



## Scrambled Control Lentifect™ Purified Lentiviral Particles •

Cat No. LPP- CSHCTR001-LVRU6MH-025, LPP- CSHCTR001-LVRU6MH-100

Ready-to-use lentiviral particles for the transduction of a variety of mammalian cells including difficult-to-transfect, primary, stem and non-dividing cells.

### Description

- Produced with a standardized protocol using highly purified plasmids and EndoFectin Lenti™ transfection and TiterBoost™ reagents
- A short non-coding stuffer instead of a specific coding insert downstream of the U6 promoter
- Hygromycin resistance marker for selection of stably transduced cells

### Contents and storage

Provided as 1 vial of 25 µl or 4 vials of 25 µl of purified lentiviral particles with titers of  $\sim 1 \times 10^8$  TU/ml.

Lentifect particles are shipped on dry ice and **must be stored at –80°C immediately upon receipt**. Avoid repeated freeze-thaw cycles as this will reduce titers.

### Quality control

The lentiviral expression construct was validated by full-length sequencing, restriction enzyme digestion and PCR-size validation using gene-specific and vector-specific primers. Product is confirmed free of bacteria, fungi and common *Mycoplasma* contamination.

### Viral titer

The transduction unit (TU or IFU) of the lentiviral particles was estimated using the formula- 1TU=100 copies of viral genomic RNA. The physical copy numbers of the viral genomic RNA was determined using qRT-PCR. The customer should test the transduction at MOI=0.3, 1, 3, 5, 10 for their specific cell lines in order to get the best transduction efficiency.

### Overview of production

The OmicsLink™ ORF lentiviral expression plasmid (GeneCopoeia Cat. No. EX-NEG-LV151) was constructed using GeneCopoeia proprietary RecJoin™ technology. This plasmid was co-transfected into 293Ta cells (GeneCopoeia Cat. No. CLv-PK-01) with the Lenti-Pac HIV Packaging Mix (GeneCopoeia Cat. No. HPK-LvTR-20). Lentivirus-containing supernatants were harvested 48 hours after transfection, the particles were purified and stored at –80°C in aliquots.

### User manual

Please contact GeneCopoeia for a copy or download at:

[http://genecopoeia.com/product/lentiviral/pdf/packageing\\_kit\\_manual.pdf](http://genecopoeia.com/product/lentiviral/pdf/packageing_kit_manual.pdf)

GeneCopoeia, Inc.  
9620 Medical Center Drive, #101  
Rockville, Maryland 20850  
Tel: 301-762-0888 Fax: 301-762-8333  
Email: [inquiry@genecopoeia.com](mailto:inquiry@genecopoeia.com)  
Web: [www.genecopoeia.com](http://www.genecopoeia.com)

GeneCopoeia Products are for Research Use Only  
Trademarks: Lentifect™, Lenti-Pac™, OmicsLink™, EndoFectin™, TiterBoost™ (GeneCopoeia Inc.)

Copyright © 2015 GeneCopoeia Inc.  
LPP CSHCTR 012815

[www.genecopoeia.com](http://www.genecopoeia.com)