation times (*i.e.* overnight incubation) will yield more colonies and are recommended.

- 6. Add 1 μ l of the Proteinase K solution to each reaction. Incubate for 10 minutes at 37°C.
- Proceed to transformation using competent *E. coli* or Transforming One Shot[®] Stbl3[™] Competent *E. coli* for transforming pLenti6/V5-DEST construct (page 87).

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

 Pick transformants and isolate plasmid DNA using PureLink[™] HQ Mini Plasmid DNA Purification Kit (page 61). Perform restriction analysis to select a clone that has incorporated the pre-miRNA expression cassette (sequence analysis is not necessary).

What You Should See

If you use *E. coli* cells with a transformation efficiency of $\ge 1 \times 10^8$ cfu/µg, the LR reaction should give > 5000 colonies if the entire LR reaction is transformed and plated. See page 88 for confirming the expression clone.

Performing the Standard LR Recombination Reaction, Continued

Use this LR recombination reaction for MultiSite Gateway[®] destination vectors. For LR recombination reaction with standard destination vectors, see previous page.

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 LR Clonase[™] Plus enzyme mix (see Step 4).

Component	Sample	Positive Control
Entry clone (60 ng/µl)	1 µl	1 µl
5' pENTR [™] vector (60 ng/µl)	1 µl	1 µl
3' pENTR [™] vector (60 ng/µl)	1 µl	1 µl
MultiSite Gateway [®] destination vector (60 ng/µl)	1 µl	1 µl
5X LR Clonase [™] Plus Buffer	3 µl	3 µl
TE Buffer, pH 8.0	to 16 µl	to 16 μl

 Remove the LR Clonase[™] Plus enzyme mix from -80°C and thaw on ice (~ 2 minutes).

- 3. Vortex the LR Clonase[™] Plus enzyme mix briefly twice (2 seconds each time).
- 4. To each sample above, add 4 μl of LR Clonase[™] Plus enzyme mix. Mix well by vortexing briefly twice (2 seconds each time).

Reminder: Return LR Clonase[™] Plus enzyme mix to -80°C immediately after use.

- 5. Incubate reactions at 25°C for 16 hours or overnight.
- 6. Add 2 μ l of the Proteinase K solution to each reaction. Incubate for 10 minutes at 37°C.
- 7. Proceed to transform a suitable *E. coli* host and select for expression clones. If you are transforming One Shot[®] TOP10 Chemically Competent *E. coli*, follow the protocol on page 138. For transforming lentiviral construct, proceed to **Transforming One Shot[®] Stb13[™] Competent** *E. coli* (page 87).

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

 Pick transformants and isolate plasmid DNA using PureLink[™] HQ Mini Plasmid DNA Purification Kit (page 61). Perform restriction analysis to select a clone that has incorporated the pre-miRNA expression cassette (sequence analysis is not necessary).

What You Should See

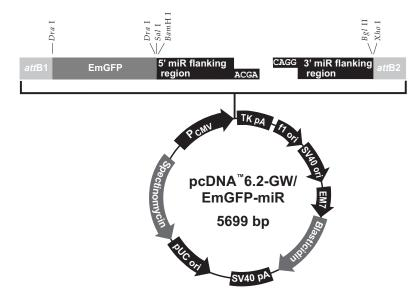
If you use *E. coli* cells with a transformation efficiency of 1×10^9 cfu/µg, the MultiSite Gateway[®] LR reaction should give approximately 2,000-8,000 colonies if the entire reaction is transformed and plated.

Map of pcDNA[™]6.2-GW/EmGFP-miR

pcDNA[™]6.2-GW/EmGFP-miR

The figure below shows the features of the pcDNA[™]6.2-GW/EmGFP-miR vector. The vector is supplied linearized between nucleotides 1518 and 1519 with 4 base pair 5' overhangs on each strand as indicated. The complete sequence of pcDNA[™]6.2-GW/EmGFP-miR is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (page 163).

Note: For an explanation of the features see page 145.



Comments for pcDNA[™] 6.2-GW/EmGFP-miR 5699 nucleotides

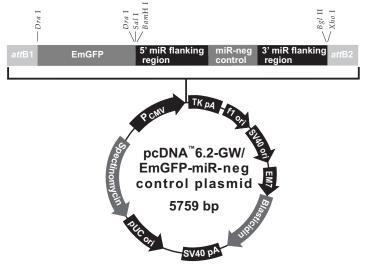
CMV promoter: bases 1-588 attB1 site: bases 680 - 704 EmGFP: bases 713-1432 EmGFP forward sequencing primer site: bases 1409-1428 5' miR flanking region: bases 1492-1518 5' overhang (C): bases 1515-1518 5'overhang: bases 1519-1522 3' miR flanking region: bases 1519-1563 attB2 site (C): bases 1592-1616 miRNA reverse sequencing primer site (C): bases 1607-1626 TK polyadenylation signal: bases 1645-1916 f1 origin: bases 2028-2456 SV40 early promoter and origin: bases 2483-2791 EM7 promoter: bases 2846-2912 Blasticidin resistance gene: bases 2913-3311 SV40 polyadenylation signal: bases 3469-3599 pUC origin (C): bases 3737-4410 Spectinomycin resistance gene (C): bases 4480-5490 Spectinomycin promoter (C): bases 5491-5624

(C) = Complementary strand

Map of pcDNA[™]6.2-GW/EmGFP-miR-neg control plasmid

pcDNA[™]6.2-GW/EmGFP-miRneg control plasmid

The figure below shows the features of the pcDNA[™]6.2-GW/EmGFP-miR-neg control plasmid. The vector contains an insert between bases 1519 and 1578 that can form a hairpin structure just as a regular pre-miRNA, but is predicted not to target any known vertebrate gene. The insert has been cloned according to the instructions in this manual. Thus, this plasmid serves as a suitable negative control for pre-miRNA experiments with pcDNA[™]6.2-GW/EmGFP-miR expression vectors. The complete sequence of pcDNA[™]6.2-GW/EmGFP-miR-neg control plasmid is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (page 163).



Comments for pcDNA[™] 6.2-GW/EmGFP-miR-neg control plasmid 5759 nucleotides

CMV promoter: bases 1-588 attB1 site: bases 680-704 EmGFP: bases 713-1432 EmGFP forward sequencing primer site: bases 1409-1428 5' miR flanking region: bases 1492-1518 miR-neg control: bases 1519-1578 3' miR flanking region: bases 1579-1623 attB2 site (C): bases 1652-1676 miRNA reverse sequencing primer site (C): bases 1667-1686 TK polyadenylation signal: bases 1705-1976 f1 origin: bases 2088-2516 SV40 early promoter and origin: bases 2543-2851 EM7 promoter: bases 2906-2972 Blasticidin resistance gene: bases 2973-3371 SV40 polyadenylation signal: bases 3529-3659 pUC origin (C): bases 3797-4470 Spectinomycin resistance gene (C): bases 4540-5550 Spectinomycin promoter (C): bases 5551-5684

(C) = Complementary strand

Features of pcDNA[™]6.2-GW/EmGFP-miR

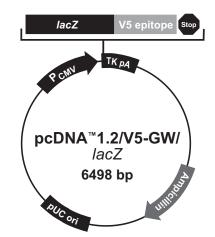
Features of the
VectorsThe pcDNA[™]6.2-GW/EmGFP-miR vector contains the following elements. All
features have been functionally tested and the vector fully sequenced.

