

Product Information

Construct:	Spike S2 (686-1212-Fc) (SARS-CoV-2)
Concentration:	0.40 mg/ml
Species:	SARS-CoV-2
Formulated In:	50 mM Tris-HCl, pH 7.5, 90 mM Glycine
Expression System:	CHO
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	MN908947
MW:	130 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

