

Product Information

Construct:	Spike S1 (P.1 Variant) (16-685-Avi-His) (SARS-CoV-2)
Mutations:	L18F, T20N, P26S, D138Y, R190S, K417T, E484K, N501Y, D614G, H655Y
Concentration:	0.28 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:	MN908947
MW:	78 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	89%

Quality Control Data

4-20% SDS-Page Coomassie Staining