Feature	Benefit
CMV promoter	Permits high-level, constitutive expression of the gene of interest (Andersson <i>et al.</i> , 1989; Boshart <i>et al.</i> , 1985; Nelson <i>et al.</i> , 1987).
miRNA forward sequencing primer	Allows sequencing of the insert (for pcDNA [™] 6.2-GW/miR vectors).
<i>att</i> B1 and <i>att</i> B2 sites	Bacteriophage λ -derived recombination sequences that allow recombinational cloning of a gene of interest in the expression construct with a Gateway [®] destination vector (Landy, 1989).
EmGFP coding sequence	Allows visual detection of transfected mammalian cells using fluorescence microscopy.
EmGFP forward sequencing primer	Allows sequencing of the insert.
5' miR flanking region	Allows formation of functional engineered pre-miRNA.
5' overhangs	Allows ligase-mediated directional cloning of the double-stranded oligonucleotide of interest.
miR-neg control	Allows formation of a pre-miRNA hairpin sequence predicted not to target any known vertebrate gene (only for pcDNA [™] 6.2-GW/EmGFP-miR-neg control). Sequence without 5' overhangs is shown below: 5' -GAAATGTACTGCGCGTGGAGACGTTTTGGCCACTGACTGA
3' miR flanking region	Allows formation of functional engineered pre-miRNA.
miRNA reverse sequencing primer	Allows sequencing of the insert.
TK polyadenylation signal	Allows transcription termination and polyadenylation of mRNA.
f1 origin	Allows rescue of single-stranded DNA.
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 (<i>bsd</i>) resistance gene	Permits selection of stably transfected mammalian cell lines (Kimura <i>et al.</i> , 1994).
SV40 polyadenylation signal	Allows transcription termination and polyadenylation of mRNA.
pUC origin	Permits high-copy replication and maintenance in <i>E. coli</i> .
Spectinomycin resistance gene (<i>aad</i> A1)	Allows selection of the plasmid in <i>E. coli</i> (Liebert <i>et al.,</i> 1999).
Spectinomycin promoter	Allows expression of the spectinomycin resistance gene in <i>E. coli</i> .

Map of pcDNA[™]1.2/V5-GW/*lacZ*

Description

pcDNA^M1.2/V5-GW/*lacZ* (6498 bp) is a control vector expressing a C-terminallytagged β -galactosidase fusion protein under the control of the human cytomegalovirus (CMV) promoter (Andersson *et al.*, 1989; Boshart *et al.*, 1985; Nelson *et al.*, 1987), and was generated using the MultiSite Gateway[®] Three-Fragment Vector Construction Kit available from Invitrogen (Catalog no. 12537-023). Briefly, a MultiSite Gateway[®] LR recombination reaction was performed with pDEST^MR4-R3 and entry clones containing the CMV promoter, *lacZ* gene, and V5 epitope and TK polyadenylation signal (Cole & Stacy, 1985) to generate the pcDNA^M1.2/V5-GW/*lacZ* vector. β -galactosidase is expressed as a C-terminal V5 fusion protein with a molecular weight of approximately 119 kDa. The complete sequence of pcDNA^M1.2/V5-GW/*lacZ* is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (page 163).



Comments for pcDNA[™]1.2/V5-GW//acZ 6498 nucleotides

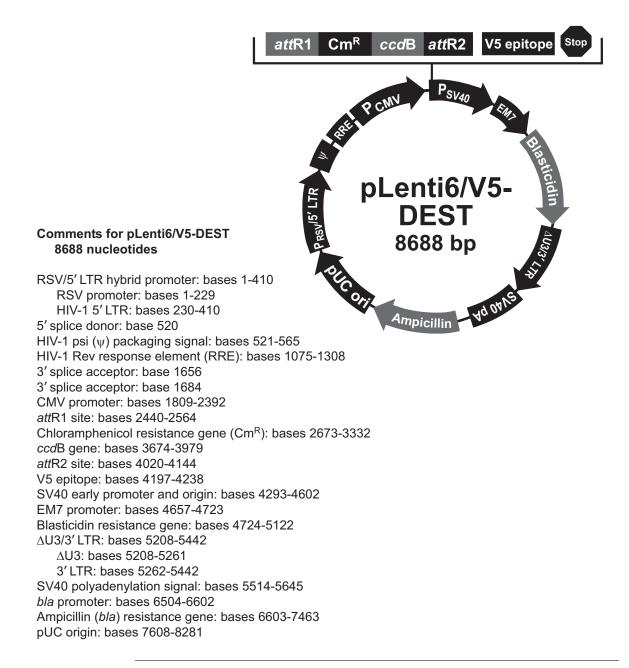
attB4: bases 5-25 CMV promoter: bases 137-724 attB1: bases 614-637 LacZ fusion protein: bases 643-3798 LacZ ORF: bases 643-3714 attB2: bases 3716-3739 V5 epitope: bases 3739-3780 lacZ forward 2 priming site: 840-859 lacZ reverse 2 priming site: 1820-1839 (C) TK polyadenylation signal: bases 3807-4078 attB3: bases 4079-4099 *bla* promoter: bases 4603-4701 Ampicillin (*bla*) resistance gene: bases 4702-5562 pUC origin: bases 5707-6380

(C) = complementary strand

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 2447 and 4130. The complete sequence for pLenti6/V5-DEST is available from our web site (www.invitrogen.com) or by contacting Technical Service (page 160).

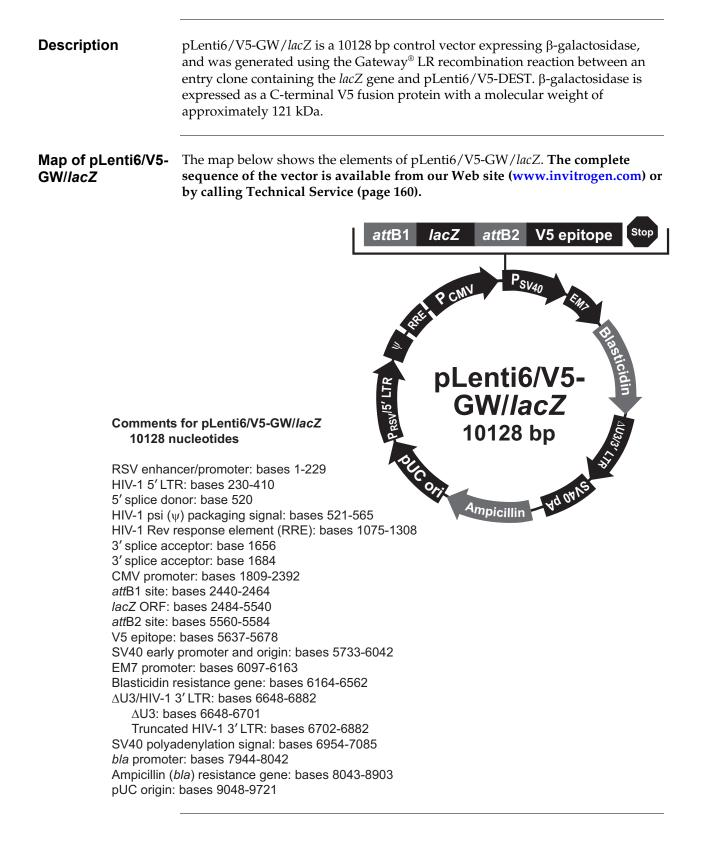


Map and Features of pLenti6/V5-DEST, Continued

Features of the	The pLenti6/V5-DEST (8688 bp) vector contains the following elements. All
Vector	features have been functionally tested and the vector is fully sequenced.

