

PRODUCT INFORMATION



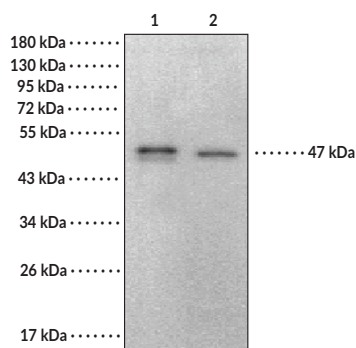
SARS-CoV-2 Nucleocapsid Protein Rabbit Monoclonal Antibody (Clone 0004)

Item No. 31801

Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	SARS-CoV-2 NP, SARS-CoV-2 Nucleoprotein, Severe Acute Respiratory Syndrome Coronavirus 2 Nucleocapsid Protein, COVID-19 NP, COVID-19 Nucleoprotein, COVID-19 Nucleocapsid Protein, 2019-nCoV NP, 2019-nCoV Nucleoprotein, 2019-nCoV Nucleocapsid Protein
Immunogen:	Recombinant SARS-CoV-2 nucleocapsid protein
Cross Reactivity:	(+) Nucleocapsid protein
Species Reactivity:	(+) SARS-CoV-2, SARS-CoV-2 Omicron (B.1.1.529), SARS-CoV-2 Omicron (BA.2), SARS-CoV-2 Omicron XE (BA.1 x BA.2), SARS-CoV-2 Omicron (BA.4), SARS-CoV, MERS-CoV; (-) HCoV-229E, HCoV-NL63, HCoV-HKU1, HCoV-OC43, Influenza A H1N1 (A/California/07/2009), Influenza A H3N2 (A/Switzerland/9715293/2013), Influenza A H2N2 (A/Ann Arbor/6/1960), Influenza A H7N9 (A/Shanghai/2/2013), Influenza B (B/Florida/4/2006))
MW:	~47 kDa
Form:	Liquid
Storage:	-80°C (as supplied)
Stability:	≥1 year
Storage Buffer:	0.2 µm filtered solution in PBS
Clone:	004
Host:	Rabbit
Isotype:	IgG
Applications:	ELISA and Western blot (WB); the recommended starting concentration is 1:5,000-1:10,000 for ELISA and 1:500-1:2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: SARS-CoV-2 Nucleocapsid Protein Rabbit Monoclonal Antibody (Clone 0004) (30 ng)
Lane 2: MERS-CoV Nucleocapsid (30 ng)

WB of SARS-CoV-2 Nucleocapsid Protein Rabbit Monoclonal Antibody (Clone 0004) at 1:10,000 dilution.

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, 11/09/2023

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 Protein Rabbit Monoclonal Antibody (Clone 0004) can be used for ELISA and Western blot (WB) applications. This recombinant antibody recognizes nucleocapsid protein at 45.6 kDa from SARS-CoV-2, several SARS-CoV-2 Omicron subvariants, SARS-CoV, and Middle East respiratory syndrome coronavirus (MERS-CoV).

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