

Product Information

Construct:	Spike S1 (B.1.351) (16-685-Avi-His) (SARS-CoV-2)
Mutations:	L18F, D80A, D215G, R246I, K417N, E484K, N501Y, D614G.
Concentration:	0.91 mg/ml
Species:	SARS-CoV-2
Formulated In:	8mM phosphate pH 7.4, 110mM NaCl, 2.2mM 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:	MN908947
MW:	78 kDa + glycans
Glycosylation:	This protein runs at higher MW by SDS-PAGE due to glycosylation.
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

