

**OCT4 Lentifect™ Purified Lentiviral Particles •**  
**Cat Nos. LPP-OCT4-Lv105-025-C, LPP-OCT4-Lv105-100-C**

Ready-to use purified lentiviral particles for the transduction of a variety of mammalian cells including difficult-to-transfect, primary, stem and non-dividing cells as well as in vivo use for transgenic animals.

**Description**

GeneCopoeia's Lentifect™ Purified Lentiviral Particles are produced from a standardized protocol using purified plasmid DNA and the proprietary reagents, EndoFectin™ Lenti (for transfection) and TiterBoost™ solution. The protocol uses a third generation self-inactivating packaging system meeting BioSafety Level 2 requirements.

The lentiviral particles include a CMV promoter for efficient expression of non-tagged, native OCT4 (variant 1) in target cells and use a puromycin 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 OCT4 lentiviral particles with titers of  $1 \times 10^7$  TU/ml.

Lentiviral particles are shipped on dry ice and must be stored at  $-80^{\circ}\text{C}$  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-  $1\text{TU}=100$  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  $\text{MOI}=0.3, 1, 3, 5, 10$  for their specific cell lines in order to get the best transduction efficiency.

**Overview of production**

The OCT4 OmicsLink™ ORF lentiviral expression plasmid (GeneCopoeia Cat. No. EX-Z0092-Lv105) was constructed using GeneCopoeia proprietary RecJoin™ technology. This plasmid was co-transfected into 293Ta cells (GeneCopoeia Cat. No. LT008) with the Lenti-Pac™ HIV packaging mix (GeneCopoeia Cat. No. LT001). Lentivirus containing supernatants were harvested 48 hours after transfection. The virus was further purified, concentrated and stored at  $-80^{\circ}\text{C}$  in aliquots.

**User manual**

Please contact GeneCopoeia for a copy or download at:  
[http://genecopoeia.com/product/lentiviral/pdf/packaging\\_kit\\_manual.pdf](http://genecopoeia.com/product/lentiviral/pdf/packaging_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)

Products are for Research Use Only  
Trademarks: Lentifect™, Lenti-Pac™, OmicsLink™, EndoFectin™, TiterBoost™ (GeneCopoeia Inc.)  
Copyright © 2019 GeneCopoeia Inc. LLPPOC-09102019