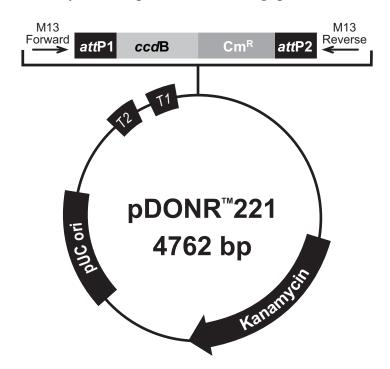
Feature	Benefit
Rous Sarcoma Virus (RSV) enhancer/promoter	Allows Tat-independent production of viral mRNA (Dull et al., 1998).
HIV-1 truncated 5' LTR	Permits viral packaging and reverse transcription of the viral mRNA (Luciw, 1996).
5' splice donor and 3' acceptors	Enhances the biosafety of the vector by facilitating removal of the Ψ 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 λ -derived DNA recombination sequences that permit recombinational cloning of the gene of interest from a Gateway [®] entry clone (Landy, 1989).
Chloramphenicol resistance gene (Cm ^R)	Allows counterselection of the plasmid.
ccdB 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). This feature is not used when expressing miRNAs.
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



Map and Features of pDONR[™]221

Map of pDONR[™]221 The map below shows the elements of pDONR[™]221 vector. **The complete** sequences of pDONR[™]221 is available for downloading from our web site (www.invitrogen.com) or by contacting Technical Service (page 160).



Comments for pDONR[™]221 4762 nucleotides

rrnB T2 transcription termination sequence: bases 268-295 (C) rrnB T1 transcription termination sequence: bases 427-470 (C) M13 Forward (-20) priming site: bases 537-552 *att*P1:bases 570-801 *ccd*B gene: bases 1197-1502 (C) Chloramphenicol resistance gene: bases 1847-2506 (C) *att*P2: bases 2754-2985 (C) M13 Reverse priming site: bases 3027-3043 Kanamycin resistance gene: bases 3156-3965 pUC origin: bases 4086-4759 (C) = complementary strand

Map and Features of pDONR[™]221, Continued

Features of pDONR[™]221

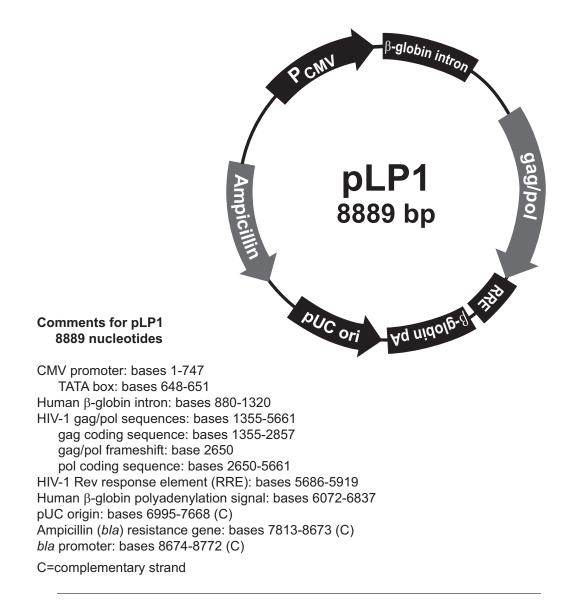
pDONR[™]221 (4762 bp) contains the following elements. All features have been functionally tested and the vectors fully sequenced.

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 λ-derived DNA recombination sequences that allow recombinational cloning of the gene of interest from a Gateway [®] expression clone or <i>att</i> B PCR product (Landy, 1989)
ccdB gene	Allows negative selection of the plasmid
Chloramphenicol resistance gene (Cm ^R)	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 then selfcleaved by the viral protease into individual Gag and Pol polyproteins. The **complete sequence of pLP1 is available for downloading from our Web site** (www.invitrogen.com) or by contacting Technical Service (see page 160).



Map and Features of pLP1, Continued

Features of pLP1

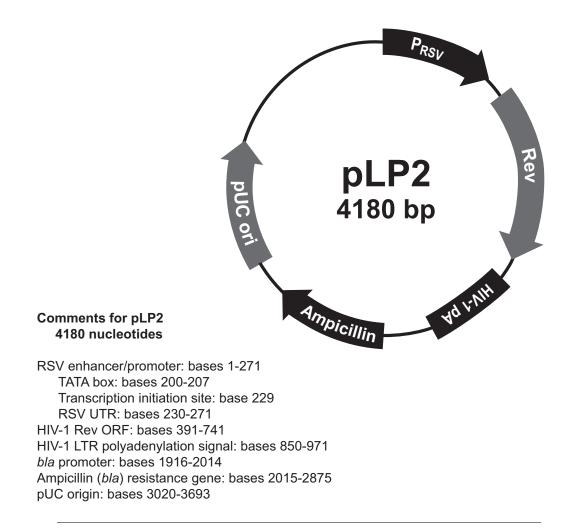
pLP1 (8889 bp) contains the following elements. All features have been functionally tested.

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 β -globin polyadenylation signal	Allows efficient transcription termination and polyadenylation of mRNA.
pUC origin of replication (<i>ori</i>)	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 complete sequence** of pLP2 is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (see page 160).



Map and Features of pLP2, Continued

Features of pLP2

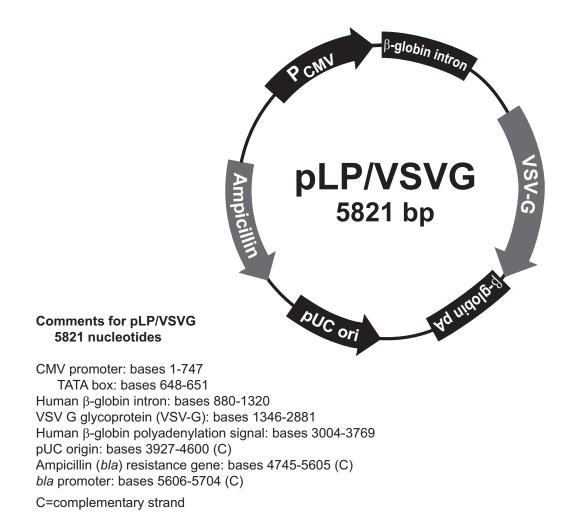
pLP2 (4180 bp) contains the following elements. All features have been functionally tested.

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 complete sequence of pLP/VSVG is available for downloading from our Web site (www.invitrogen.com) or by contacting Technical Service (see page 160).



Map and Features of pLP/VSVG, Continued

Features of pLP/VSVG

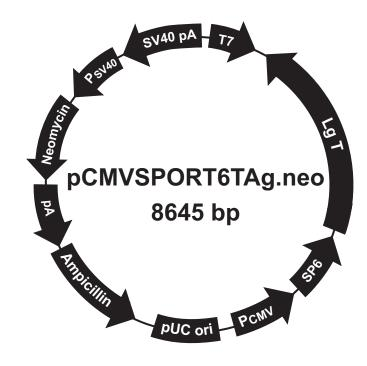
 $\rm pLP/VSVG$ (5821 bp) contains the following elements. All features have been functionally tested.

