

PRODUCT INFORMATION

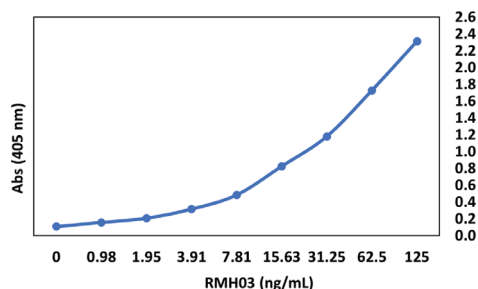
SARS-CoV-2 Nucleocapsid Chimeric Human Monoclonal Antibody (Clone RMH03)

Item No. 42860

Overview and Properties

Contents:	This vial contains 100 µg of protein A-affinity purified recombinant monoclonal antibody
Synonyms:	2019-nCoV NP, 2019-nCoV Nucleocapsid Protein, 2019-nCoV Nucleoprotein, COVID-19 NP, COVID-19 Nucleocapsid Protein, COVID-19 Nucleoprotein, SARS-CoV-2 NP, SARS-CoV-2 Nucleoprotein, Severe Acute Respiratory Syndrome Coronavirus 2 Nucleocapsid Protein
Immunogen:	Recombinant SARS-CoV-2 nucleocapsid protein
Cross Reactivity:	(+) SARS-CoV-2 nucleocapsid protein; (-) SARS-CoV nucleocapsid protein
Species Reactivity:	(+) SARS-CoV-2
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RMH03
Host:	<i>E. coli</i>
Isotype:	IgG
Applications:	Sandwich ELISA with SARS-CoV-2 Nucleocapsid Rabbit Monoclonal Antibody (Clone RM420). Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Binding curve of SARS-CoV-2 Nucleocapsid Chimeric Human Monoclonal Antibody (Clone RMH03) and recombinant 2019-nCoV Nucleocapsid protein. ELISA plate was coated with 50 ng recombinant 2019-nCoV Nucleocapsid protein at a concentration of 1 µg/ml. A two-fold serial dilution from 125 ng/ml was performed using RMH03. For detection, AP-conjugated goat anti-human IgG secondary antibody was used.

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, 03/24/2025

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) nucleocapsid protein is a viral protein encoded by the *N* gene in SARS-CoV-2 RNA.¹ SARS-CoV-2 is a member of the *Betacoronavirus* genus of viruses and has 88% sequence identity with two bat-derived SARS-like CoVs.² SARS-CoV-2 nucleocapsid protein is composed of an N-terminal domain (NTD) that binds RNA, a disordered linker region that contains regulatory phosphorylation sites, a C-terminal domain (CTD) that mediates dimerization, and a C-terminal tail.³ In a similar virus, SARS-CoV, the nucleocapsid protein packages the viral RNA into a helical ribonucleoprotein (RNP) complex that is a template for viral replication.⁴ The SARS-CoV nucleocapsid protein is integral for viral self-assembly and is involved in cell cycle regulation. The SARS-CoV-2 nucleocapsid protein gene sequence is greater than 90% similar to the SARS-CoV nucleocapsid protein, and it contains 27 T cell epitopes that are identical to SARS-CoV T cell epitopes.⁵ SARS-CoV-2 nucleocapsid protein inhibits the formation of stress granules, which sequester viral factors and are involved in the host antiviral response.⁶ SARS-CoV-2 is 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.⁷⁻⁹ Cayman's SARS-CoV-2 Nucleocapsid Chimeric Human Monoclonal Antibody (Clone RMH03) is composed of human IgG1 constant domains fused to the antigen-binding domain of a rabbit anti-SARS-CoV-2 nucleocapsid monoclonal antibody. It can be used as a detection antibody paired with the capture antibody SARS-CoV-2 Nucleocapsid Rabbit Monoclonal Antibody (Clone RM420) (Item No. 42859) in sandwich ELISAs.

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. Zhang, B., Tian, J., Zhang, Q., *et al.* Comparing the nucleocapsid proteins of human coronaviruses: Structure, immunoregulation, vaccine, and targeted drug. *Front. Mol. Biosci.* **9**, 761173 (2022).
4. Chang, C.-K., Hou, M.-H., Chang, C.-F., *et al.* The SARS coronavirus nucleocapsid protein – Forms and functions. *Antiviral Res.* **103**, 39-50 (2014).
5. Ahmed, S.F., Quadeer, A.A., and McKay, M.R. Preliminary identification of potential vaccine targets for the COVID-19 coronavirus (SARS-CoV-2) based on SARS-CoV immunological studies. *Viruses* **12**(3), E254 (2020).
6. Zheng, Z.-Q., Wang, S.-Y., Xu, Z.-S., *et al.* SARS-CoV-2 nucleocapsid protein impairs stress granule formation to promote viral replication. *Cell Discov.* **7**(1), 38 (2021).
7. 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).
8. Klok, F.A., Kruip, 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).
9. Yang, F., Shi, S., Zhu, J., *et al.* Analysis of 92 deceased patients with COVID-19. *J. Med. Virol.* **92**(11), 2511-2515 (2020).

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