



Lentifect™ Pre-made Lentiviral Particles

GeneCopoeia offers Lentifect™ pre-made purified and ultra-purified lentivirus for *in vitro* and *in vivo* research, includes reporter gene lentivirus, cell immortalization lentivirus, Cas9 nuclease lentivirus, iPSc reprogramming lentivirus, exosome labeling lentivirus and other commonly used lentivirus.

Advantages

- Low cost and ideal for cell-line specific MOI testing or as transduction controls.
- Hundreds of pre-made viral particles – ship next day.
- Including – fluorescent proteins, luciferase reporters, empty vectors controls, cell immortalization genes and more.

Fluorescent proteins, luciferase reporters lentiviral particles

Promoter	Selection	Available genes
CMV	Puromycin / Hygromycin / Neomycin / Blasticidin	eGFP
CMV	Puromycin	eYFP
CMV	Puromycin	eCFP
CMV	Puromycin	mCherry
CMV	Puromycin	PLUM
CMV	Puromycin	Firefly luciferase
CMV	Puromycin	Renilla luciferase
CMV	Puromycin	Gaussia luciferase
CMV	Puromycin / N/A	Firefly luciferase + eGFP/mCherry
CMV	Puromycin / N/A	Renilla luciferase + eGFP/mCherry

Visit our website, <https://www.genecopoeia.com/product/lentivirus-system/premade-lentiviral-particles/>, to see a complete list.

Transduction of H1299 cells with GeneCopoeia's Lentifect™ lentiviral particles expressing eGFP

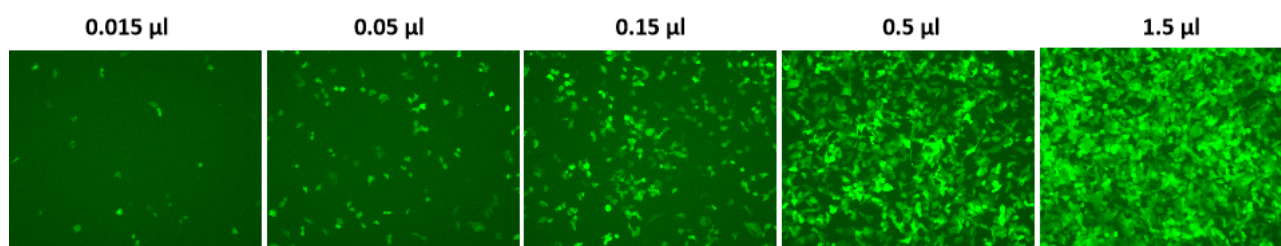


Figure 1. H1299 cells in 24-well plates were transduced with the indicated amounts of eGFP purified particles, (CAT#.LP303) in the presence of 5 µg/ml of polybrene. The expression of eGFP was visualized with a fluorescence microscope 72 hours post-transduction.

Lentifect™ Pre-made Lentiviral Particles

Cell immortalization lentiviral particles

Promoter	Selection	Available genes
CMV / EF1α	Puromycin	SV40 small T & large T antigen
CMV / EF1α	Puromycin / Hygromycin / N/A	SV40 large T antigen
CMV / EF1α	Puromycin / Hygromycin / N/A	hTERT
CMV	Puromycin / N/A	CDK4
CMV	Puromycin / N/A	HOXB8
CMV	Puromycin / N/A	HOXA9
CMV	Puromycin / N/A	HOXA10
CMV	Puromycin	c-Myc
CMV	Puromycin / N/A	Bmi1
CMV	Puromycin / N/A	HPV-16 E6-E7

Visit our website, <https://www.genecopoeia.com/product/lentivirus-system/cell-immortalization-reagents/>, to see a complete list.

Related products:

- Custom lentivirus including ORF cDNA, shRNA, sgRNA, miRNA and promoters
- Cas9 nuclease lentivirus and control sgRNA lentivirus
- iPSc reprogramming lentivirus
- Exosome labeling lentivirus
- SARS-CoV-2 spike-pseudotyped lentivirus

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

Lenti-101823