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 β -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 pCMVSPORT6TAg.neo

Description The pCMVSPORT6TAg.neo plasmid is derived from pCMVSPORT6, which has been modified to include the following features:

- The neomycin resistance gene for stable selection in mammalian cells (Southern & Berg, 1982). Expression of the neomycin resistance gene is controlled by the SV40 early enhancer/promoter from which the SV40 origin of replication has been removed.
- The gene encoding the SV40 large T antigen to facilitate optimal virus production (*e.g.* Invitrogen's ViraPower[™] Lentiviral Expression System) and to permit episomal replication of plasmids containing the SV40 early promoter and origin. Expression of the SV40 large T antigen is controlled by the human cytomegalovirus (CMV) promoter.



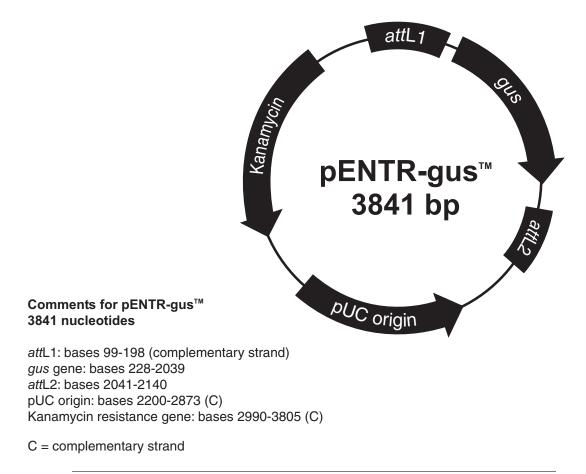
Map of pENTR[™]-gus

Description

pENTRTM-gus is a 3841 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 *att*B recombination sites. The amplified PCR product was then used in a BP recombination reaction with pDONR201TM to generate the entry clone. For more information about the BP recombination reaction, refer to the Gateway[®] Technology with ClonaseTM II manual which is available for downloading from our Web site or by contacting Technical Service.

Map of Control Vector

The figure below summarizes the features of the pENTR[™]-gus vector. The complete sequence for pENTR[™]-gus is available from our Web site (www.invitrogen.com) or by contacting Technical Service (see page 160).



Appendix II

Accessory Products

BioModule[™] Units

Additional BioModule[™] Units that can be used for validation experiments or gene expression profiling are available from Invitrogen. Ordering information is provided below. For more information, visit www.invitrogen.com or call Technical Service (page 163).

Product	Amount	Catalog no.
BioModule [™] Western Analysis Unit		
for chromogenic detection	20 transfers	WFGE09
for chemiluminescent detection	20 transfers	WFGE10
BioModule [™] qRT-PCR Unit	100 reactions	WFGE01
	1000 reactions	WFGE02
BioModule [™] Immunohistochemical (IHC) Staining for Tissue	150 slides	WFGE11
BioModule [™] Transfection and Control Unit with BLOCK-iT [™] Technology	1 unit	WFGE06
BioModule [™] Microarray Analysis		
with indirect labeling	15 arrays	WFGE03
with direct labeling	15 arrays	WFGE04
BioModule [™] Lentiviral 293 Unit (for contents, see below)	1 Unit	WFGE08-S

BioModule[™] Lentiviral 293 Unit

The BioModule[™] Lentiviral 293 Unit contains all necessary media, antibiotics, serum, and Trypan Blue for growth, maintenance, and checking cell viability of 293 FT cells.

Product	Amount
Fetal Bovine Serum	2 x 100 ml
200 mM L-Glutamine (100X)	100 ml
Penicillin-Streptomycin, liquid (5,000 units of penicillin and 5,000 μ g of streptomycin)	100 ml
Trypsin-EDTA	100 ml
Dulbecco's Modified Eagle Medium	2 x 1000 ml
10 mM MEM Non-Essential Amino Acids Solution (100X)	100 ml
MEM Sodium Pyruvate Solution (100X)	100 ml
Phosphate-Buffered Saline, pH 7.4	500 ml
Opti-MEM® I Reduced Serum Medium	500 ml
Geneticin [®] Selective Antibiotic (50 mg/ml)	20 ml
Trypan Blue Stain	100 ml

Accessory Products, Continued

AccessoryMany of the reagents supplied in the BioModule™ BLOCK-iT™ RNAi Units as
well as other products suitable for use with the units are available separately
from Invitrogen. Ordering information is provided below.

Product	Amount	Catalog no.
Lipofectamine [™] 2000 Reagent	0.75 ml	11668-027
	1.5 ml	11668-019
pDONR [™] 221 Vector	6 µg	12536-017
Opti-MEM [®] I Reduced Serum Medium	100 ml	31985-062
Gateway® BP Clonase™ II Enzyme Mix	20 reactions	11789-020
	100 reactions	11789-100
Gateway [®] LR Clonase [™] II Enzyme Mix	20 reactions	11791-020
	100 reactions	11791-100
Phosphate-Buffered Saline (PBS), pH 7.4	500 ml	10010-023
One Shot [®] Stbl3 [™] Chemically Competent <i>E. coli</i>	20 x 50 μl	C7373-03
One Shot® TOP10 Chemically Competent E. coli	20 x 50 μl	C4040-03
Library Efficiency [®] DB3.1 [™] Competent Cells	5 x 200 µl	11782-018
One Shot [®] ccdB Survival [™] T1 ^R Chemically Competent Cells	10 reactions	C7510-03
ViraPower [™] Lentiviral Packaging Mix	60 reactions	K4975-00
Blasticidin	50 mg	R210-01
Kanamycin Sulfate (100X)	100 ml	15160-054
Ampicillin	5 g	Q100-16
Geneticin [®] Selective Antibiotic (50 mg/ml)	20 ml	10131-035
LB Broth (1X)	500 ml	10855-021
LB Agar (Lennox L Agar) [®]	500 g	22700-025
PureLink [™] HQ Mini Plasmid Purification Kit	100 preps	K2100-01
PureLink [™] Quick Gel Extraction Kit	50 preps	K2100-12
BlueJuice [™] Gel Loading Buffer	3 x 1 ml	10816-015
TrackIt [™] Cyan/Orange Loading Buffer	3 x 0.5 ml	10482-028
TrackIt [™] Cyan/Yellow Loading Buffer	3 x 0.5 ml	10482-035
TrackIt [™] 10 bp DNA Ladder	20 applications	10488-019
TrackIt [™] 1 Kb Plus DNA Ladder	100 applications	10488-085
Quant-iT [™] DNA Assay Kit, Broad-Range	1000 assays	Q33130

Accessory Products, Continued

BLOCK-iT [™] RNAi Designer	The BLOCK-iT [™] RNAi Designer is an online tool (<u>www.invitrogen.com/rnaidesigner</u>) to help you design and order microRNA sequences for any target gene of interest. The RNAi Designer incorporates published rules on RNAi design into a proprietary algorithm to design most effective microRNA sequences to obtain high level gene knockdown.
BLOCK-iT [™] RNAi Products	A large variety of BLOCK-iT [™] RNAi products are available from Invitrogen to facilitate RNAi analysis including Stealth [™] RNAi, the Validated Stealth [™] RNAi, Stealth [™] RNAi Collection, and a large selection of RNAi vectors.
	For details, visit the RNAi Central portal at www.invitrogen.com/rnai or contact Technical Service (page 163).
Gateway [®] Destination Vectors	A large selection of Gateway [®] destination vectors are available from Invitrogen to facilitate the transfer of the pre-miRNA sequence into a suitable destination vector to allow miRNA expression in multiple systems including viral expression systems and tissue-specific expression. See page 77 for a list of destination vectors compatible with the pcDNA [™] 6.2-GW/EmGFP-miR Vector.
Antibodies	A large variety of high-quality antibodies including the Zymed [®] Antibodies is available from Invitrogen for use in Western immunodetection, immunohistochemistry, or ELISA assays. For details, visit <u>www.invitrogen.com</u> or contact Technical Service (page 163).

