

---

# Recombinant SARS-CoV-2 Envelope Protein with His and Avi Tag

(Catalog # E80005)

---

## Description

Recombinant SARS-CoV-2 Envelope Protein is produced by E. coli expression system. The target protein is expressed with sequence (Met1-Val175) of 2019-ncov envelope (Accession #QHD43418.1) fused with an initial Met, a 6xHis, Avi tag at the N-terminus.

## Bio-Activity

Measured by its binding ability in a functional ELISA. Immobilized Recombinant SARS-COV-2 Envelope at 2ug/mL (100 uL/well) can bind Recombinant SARS-COV-2 Nucleocapsid with a linear range of 1.22-166.1 ng/mL.

## Purity

> 95% by SDS-PAGE.

## Endotoxin

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

## Formulation

Supplied as a 0.22 um filtered solution in 20mM Tris, 250mM NaCl, 0.5%TritonX-100, pH 8.0.

## Species

SARS-COV-2

## Calculated MW

11.4kDa

## Expressed Host

*E. coli*

## Tag

6xHis, Avi tag at the N-terminus

## Gene ID

43740570

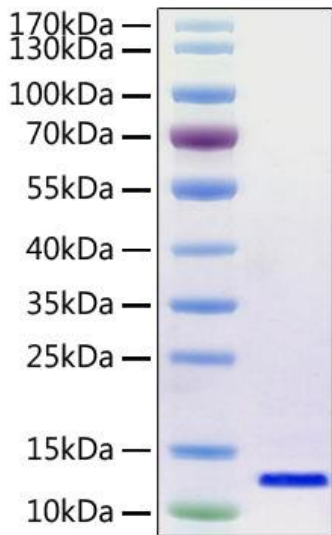
## Storage

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.

## Alternative Names

2019-nCoV E protein; 2019-nCoV sM protein; Envelope protein; Env polyprotein; Envelope glycoprotein; env; COVID-19; SARS-CoV-2; 2019-nCoV

---



Recombinant SARS-CoV-2 envelope Protein with His and Avi Tag was determined by SDS-PAGE with Coomassie Blue, showing a band at 12 kDa.