



## **CRE-excisable mouse OSKM Lentifect™ Purified Lentiviral Particles**

**Cat. No. LP810-025, LP810-100**

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: mouse Oct4-Sox2-Klf4-Myc

Promoter: EF1 $\alpha$

Tag: N/A

Reporter: N/A

Resistance marker: N/A

Additional note: N/A

LoxP sequences were inserted to the 5'LTR and 3'LTR region. The viral transgene insertion can be excised by CRE recombinase.

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  $\mu$ l or 4 vials of 25  $\mu$ l of purified mouse OSKM lentiviral particles with titers of  $\sim 1 \times 10^8$  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=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

The mouse OSKM OmicsLink™ ORF lentiviral expression plasmid (GeneCopoeia Cat. No. CS-Mm24188-Lv195-01). 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:

<https://www.genecopoeia.com/wp-content/uploads/2018/03/Lentivirus-protocol-GeneCopoeia.pdf>

GeneCopoeia, Inc.

9620 Medical Center Drive, Suite 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 Copyright © 2020 GeneCopoeia Inc. Trademarks: Lentifect™, Lenti-Pac™, OmicsLink™, EndoFectin™, TiterBoost™ (GeneCopoeia Inc.)

LP810-030320