

Overview

OmicsLink™ shRNA clone collections consist of lentiviral and other mammalian expression vector-based small hairpin RNA (shRNA) clones against genome-wide target genes from human, mouse and rat. A set of four expression constructs for every target gene ensures high knockdown efficiency with minimal off-target effects.

GeneCopoeia provides four shRNA constructs for every target gene and guarantees that at least one of the four will have a knockdown effect of 70% on corresponding gene expression as determined by qRT-PCR.

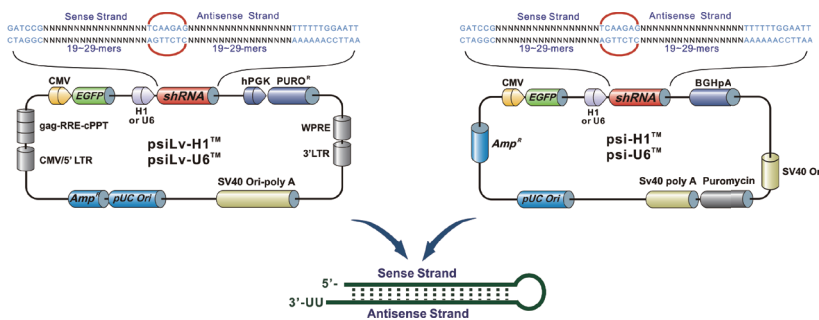


Figure 1. Lentiviral expression vector-based shRNA clones with H1 or U6 promoter.

Selected publications

1. Wong, L. C. et al. (2010) Fully-automated image processing software to analyze calcium traces in populations of single cells. *Cell Calcium*, 48 (5): 270-274
2. St John, MA et al (2009) Proinflammatory mediators upregulate snail in head and neck squamous cell carcinoma. *Clin Cancer Res* 15: 6018
3. Kim, YD et al (2011) NSrp70 is a novel nuclear speckle-related protein that modulates alternative pre-mRNA splicing in vivo. *Nucl. Acids Res.* (2011) doi: 10.1093/nar/gkq1267
4. Gupta, S. et al. (2010) HSP72 Protects Cells from ER Stress-induced Apoptosis via Enhancement of IRE1α-XBP1 Signaling through a Physical Interaction. *PLoS Biol* 8 (7): e1000410

Advantages

Guaranteed knockdown

- Four shRNA sequences are selected through a proprietary algorithm
- Guarantee at least one of the four will have a knockdown effect of 70% on corresponding gene expression as determined by qRT-PCR

Versatile deliveries

- Available in expression plasmid or lentiviral particles for gene silencing in virtually all cell types including difficult-to-transfect and non-dividing cells

Markers and reporters

- Enable stable cell line selection and expression verification

Fully sequenced

- The expression cassettes of all shRNA clones are fully sequenced

Vector list

Vector	Promoter	Selection marker	Reporter gene	Viral type
psiLv-H1	H1	Puromycin	eGFP	Lenti
psiLv-U6	U6	Puromycin	eGFP	Lenti
psiLv-mH1	H1	Puromycin	mCherry	Lenti
psiLv-mU6	U6	Puromycin	mCherry	Lenti
psiLV-nH1	H1	Puromycin	N/A	Lenti
psiLV-nU6	U6	Puromycin	N/A	Lenti
psi-H1	H1	Puromycin	eGFP	N/A
psi-U6	U6	Puromycin	eGFP	N/A
psi-mH1	H1	Puromycin	mCherry	N/A
psi-mU6	U6	Puromycin	mCherry	N/A
psi-nH1	H1	Puromycin	N/A	N/A
psi-nU6	U6	Puromycin	N/A	N/A

Related products

Product	Description
Lentifect™ Custom Lentivirus Production Services	Up to 10 ¹⁰ copies/ml, crude or purified lentiviral particles
Lenti-Pac™ HIV or FIV Expression Packaging Kits	HIV or FIV-based lentiviral packaging plasmids, EndoFectin™ Lenti Transfection Reagent, eGFP control clone, TiterBoost™ Reagent
293Ta Lentiviral Packaging Cell Line	1.5 x 10 ⁶ cells
GCI-L3 Chemically Competent E. coli Cells	Competent cells optimized for lentiviral expression clones
OmicsLink™ Lentiviral ORF Expression Clones	45,000 human and mouse clones available in lentiviral and non-viral expression vectors
miExpress™ Precursor miRNA Expression Clones	1048 human, 672 mouse, and 408 rat clones available in lentiviral and non-viral expression vectors

To order

To search and order shRNA clones, please visit www.genecopoeia.com.

GeneCopoeia, Inc.
 9620 Medical Center Drive, Suite 101
 Rockville, MD 20850, USA
 Email inquiry@genecopoeia.com
 Tel +1 (301) 762-0888
 Toll free +1 (866) 360-9531
 Fax +1 (301) 762-3888
 Website www.genecopoeia.com