

Product Information

Construct:	Spike Trimer (S1+S2) (B.1.429 Variant) (16-1213(end)-His) (SARS-CoV-2)
Mutations:	W152C, L452R, D614G
Concentration:	0.26 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:	139 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	85%

Quality Control Data

4-20% SDS-Page Coomassie Staining

