



Infect All Cells

LentiFect™ Lentiviral Particles

For *in vitro* and *in vivo* applications

Features

- Titers ranging from 10^8 to 10^9 TU/ml
- Purified and ultra-purified for *in vitro* and *in vivo* studies, respectively
- Largest collection of sequence-verified ORF cDNA, miRNA, shRNA, promoter, and sgRNA lentiviral clones available for packaging
- Collection of 100+ premade lentiviral particles available for *Next-day*™ shipment, including GFP, mCherry, YFP, CFP and luciferase

Purified particles for *in vitro* (cell culture) studies

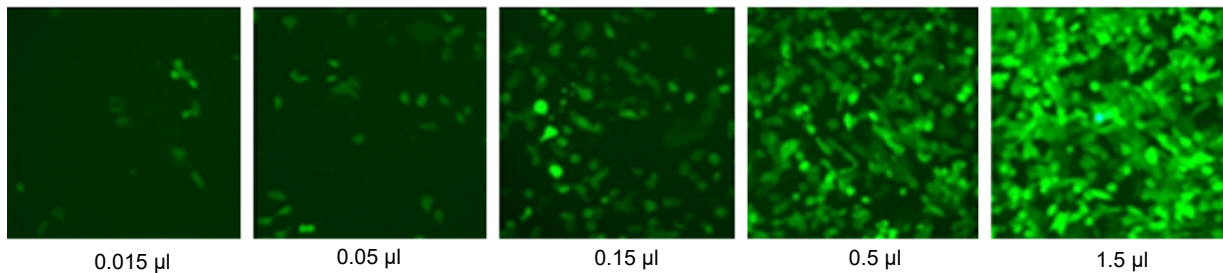


Figure 1. H1299 cells in 24-well plates were transduced with the indicated amounts of eGFP purified particles, (catalog number LPP-EGFP-LV105) in the presence of 5 µg/ml of polybrene. The expression of eGFP was visualized with a fluorescence microscope 72 hours post-transduction.

Ultra-purified particles for *in vivo* (animal) studies

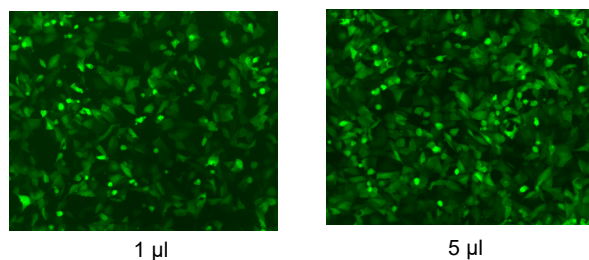


Figure 2. H1299 cells in 24-well plates were transduced with the indicated amounts of eGFP ultra purified particles. The expression of eGFP was checked with a fluorescence microscope 48 hours post-transduction. Exposure time was 0.2 s.

LentiFect™ Lentiviral Particles

High quality • Quick delivery • Low price

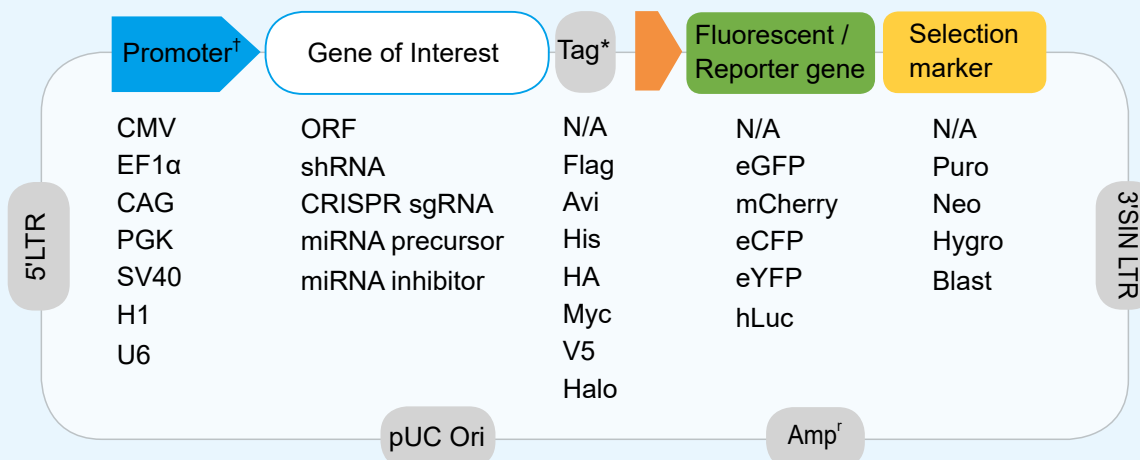
Lentiviral Particles for *in vivo* and *in vitro* Studies

Custom lentiviral particles Titer of 10 ⁸ TU/mL, purified and concentrated	Features
Any ORF, shRNA, sgRNA, precursor miRNA, miRNA inhibitor, or promoter reporter	Third generation packaging system Titer determined by qPCR
Premade particles Titer over 10 ⁸ TU/mL, purified and concentrated	Features
Positive and negative controls	Fluorescent tags: eGFP, eYFP, eCFP, mCherry, PLUMLuciferase reporters: Firefly, Renilla, Gaussia
iPSC ORF clones	OSKM transcription factors: Oct4, Sox2, Klf4, c-Myc, Nanog and Lin28
Cell immortalization	SV40 large T antigen, hTERT, c-Myc, CDK4, HOX, Bmi1, and more
Utility ORF clones	CRE recombinase, FLP recombinase, HIV tat, Cas9 nuclease*

*Cas9 nuclease titer ranges from 10⁶ to 10⁸ TU/mL

Premade particles: Next-day shipment and clone price included!

GeneCopoeia lentiviral clone features



*Not all tags, fluorescent/reporter genes and selection markers might be available for your product/vector of choice. Please refer to our website for detailed information.

†You can also choose the promoter of a specific gene to study its transcription and regulation. Please refer to "promoter reporter" product on our website for more details.

Contact us

GeneCopoeia, Inc.

9620 Medical Center Dr. Suite 101,
Rockville, MD 20850 USA



Email inquiry@genecopoeia.com
Tel +1 (301) 762-0888
Fax +1 (301) 762-3888
Website www.genecopoeia.com

LPP-030722