Technical Service

Web Resources	Technica applicatiCompletAccess to	trogen Web site at <u>www.invitrog</u> l resources, including manuals, v on notes, MSDSs, FAQs, formula e technical service contact inform o the Invitrogen Online Catalog nal product information and spec	vector maps and sequences, ations, citations, handbooks, etc. nation
Contact Us		ormation or technical assistance, l offices are listed on our Web pa	call, write, fax, or email. Additional ge (<u>www.invitrogen.com</u>).
Corporate Headquarter Invitrogen Corporation 1600 Faraday Avenue Carlsbad, CA 92008 US Tel: 1 760 603 7200 Tel (Toll Free): 1 800 95 Fax: 1 760 602 6500 E-mail: tech_service@invitroge	n 5A 55 6288	Japanese Headquarters: Invitrogen Japan LOOP-X Bldg. 6F 3-9-15, Kaigan Minato-ku, Tokyo 108-0022 Tel: 81 3 5730 6509 Fax: 81 3 5730 6519 E-mail: jpinfo@invitrogen.com	European Headquarters: Invitrogen Ltd Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF, UK Tel: +44 (0) 141 814 6100 Tech Fax: +44 (0) 141 814 6117 E-mail: <u>eurotech@invitrogen.com</u>
Material Data Safety Sheets (MSDSs)	page, click o	vailable on our Web site at <u>www</u> n Technical Resources and follov le MSDS for your product.	U
Limited Warranty	Limited Warranty Invitrogen is committed to providing our customers with high-quality goods and service Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about an Invitrogen product or service, contact our Technical Service Representatives. Invitrogen warrants that all of its products will perform according to specifications stated on the certificate of analysis. The company will replace, free of charge, any product that does not meet those specifications. This warranty limits Invitrogen Corporation's liability only to the cost of the product. No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. Invitrogen reserves the right to select the method(s) used to analyze a product unless Invitrogen agrees to a specified method in writing prior to acceptance of the order. Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore Invitrogen makes no warranty of any kind regarding the contents of any publications or documentation. If you		
	discover an er Representative Invitrogen ass consequential exclusive. No	ror in any of our publications, please	e report it to our Technical Service or any special, incidental, indirect or ove limited warranty is sole and pressed or implied, including any

