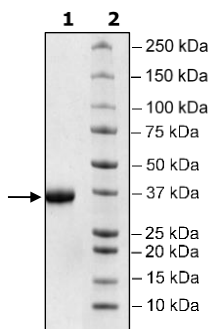


## Product Information

<b>Construct:</b>	Spike S1 RBD (V367F) (319-541-Avi-His) (SARS-CoV-2)
<b>Mutation:</b>	V367F
<b>Concentration:</b>	2.94 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>	28 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Purity:</b>	≥90%
<b>Aggregation:</b>	<10%

## Quality Control Data

### 4-20% SDS-Page Coomassie Staining



## Gel Filtration Trace

