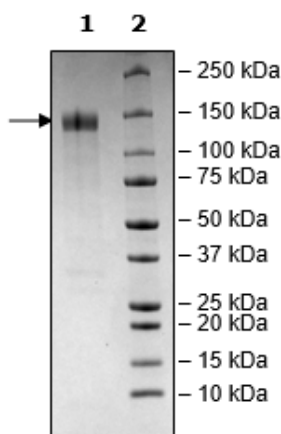


Product Information

Construct:	Spike S1 ((13-665)-Fc(IgG1)-Avi) (SARS-CoV-2)
Concentration:	0.88 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:	102 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	≥90%
Assay Conditions:	Assay was done according to the Spike S1 (SARS-CoV-2): ACE2 Inhibitor Screening Colorimetric Assay Kit (BPS Bioscience #79954). 100 ng of Spike S1 #100678 was coated overnight. Various amounts of ACE2 [Biotinylated] (BPS Bioscience #100665) added and incubated for 1 hour, followed by 1 hour incubation with streptavidin-HRP, and addition of colorimetric HRP substrate and 1 N HCl. Absorbance measure at 450 nm.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Spike S1 Fc:ACE2 [Biotinylated] Binding

