

CRE Recombinase Lentifect™ Purified Lentiviral Particles •

Cat Nos. LPP-CRE-Lv201-025-C, LPP-CRE-Lv201-100-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 CRE recombinase (NC_005856.1)
- SV40 promoter for the expression of GFP
- Puromycin resistance marker (Pac gene) for selection of stably transduced cells
- Pac gene bicistronically coexpressed with GFP

Contents and storage

Provided as 1 vial of 25 µl or 4 vials of 25 µl of purified lentiviral particles with titers of 1×10^8 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- $1\text{TU}=100$ 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 MOI=0.3, 1, 3, 5, 10 for their specific cell lines in order to get the best transduction efficiency.

Overview of production

The CRE Recombinase OmicsLink™ ORF lentiviral expression plasmid (GeneCopoeia Cat. No. EX-CRE-Lv201) 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 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. LPPCRE082219