

EF1α-SV40T (Puro) Lentifect™ Purified Lentiviral Particles

Cat. No. LP722-025, LP722-100

(Old Cat. No. LPP-SV40T-Lv195-025-C, LPP-SV40T-Lv195-100-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.

Description

Gene: SV40 small T and large T antigen Promoter: EF1α Tag: N/A Reporter: N/A Resistance marker: Puromycin Additional note: N/A

GeneCopoeia Lentifect[™] 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.

Contents and storage

Provided as 1 vial of 25 μ l or 4 vials of 25 μ l of purified SV40 small T and large T antigen lentiviral particles with titers of ~1 x 10⁸ 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=30 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 SV40 Small T and Large T Antigen OmicsLink[™] ORF lentiviral expression plasmid (GeneCopoeia Cat. No. EX-SV40T-Lv195). 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: <u>https://www.genecopoeia.com/wp-content/uploads/2018/03/Lentivirus-protocol-GeneCopoeia.pdf</u>

GeneCopoeia, Inc. 9620 Medical Center Drive, Suite 101 Rockville, Maryland 20850 Tel: 301-762-0888 Fax: 301-762-8333 Email: <u>inquiry@genecopoeia.com</u> Web: <u>www.genecopoeia.com</u>

Copyright © 2023 GeneCopoeia Inc.