

# PRODUCT INFORMATION



## SARS-CoV-2 nsp16 Methyltransferase (recombinant)

Item No. 31816

### Overview and Properties

**Synonyms:** 2019-nCoV nsp16, 2'-O-MTase, ME, nsp16, 2'-O-methyltransferase, SARS-CoV-2 nsp16, SARS-CoV-2 nsp16, Severe Acute Respiratory Syndrome Coronavirus 2 nsp16 Methyltransferase

**Source:** Recombinant SARS-CoV-2 C-terminal His-tagged nsp16 methyltransferase expressed in *E. coli*

**Uniprot No.:** P0DTC2

**Amino Acids:** 6,799-7,096

**Molecular Weight:** 33.5 kDa

**Storage:** -20°C (as supplied)

**Stability:** ≥1 year

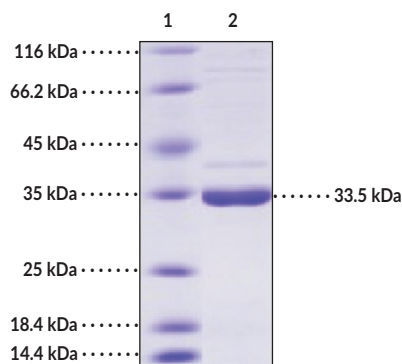
**Purity:** ≥85% estimated by SDS-PAGE

**Supplied in:** Lyophilized from sterile 20mM Tris 150mM sodium chloride, pH 8.3, with 5% trehalose, 5% mannitol, and 0.01% tween-80

**Reconstitution:** Add 670 µl of sterile water to the vial to prepare a stock solution of 0.15 mg/ml

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

### Image



Lane 1: MW Markers

Lane 2: SARS-CoV-2 nsp16 Methyltransferase (recombinant)

**SDS-PAGE Analysis of SARS-CoV-2 nsp16 Methyltransferase (recombinant).** This protein has a calculated molecular weight of 33.5 kDa.

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, 03/27/2024

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

# PRODUCT INFORMATION



## Description

---

Severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) (recombinant) 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 nsp16 methyltransferase (recombinant) is a 2'-O-methyltransferase (2'-O-MTase) and has a role in 5'-end capping of viral mRNAs.<sup>7,8</sup> In coronaviruses, nsp16 forms a complex with nsp10 to methylate nascent mRNAs at the ribose 2'-O position, creating a Cap-1 structure that facilitates increased translation of viral mRNAs and reduced innate immune recognition by the host cell.<sup>8,9</sup> Deficiency of nsp16 in the related virus SARS-CoV reduces viral RNA synthesis by approximately 10-fold *in vitro*, and mutations in the nsp16 KDKE catalytic tetrad in mouse-adapted SARS-CoV attenuate the virus *in vivo*.<sup>8,10</sup> An nsp10-derived peptide inhibitor of the nsp16/nsp10 complex increases survival in a mouse model of infection with the coronavirus mouse hepatitis virus (MHV).<sup>9</sup> Cayman's SARS-CoV-2 nsp16 Methyltransferase (recombinant) protein consists of 299 amino acids and has a calculated molecular weight of 33.5 kDa.

## References

---

1. Kandeel, M., Ibrahim, A., Fayez, M., *et al.* From SARS and MERS CoVs to SARS-CoV-2: Moving toward more biased codon usage in viral structural and nonstructural genes. *J. Med. Virol.* **92(6)**, 660-666 (2020).
2. Lu, R., Zhao, X., Li, J., *et al.* Genomic characterisation and epidemiology of 2019 novel coronavirus: Implications for virus origins and receptor binding. *Lancet* **395(10224)**, 565-574 (2020).
3. Meo, S.A., Alhowikan, A.M., Al-Khlaiwi, T., *et al.* Novel coronavirus 2019-nCoV: Prevalence, biological and clinical characteristics comparison with SARS-CoV and MERS-CoV. *Eur. Rev. Med. Pharmacol. Sci.* **24(4)**, 2012-2019 (2020).
4. Klok, F.A., Kruijff, M.J.H.A., van der Meer, N.J.M., *et al.* Incidence of thrombotic complications in critically ill ICU patients with COVID-19. *Thromb. Res.* **191**, 145-147 (2020).
5. Yang, F., Shi, S., Zhu, J., *et al.* Analysis of 92 deceased patients with COVID-19. *J. Med. Virol.* (2020).
6. Romano, M., Ruggiero, A., Squeglia, F., *et al.* A structural view of SARS-CoV-2 RNA replication machinery: RNA synthesis, proofreading and final capping. *Cells* **9(5)**, 1267 (2020).
7. Krafcikova, P., Silhan, J., Nancka, R., *et al.* Structural analysis of the SARS-CoV-2 methyltransferase complex involved in RNA cap creation bound to sinefungin. *Nat. Commun.* **11(1)**, 3717 (2020).
8. Viswanathan, T., Arya, S., Chan, S.-H., *et al.* Structural basis of RNA cap modification by SARS-CoV-2. *Nat. Commun.* **11(1)**, 3718 (2020).
9. Wang, Y., Sun, Y., Wu, A., *et al.* Coronavirus nsp10/nsp16 methyltransferase can be targeted by nsp10-derived peptide *in vitro* and *in vivo* to reduce replication and pathogenesis. *J. Virol.* **89(16)**, 8416-8427 (2015).
10. Menachery, V.D., Yount, B.L.J., Josset, L., *et al.* Attenuation and restoration of severe acute respiratory syndrome coronavirus mutant lacking 2'-O-methyltransferase activity. *J. Virol.* **88(8)**, 4251-4264 (2014).

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