

Product Information

Construct:	Spike S1 RBD (V483A) (319-541-Avi-His) (SARS-CoV-2)
Concentration:	0.83 mg/ml
Species:	SARS-CoV-2
Formulated In:	8 mM phosphate, 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:	28 kDa +glycans
Purity:	>90%
Notes:	This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

Quality Control Data

4-20% SDS-Page Coomassie Staining

