

Product Information

MISSION® Lenti-miRs – Lentiviral Constructs for miRNA Expression

Catalog Numbers **HLMIR0001 – HLMIR2781, MLMIR0001 – MLMIR2049**

Storage Temperature -70 °C

Product Description

MicroRNAs (miRNAs) are a class of genome-encoded nucleic acids that have been shown to regulate gene expression in a variety of organisms. These small, noncoding RNA molecules function by mediating transcript degradation, inhibiting translation, or a combination of these mechanisms. This type of regulation generally occurs by base pairing of the miRNA to a target sequence within the 3' UTR of a transcript. Thousands of miRNAs have been identified and classified via sequencing or bioinformatics approaches that are based on strongly conserved sequence motifs. The University of Manchester operates the publicly available *miRBase Sequence Database*, where microRNA data are managed and annotated.

Sigma's Mission Lenti-miRs express miRNAs from a common backbone, whose structure meets requirements for accurate Dicer processing.¹ Mature microRNA sequences are obtained from miRBase, and a partially complementary strand is designed to mimic the base pairing pattern in the backbone structure using a proprietary algorithm. Oligos containing the microRNA sequences are cloned into the TRC2-pLKO-puro vector (see figure 1). Each miRNA construct has been cloned and sequence verified.

Lentiviral transduction particles are produced from sequence-verified lentiviral plasmid vectors. Oligos containing the microRNA sequences are cloned into the TRC2-pLKO-puro vector (see figure 1). Co-transfection of this vector into the appropriate cell line with compatible packaging plasmids produces viral particles that can be used to transduce mammalian cells. The polymerase II promoter, elongation factor 1 alpha (EF1A), was chosen to drive miRNA expression.² The vector also contains elements needed for reverse transcription of viral RNA and integration of viral DNA into the host cell genome. Additionally, the

Woodchuck Hepatitis Post-Transcriptional Regulatory element³ (WPRE) is included, allowing for enhanced expression of transgenes delivered by lentiviral vectors.⁴ This lentiviral vector also carries a puromycin resistance gene for selection of cells. Unlike murine-based MMLV or MSCV retroviral systems, lentiviral-based particles permit efficient infection and integration of the specific miRNA construct into differentiated and non-dividing cells, such as neurons and dendritic cells,⁵ overcoming low transfection and integration difficulties when using these cell lines. Self-inactivating replication incompetent viral particles are produced in packaging cells (HEK293T) by co-transfection with compatible packaging plasmids.⁶⁻⁷

In addition, the lentiviral transduction particles are pseudotyped with an envelope G glycoprotein from Vesicular Stomatitis Virus (VSV-G), allowing transduction of a wide variety of mammalian cells including primary and embryonic stem cells.⁸ The lentiviral transduction particles are titered via a p24 antigen ELISA assay and pg/ml of p24 are then converted to transducing units per ml using a conversion factor.

To use Lenti-miRs, one needs to first select a cell type in which the miRNA of interest is expressed at low or undetectable levels. If unsure about the levels of miRNA expression in a particular cell line, it is critical that you determine those levels by conducting qRT-PCR or similar assay. We recommend using the MystiCq™ microRNA cDNA synthesis Mix. Starting with total RNA or RNA preparations pre-enriched for microRNAs, this kit provides all the components necessary to convert mature miRNAs into cDNA templates for qPCR. Once miRNAs have been chosen and quantified in the cell line of choice, cells can be transduced with lentiviral particles containing the miRNAs.

MISSION Lenti-miR activity may be assessed by the Dual-Luciferase Assay System which requires co-transfecting a Dual-Luciferase[®] Reporter (DLR[™]) plasmid with a complementary miRNA binding sequence (see Figure 2 in Appendix). Levels of target mRNA or protein can be assessed using a qRT-PCR assay or western Blotting, respectively.

Components/Reagents

The individual constructs are provided in Dulbecco's Modified Eagle's Medium with 10% heat-inactivated fetal bovine serum and penicillin-streptomycin.

Standard volume: 0.2 mL

Precautions and Disclaimer

These products are for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices. Though the lentiviral transduction particles produced are replication incompetent, it is highly recommended that they be treated as **Risk Group Level 2 (RGL-2)** organisms.⁹ Follow all published RGL-2 guidelines for handling and waste decontamination.

Storage/Stability

All components are stable for at least six months after receipt when stored at -70 °C. Avoid repeated freeze/thaw cycles, which will severely reduce functional viral titer.