Product Qualification

described in this manual and the reaction is transformed into One Shot® TOP10 Chemically Competent E. coli. Sequence analysis is performed on 20 colonies using the appropriate forward and reverse sequencing primers. The qualification must meet the following criteria: • Cloning efficiency of ≥95% • Sequencing analysis must show ≥95% clones contain the insert in the correct orientation and ≥80% clones must have the correct sequence in both directions (forward and reverse) The attB1 and attB2 sequences should be correct pLenti6/V5-DEST vector The pLenti6/V5-DEST vector is qualified in a recombination assay using Gateway® LR Clonase [™] II Enzyme Mix. The ccdB gene is assayed by transformation using an appropriate E. coli strain. pDONR [™] 221 Vector pDONR [™] 221 vector is qualified in a BP recombination reaction using the Gateway® BP Clonase [™] II. The ccdB gene is assayed by transformation using an appropriate E. coli strain. gboot Lipofectamine [™] 2000 is tested for the absence of microbial contamination using blood agar plates, Sabaraud dextrose agar plates, and fluid thioglycolate medium, and functionally by transfection with a luciferase reporter-containing plasmid. Antibiotics Kanamycin, Ampicillin, and Spectinomycin Competent E. coli cells are transformed with an appropriate antibiotic-resistant plasmid and appropriate antibiotic. Colony growth should only be observed on LB plates plated with antibiotic-resistant plasmid donly be observed on LB plates plated with antibiotic-resistant plasmid transformation mixture. Blasticidin Blasticidin-sens	Introduction	This section describes the criteria used to qualify the components of the BioModule [™] BLOCK-iT [™] RNAi Units.				
The miR-lacZ positive ds control oligo is annealed and cloned into each vector as described in this manual and the reaction is transformed into One Shot [®] TOP10 Chemically Competent <i>E. coli</i> . Sequence analysis is performed on 20 colonies using the appropriate forward and reverse sequencing primers. The qualification must meet the following criteria: • Cloning efficiency of 295% • Sequencing analysis must show ≥95% clones contain the insert in the correct orientation and ≥80% clones must have the correct sequence in both directions (forward and reverse) The attB1 and attB2 sequences should be correct pLenti6/V5-DEST vector The pLenti6/V5-DEST vector is qualified in a recombination assay using Gateway [®] LR Clonase [®] II Enzyme Mix. The <i>ccdB</i> gene is assayed by transformation using an appropriate <i>E. coli</i> strain. pDONR [™] 221 vector pDONR [™] 221 vector is qualified in a BP recombination reaction using the Gateway [®] BP Clonase [™] II. The <i>ccdB</i> gene is assayed by transformation using an appropriate <i>E. coli</i> strain. 2000 Lipofectamine [™] 2000 Kanamycin, Ampicillin, and Spectinomycin Competent <i>E. coli</i> cells are transformed with an appropriate antibiotic-resistant plasmid and appropriate antibiotic-resistant plasmid and appropriate antibiotic-resistant plasmid and appropriate antibiotic-resistant plasmid and appropriate antibiotic-resistant plasmid transformation mixture. Blasticidin Blasticidin is lot-qualified by performing a kill curve on Blasticdin-sensitive and resistant mammalian cell lines. Blast	Vectors	, , ,				
described in this manual and the reaction is transformed into One Shot® TOP10 Chemically Competent E. coli. Sequence analysis is performed on 20 colonies using the appropriate forward and reverse sequencing primers. The qualification must meet the following criteria: • Cloning efficiency of ≥95% • Sequencing analysis must show ≥95% clones contain the insert in the correct orientation and ≥80% clones must have the correct sequence in both directions (forward and reverse) The attB1 and attB2 sequences should be correct pLenti6/V5-DEST vector The pLenti6/V5-DEST vector is qualified in a recombination assay using Gateway® LR Clonase [™] II Enzyme Mix. The ccdB gene is assayed by transformation using an appropriate E. coli strain. pDONR [™] 221 Vector pDONR [™] 221 vector is qualified in a BP recombination reaction using the Gateway® BP Clonase [™] II. The ccdB gene is assayed by transformation using an appropriate E. coli strain. gboot Lipofectamine [™] 2000 is tested for the absence of microbial contamination using blood agar plates, Sabaraud dextrose agar plates, and fluid thioglycolate medium, and functionally by transfection with a luciferase reporter-containing plasmid. Antibiotics Kanamycin, Ampicillin, and Spectinomycin Competent E. coli cells are transformed with an appropriate antibiotic-resistant plasmid and appropriate antibiotic. Colony growth should only be observed on LB plates plated with antibiotic-resistant plasmid donly be observed on LB plates plated with antibiotic-resistant plasmid transformation mixture. Blasticidin Blasticidin-sens		pcDNA [™] 6.2-GW/EmGFP-miR Vector				
 Sequencing analysis must show ≥95% clones contain the insert in the correct orientation and ≥80% clones must have the correct sequence in both directions (forward and reverse) The <i>att</i>B1 and <i>att</i>B2 sequences should be correct pLenti6/V5-DEST vector The <i>pLenti6/V5-DEST vector</i> is qualified in a recombination assay using Gateway® LR Clonase[™] II Enzyme Mix. The <i>ccdB</i> gene is assayed by transformation using an appropriate <i>E. coli</i> strain. pDONR[™]221 Vector pDONR^{™221} vector is qualified in a BP recombination reaction using the Gateway® BP Clonase[™] II. The <i>ccdB</i> gene is assayed by transformation using an appropriate <i>E. coli</i> strain. Dipofectamine[™] Lipofectamine[™] 2000 is tested for the absence of microbial contamination using an appropriate <i>E. coli</i> strain. Lipofectamine[™] Lipofectamine[™] 2000 is tested for the absence of microbial contamination using blood agar plates, Sabaraud dextrose agar plates, and fluid thioglycolate medium, and functionally by transfection with a luciferase reporter-containing plasmid. Antibiotics Kanamycin, Ampicillin, and Spectinomycin Competent <i>E. coli</i> cells are transformed with an appropriate antibiotic-resistant plasmid and appropriate antibiotic-sensitive plasmid. Cells are plated on LB medium containing the appropriate antibiotic. Colony growth should only be observed on LB plates plated with antibiotic-resistant plasmid transformation mixture. Blasticidin Blasticidin is lot-qualified by performing a kill curve on Blasticidin-sensitive and resistant mammalian cell lines. Blasticidin-sensitive cells should be killed at all		Chemically Competent <i>E. coli</i> . Sequence analysis is performed on 20 colonies using the appropriate forward and reverse sequencing primers. The qualification				
pLenti6/V5-DEST vector The pLenti6/V5-DEST vector is qualified in a recombination assay using Gateway [®] LR Clonase [™] II Enzyme Mix. The <i>ccdB</i> gene is assayed by transformation using an appropriate <i>E. coli</i> strain. pDONR [™] 221 Vector pDONR [™] 221 vector is qualified in a BP recombination reaction using the Gateway [®] BP Clonase [™] II. The <i>ccdB</i> gene is assayed by transformation using an appropriate <i>E. coli</i> strain. Lipofectamine [™] 2000 Lipofectamine [™] Lipofectamine [™] Lipofectamine [™] Competent <i>E. coli</i> strain. Lipofectamine [™] Conse [™] Competent <i>E. coli</i> strain. Conse [™] Competent <i>E. coli</i> strain. Antibiotics Kanamycin, Ampicillin, and Spectinomycin Competent <i>E. coli</i> cells are transformed with an appropriate antibiotic-resistant plasmid and appropriate antibiotic-sensitive plasmid. Cells are plated on LB medium containing the appropriate antibiotic. Colony growth should only be observed on LB plates plated with antibiotic-resistant plasmid transformation mixture. Blasticidin Blasticidin is lot-qualified by performing a kill curve on Blasticidin-sensitive and resistant mammalian cell lines. Blasticidin-sensitive cells should be killed at all		• Sequencing analysis must show ≥95% clones contain the insert in the correct orientation and ≥80% clones must have the correct sequence in both				
The pLenti6/V5-DEST vector is qualified in a recombination assay using Gateway [®] LR Clonase [™] II Enzyme Mix. The <i>ccdB</i> gene is assayed by transformation using an appropriate <i>E. coli</i> strain. pDONR [™] 221 Vector pDONR [™] 221 vector is qualified in a BP recombination reaction using the Gateway [®] BP Clonase [™] II. The <i>ccdB</i> gene is assayed by transformation using an appropriate <i>E. coli</i> strain. Lipofectamine [™] Lipofectamine [™] 2000 is tested for the absence of microbial contamination using blood agar plates, Sabaraud dextrose agar plates, and fluid thioglycolate medium, and functionally by transfection with a luciferase reporter-containing plasmid. Antibiotics Kanamycin, Ampicillin, and Spectinomycin Competent <i>E. coli</i> cells are transformed with an appropriate antibiotic-resistant plasmid and appropriate antibiotic. Colony growth should only be observed on LB plates plated with antibiotic. Colony growth should only be observed on LB plates plated with antibiotic. Colony growth should only be observed on LB plates plated with antibiotic cells and transformation mixture. Blasticidin Blasticidin is lot-qualified by performing a kill curve on Blasticidin-sensitive and resistant mammalian cell lines. Blasticidin-sensitive cells should be killed at all		The <i>att</i> B1 and <i>att</i> B2 sequences should be correct				
Gateway [®] LR Clonase [™] II Enzyme Mix. The <i>ccdB</i> gene is assayed by transformation using an appropriate <i>E. coli</i> strain. pDONR™221 Vector pDONR™221 vector is qualified in a BP recombination reaction using the Gateway [®] BP Clonase [™] II. The <i>ccdB</i> gene is assayed by transformation using an appropriate <i>E. coli</i> strain. Lipofectamine [™] Coold agar plates, Sabaraud dextrose agar plates, and fluid thioglycolate medium, and functionally by transfection with a luciferase reporter-containing plasmid. Antibiotics Kanamycin, Ampicillin, and Spectinomycin Competent <i>E. coli</i> cells are transformed with an appropriate antibiotic-resistant plasmid and appropriate antibiotic-sensitive plasmid. Cells are plated on LB medium containing the appropriate antibiotic. Colony growth should only be observed on LB plates plated with antibiotic-resistant plasmid transformation mixture. Blasticidin Blasticidin is lot-qualified by performing a kill curve on Blasticidin-sensitive and resistant mammalian cell lines. Blasticidin-sensitive cells should be killed at all		-				
pDONR [™] 221 vector is qualified in a BP recombination reaction using the Gateway [®] BP Clonase [™] II. The <i>ccd</i> B gene is assayed by transformation using an appropriate <i>E. coli</i> strain. Lipofectamine [™] Lipofectamine [™] 2000 is tested for the absence of microbial contamination using blood agar plates, Sabaraud dextrose agar plates, and fluid thioglycolate medium, and functionally by transfection with a luciferase reporter-containing plasmid. Antibiotics Kanamycin, Ampicillin, and Spectinomycin Competent <i>E. coli</i> cells are transformed with an appropriate antibiotic-resistant plasmid and appropriate antibiotic-sensitive plasmid. Cells are plated on LB medium containing the appropriate antibiotic. Colony growth should only be observed on LB plates plated with antibiotic-resistant plasmid transformation mixture. Blasticidin Blasticidin is lot-qualified by performing a kill curve on Blasticidin-sensitive and resistant mammalian cell lines. Blasticidin-sensitive cells should be killed at all		Gateway [®] LR Clonase [™] II Enzyme Mix. The <i>ccd</i> B gene is assayed by				
Gateway® BP Clonase™ II. The ccdB gene is assayed by transformation using an appropriate E. coli strain. Lipofectamine™ 2000 Lipofectamine™ Lipofectamine™ Lipofectamine™ Lipofectamine™ Lipofectamine™ Lipofectamine™ 2000 Lipofectamine™ Lipofectamine™ 2000 <		pDONR [™] 221 Vector				
 2000 blood agar plates, Sabaraud dextrose agar plates, and fluid thioglycolate medium, and functionally by transfection with a luciferase reporter-containing plasmid. Antibiotics Kanamycin, Ampicillin, and Spectinomycin Competent <i>E. coli</i> cells are transformed with an appropriate antibiotic-resistant plasmid and appropriate antibiotic-sensitive plasmid. Cells are plated on LB medium containing the appropriate antibiotic. Colony growth should only be observed on LB plates plated with antibiotic-resistant plasmid transformation mixture. Blasticidin Blasticidin is lot-qualified by performing a kill curve on Blasticidin-sensitive and resistant mammalian cell lines. Blasticidin-sensitive cells should be killed at all 		Gateway [®] BP Clonase [™] II. The <i>ccd</i> B gene is assayed by transformation using an				
 Competent <i>E. coli</i> cells are transformed with an appropriate antibiotic-resistant plasmid and appropriate antibiotic-sensitive plasmid. Cells are plated on LB medium containing the appropriate antibiotic. Colony growth should only be observed on LB plates plated with antibiotic-resistant plasmid transformation mixture. Blasticidin Blasticidin is lot-qualified by performing a kill curve on Blasticidin-sensitive and resistant mammalian cell lines. Blasticidin-sensitive cells should be killed at all 	-	blood agar plates, Sabaraud dextrose agar plates, and fluid thioglycolate medium, and functionally by transfection with a luciferase reporter-containing				
plasmid and appropriate antibiotic-sensitive plasmid. Cells are plated on LB medium containing the appropriate antibiotic. Colony growth should only be observed on LB plates plated with antibiotic-resistant plasmid transformation mixture. Blasticidin Blasticidin is lot-qualified by performing a kill curve on Blasticidin-sensitive and resistant mammalian cell lines. Blasticidin-sensitive cells should be killed at all	Antibiotics	Kanamycin, Ampicillin, and Spectinomycin				
Blasticidin is lot-qualified by performing a kill curve on Blasticidin-sensitive and resistant mammalian cell lines. Blasticidin-sensitive cells should be killed at all		plasmid and appropriate antibiotic-sensitive plasmid. Cells are plated on LB medium containing the appropriate antibiotic. Colony growth should only be observed on LB plates plated with antibiotic-resistant plasmid transformation				
resistant mammalian cell lines. Blasticidin-sensitive cells should be killed at all		Blasticidin				
concentrations tested (2.5-10 μ g/ml) within 10 days after addition of Blasticidin.		Blasticidin is lot-qualified by performing a kill curve on Blasticidin-sensitive and resistant mammalian cell lines. Blasticidin-sensitive cells should be killed at all concentrations tested (2.5-10 μ g/ml) within 10 days after addition of Blasticidin.				

