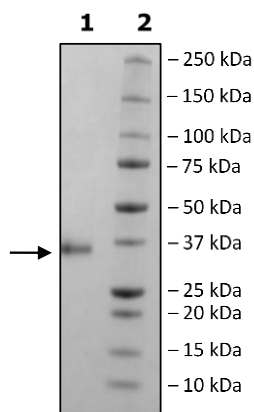


Product Information

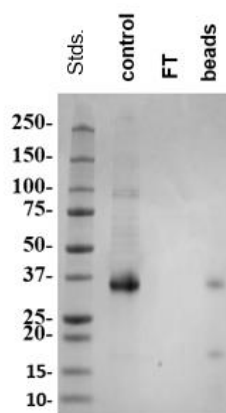
Construct:	Spike S1 RBD (B.1.617.2.1 Variant) (319-541-Avi-His)-(Biotin) (SARS-CoV-2)
Mutations:	K417N, L452R, T478K
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
Concentration:	0.75 mg/ml
Species:	SARS-CoV-2
Formulated In:	8 mM phosphate 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:	28 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	≥90%
Aggregation:	<10%

Quality Control Data

4-20% SDS-Page Coomassie Staining



Biotin-Avidin Pulldown



Gel Filtration Trace