Related Products

- Hexadimethrine Bromide, Catalog Number H9268
- Puromycin dihydrochloride, Ready Made solution, 10 mg/ml in H₂O, Catalog Number P9620
- Minimum Essential Medium containing 10% fetal calf serum or growth medium optimized for the specific cell line
- MISSION ExpressMag[®] 96-Well Magnetic Kit, Catalog Number SHM02
- MystiCq[®] microRNA cDNA Synthesis Mix, Catalog Number MIRRT
- KiCqStart[®] SYBR[®] Green qPCR ReadyMix[™], Catalog Number KCQS00
- mirPremier[®] microRNA Isolation Kit, Catalog Number SNC10, SNC50

- MISSION microRNA Mimics
- MISSION Lenti microRNA Inhibitors
- MISSION Synthetic microRNA Inhibitors
- MISSION TRC2 pLKO.5-puro-CMV-TurboGFP[™] Positive Control Transduction Particles, Catalog Number SHC203V
- MISSION pLKO.1-puro Non-Mammalian shRNA Negative Control Transduction Particles, Catalog Number SHC002V
- MISSION Lenti microRNA, ath-miR416, Negative Control 1 Transduction Particles, Catalog Number NCLMIR001
- MISSION Lenti microRNA, cel-miR-243-3p, Negative Control 2 Transduction Particles, Catalog Number NCLMIR002

Procedure for the Use of MISSION Lentiviral microRNA Transduction Particles:

Puromycin titration (kill curve) should be performed before starting transduction when working with a new cell type.

1. Plate 1.6×10^4 cells into wells of a 96-well plate with 120 μ l fresh media.
2. The next day add 0.5-10 μ g/ml of puromycin to selected wells.
3. Examine viability every 2 days.
4. Culture for 3 – 14 days depending on the growth rate of the cell type and the length of time that cells would typically be under selection during a normal experimental protocol. Replace the media containing puromycin every 3 days. The minimum concentration of puromycin that causes complete cell death after the desired time should be used for that cell type and experiment.

Day 1

Seed the mammalian cell line of choice in complete medium into culture dishes or multiwell plates 24 hours prior to transduction. Plating density should be such that cells are 50-80% confluent at the time of transduction.

Day 2

Thaw the lentiviral stock slowly on ice. Gently spin down material in tubes before opening. Add hexadimethrine bromide (the chemical equivalent of Polybrene) to the cells at a final concentration of 8 µg/ml.

Note: Hexadimethrine bromide enhances transduction of most cell types. However, some cells, such as primary neurons, are sensitive to hexadimethrine bromide. When using sensitive cells, do not add the hexadimethrine bromide and the cells should still be transduced.

Following addition of hexadimethrine bromide, gently swirl the plate to mix. Add the appropriate amount of viral particles at a suitable multiplicity of infection (MOI) and swirl the plate gently to mix. Incubate the cell-viral particle mixture at 37 °C overnight.

Multiplicity of Infection (MOI) is the number of transducing lentiviral particles per cell. It is highly recommended that for each new cell type to be transduced, a range of MOI be tested. To calculate MOI: (total number of cells per well) x (desired MOI) = total transducing units needed (TU)

(total TU needed) / (TU/ml reported on C of A) = total ml of lentiviral particles to add to each well

Notes

- When transducing a lentiviral construct into a cell line for the first time, it is recommended that a range of MOIs (0.5-20) be used to find the optimum degree of transduction efficiency.
- When overnight incubation presents a toxicity concern, cells may be incubated for as little as 4 hours before changing the medium. Cells can be transduced in reduced medium volumes to increase transduction efficiencies.

Day 3

Remove the viral particle-containing medium and replace it with fresh, pre-warmed complete culture medium.

Day 4

For puromycin selection, remove the medium and replace it with fresh, complete medium that contains the appropriate amount of puromycin for selection of transduced cells. Proceed to day 5.

Note: When the appropriate concentration of puromycin for a specific cell type is unknown, perform a kill curve experiment, as described above. Typically, puromycin concentrations ranging from 0.5-10 µg/ml are sufficient to kill most untransduced mammalian cell lines. Excess puromycin will cause undesired phenotypic responses in most cell types.

Day 5 and forward

Replace medium with fresh, puromycin-containing medium every 3-4 days until resistant colonies can be identified (generally, 10-12 days after selection). If isolated colonies are desired, pick a minimum of 5 puromycin-resistant colonies and expand each clone to assay. Alternatively, miRNA expression and function may be evaluated with a puro-selected culture, containing all transduced cells.

