

# PRODUCT INFORMATION



## SARS-CoV-2 nsp1 (recombinant)

Item No. 40879

### Overview and Properties

<b>Synonyms:</b>	SARS-CoV-2 Host Translation Inhibitor nsp1, SARS-CoV-2 Non-structural Protein 1, Severe Acute Respiratory Syndrome Coronavirus 2 nsp1
<b>Source:</b>	Recombinant SARS-CoV-2 N-terminal His-tagged nsp1 expressed in <i>E. coli</i>
<b>Amino Acids:</b>	13-128
<b>Uniprot No.:</b>	PODTD1
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Purity:</b>	≥90%
<b>Supplied in:</b>	50 mM Tris-HCl, pH 7.5, 200 mM sodium chloride, and 20% glycerol
<b>Endotoxin Testing:</b>	<1.0 EU/μg, determined by the LAL endotoxin assay
<b>Protein Concentration:</b>	<i>batch specific</i> mg/ml
<b>Special Conditions:</b>	Avoid repeated freeze/thaw cycles.

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description

Severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) is an enveloped positive-stranded RNA virus and the causative agent of COVID-19, a primarily respiratory illness characterized by fever, cough, and shortness of breath that can lead to life-threatening complications.<sup>1-5</sup> The SARS-CoV-2 genome contains approximately 30 kilobases and 14 open reading frames (ORFs) that encode four structural proteins: spike, envelope, membrane, and nucleocapsid, as well as 16 non-structural proteins and 9 accessory factors.<sup>6</sup> SARS-CoV-2 non-structural protein 1 (nsp1) is a multifunctional protein and virulence factor that inhibits host mRNA translation, including translation of antiviral proteins such as IFN-β and IFN-λ1.<sup>7-9</sup> It is composed of an N-terminal domain and a C-terminal domain connected *via* an unstructured 20-amino acid residue linker.<sup>9,10</sup> The SARS-CoV-2 nsp1 C-terminal domain binds to the mRNA entry site of the 40S subunit of the host ribosome, which prevents host mRNA from entering, while the N-terminal binds to the 40S decoding center preventing accurate alignment of tRNA and mRNA.<sup>8-10</sup> SARS-CoV-2 nsp1 also induces host mRNA degradation, prevents mRNA export from the nucleus to the cytoplasm, and binds to pre-rRNA in the nucleolus to disrupt rRNA biogenesis.<sup>7,9,11,12</sup>

### References

1. Kandeel, M., Ibrahim, A., Fayed, M., *et al. J. Med. Virol.* **92(6)**, 660-666 (2020).
2. Lu, R., Zhao, X., Li, J., *et al. Lancet* **395(10224)**, 565-574 (2020).
3. Meo, S.A., Alhowikan, A.M., Al-Khlaiwi, T., *et al. Eur. Rev. Med. Pharmacol. Sci.* **24(4)**, 2012-2019 (2020).
4. Klok, F.A., Kruip, M.J.H.A., van der Meer, N.J.M., *et al. Thromb. Res.* **191**, 145-147 (2020).
5. Yang, F., Shi, S., Zhu, J., *et al. J. Med. Virol.* **92(11)**, 2511-2515 (2020).
6. Romano, M., Ruggiero, A., Squeglia, F., *et al. Cells* **9(5)**, 1267 (2020).
7. Yerlici, V.T., Astori, A., Kejiou, N.S., *et al. Cell Rep.* **43(3)**, 113891 (2024).
8. Thoms, M., Buschauer, R., Ameisemeier, M., *et al. Science* **369(6508)**, 1249-1255 (2020).
9. Karousis, E.D. *Biochem. Soc. Trans.* **52(1)**, 481-490 (2024).
10. Schubert, K., Karousis, E.D., Jomaa, A., *et al. Nat. Struct. Mol. Biol.* **27(10)**, 959-966 (2020).
11. Mendez, A.S., Ly, M., González-Sánchez, A.M., *et al. Cell Rep.* **37(3)**, 109841 (2021).
12. Zhang, K., Miorin, L., Makio, T., *et al. Sci. Adv.* **7(6)**, eabe7386 (2021).

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/01/2024

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM