

## **CD63-mCherry-Flag Lentifact™ Purified Lentiviral Particles • Catalog #: LP781-025, LP781-100**

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: CD63-mCherry-Flag
- Promoter: CMV
- Tag: 3xFLAG
- Reporter: mCherry
- Resistance marker: Blastidicin 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 10<sup>8</sup> 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=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: pEZ-Lv197. 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](mailto:inquiry@genecopoeia.com) Web: [www.genecopoeia.com](http://www.genecopoeia.com)

GeneCopoeia Products are for Research Use Only.

Trademarks: EndoFectin™, Lenti-Pac™, TiterBoost™ (GeneCopoeia Inc.) Copyright© 2020 GeneCopoeia Inc.

LP-20200206