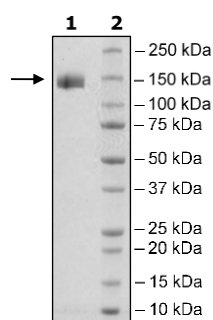


Product Information

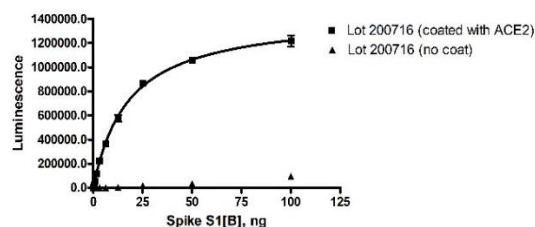
Construct:	Spike S1 (13-665-Fc(IgG1)-Avi)-(Biotin) (SARS-CoV-2)
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%.
Concentration:	0.65 mg/ml
Species:	Human
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
Expression 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 molecular weight due to glycosylation.
Purity:	≥90%
Assay Conditions:	Reaction done according to ACE2:SARS-CoV-2 Spike S1 Inhibitor Screening Assay Kit (BPS Bioscience #79945) protocol. Buffer used is PBS + 0.1% BSA. Blocking buffer is Superblock PBS + 0.05% Tween-20. Coat protein (ACE2, His-tag) was added overnight at 4° 50 ng/well. Various amounts of biotinylated ligand (Spike S1[Biotin-labeled]) (0-400 ng) added for 1h at room temperature. Strep-HRP added for 1h at room temperature. Luminescence developed with ELISA ECL substrates.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



ACE2-His: Spike S1 (SARS-CoV-2) Biotinylated Binding



Biotin-Avidin Pulldown

