

# Spike Trimer (S1+S2) (D614G), His-Tag (SARS-CoV-2) Recombinant

Catalog: 100810

Lot: 210325

## Product Information

<b>Description:</b>	Recombinant SARS-CoV-2 Spike protein in its homotrimeric form, containing S1+S2 subunits and encompassing amino acids 16-1213. This protein contains mutation of interest D614G. The construct also contains mutations 682RRAR685>A, K986P and V987P, and a T4 trimerization domain followed by a His-tag (6xHis) in C-terminal. The recombinant protein was affinity purified.
<b>Species</b>	SARS-CoV-2
<b>Construct:</b>	Spike Trimer (S1+S2) (D614G) (16-1213-His) (SARS-CoV-2)
<b>Mutation:</b>	D614G
<b>Concentration:</b>	0.94 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>	139 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

