

# **A-Tubulin RFP fusion Lentifect<sup>™</sup> Lentiviral Particles** Cat No. LP506, LP507

Ready-to-use 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 Lentifect<sup>™</sup> Lentiviral Particles are produced from a standardized protocol using purified plasmid DNA and the proprietary reagents, EndoFectin<sup>™</sup> Lenti (for transfection) and TiterBoost <sup>™</sup> solution. The protocol uses a third generation self-inactivating packaging system meeting BioSafety Level 2 requirements.

The Lentifect particles include a CMV promoter for efficient expression of RFP-tagged, A-Tubulin protein 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 or purified A-Tubulin lentiviral particles with titers of ~1 x 10<sup>7</sup> 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 A-Tubulin OmicsLink<sup>™</sup> ORF lentiviral expression plasmid (GeneCopoeia Cat. No. EX-CS-RFP-A4294-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 and stored at -80°C.

#### **User manual**

Please contact GeneCopoeia for a copy or download at: 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

Web: www.genecopoeia.com