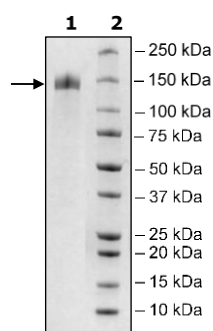


Product Information

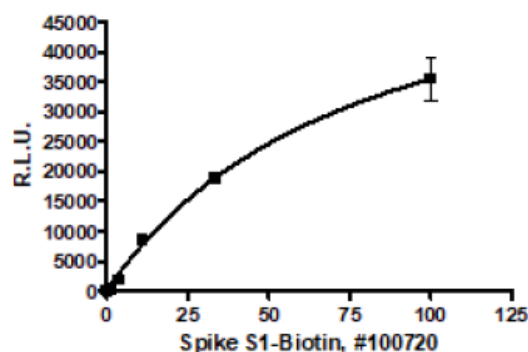
Construct:	Spike S1 (16-685-Fc (IgG1)-Avi)-(Biotin) (SARS-CoV-2)
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%.
Concentration:	1.02 mg/ml
Species:	SARS-CoV-2
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
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 MW by SDS-PAGE due to glycosylation.
Purity:	≥90%
Assay Conditions:	SARS-CoV-2 S1 (BPS Bioscience #100720) 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 #100720) protein and Streptavidin-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

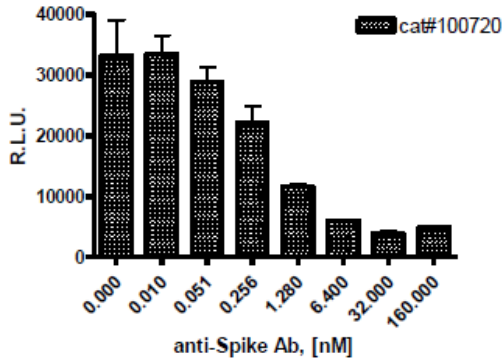
4-20% SDS-Page Coomassie Staining



ACE2: SARS-CoV-2 Inhibition



Biotin-Avidin Pulldown



SARS-CoV-2 S1(BPS Bioscience #100720, 20 nM) was incubated with various concentrations of the anti-Spike Ab for 1 hour at room temperature. After then, S1/Ab mixture was added to the plate coated with ACE2 as described in the assay kit (BPS Bioscience #79936). Streptavidin-HRP were used to measure SARS-CoV-2-S1/ACE2 binding.

