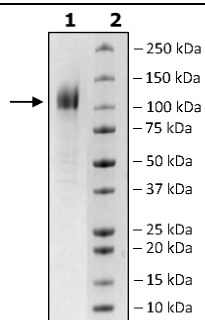


Product Information

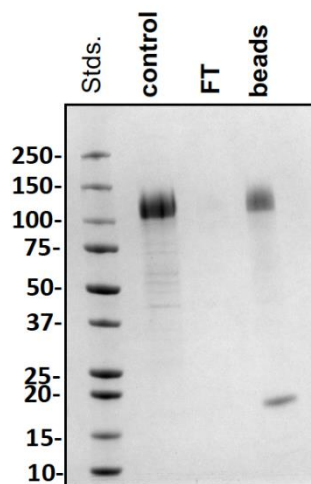
Construct:	Spike S1 protein (SARS-CoV-2) (16-685-Avi-His)-(Biotin)
Concentration:	0.21 mg/ml
Species:	SARS-CoV-2
Formulated In:	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2mM KCl, and 20% glycerol.
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	MN908947
MW:	78 kDa +glycans
Purity:	85%
Notes:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Biotin-Avidin Pulldown



This protein is enzymatically biotinylated using
Avi-Tag™ technology.
Biotinylation is confirmed to be $\geq 90\%$.