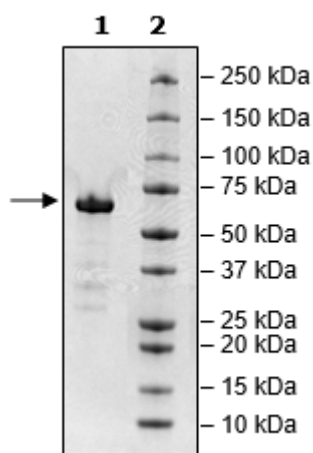


Product Information

Construct:	Spike S1 RBD (319-541-Fc (IgG1)-Avi)-(Biotin) (SARS-CoV-2)
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%.
Concentration:	0.61 mg/ml
Species:	SARS-CoV-2
Formulated In:	8 mM Phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 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:	MN_908947
MW:	54 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining



Biotin-Avidin Pulldown

