

Product Information

MISSION® Lentiviral Transduction Particles – Lentiviral MicroRNA Inhibitors

Catalog Numbers **HLTUD0001-HLTUD2235,**
MLTUD0001-MLTUD1405, CSTTUD,
HLTUD001C, HLTUD002C

Storage Temperature –70 °C

TECHNICAL BULLETIN

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.

Individual microRNA inhibitors are designed using a proprietary algorithm, which is based on the work of Haraguchi, T, *et al.*¹, and in collaboration with Dr. Hideo Iba, University of Tokyo. This algorithm utilizes the tough decoy (TuD) design. Each miRNA inhibitor construct has been cloned and sequence verified to ensure a match to the target miRNA.

The Lentiviral Transduction Particles are produced from sequence-verified lentiviral plasmid vectors. The Lentiviral microRNA Inhibitors 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 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 Element² (WPRE) is included, allowing for enhanced expression of transgenes delivered by lentiviral vectors.³ 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 inhibitor construct into differentiated and non-dividing cells, such as neurons and dendritic cells,⁴ 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.⁵⁻⁶

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.⁷ 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 the lentiviral microRNA inhibitors one needs to first select the miRNA targets to be inhibited/ investigated. 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 microRNAs into cDNA templates for qPCR. Once miRNA targets have been chosen and quantified in the cell line of choice, cells can be transduced with lentiviral particles containing the miRNA inhibitors. Levels of target mRNA or protein can then be assessed using a qRT-PCR assay (optional) or Western Blotting, respectively. The lentiviral miRNA inhibitors can also be used to investigate the functional role of the miRNA being studied.

Components/Reagents

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

Standard Volume:

- 0.2 mL

Custom Volumes:

- 0.1 mL
- 1.0 mL
- 2.0 mL
- 5.0 mL
- 10.0 mL

Orders of 25 or fewer lenti microRNA inhibitors are provided in individual vials. Orders of >25 lenti microRNA inhibitors are provided in a 96-well plate. 96-well plates are provided with a CD containing plate map positions.

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.⁸ 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 Inhibitor, ath-miR416, Negative Control 1, Sequence from *Arabidopsis thaliana* with no homology to human and mouse gene sequences, Catalog Number [HLTUD001C](#)
- MISSION Lenti microRNA Inhibitor, cel-miR-243-3p, Negative Control 2, Sequence from *Caenorhabditis elegans* with no homology to human and mouse gene sequences, Catalog Number [HLTUD002C](#)
- MISSION TRC2 pLKO.5-puro-CMV-TurboGFP[™] Positive Control Transduction Particles, Catalog Number [SHC203V](#) or MISSION[®] TRC2 pLKO.5-puro Empty Vector Control Transduction Particles puro, Catalog Number [SHC201V](#)
- qRT-PCR Reagents, please visit <http://www.sigma-aldrich.com/pcr>
- Prestige Antibodies, please visit <http://www.sigma-aldrich.com/prestige>

Procedure for the Use of MISSION Lentiviral microRNA Inhibitor Transduction Particles

Day 1

Plate the mammalian cell line of choice in complete medium 24 hours prior to transduction. Take into account the length of time that the cells will be cultured prior to performing miRNA target inhibition analysis when determining plating density. Typically cells are transduced at 50-80% confluency.

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 $\mu\text{g/ml}$.

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

If placing cells under 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. Typically, puromycin concentrations ranging from 0.5-10 $\mu\text{g/ml}$ are sufficient to kill most untransduced mammalian cell lines.

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

- Plate 1.6×10^4 cells into wells of a 96-well plate with 120 μL fresh media.
- The next day add 0.5-10 $\mu\text{g/ml}$ of puromycin to selected wells.
- Examine viability every 2 days.
- 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.

Note: Excess puromycin can cause many 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). Pick a minimum of 5 puromycin-resistant colonies and expand each clone to assay for inhibition of the target miRNA.

