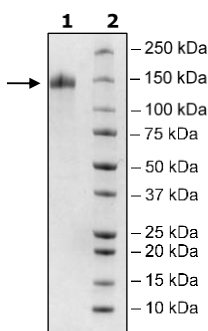


Product Information

Construct:	Spike S1 (16-685-Fc(IgG1)-Avi) (SARS-CoV-2)
Concentration:	0.90 mg/ml
Species:	SARS-CoV-2
Formulated In:	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	MN908947
MW:	104 kDa + glycans
Glycosylation:	This protein runs at a higher molecular weight due to glycosylation.
Purity:	≥90%
Assay Conditions:	SARS-CoV-2 S1 (BPS Bioscience #100719) and ACE2 binding assay was done using BPS ACE2: SARS-CoV-2 Spike Inhibitor Screening Assay Kit (BPS Bioscience #79936). SARS-CoV-2 S1 (BPS Bioscience #100719) protein and antihuman Fc-HRP were used instead of SARS-CoV-2 Spike protein (RBD) and anti-mouse Fc-HRP respectively.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



ACE2: SARS-CoV-2 Inhibition