Product Qualification, Continued

Gateway [®] Clonase [™] II Enzyme Mix	Gateway [®] BP and LR Clonase [™] II Enzyme Mixes are functionally tested in a one hour recombination reaction followed by a transformation assay.	
One Shot [®] Competent Cells	Each lot of One Shot [®] Competent cells is tested for transformation efficiency using the pUC19 control plasmid included in the kit and following the procedure described in this manual. Test transformations are performed on 3 to 20 vials per lot, depending on batch size. Transformed cultures are plated on LB plates containing 100 μ g/ml ampicillin and incubated overnight. Transformation efficiency should be greater than 1 x 10 ⁸ cfu/ μ g plasmid DNA. In addition, untransformed cells are tested for the appropriate antibiotic sensitivity and the absence of phage contamination.	
293FT Cell Line	Each lot of cells is tested for cell growth and viability post-recovery from cryo- preservation. Master Cell Banks are screened for viruses, mycoplasma, and sterility, and expression of the SV40 large T antigen is confirmed by western blot.	
Lentivirus Production	Using the reagents provided in the kit, the ViraPower [™] Packaging Mix and control lentiviral construct are cotransfected into 293FT cells using the protocol on page 105. Lentiviral supernatants are harvested 48 hours post-transfection, and the titer is determined using HT1080 cells. The control lentiviral construct must demonstrate a titer of greater than 1 x 10 ⁵ TU/ml.	
miR- <i>lacZ</i> ds oligo	The miR- <i>lacZ</i> positive ds control Oligo is functionally qualified by use in a ligation reaction as described in this manual.	
Sequencing Primers	Primers are lot-qualified by DNA sequencing experiments using the dideoxy chain termination technique.	
T4 DNA Ligase and 5X DNA Ligase Buffer	 T4 DNA Ligase is lot-qualified and must pass the following quality control assays: Functional absence of endonuclease and exonuclease activities Ligation/recut Ligation efficiency 5X DNA Ligase Buffer is functionally qualified with the enzyme and meets quality control specifications. 	

Product Qualification, Continued

PureLink [™] HQ Mini Plasmid Purification Kit	The PureLink TM HQ Mini Plasmid Purification Kit is functionally qualified by isolating a high copy number plasmid DNA from 1–2 $\times 10^9 E$. <i>coli</i> cells as described in this manual. The kit must produce the following results: • $A_{260}/A_{280} \ge 1.80$		
	• No detectable genomic DNA or RNA contamination on a 0.8% agarose gel		
	In addition, each kit component is sterile and is lot qualified for optimal performance.		
S.N.A.P. [™] MidiPrep Kit	Each component of the S.N.A.P. [™] MidiPrep is lot qualified for maximum performance in accordance with the S.N.A.P. [™] MidiPrep protocol. A sample of 4 columns from each lot is qualified in the following manner:		
	Binding Capacity $300 \ \mu g$ of pre-purified control plasmid is applied to the column and eluted according to the protocol. The quantity of plasmid DNA eluted from the column must be >200 μg .		
	Kit Performance		
	50 ml of control plasmid is grown in TOP10F´ cells and plasmid DNA is isolated using the S.N.A.P. [™] MidiPrep Kit.		
	Plasmid DNA is qualified in the following manner:		
	Visual inspection		
	Four samples are run on 0.8% agarose gel for a visual inspection. Each isolated plasmid must show supercoiled plasmid DNA with no RNA contamination.		
	Restriction Digest		
	Four 500 ng plasmid DNA samples are digested with <i>Apa</i> I restriction enzyme for 90 minutes. Restriction digest is run on a 0.8% agarose gel and must be >99% complete.		
	Endonuclease Activity		
	Four 500 ng plasmid DNA samples are incubated with 10 mM Mg ²⁺ for 4 hours and then run on a 0.8% agarose gel. Each sample must show no endonuclease contamination of supercoiled DNA present.		
DNA Ladders	Agarose gel analysis must show that bands are distinguishable after ethidium bromide staining.		
E-Gel [®] Agarose Gels	E-Gel [®] agarose gels are tested by running a mixture of DNA markers under standard running conditions described in this manual. Gels are visualized under UV light for proper staining, resolution, and migration of bands.		

