

Spike Trimer (S1+S2), His-Tag (SARS-CoV-2) Recombinant

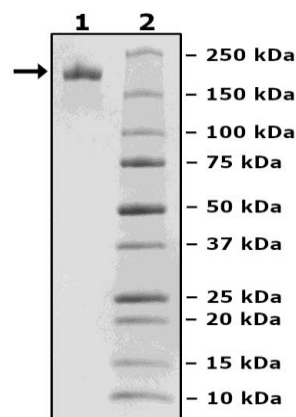
Catalog: 100728
Lot: 230228-2

Product Information

Description:	Recombinant SARS-CoV-2 Spike protein in its homotrimeric form, containing S1+S2 subunits and encompassing amino acids 16-1213. This protein corresponds to the wild-type (strain of reference). The construct contains mutations 682RRAR685>A, K986P and V987P, and a T4 trimerization domain followed by a His-tag (6xHis) in C-terminal. The recombinant protein was affinity purified.
Species	SARS-CoV-2
Construct:	Spike Trimer (S1+S2) (16-1213-His) (SARS-CoV-2)
Concentration:	0.16 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	137 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	MN908947
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Assay Conditions:	Assay was done according to SARS-CoV-2 Spike Trimer (S1+S2):ACE2 Inhibitor Screening Colorimetric Assay Kit (BPS Bioscience #79999) with ACE2-Biotin titrated at various concentrations.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Spike Trimer (S1+S2): ACE2-Biotin Activity

