

**Firefly Luciferase + IRES-mCherry Lentifect™ Purified Lentiviral Particles •
Cat Nos. LPP-HLUC-Lv214-025-C, LPP-HLUC-Lv214-100-C**

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

- Produced with a standardized protocol using highly purified plasmids and EndoFectin-Lenti™ transfection and TiterBoost™ reagents. The protocol uses a third generation self-inactivating packaging system meeting BioSafety Level 2 requirements.
- CMV promoter for the expression of HLUC (humanized firefly luciferase) in target cells
- mCherry bicistronically co-expressed with FLUC
- Drug selection marker: none

Contents and storage

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

Lentiviral 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 Firefly Luciferase + IRES-mCherry OmicsLink™ ORF lentiviral expression plasmid (GeneCopoeia Cat. No. EX-HLUC-Lv214) was constructed using GeneCopoeia proprietary RecJoin™ technology. This plasmid was cotransfected 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 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

GeneCopoeia Products are for Research Use Only
Trademarks: Lentifect™, Lenti-Pac™, OmicsLink™, EndoFectin™, TiterBoost™ (GeneCopoeia Inc.)
Copyright © 2019 GeneCopoeia Inc. LPPHLUCM082219