

Data Sheet ***Spike (SARS-CoV-2) Lentivirus*** **Catalog #: 78010**

Product Description

Cell entry of SARS-CoV-2 depends on the binding of viral spike protein to cellular receptor ACE2. The SARS-CoV-2 Spike Lentivirus are replication incompetent, HIV-based, VSV-G pseudotyped lentiviral particles that are ready to be transduced into almost all types mammalian cells, including primary and non-dividing cells. The particles contain the full length SARS-CoV-2 spike gene (QHD43416.1) driven by an EF1A promoter (Figure 1).

Application

1. Transient expression of full length SARS-CoV-2 spike in target cells.
2. Generation of stable cell line expressing SARS-CoV-2 spike with puromycin selection.

Formulation

The lentiviruses were produced from HEK293T cells in medium containing 90% DMEM + 10% FBS.

Titer

The titer will vary with each lot; the exact value is provided with each shipment.

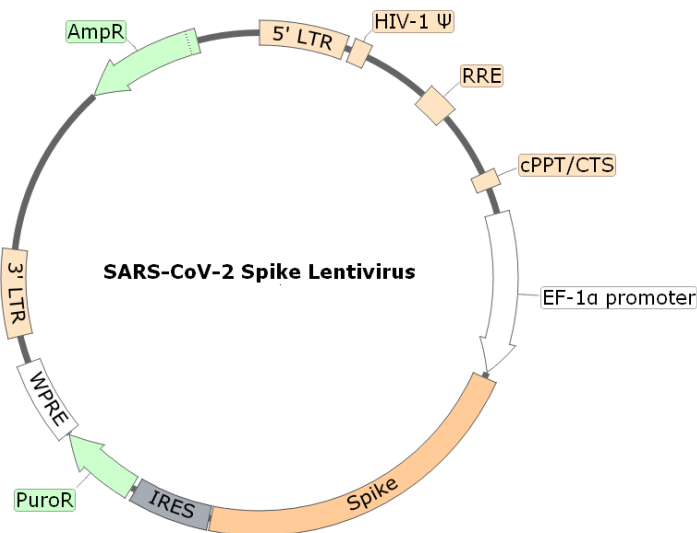


Figure 1. Schematic of the lenti-vector used to generate the SARS-CoV-2 spike lentivirus

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

Storage

Lentiviruses are shipped with dry ice. For long term storage, it is recommended to store the virus at -80°C . Avoid repeated freeze-thaw cycles. Titers can drop significantly with each freeze-thaw cycle.

Biosafety

None of the HIV genes (gag, pol, rev) will be expressed in the transduced cells, as they are expressed from packaging plasmids lacking the packing signal. Although the pseudotyped lentiviruses are replication-incompetent, they require the use of a Biosafety Level 2 facility. BPS recommends following all local federal, state, and institutional regulations and using all appropriate safety precautions.

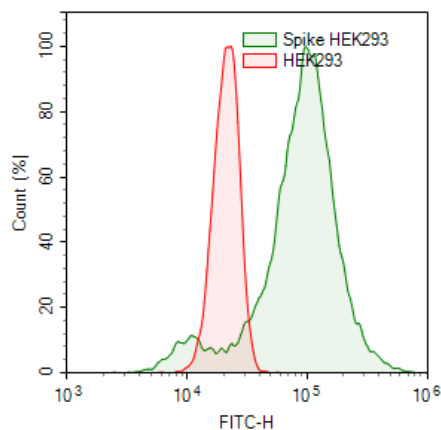


Figure 2. The expression of SARS-CoV-2 Spike protein in HEK293 cells transduced with SARS-CoV-2 spike lentivirus. Appropriate 500,000 cells/well (6-well culture plate) were transduced with 2,000,000 TU/well SARS-CoV-2 spike lentivirus in the presence of $5\ \mu\text{g/ml}$ of polybrene. After 52 hours of transduction, the transduced cells were selected in Growth Medium 1N (BPS Bioscience, #79801) which contains $0.5\ \mu\text{g/ml}$ puromycin. The spike expression in the puromycin-resistant cell pool was analyzed by intracellular flow cytometry staining using anti-Spike S1 primary antibody (BPS Bioscience, #100715-1) and FITC-conjugated goat anti-human IgG secondary antibody (ThermoFisher, #H10101C). Red, HEK293 parental cells; Green, HEK293 cells transduced with SARS-CoV-2 spike lentivirus.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence

