

Lentifect™ SARS-CoV-2 Spike Protein-pseudotyped Lentiviral Particle Kit, Full Length (*in vitro* grade)

Catalog #: LT014

The SARS-CoV-2 spike protein-pseudotyped lentiviral particle kit contains one pseudotyped lentiviral particle, two controls and an ACE2/HEK293T stable cell line, which can be used to test the ability of the pseudovirus to infect the cells via the human ACE2 receptor on the cell surface.

Description

Lentifect™ SARS-CoV-2 Spike-Pseudotyped Lentiviral Particle Kit, **full length** (Cat No.LT014)

Reagent Cat No.	Contents	Viral titer
SP101-100	SARS-CoV-2 spike-pseudotyped lentivirus, full length (100 µl)	> 1×10 ⁵ TU/ml or > 1×10 ⁷ RLU/ml ¹
SP401-025	Lentivirus without envelope protein negative control (25 µl, 1 vial)	> 1 × 10 ¹¹ copies/ml ²
LP464-025	Standard VSV-G lentivirus (25 µl, 1 vial)	> 1×10 ⁸ TU/ml ³
SL221	ACE2-expressing HEK293T cells (2 x 10 ⁶ cells x 1 vial)	-

¹The functional titers of SARS-CoV-2 Spike pseudotyped lentivirus are determined by FACS sorting of GFP positive ACE2-expressing ACE2/HEK293T cells (GeneCopoeia Cat.No. SL221) and by luciferase analysis (GeneCopoeia Cat.No. LF007, LF008, LF009) of transduced ACE2/HEK293T.

²Titers of standard VSV-G pseudotyped lentivirus are determined by RT-qPCR to determine physical titer (copies/ml) and extrapolated to functional titer (TU/ml).

³The titer of the lentiviral particles without envelope protein is determined by RT-qPCR.

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.

Datasheet

Please contact GeneCopoeia for a copy or download at:

1) SARS-CoV-2 spike-pseudotyped lentivirus (SP101-100)

<https://www.genecopoeia.com/wp-content/uploads/2020/08/SP101-025-G107142K2009.pdf>

For the lot# specific datasheet, please download a copy at: <https://www.genecopoeia.com/tech/datasheet/>

2) Lentivirus without envelope protein negative control (SP401-025)

<https://www.genecopoeia.com/wp-content/uploads/2020/08/SP401.G.pdf>

3) Standard VSV-G lentivirus (LP464-025)

https://www.genecopoeia.com/wp-content/uploads/2017/01/hluc_lv201.pdf

4) ACE2-expressing HEK293T cells (SL221)

<https://www.genecopoeia.com/wp-content/uploads/2020/04/Datasheet-SL221-HEK293T-ACE2.pdf>

GeneCopoeia, Inc.

9620 Medical Center Drive, #101

Rockville, Maryland 20850

Tel: 301-762-0888 Fax: 301-762-8333

Email: inquiry@genecopoeia.com

Web: www.genecopoeia.com