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

Assessing loss of miRNA function

miRNA targets can be validated by quantitating target protein and/ or messenger RNA levels in response to miRNA downregulation. 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. Many researchers find it is necessary to use a combination of assays to assess miRNA target inhibition.

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

Name	Description
hU6	U6 Promoter
cppt	Central polypurine tract
hPGK	Human phosphoglycerate kinase eukaryotic promoter
PuroR	Puromycin resistance gene for mammalian selection
WPRE	Woodchuck Hepatitis Post-Transcriptional 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
RRE	Rev response element
miRNA inhibitor	miRNA inhibitor sequence (clone specific)

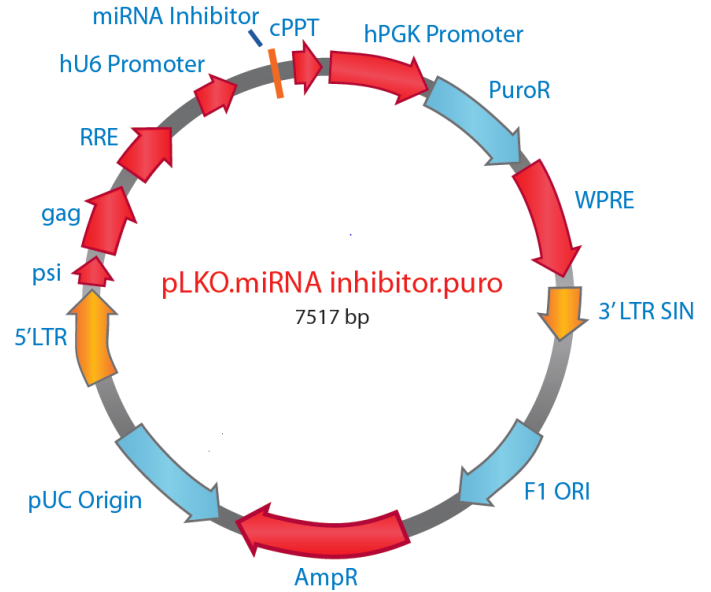
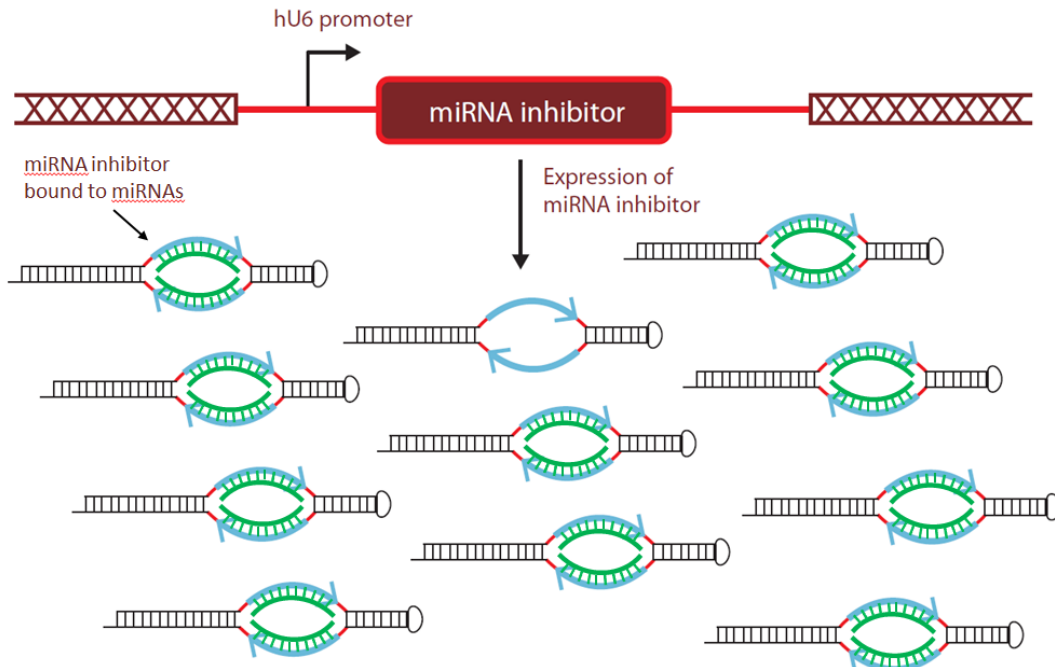


