

## Recombinant SARS-CoV-2 NSP9 protein

**Catalog No:** 81322, 81622

**Expressed In:** *E. coli*

**Quantity:** 50, 1000 µg

**Concentration:** 1.5 µg/µl

**Source:** SARS-CoV-2

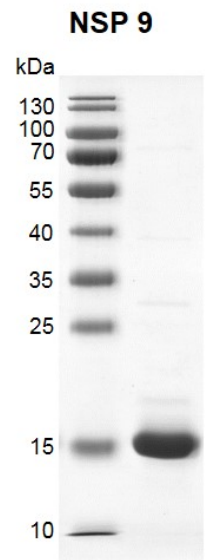
**Buffer Contents:** Recombinant SARS-CoV-2 NSP9 protein is supplied in 25 mM Tris pH 7.4, 300 mM NaCl, 10% glycerol, and 0.5 mM TCEP.

**Background:** SARS-CoV-2 NSP9 (Nonstructural Protein 9) is one of the nonstructural proteins encoded by SARS-CoV-2 orf1ab. The polyproteins of CoVs are cleaved by virus-encoded cysteine proteinases comprise papain- and chymotrypsinlike proteases into 16 nonstructural proteins including the expression of NSP1 to NSP11 by orf1a and encoding NSP12 to NSP16 by orf1b. According to BLAST analysis, the sequence identity of ORF1ab protein between SARS-CoV-2 and SARS-CoV is more than 90% with the query cover of about 100%, the sequence identity of NSP9 between these two viruses is about 100%.

**Protein Details:** Recombinant SARS-CoV-2 NSP9 protein was expressed in *E. coli* cells as the full length protein (accession number YP\_009725305.1) with a C-terminal Avi and a C-terminal 6×His tag. The molecular weight of the protein is 12.38 kDa.

**Application Notes:** Recombinant SARS-CoV-2 NSP9 protein is suitable for use in the study of SARS-CoV-2. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for this product is shown.

**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of arrival.



**Recombinant SARS-CoV-2 NSP9 protein gel**  
12.5% SDS-PAGE with Coomassie blue staining

MW: 12.38 kDa

Purity: >90%