

# Spike S1, Avi-His-Tag, iFluor-488-labeled (SARS-CoV-2) Recombinant

Catalog: 100936

Lot: 230109

## Product Information

<b>Description:</b>	Recombinant SARS-CoV-2 Spike protein, S1 subunit encompassing amino acids 16-685. It contains a C-terminal Avi-Tag™ followed by a His-tag (6xHis). The recombinant protein has been conjugated with iFluor-488™ and affinity purified.
<b>Species</b>	SARS-CoV-2
<b>Construct:</b>	Spike-S1 (16-685-Avi-His)-(iFluor-488) (SARS-CoV-2)
<b>Concentration:</b>	0.74 mg/ml
<b>Expression System:</b>	HEK293
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>MW:</b>	78 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Genbank Accession:</b>	MN908947
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	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