Figure 2. Lentiviral miRNA inhibitor expression regulates miRNA function

Expression of the miRNA inhibitor is driven by the U6 promoter upon genomic integration of the lentiviral transfer vector into the host cell post-transduction. miRNA inhibitors are able to competitively bind specific miRNAs and prevent them from regulating their endogenous targets¹.



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 inhibition 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 ExpressMag.
	MOI is too low	Transduce at a higher MOI, or evaluate transduction enhancement reagents, such as ExpressMag.
	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 [®] Lenti microRNA Inhibitor, ath-miR416, Negative Control 1, Catalog Number HLTUD001C, MISSION [®] TRC2 pLKO.5-puro-CMV-TurboGFP [™] Positive Control Transduction Particles (SHC203V), or MISSION [®] TRC2 pLKO.5-puro Empty Vector Control Transduction Particles puro (SHC201V) 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 inhibition detected due to target choice or cell line variabilities in expression of the miRNA	microRNA 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. 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	Transfect cells with ready-to-use MISSION miRNA mimics along with reporter plasmid. These small, double-stranded RNA molecules, designed to mimic endogenous mature miRNA molecules when introduced into cells, will aid in assessing if the assay is working.
Cytotoxic effects observed after transduction	Target miRNA is essential for cell viability	Be sure that target miRNA is not essential for cell growth or viability.
	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 miRNA inhibitor sequence from <i>Arabidopsis</i>	MISSION Lenti microRNA Inhibitor, ath-miR416, Negative Control 1, Catalog Number HLTUD001C These viral particles contain a sequence designed to inhibit a miRNA found in <i>Arabidopsis</i> that will not target any mammalian miRNAs. Side-by-side transductions using these control particles and experimental lenti microRNA inhibitors will allow for observation of cellular effects of the transduction process, the puromycin selection process, and will provide control cells for downstream applications such as qRT-PCR and Western blot analysis.
Negative Control: Transduction with lentiviral particles containing a miRNA inhibitor sequence from <i>C. elegans</i>	MISSION Lenti microRNA Inhibitor, cel-miR-243-3p, Negative Control 2, Catalog Number HLTUD002C These viral particles contain a sequence designed to inhibit a miRNA found in <i>Caenorhabditis elegans</i> that will not target any mammalian miRNAs. Side-by-side transductions using these control particles and experimental lenti microRNA inhibitors will allow for observation of cellular effects of the transduction process, the puromycin selection process, and will provide control cells for downstream applications such as qRT-PCR and Western blot analysis.

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.sigmaldrich.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 epithelioid carcinoma	Primary cells, human	Cell Type	cardiomyocytes	rat neonatal cardiomyocytes
PC3	prostate carcinoma	astrocytes	normal		
DU145	prostate carcinoma	C3H10T1/2	mesenchymal		

References

1. Haraguchi, T., *et al.*, Vectors expressing efficient RNA decoys achieve the long-term suppression of specific microRNA activity in mammalian cells. *Nucleic Acids Res.*, 37, (2009).
2. Donello, J.E., *et al.*, Woodchuck hepatitis virus contains a tripartite posttranscriptional regulatory element. *J Virol.*, 72, 5085-5092 (1998).
3. Zufferey, R., *et al.*, Woodchuck hepatitis virus posttranscriptional regulatory element enhances expression of transgenes delivered by retroviral vectors. *J Virol.*, 73, 2886-2892 (1999).
4. Stewart, S.A., *et al.*, Lentivirus-delivered stable gene silencing by RNAi in primary cells. *RNA*, 9, 493-501 (2003).
5. Zufferey, R., *et al.*, Multiply attenuated lentiviral vector achieves efficient gene delivery *in vivo*. *Nat. Biotechnol.* 15, 871-885 (1997).
6. Zufferey, R., *et al.*, Self-inactivating lentivirus vector for safe and efficient *in vivo* gene delivery. *J Virol.*, 72, 9873-9880 (1998).
7. Burns, J.C., *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).
8. NIH Guidelines for Research Involving Recombinant DNA Molecules (NIH Guidelines) 2002 (<http://www4.od.nih.gov/oba>).

