

Lentiviral particles for HUWE

Catalog #: LPP-HUWE-LvSG03-A00

Lot #: GC08132K1802

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: HUWE
- Promoter: U6
- Tag: N/A
- Reporter: mcherry
- Resistance marker: Puromycin resistance gene
- Additional note: N/A

Contents and storage

Provided as 50 µl×15 of purified lentiviral particles

Titer: 2.53×10^8 TU/ml

Lentivirus 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=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

Identity of lentiviral transfer vector: pCRISPR-LvSG03

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 lentivirus 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:

http://www.genecopoeia.com/product/lentiviral/pdf/packaging_kit_manual.pdf (HIV)

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.)

Copyright© 2018 GeneCopoeia Inc.