



Renilla Luciferase + IRES-mCherry Lentifact™ Purified Lentiviral Particles • Cat No. LP-RLUC-LV177-025

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 RLUC in target cells
- mCherry bicistronically coexpressed with RLUC
- Drug selection marker: none

Contents and storage

Provided as 1 vial of 25 µl of purified lentiviral particles with titers of ~6x 10⁹ 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 Renilla Luciferase + IRES-mCherry OmicsLink™ ORF lentiviral expression plasmid (GeneCopeia Cat. No. EX-RLUC-LV177) was constructed using GeneCopeia proprietary RecJoin™ technology. This plasmid was co-transfected into 293Ta cells (GeneCopeia Cat. No. CLV-PK-01) with the Lenti-Pac HIV Packaging Mix (GeneCopeia Cat. No. HPK-LvTR-20). Lentivirus-containing supernatants were harvested 48 hours after transfection and stored at –80°C.

User manual

Please contact GeneCopeia for a copy or download at:
http://genecopeia.com/product/lentiviral/pdf/packaging_kit_manual.pdf

GeneCopeia, Inc.
9620 Medical Center Drive, #101
Rockville, Maryland 20850
Tel: 301-762-0888 Fax: 301-762-8333
Email: inquiry@genecopeia.com
Web: www.genecopeia.com

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

Copyright © 2011 GeneCopeia Inc.
LPPRLUCM061311