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

Limited Use Licenses

Use of this product for Commercial Purposes requires a license from Sigma-Aldrich Corporation. The purchase of this product conveys to the buyer the nontransferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The 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. Commercial Purposes means any activity by a party for consideration, but excludes not-for-profit core facilities providing services within their own research institutions at cost. Core facilities are invited to join Sigma-Aldrich's RNAi Partnership Program. Details of Sigma-Aldrich's RNAi Partnership Program can be found at www.sigma.com/rpp.

This product is licensed under U.S. Pat. Nos. 5,817,491; 5,591,624; 5,716,832; 6,312,682; 6,669,936; 6,235,522; 6,924,123 and foreign equivalents from Oxford BioMedica (UK) Ltd., Oxford, UK, and is provided for use in academic and commercial *in vitro* and *in vivo* research for elucidating gene function, and for validating potential gene products and pathways for drug discovery and development, but excludes any use of LentiVector® technology for: creating transgenic birds for the purpose of producing useful or valuable proteins in the eggs of such transgenic birds, the delivery of gene therapies, and for commercial production of therapeutic, diagnostic or other commercial products not intended for research use where such products do not consist of or incorporate a lentiviral vector. Information about licenses for commercial uses excluded under this license is available from Oxford BioMedica (UK), Ltd., Medawar Center, Oxford Science Park, Oxford OX4 4GA UK enquiries@oxfordbiomedica.co.uk or BioMedica Inc., 11622 El Camino Real #100, San Diego CA 92130-2049 USA. LentiVector is a registered US and European Community trademark of Oxford BioMedica plc.

This product (based upon the lentikat system) is sub-licensed from Invitrogen Corporation under U.S. Patent Nos. 5,686,279, 5,834,256, 5,858,740; 5,994,136; 6,013,516; 6,051,427, 6,165,782, and 6,218,187 and corresponding patents and applications in other countries for internal research purposes only. Use of this technology for gene therapy applications or bioprocessing other than for nonhuman research use requires a license from Cell Genesys, Inc. Please contact Cell Genesys, Inc. at 342 Lakeside Drive, Foster City, California 94404. Use of this technology to make or sell products or offer services for consideration in the research market requires a license from Invitrogen Corporation, 1600 Faraday Ave., Carlsbad, CA 92008.

This product is for non-clinical research use only. It is not to be used for commercial purposes. Use of this product to produce products for sale or for diagnostic, therapeutic or high throughput drug discovery purposes (the screening of more than 10,000 compounds per day) is prohibited. This product is sold under license from Invitrogen Corporation. In order to obtain a license to use this product for these commercial purposes, contact The Regents of the University of California. This product or the use of this product is covered by U.S. Patent No. 5,624,803 owned by The Regents of the University of California.

All Mission TRC II Lentiviral backbone-containing products contain a specific genetic component (WPRE), which is licensed from the Salk Institute for Biological Studies and covered under the following patents:

U.S. Patent No. 6,136,597, U.S. Patent No. 6,284,469, U.S. Patent No. 6,312,912, U.S. Patent No. 6,287,814.

Purchaser Notification:

Licensee has a license to sell the Product containing WPRE, under the terms described below. Any use of WPRE outside of Licensee's Product or the Product's intended use, requires a license as detailed below. Before using the Product containing WPRE, please read the following license agreement. If you do not agree to be bound by its terms, contact Licensee within 10 days for authorization to return the unused Product containing WPRE and to receive a full credit.

