

Product Information

Construct:	Spike Trimer (S1+S2) (B.1.617.1 Variant) (16-1213-His) (SARS-CoV-2)
Mutations:	G142D, E154K, L452R, E484Q, D614G, P681R, Q1071H
Concentration:	1.17 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 + 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