MFVFLVLLPLVSSQCVNLTTRTQLPPAYTNSFTRGVVYYPDKVFRSSVLHSTQDLFLPFFSNVTW
FHAIHVSNGTNGTKRFDNPVLPFNDGVYFASTEKSNIRGWIFGTTLDSTQSLNATNVVIVK
CEFQFCNDPFLGVYHKNKSWMESEFRVYSSANNCTFEYVSQPFLMDLEGKQGNFKNLREF
VFKNIDGYFKIYSKHTPINLVRDLPQGFSALEPLVDLPIGINITRFQTLALHRSYLTGDSSSGW
TAGAAAYYVGYLQPRTFLLKYNENGTITDAVDCALDPLSETKCTLKSFTVEKGIYQTSNFRVQP
TESIVRFPNITNLCPFGEVFNATRFASVYAWNRKRISNCVADYSVLYNSASFSTFKCYGVSPTKL
NDLCFTNVYADSFVIRGDEVQRQIAPGQTGKIADYNYKLPDDFTGCVIAWNSNLDKSVGGNYN
YLYRFRKSNLKPFRDISTEIQAGSTPCNGVEGFNCYFPLQSYGFQPTNGVGYQPYRVVVL
SFELLHAPATVCGPKKSTNLVKNKCVNFNGLTGTGVLTESNKKFLPFQFGRDIADTTDAVR
DPQTLEILDITPCSFGGVSVITPGTNTSNQVAVLYQDVNCTEVPVAIHADQLTPTWRVYSTGSN
VFQTRAGCLIGAEHVNSYECDIPIGAGICASYQTQTNPRRARSVASQSIIAYTMSLGAENSV
YSNSIAIPTNFTISVTTEILPVSMKTSVDCTMYICGDSTECSNLLQYGSFCTQLNRALTGI
EQDKNTQEVFAQVKQIYKTPPIKDFGGFNFSQILPDPSKPSKRSFIEDLLFNKVTLADAGFIKQY
GDCLGDIARDLICAQKFNGLTVLPPLTDEMIQYTSALLAGTITSGWTFGAGAALQIPFAMQM
AYRFNGIGVTQNVLYENQKLIANQFNVAIGKIQDLSSTASALGKLQDVVNQNAQALNTLVKQL
SSNFGAISSVLNDILSRDKVEAEVQIDRLITGRLQSLQTYVTQQLIRAAEIRASANLAATKMSEC
VLGQSKRVDFCGKGYHLSFPQSAPHGVVFLHVTVYVPAQEKNTTAPAICHDKAHFPREGV
FVSNGTHWFVTQRNFYEPQIITDNTFVSGNCDVVIGIVNNTVYDPLQPELDSFKEELDKYFKN
HTSPDVDLGDIGINASVNIQKEIDRLNEVAKNLNESLIDLQELGKYEYIKWPWYIWLGFIAGL
IAIVMVTIMLCCMTSCCCLKGCCSCGSCCKFDEDDSEPVKGVKLHYT

Related Products

Product	Cat. #	Size
Spike Pseudotyped Lentivirus (Luciferase Reporter)	79942	500 µl x2
Bald Lentiviral Pseudovirion (Luciferase Reporter)	79943	500 µl x2
Spike Pseudotyped Lentivirus (eGFP Reporter)	79981	500 µl x2
Bald Lentiviral Pseudovirion (eGFP Reporter)	79987	500 µl x2
Spike Pseudotyped Lentivirus (Luciferase-eGFP Dual Reporter)	79982	500 µl x2
Bald Lentiviral Pseudovirion (Luciferase-eGFP Dual Reporter)	79988	500 µl x2
eGFP Lentivirus	79979	500 µl x2
Firefly Luciferase-eGFP Lentivirus	79980	500 µl x2
Negative Control Lentivirus	79578	500 µl x2
Renilla Luciferase (Rluc) Lentivirus	79565	500 µl x2
Firefly Luciferase (Fluc) Lentivirus	79692	500 µl x2
ACE2 Lentivirus	79944	500 µl x2

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com