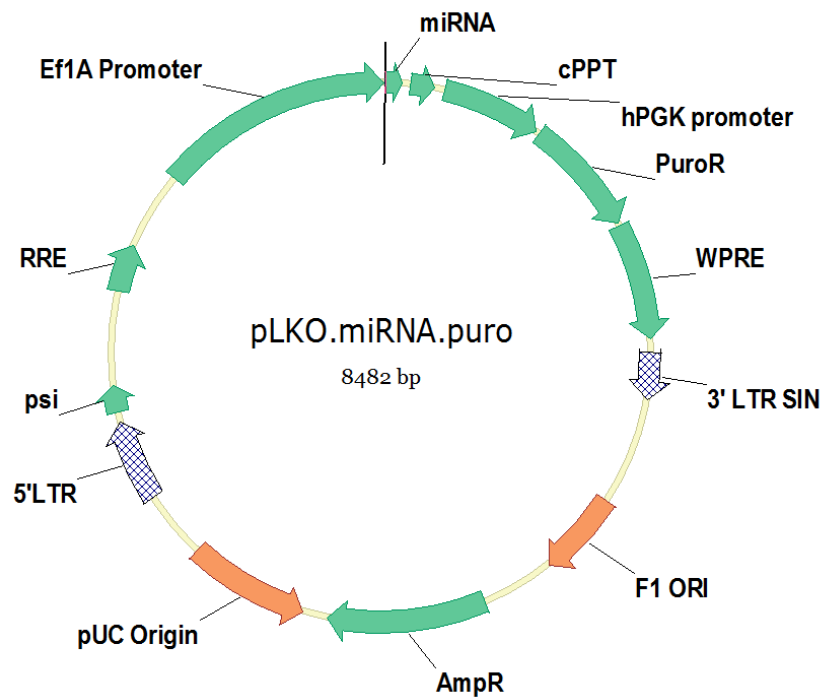
Note: Due to the random integration of the lentivirus into the genome, varying levels of miRNA may be seen from different puromycin-resistant clones. Testing a number of puromycin-resistant clones will allow a determination of which clone is optimal.

Assessing miRNA expression and function

miRNA expression from Lenti-miRs can be evaluated by RT-qPCR. We recommend the MystiCq® microRNA cDNA Synthesis Mix and MystiCq® microRNA qPCR assay primers. To evaluate function, reporter assays or levels of target protein and/or messenger RNA levels may be evaluated. miRNA functional studies may require simultaneous analyses of both mRNA and protein expression. While qRT-PCR can be used to assess levels of target transcript, Western analysis or other validated immunoassays are used to investigate the impact on protein quantity. Reporter assays, such as a dual luciferase reporter assay, are used to study the interaction between miRNAs and their target sites.

Figure 1. TRC2 Lentiviral Plasmid vector TRC2-pLKO-puro Features

Name	Description
EF-1alpha	Elongation factor-1alpha Promoter
cppt	Central polypurine tract
hPGK	Human phosphoglycerate kinase eukaryotic promoter
PuroR	Puromycin resistance gene for mammalian selection
WPRE	Woodchuck Hepatitis PostTranscriptional Regulatory element
SIN/3' LTR	3' self inactivating long terminal repeat
f1 ori	f1 origin of replication
AmpR	Ampicillin resistance gene for bacterial selection
pUC ori	pUC origin of replication
5' LTR	5' long terminal repeat
Psi	RNA packaging signal
RPE	Rev response element
miRNA	miRNA sequence (clone specific)



Troubleshooting Guide

Problem	Cause	Solution
No Transduction of cells	Viral stock stored incorrectly	Store stocks at -70 °C. Do not freeze/thaw more than 2 times.
Low miRNA expression detected due to low transduction efficiency	Population of cells not transduced efficiently	Select for transduced cells with puromycin, and allow selected population to become mitotic.
	Hexadimethrine bromide not included during transduction	Transduce in the presence of hexadimethrine bromide.
	Non-dividing cell type used	Transduce at a higher MOI, or evaluate transduction enhancement reagents, such as Express Mag.
	It is unknown how efficiently cells can be transduced with VSV-G pseudotyped lentiviral particles	Try transducing cells with control lentiviral particles such as MISSION® TRC2 pLKO.5-puro-CMV-TurboGFP™ Positive Control Transduction Particles (SHC203V) to establish experimental parameters for transductions.
	Cells were harvested and assayed too soon after transduction	Harvest cells 72 hours after transduction, and not earlier. Alternatively, results may be improved by placing cells under puromycin selection because untransduced cells will be killed.
Low miRNA impact due to target choice or cell line variability in expression of the miRNA	miRNA is not expressed at a high enough level for analysis in cell line	Evaluate miRNA levels in cell type of choice via qRT-PCR using the MystiCq™ microRNA cDNA Synthesis Mix kit. Consider alternative cell lines if endogenous miRNA levels are high. Use a combination of assays to test the miRNA target, such as qRT-PCR, Western analysis, and reporter assays such as a dual luciferase assay
No signal from reporter assay (such as a dual luciferase assay)	Reporter assay might not be working correctly	Double check to make sure reporter plasmid target sequence and orientation is correct.
Cytotoxic effects observed after transduction	Expressed miRNA causes cell cycle arrest or apoptosis	Be sure expressed miRNA does not have an adverse phenotype on your cell line of interest.