Product Qualification, Continued

Opti-MEM [®] I Reduced Serum Medium	Opti-MEM [®] I Reduced Serum Medium is subjected to pH, osmolality, endotoxin, bacterial, fungal, and mycoplasma testing. The endotoxin level must be less than 1.0 EU/ml.
	Each lot of Opti-MEM [®] I is evaluated utilizing sensitive quantitative assays for its ability to support cloning efficiency of a murine myeloma cell line, and growth over multiple subcultures of an adherent cell line. Test lots of Opti- MEM [®] I Reduced Serum Medium at 2% (CHO growth) and 4% (Sp2 cloning) serum supplementation are compared to a previously approved Opti-MEM [®] I Reduced Serum Medium control.
	GIBCO [®] cell culture liquid products are prepared by an aseptic process for which each step has been validated to ensure that all products meet the industry standard sterility assurance level of 10 ⁻³ ; i.e., product that demonstrates a contamination level of no more than 1 of 1000 units during the manufacturing process. The highest level of sterility assurance (equal to or greater than 10 ⁻⁶) cannot be achieved without terminal sterilization which is harmful to the performance of cell culture products.

Purchaser Notification

Introduction	Use of the BioModule TM BLOCK-iT TM RNAi Units is covered under the licenses detailed below.
Information for European Customers	The 293FT Cell Line is genetically modified and carries the pUC-derived plasmid, pCMVSPORT6TAg.neo. As a condition of sale, use of this product must be in accordance with all applicable local legislation and guidelines including EC Directive 90/219/EEC on the contained use of genetically modified organisms.
Limited Use Label License No. 19: Gateway [®] Cloning Products	This product and its use is the subject of U.S. patents and/or other pending U.S. and foreign patent applications owned by Invitrogen Corporation. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for profit entity). The purchase of this product does not convey a license under any method claims in the foregoing patents or patent applications, or to use this product with any recombination sites other than those purchased from Invitrogen Corporation or its authorized distributors. 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 ¹⁰ purchased from Invitrogen Corporation or its authorized distributors. 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 or materials made by the employment of this product or its components or materials made by the employment of this product or its components for a scientific collaborator, provided that such transfer is not for any Commercial Purposes. The buyer may transfer information or materials made in thransfer is not for any Commercial Purposes, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Notwithstanding the preceding, any buyer who is employed in an academic or government institution may transfer materials made with this product to a third party who has a license from Invitrogen under the patents identified above to distributes such materials to rus of the product or its components to provide a service, information, or data; (3) use of the product or its components to provide a service, information, or data; (3) use of

Gateway [®] Clone Distribution Policy	For additional information about Invitrogen's policy for the use and distribution of Gateway [®] clones, see the section entitled Gateway[®] Clone Distribution Policy , page 176.
Limited Use Label License No. 23: GUS Control Vector	The GUS positive control vector in these products is claimed in patents and patent applic- ations licensed to Invitrogen 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: Lipofectamine™ 2000	The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components should determine if any licenses are required. Invitrogen Corporation will not assert a claim against the buyer of infringement of patents owned by Invitrogen and claiming this product based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic purposes is write product nor any of its components was used in the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product dueloped in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product dueloped in research

Limited Use Label License No. 51: Blasticidin and the Blasticidin Selection Marker Blasticidin and the blasticidin resistance gene (*bsd*) are the subject of U.S. patents 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, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500.

Limited Use Label License No. 61: E-Gel[®] Agarose Gels This product is the subject of U.S. and foreign patents owned by Invitrogen Corporation. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of 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 neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500.

Limited Use Label License No. 108: Lentiviral Technology The Lentiviral Technology (based upon the lentikat[™] system) is exclusively licensed from Cell Genesys, Inc., under U.S. patents and 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, including non-gene therapy research and target validation applications in laboratory animals (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of 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 neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500.

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 patents and applications in other countries for internal research purposes only. Use of these cell lines for Commercial Purposes requires a license from Invitrogen. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of 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 neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500.

Limited Use Label License No. 127: GFP with Heterologous Promoter

This product and its use is the subject of U.S. and foreign patents. 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. 181: Gateway[®]-adapted Viral Vectors The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of patents owned by Invitrogen and claiming this product based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500.

Limited Use Label License No. 198: Fluorescent Protein Products This product and its use is the subject of U.S. and foreign patents. Any use of this product by a commercial (for-profit) entity requires a separate license from either GE Healthcare or Invitrogen Corporation. For information on obtaining a commercial license to use this product, please refer to the contact information located at the bottom of this statement. 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). No rights are conveyed to modify or clone the gene encoding GFP contained in this product. 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. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that none of this product, or any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to use this product for purposes other than those permitted above, contact Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200.

Limited Use Label License No. 267: Mutant GFP Products This product and its use is the subject of U.S. and foreign patents.

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. The purchase or transfer of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of 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 neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500.

Limited Use Label License No. 272: Humanized GFP

This product is the subject of one or more of U.S. and foreign patents licensed by Invitrogen Corporation. This product is sold for research use only. Not for therapeutic or diagnostic use in humans.

Gateway[®] Clone Distribution Policy

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

References

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
- Ausubel, F. M., Brent, R., Kingston, R. E., Moore, D. D., Seidman, J. G., Smith, J. A., and Struhl, K. (1994) *Current Protocols in Molecular Biology*, Greene Publishing Associates and Wiley-Interscience, New York
- 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[™] 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
- Cole, C. N., and Stacy, T. P. (1985) Identification of Sequences in the Herpes Simplex Virus Thymidine Kinase Gene Required for Efficient Processing and Polyadenylation. Mol. Cell. Biol. 5, 2104-2113
- 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

References, Continued

- 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
- Gitlin, L., Karelsky, S., and Andino, R. (2002) Short Interfering RNA Confers Intracellular Antiviral Immunity in Human Cells. Nature 418, 430-434
- 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
- Graham, F. L., Smiley, J., Russell, W. C., and Nairn, R. (1977) Characteristics of a Human Cell Line Transformed by DNA from Human Adenovirus Type 5. J. Gen. Virol. *36*, 59-74
- 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
- Harrison, T., Graham, F., and Williams, J. (1977) Host-range Mutants of Adenovirus Type 5 Defective for Growth in HeLa Cells. Virology 77, 319-329
- 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 b-glucuronidase 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
- Lagos-Quintana, M., Rauhut, R., Yalcin, A., Meyer, J., Lendeckel, W., and Tuschl, T. (2002) Identification of tissue-specific microRNAs from mouse. Curr Biol 12, 735-739.
- 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

References, Continued

Li, W. X., and Ding, S. W. (2001) Viral Suppressors of RNA Silencing. Curr. Opin. Biotechnol. 12, 150-154

- Liebert, C. A., Watson, A. L., and Summers, A. O. (1999) Transposon Tn21, Flagship of the Floating Genome. Microbiol. Mol. Biol. Rev. 63, 507-522
- 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
- Miller, J. H. (1972) *Experiments in Molecular Genetics*, Cold Spring Harbor Laboratory, Cold Spring Harbor, New York
- 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
- Southern, P. J., and Berg, P. (1982) Transformation of Mammalian Cells to Antibiotic Resistance with a Bacterial Gene Under Control of the SV40 Early Region Promoter. J. Molec. Appl. Gen. 1, 327-339
- 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
- Weiss, B., Jacquemin-Sablon, A., Live, T. R., Fareed, G. C., and Richardson, C. C. (1968) Enzymatic Breakage and Joining of Deoxyribonucleic Acid. VI. Further Purification and Properties of Polynucleotide Ligase from *Escherichia coli* Infected with Bacteriophage T4. J. Biol. Chem. 243, 4543-4555
- 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 immunodeficiency 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

References, Continued

- 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

©2005, 2010 Invitrogen Corporation. All rights reserved.

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

Polybrene[®] is a registered trademark of Abbott Laboratories BigDye[®] is a registered trademark of Applied Biosystems