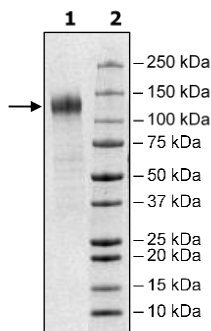


## Product Information

<b>Construct:</b>	Spike S1 (16-685-Avi-His)(SARS-CoV-2)
<b>Concentration:</b>	2.13 mg/ml
<b>Species:</b>	SARS-CoV-2
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
<b>Expression System:</b>	HEK293
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	MN908947
<b>MW:</b>	78 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher molecular weight due to glycosylation.
<b>Purity:</b>	≥90%
<b>Assay Conditions:</b>	SARS-CoV-2 Spike S1 protein (BPS Bioscience #100730) was coated on the plate for overnight at 4°C, and tested for the binding to ACE2 using ACE2-His, Avi-Biotin labeled (BPS Bioscience #100665) at various concentrations. Streptavidin-HRP was used to develop the luminescence signal.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

## 4-20% SDS-Page Coomassie Staining



## SARS-CoV-2 Spike S1:ACE2 Binding

