
Recombinant SARS-CoV-2 S1+S2 ECD (S-ECD) Protein with His-Tag

(Catalog # E80002)

Background

The spike protein (S) of coronavirus (CoV) attaches the virus to its cellular receptor, angiotensin-converting enzyme 2 (ACE2). A defined receptor-binding domain (RBD) on S mediates this interaction. The S protein plays key parts in the induction of neutralizing-antibody and T-cell responses, as well as protective immunity.

Description

Recombinant SARS-CoV-2 S1+S2 ECD (S-ECD) Protein is produced by HEK293 cells expression system. The target protein is expressed with sequence (Val11-Gln1208) of 2019-ncov S1+S2 ECD (S-ECD) (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+S2 ECD-His at 2ug/mL (100 uL/well) can bind Recombinant Human ACE2 with a linear range of 0.15-9.92 ng/mL

Purity

>97% by SDS-PAGE.

Endotoxin

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

Formulation

Supplied as a 0.22 um filtered solution in PBS, pH 7.4.

Species

SARS-COV-2

Calculated MW

133.8kDa

Expressed Host

HEK293 cells

Tag

6xHis tag at the C-terminus

Gene ID

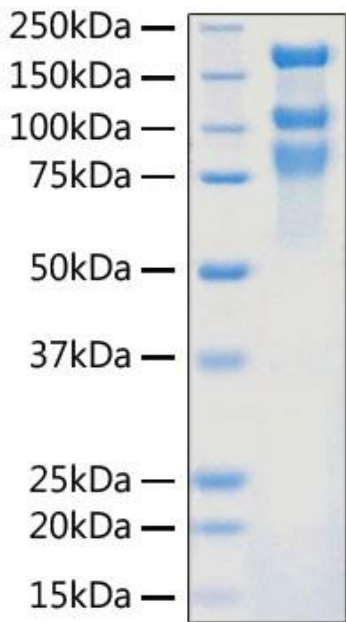
43740568

Storage

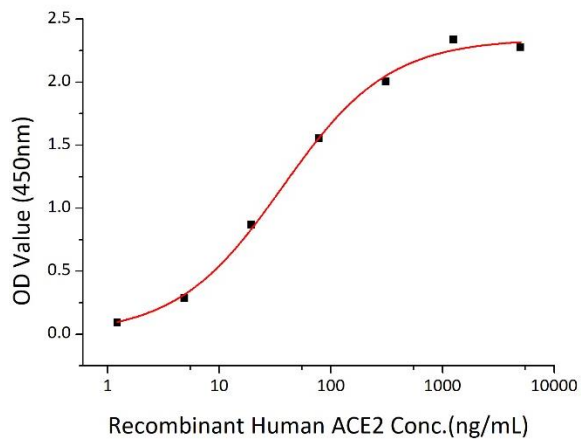
This product is stable at $\leq -70^{\circ}\text{C}$ for up to 6 months from the date of receipt. For optimal storage, aliquot into smaller quantities after centrifugation and store at the recommended temperature. Avoid repeated freeze-thaw cycles.

Alternative Names

Cov spike Protein; 2019-nCoV; ncov spike Protein; 2019-nCoV; coronavirus spike Protein; 2019-nCoV; S glycoprotein; COVID-19; S protein; Spike glycoprotein; SARS-CoV-2



Recombinant SARS-CoV-2 S1+S2 ECD (S-ECD) Protein with His-Tag was determined by SDS-PAGE with Coomassie Blue, showing bands around 80,110,180 kDa.



Immobilized Recombinant SARS-CoV-2 Spike S1+S2 ECD-His at 2 μ g/mL (100 μ L/well) can bind Recombinant Human ACE2 with a linear range of 1.5-36.9 ng/mL.