
Recombinant SARS-CoV-2 Spike S2 ECD Protein with His-Tag

(Catalog # E80009)

Description

Recombinant SARS-CoV-2 Spike S2 ECD Protein is produced by the Baculovirus-Insect Cells expression system. The target protein is expressed with sequence (Ser686-Pro1213) of 2019-nCoV Spike S2 ECD (Accession #YP_009724390.1) fused with a His-tag at the C-terminus.

Purity

> 90% by SDS-PAGE.

Endotoxin

< 1.0 EU/ug of the protein by LAL method.

Formulation

Lyophilized from a 0.22 um filtered solution of sterile 20 mM PB, 300 mM NaCl, pH 7.0. or Supplied as a 0.22 um filtered solution in 20 mM PB, 300 mM NaCl, pH 7.0

Species

Human

Calculated MW

58.8kDa

Expressed Host

Baculovirus-Insect Cells

Tag

His tag at the C-terminus

Gene ID

43740568

Storage

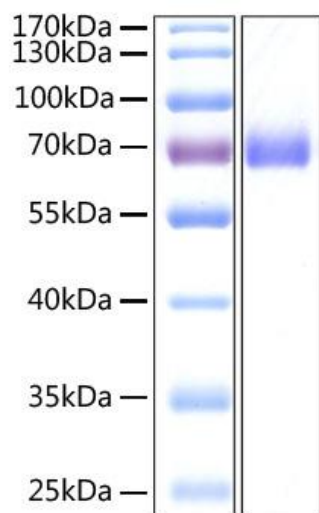
Store the lyophilized protein at -20°C to -80°C for long term. After reconstitution, the protein solution is stable at -20°C for 3 months, at 2-8°C for up to 1 week. or This product is stable at ≤ -70°C for up to 6 months from the date of receipt. For optimal storage, aliquot into smaller quantities after centrifugation and store at the recommended temperature.

Reconstitution

For lyophilized protein: Reconstitute to a concentration of 0.1-0.5 mg/mL in sterile distilled water.

Alternative Names

S2 protein; Spike glycoprotein Subunit2; S glycoprotein Subunit2; Spike protein S2; COVID-19; SARS-CoV-2; 2019-nCoV



Recombinant SARS-CoV-2 Spike S2 ECD Protein with His-Tag was determined by SDS-PAGE with Coomassie Blue, showing a band at 70 kDa.