

Spike S1 RBD, Avi-His-Tag, Biotin-labeled (SARS-CoV-2) Recombinant

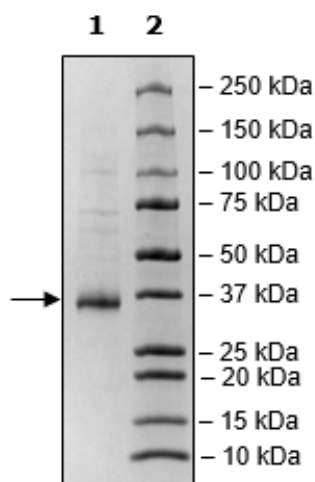
Catalog: 100697
Lot: 201105-1

Product Information

Description:	Recombinant SARS-CoV-2 Spike protein, RBD (Receptor Binding Domain) encompassing amino acids 319-541. The construct contains a C-terminal Avi-Tag™ followed by a His-tag (6xHis). The protein was enzymatically biotinylated using the Avi-Tag™ and affinity purified.
Species	Human
Construct:	Spike S1 RBD (319-541-Avi-His)-(Biotin)
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
Concentration:	2.41 mg/ml
Expression System:	HEK293
Purity:	86%
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	28 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	MN908947
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown

