



## eGFP Lentifact™ Purified Lentiviral Particles • Cat No. LPP-EGFP-LV105-025

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 Lentifact™ 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 Lentifact particles include a CMV promoter for efficient expression of non-tagged eGFP 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 of purified eGFP lentiviral particles with titers of  $\sim 1 \times 10^{11}$  copies/ml.

Lentifact 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

Lentivirus products were titrated using qRT-PCR, which determines the copy number of viral genomic RNA.

### Overview of production

The eGFP OmicsLink™ ORF lentiviral expression plasmid (GeneCopoeia Cat. No. EX-EGFP-Lv105) 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 virus was further purified, concentrated and stored at –80°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)

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

Copyright © 2011 GeneCopoeia Inc.  
LLPPEG-061311