Cytotoxic effects observed after transduction (continued)	Hexadimethrine bromide was used during transduction	Be sure that cells are not sensitive to hexadimethrine bromide. Omit the hexadimethrine bromide during the transduction.
	Too much puromycin was used for selection	Determine the puromycin sensitivity of the cells by performing a kill curve and use the minimum concentration required to kill the untransduced cells.

Control Selection Table

Recommended Control	Objective
Negative Control: Untreated Cells	Untreated cells will provide a reference point for comparing all other samples.
Negative control: Transduction with lentiviral particles containing a sequence targeting no known mammalian genes (designed against Turbo GFP sequence), Catalog Number SHC002V	MISSION® pLKO.1-puro Non-Mammalian shRNA Control Transduction Particles provides a reference point to compare lentiviral transduced samples.
Negative control: Transduction with lentiviral particles containing a sequence targeting no known mammalian genes, Catalog Number SHC016V	MISSION® pLKO.1-puro Non-Mammalian shRNA Control Transduction Particles provides a reference point to compare lentiviral transduced samples.
Negative control: Transduction with MISSION Lenti microRNA, ath-miR416, Negative Control 1, Sequence from <i>Arabidopsis thaliana</i> with no homology to human and mouse gene sequences, Catalog Number NCLMIR001	MISSION® pLKO.miRNA-puro Non-Mammalian microRNA Control Transduction Particles provides a reference point to compare lentiviral transduced samples.
Negative control: Transduction with MISSION Lenti microRNA, cel-miR-243-3p, Negative Control 2, Sequence from <i>Caenorhabditis elegans</i> with no homology to human and mouse gene sequences, Catalog Number NCLMIR2	MISSION® pLKO.miRNA-puro Non-Mammalian microRNA Control Transduction Particles provides a reference point to compare lentiviral transduced samples.
Positive control: Transduction with lentiviral particles containing Turbo GFP, Catalog Number SHC203V	MISSION® TRC2 pLKO.5-puro-CMV-TurboGFP™ Positive Control Transduction Particles can be used to monitor transduction efficiency.

Cell Type Table

The cell types listed below have been successfully infected by pLKO.2-puro based lentiviral particles. Optimal conditions will need to be determined for your experimental needs. For the most updated cell line list, and some guidelines for conditions, please visit:

<http://www.sigmaaldrich.com/life-science/functional-genomics-and-rnai/shrna/learning-center/getting-started.html>

Cell lines, human	Cell Type	Cell lines, human	Cell Type	Primary cells human	Cell Type
HEK293	embryonic kidney cells	A431	epidermal carcinoma	dendritic	immature dendritic
HeLa	cervical adenocarcinoma	THP1	monocytic	T-cells	lymphocytes
A549	lung adenocarcinoma	RAW264.7	macrophage	epithelial	prostate
H1299	lung carcinoma	SH-SY5Y	brain neuroblastoma	fibroblasts	primary mammary
HT29-D4	colon carcinoma	HCN-1A	brain cortical neuron	Primary cells, other species	Cell Type
HepG2	hepatocellular carcinoma	SupT1	T-cells	ECS	mouse embryonic stem cells
HCT116	colon carcinoma	BJ-TERT	diploid fibroblasts	fibroblasts	Mouse embryonic fibroblasts
MCF7	breast carcinoma	Cell lines, mouse	Cell Type	MC3T3-E1	mouse bone marrow derived
MCF10A	breast carcinoma	NIH3T3	fibroblast	molar mesenchymal	mouse embryonic mesenchymal
Panc-1	pancreatic epitheliod carcinoma	Primary cells, human	Cell Type	cardiomyocytes	rat neonatal cardiomyocytes
PC3	prostate carcinoma	astrocytes	normal		
DU145	prostate carcinoma	C3H10T1/2	mesenchymal		

