

Infect All Cells

LentiFect™ Lentiviral Particles

Pre-made particles and custom services available

Main Features

- Titer above 10^8 TU/mL for premade particles
- All lentiviral products purified and ready to transduce
- High transduction efficiency and long-term expression
- Third generation packaging system ensures safety
- Largest collection of ORF, miRNA, shRNA, promoter and sgRNA lentiviral clones to choose from
- Cost-effective, high quality and fast delivery
- All lentiviral clones are sequence verified

Completely customizable with modular components.

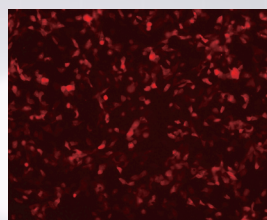
Type of clones	ORF	Precursor miRNA	miRNA inhibitor	shRNA	Promoter	sgRNA
Promoter	CMV, CMV5, EF1 α , CAG, PDK, SV40, Inducible	H1, CMV, EF1 α	H1, U6, CMV	H1, U6, Inducible	Various	U6
Reporter gene	eGFP, mCherry, Gluc, eCFP, eYFP	eGFP, mCherry		eGFP, mCherry, Gluc	mCherry	
Selection marker	Puromycin, Neomycin, Hygromycin, Zeomycin	Puromycin, Neomycin, Hygromycin		Puromycin		

LentiFect™ Lentiviral Particles

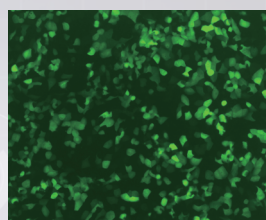
Quick delivery and low price

Custom lentiviral particles	Titer of 10^7 - 10^9 TU/mL, purified and concentrated
Any ORF, shRNA, sgRNA, precursor miRNA or miRNA inhibitors	Third generation packaging system Titer determined by qPCR
Premade particles	Titer over 10^8 TU/mL, purified and concentrated
Positive and negative controls	Fluorescent tags: eGFP, eYFP, eCFP, mCherry, PLUM Luciferase reporters: Firefly, Renilla, Gaussia
iPSC ORF clones	OSKM transcription factors: Oct4, Sox2, Klf4, c-Myc, Nanog and Lin28
Utility ORF clones	CRE recombinase, FLP recombinase, SV40 large T antigen, HIV tat, Cas9 nuclease

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

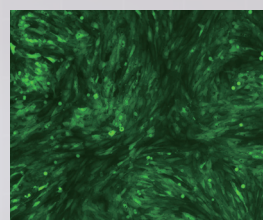


mCherry

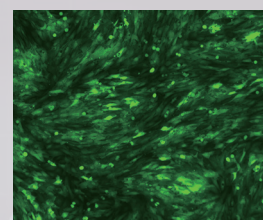


eGFP

Figure 1. H1299 cells were transduced in a 24-well plate with mCherry (left panel: LPP-MCHR-Lv105) or eGFP (right panel: LPP-EGFP-Lv105). Exposure time was ≤ 1 s with an MOI=1.

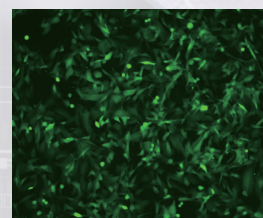


MOI=10

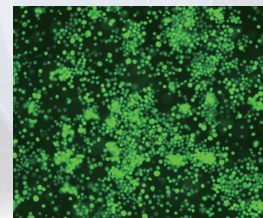


MOI=30

Figure 2. NIH/3T3 cells were transduced in a 24-well plate at two different multiplicities of infection (MOI) using LPP-NEG-Lv201 particles.



HOS



K562

Figure 3. Two cancer cell lines (Left: HOS, Right: K562) were transduced and puromycin selection was applied after 48 h. Inquire today for custom stable cell lines.

Contact us:

GeneCopoeia, Inc.

9620 Medical Center Drive, Suite 101
Rockville, MD 20850, USA

Email inquiry@genecopoeia.com
Tel +1 (301) 762-0888
Toll free +1 (866) 360-9531
Fax +1 (301) 762-3888
Website www.genecopoeia.com