

Product Information

Construct:	Spike S1 RBD (319-541-Fc(IgG1)) (SARS-CoV-2)
Concentration:	0.20 mg/ml
Species:	SARS-CoV-2
Formulated In:	50 mM Tris-HCl, pH 7.5, 90 mM Glycine, 01 mM PMSF
Expression System:	CHO
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	MN908947
MW:	50 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

