

Product Information

Construct:	Spike Trimer (B.1.1.7 Variant) (SARS-CoV-2) (16-1213-His)
Mutations:	Deletions: del69–70 HV, del144 Y Substitutions: N501Y, A570D, D614G, P681H, T716I, S982A, D1118H
Concentration:	0.59 mg/ml
Species:	SARS-CoV-2
Formulated In:	8 mM phosphate pH 7.4, 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:	137 kDa
Glycosylation	This protein runs at higher M.W. by SDS-PAGE due to glycosylation.
Purity:	>90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

