

Function Validated Lentiviral sgRNA for human PD-1 Gene Knock Out

Catalog #: LPP-HCP001055-LvSG03-15-10-h-50 Lot #: GC07262K1711

Ready-to-use purified lentiviral particles, used together with Cas9 nuclease lentiviral particles (See Cat No. followed) for effective human PD-1 gene knock out in a variety of mammalian cells including difficult-to-transfect, primary, stem and non-dividing cells as well as in vivo use for transgenic animals.

PD-1 gene knock out validated in HEK293 cells

Description

- . Gene: sgRNA (target in human PD-1 gene)
- . Promoter: U6
- . Tag: N/A
- . Reporter: mcherry
- . Resistance marker: Puromycin resistance gene
- . Additional note: sgRNA sequence available after order placed

Suggested Cas9 Nuclease lentiviral Particles from GeneCopia

The product can be combined with any of the following Ready-to-use Cas9 nuclease lentiviral particles for co-transduction or Cas9: sgRNA sequential transduction:

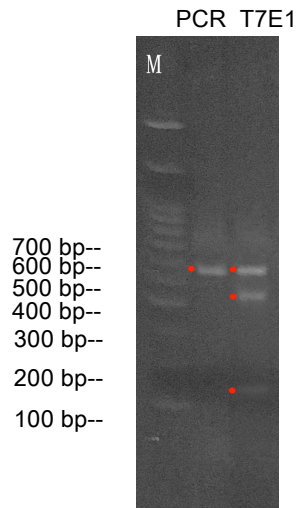
- . LPP-CP-LvC9NU-01-100-C
- . LPP-CP-LvC9NU-02-100-C
- . LPP-CP-LvC9NU-09-100-C
- . LPP-CP-LvC9NU-10-100-C

Product QC:

T7 Endonuclease I (T7 E1) Assay

PD-1 Site T7 E1 Assay

sgRNA targeting to PD-1 gene was transduced into HEK293/Cas9 Stable Cell Line by transduction. PD-1 gene was cut by CAS9 expressed inside the cells and repaired through NHEJ with mutation. A 577 bp PD-1 gene fragment from PCR was then tested by T7 Endonuclease I (T7 E1) Assay. The T7 E1 cleavage will result in two additional bands: one ~441 bp and the other ~136 bp.



Contents and storage

Provided as 50 μ l \times 1 of purified lentiviral particles

Titer: 1.9×10^8 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 .

GeneCopia, Inc.

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

GeneCopia Products are for Research Use Only
Trademarks: EndoFectinTM, Lenti-PacTM, TiterBoostTM Genome-CRISPTM (GeneCopia Inc.)
Copyright© 2017 GeneCopia Inc.

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 #: CLv-PK-01) with Lenti-Pac™ HIV packaging mix (Cat #: HPK-LvTR-20) . The lentivirus particles were purified and stored at -80°C in aliquots (purified particles).

User protocol

The large insertion size of cas9 expression cassette leads to the sharp titer decline of Cas9 lentiviral particles. To get effective desired genome editing, we strongly suggest antibiotic selection with indicate resistance gene to collect cells genomically integrated cas9 after transduction.

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™ Genome-CRISP™ (GeneCopoeia Inc.)
Copyright© 2017 GeneCopoeia Inc.