

## Lentiviral Particles for Firefly Luciferase

### Catalog #: SP105-100

Lentiviral particles pseudotyped with Full Length SARS-CoV-2 spike glycoprotein (D614G & N439K) for transferring genes to a variety of mammalian cells expressing human ACE2 receptor on the cell surface.

#### Description

- **Gene:** Firefly Luciferase
- **Promoter:** CMV
- **Tag:** N/A
- **Reporter:** EGFP
- **Resistance marker:** Puromycin resistance gene
- **Additional note:** N/A

#### Contents and storage

Provided as 50 µl x 2 of purified lentiviral particles

**Titer 1:** > 1×10<sup>6</sup> TU/ml      **Titer 2:** > 1×10<sup>8</sup> RLU/ml

For lot-specific datasheet and titers, please visit <https://www.genecopia.com/tech/datasheet/>

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 titers of Transduction Unit/ml (TU/ml) and Relative Fluorescent Unit/ml (RLU/ml) of the spike protein-pseudotyped lentiviral particles were provided for each particle, based on the transduction efficiency using GeneCopia's ACE2/HEK293T stable cell line (Cat.# SL221).

The luciferase titer (RLU/ml) was based on the viral particle's firefly luciferase signal after transducing the ACE2/HEK293T cells, and was determined using GeneCopia's Luc-Pair<sup>TM</sup> Firefly Luciferase HS Assay Kit (Cat.# LF007). The eGFP transduction titer (TU/ml) was determined by the FACS sorting of the eGFP expressing ACE2/HEK293T cells after pseudotyped viral transduction.

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-Lv201

The lentiviral particles were generated by following a standardized protocol using highly purified plasmids and EndoFectin-Lenti<sup>TM</sup> and TiterBoost<sup>TM</sup> reagents.

The lentiviral transfer vector was co-transfected into 293Ta cells (Cat.#: LT008) with Lenti-Pac SARS-CoV-2 Full Length Spike protein (D614G & N439K) packaging mix (Cat.#: LT031). 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 GeneCopia for a copy or download at:

[Lenti-Pac<sup>TM</sup> SARS-CoV-2 Full Length Spike protein-pseudotyped \(D614G & N439K\) Lentivirus Packaging Kit](#)

GeneCopia, Inc.  
9620 Medical Center Drive, #101, Rockville, Maryland 20850  
Tel: 301-762-0888 Fax: 301-762-8333  
Email: [inquiry@genecopia.com](mailto:inquiry@genecopia.com) Web: [www.genecopia.com](http://www.genecopia.com)