
Recombinant 2019-nCoV Spike S1 Protein with His-Tag

(Catalog # E80007)

Description

Recombinant 2019-nCoV Spike S1 Protein is produced by HEK293 cells expression system. The target protein is expressed with sequence (Val11-Arg682) of 2019-ncov Spike S1 (Accession #YP_009724390.1) fused with a 6xHis tag at the C-terminus.

Bio-Activity

Measured by its binding ability in a functional ELISA. Immobilized Recombinant SARS-COV-2 Spike S1-His at 2ug/mL (100 uL/well) can bind Recombinant Human ACE2 with a linear range of 1.5-36.5 ng/mL. Immobilized Human ACE2 on COOH Chip, can bind SARS-COV-2 Spike S1 with an affinity constant of 27.3 nM as determined in a SPR assay (Nicoya OpenSPR).

Purity

> 90% by SDS-PAGE.

Endotoxin

< 0.1 EU/ug of the protein by LAL method.

Formulation

Lyophilized from a 0.22 um filtered solution of PBS, pH 7.4. or Supplied as a 0.22 um filtered solution in PBS, pH 7.4.

Species

SARS-COV-2

Calculated MW

76kDa

Expressed Host

HEK293 cells

Tag

6xHis 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.

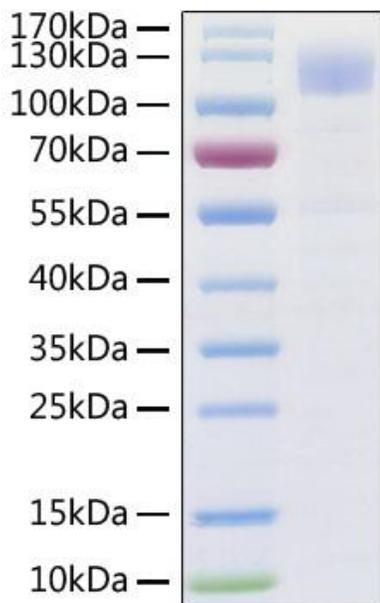
or This product is stable at ≤ -70°C for up to 6 months from the date of receipt. For optimal storage, aliquot into smaller quantities after centrifugation and store at recommended temperature.

Reconstitution

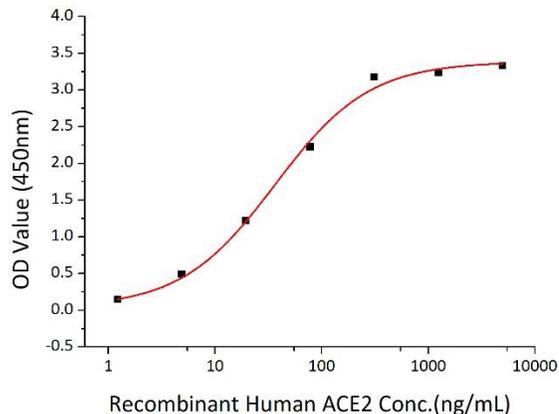
For lyophilized protein: Reconstitute to a concentration of 0.1-0.5 mg/mL in sterile distilled water.

Alternative Names

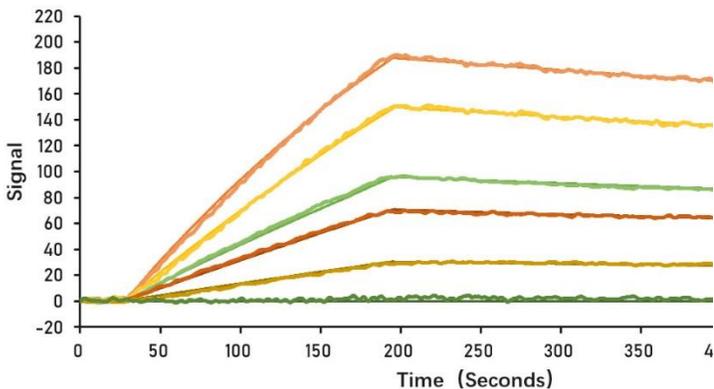
S1 protein; Spike glycoprotein Subunit1; S glycoprotein Subunit1; Spike protein S1; novel coronavirus s1 Protein



Recombinant 2019-nCoV Spike S1 Protein with His-Tag was determined by SDS-PAGE with Coomassie Blue, showing a band at 110-130 kDa.



Immobilized Recombinant 2019-nCoV Spike S1-His at 2µg/mL (100 µL/well) can bind Recombinant Human ACE2 with a linear range of 1.5-36.5 ng/mL.



Immobilized Human ACE2 on COOH Chip, can bind 2019-nCoV Spike S1 with an affinity constant of 27.3 nM as determined in a SPR assay (Nicoya OpenSPR).