References

1. Gu, S., et al., The Loop Position of shRNAs and Pre-miRNAs is Critical for the Accuracy of Dicer Processing in Vivo. *Cell*, 151(4), 900-911 (2012).
2. Lebbink, R. J., et al., Polymerase II Promoter Strength Determines Efficacy of microRNA Adapted shRNAs. *Plos One*, 6(10), e260213 (2011).
3. Donello, J.E., et al., Woodchuck hepatitis virus contains a tripartite posttranscriptional regulatory element. *J Virol.*, 72, 5085-5092 (1998).
4. Zufferey, R., et al., Woodchuck hepatitis virus posttranscriptional regulatory element enhances expression of transgenes delivered by retroviral vectors. *J Virol.*, 73, 2886-2892 (1999).
5. Stewart, S.A., et al., Lentivirus-delivered stable gene silencing by RNAi in primary cells. *RNA*, 9, 493-501 (2003).
6. Zufferey, R., et al., Multiply attenuated lentiviral vector achieves efficient gene delivery *in vivo*. *Nat. Biotechnol.* 15, 871-885 (1997).
7. Zufferey, R., et al., Self-inactivating lentivirus vector for safe and efficient *in vivo* gene delivery. *J Virol.*, 72, 9873-9880 (1998).
8. Burns, JC., et al., 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 (1993).
9. NIH Guidelines for Research Involving Recombinant DNA Molecules (NIH Guidelines) 2002 (<http://11www4.od.nih.gov/oba>).

KiCqStart and MystiCq are registered trademarks of Quanta BioSciences Inc. ExpressMag, MISSION and mirPremier are registered trademarks, and ReadyMix is a trademark, of Sigma-Aldrich Co. LLC SYBR is a registered trademark of Life Technologies TurboGFP is a trademark of Evrogen Co.

IDC,SS,EE,PHC 05/14-1

Appendix

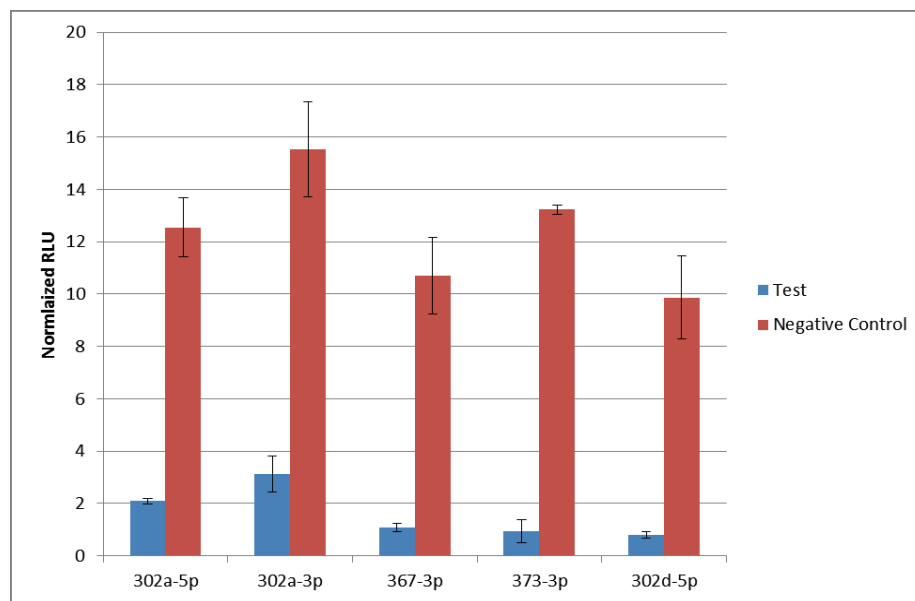


Figure 2. Lenti-miRs were transduced into HeLa cells. After two weeks under puromycin selection, the Dual-Luciferase® Reporter (DLR™) plasmid, with a complementary miRNA binding sequence, was transfected into the cells. miRNA activity is shown for 5 miRNAs, as compared to a negative control, using the Dual-Luciferase Assay System.

©2014 Sigma-Aldrich Co. LLC. All rights reserved. SIGMA-ALDRICH is a trademark of Sigma-Aldrich Co. LLC, registered in the US and other countries. Sigma brand products are sold through Sigma-Aldrich, Inc. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see product information on the Sigma-Aldrich website at www.sigmaaldrich.com and/or on the reverse side of the invoice or packing slip.