Licensee grants you a non-exclusive license to use the enclosed Product containing WPRE in its entirety for its intended research use. The Product containing WPRE is being transferred to you in furtherance of, and reliance on, such license. Any use of WPRE outside of Licensee's Product or the Product's intended use including for Commercial Purposes, requires a license from the Salk Institute for Biological Studies. Commercial Purposes means any activity by a party for consideration, but excludes not-for-profit core facilities providing services within their own research institutions at cost. This license agreement is effective until terminated. You may terminate it at any time by destroying all Products containing WPRE in your control. It will also terminate automatically if you fail to comply with the terms and conditions of the license agreement. You shall, upon termination of the license agreement, destroy all Products containing WPRE in your control, and so notify Licensee in writing.

This License shall be governed in its interpretation and enforcement by the laws of the State of California.

Contact for WPRE Licensing:

The Salk Institute for Biological Studies
10010 North Torrey Pines Road
La Jolla, CA 92037
Attn.: Office of Technology Management
Phone: (858) 453-4100 extension 1703
Fax: (858) 546-8093

This Product is covered by US and foreign patent applications or patents and other proprietary intellectual property rights owned by CSHL ("CSHL shRNA IP Rights"), including U.S. Patent Nos. 8,153,776, 8,202,846, 8,383,599, 8,829,264, and EP1546174.

Subject to acceptance and all terms and conditions of this License, sale of the Product to Buyer by Sigma-Aldrich, Co. (acting under its license, an "Authorized Sale") conveys to Buyer only the nonexclusive, nontransferable right (with no right to sublicense) under the shRNA IP Rights to use the Product solely for Customer's internal research purposes, and only at its facility where the Product is delivered by Sigma-Aldrich, Co.

The Product is for research use only and may not be used *in vitro* or *in vivo* for any diagnostic, preventative, therapeutic or vaccine application, or used (directly or indirectly) in humans for any purpose.

Non-Profit Buyers. If Buyer is a Non-Profit Entity, then the following additional restrictions will apply:

Customer obtains no right to use, develop or otherwise exploit the product for any commercial purpose.

Commercial Buyers. If Buyer is a Commercial Entity, then the following additional restrictions will apply:

A Product sale is an Authorized Sale only if Buyer has already entered into a separate written agreement that has been executed by CSHL or Hairpin Technologies, that covers the CSHL shRNA IP Rights, and that is then currently in effect. Any delivery or transfer of Product to Customer outside of an Authorized Sale is void, conveys no implied or express right under this license and Customer will immediately return Product to Sigma-Aldrich for a refund.

"Commercial Entity" means any entity or organization other than a Non-Profit Entity.

"CSHL" means Cold Spring Harbor Laboratory.

"Hairpin Technologies" means Hairpin Technologies, Inc. located at 2200 Smithtown Avenue, Ronkonkoma, NY 11779, www.hairpintechnologies.com.

"Non-Profit Entity" means any college, university or governmental entity (including without limitation, governmental and quasi-governmental institutes and research laboratories), or any non-profit scientific, research or educational organization of the type described in section 501(c)(3) of the Internal Revenue Code or qualified under a state non-profit organization statute.

"Product" means a product (including, without limitation, expression vectors encoding a shRNA, the design, manufacture or use of which (in whole or in part) is the subject of the shRNA IP Rights, and is deemed to include all components, progeny, reproductions, modified versions and other derivatives thereof.

This license is subject to a license from CSHL or Hairpin Technologies, and CSHL and Hairpin Technologies reserves all other rights under its license. For information on licensing rights for Commercial Entities, including use of this product for purposes other than research and trial licenses, please contact Hairpin Technologies, Inc. at info@hairpintechnologies.com or call (631) 881-0844.

KT,RC,PHC,MAM 08/17-1