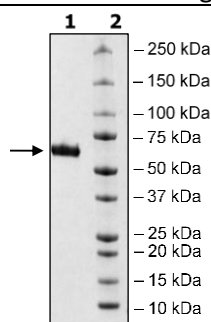


## Product Information

<b>Construct:</b>	Spike S1 RBD, Mouse Fc-Fusion (SARS-CoV-2) (319-541-mouse Fc)
<b>Concentration:</b>	0.7 mg/ml
<b>Species:</b>	SARS-CoV-2
<b>Formulated In:</b>	8 mM phosphate, 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>	QHD43416.1
<b>MW:</b>	50 kDa
<b>Purity:</b>	≥90%
<b>Assay Conditions:</b>	Assay was performed according to the ACE2: Spike RBD (SARS-CoV-2) Inhibitor Screening Assay Kit (BPS Bioscience, #79936) with SARS-CoV-2 Spike titrated from 80 ng- 0 ng.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.
<b>Notes:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.

## Quality Control Data

## 4-20% SDS-Page Coomassie Staining



## ACE2:SARS-CoV-2 Spike Interaction

