

Product Information

Construct:	Spike Trimer (K417T, E484K, N501Y)(16-1213-His) (SARS-CoV-2)
Mutation:	K417T, E484K, N501Y
Concentration:	0.29 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 higher M.W. by SDS-PAGE due to glycosylation.
Purity:	>90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

