

Halo-CD63 Lentifect[™] Purified Lentiviral Particles • Catalog #: LP777-025, LP777-100

Lentiviral particles pseudotyped with VSV-G protein for transferring genes to a variety of mammalian cells including primary, stem and non-dividing cells.

Description

Gene: Halo-CD63Promoter: EF1aTag: N-terminal HaloReporter: N/A

• Resistance marker: Puromycin resistance gene

· Additional note: N/A

Contents and storage

Provided as 1 vial of 25 µl or 4 vials of 25 µl of purified CD63 lentiviral particles with titers of ~1 x 108 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. Divide into useful aliquots if necessary and store at -80°C.

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

Identity of lentiviral transfer vector: pEZ-Lv195. The lentiviral particles were generated by following a standardized protocol using highly purified plasmids and EndoFectin-Lenti[™] and TiterBoost[™] reagents. The lentiviral transfer vector was co-transfected into 293Ta cells (Cat #: LT008) with Lenti-Pac[™] HIV packaging mix (Cat #: LT001). The lentiviral particles were purified and stored at -80°C in aliquots (purified particles).

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.

User manual

Please contact GeneCopoeia for a copy or download at: https://www.genecopoeia.com/wp-content/uploads/2018/03/Lentivirus-protocol-GeneCopoeia.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: EndoFectin™, Lenti-Pac™, TiterBoost™ (GeneCopoeia Inc.)