

# PRODUCT INFORMATION

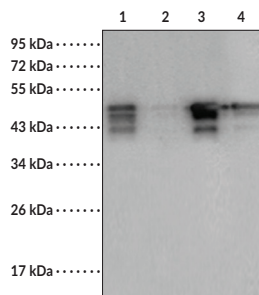
## SARS-CoV/SARS-CoV-2 Nucleocapsid Protein Rabbit Monoclonal Antibody (Clone 004)

Item No. 31986

### Overview and Properties

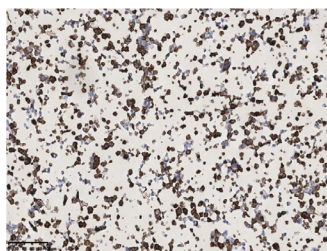
<b>Contents:</b>	This vial contains 50 or 100 µl of protein A-affinity purified recombinant monoclonal antibody.
<b>Synonyms:</b>	COVID-19 NP, COVID-19 Nucleocapsid Protein, COVID-19 Nucleoprotein, 2019-nCoV NP, 2019-nCoV Nucleocapsid Protein, 2019-nCoV Nucleoprotein, SARS-CoV-2 NP, SARS-CoV-2 Nucleoprotein, Severe Acute Respiratory Syndrome Coronavirus 2 Nucleocapsid Protein
<b>Immunogen:</b>	Recombinant SARS-CoV nucleocapsid protein
<b>Cross Reactivity:</b>	See page 2
<b>Species Reactivity:</b>	See page 2
<b>Molecular Weight:</b>	45.6 kDa
<b>Uniprot No.:</b>	PODTC9
<b>Form:</b>	Liquid
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	0.2 µm filtered solution in PBS
<b>Clone:</b>	004
<b>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Applications:</b>	ELISA, Flow cytometry (FC), Immunocytochemistry (ICC), Immunofluorescence (IF), Immunohistochemistry-paraffin (IHC-P), and Western blot (WB); the recommended starting dilution is 1:5,000-1:10,000 for ELISA, 1:25-1:100 for FC, 1:20-1:100 for IF, 1:100-1:500 for IHC-P, and 1:1,000-1:10,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images

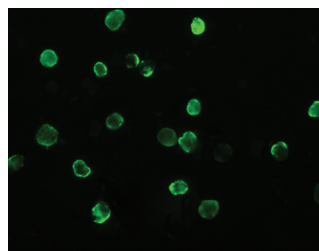


Lane 1: SARS-CoV NP Protein (30 ng)  
 Lane 2: SARS-CoV NP Protein (5 ng)  
 Lane 3: SARS-CoV-2 (2019-nCoV) NP Protein (30 ng)  
 Lane 4: SARS-CoV-2 (2019-nCoV) NP Protein (5 ng)

WB of SARS-CoV/SARS-CoV-2 Nucleocapsid Protein Rabbit Monoclonal Antibody (Clone 004) at 1:1,000 dilution.



Immunohistochemical analysis of nucleocapsid protein overexpressed in HEK293 cells. Cells were stained with Cayman's SARS-CoV/SARS-CoV-2 Nucleocapsid Protein Rabbit Monoclonal Antibody (Clone 004), then a HRP-conjugated second step antibody.



Immunofluorescence analysis of nucleocapsid protein overexpressed in HEK293 cells. Cells were stained with Cayman's SARS-CoV/SARS-CoV-2 Nucleocapsid Protein Rabbit Monoclonal Antibody (Clone 004), then an AlexaFluor®488-conjugated second step antibody.

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

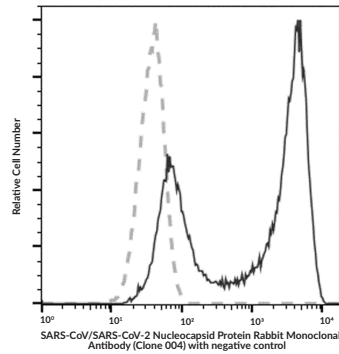
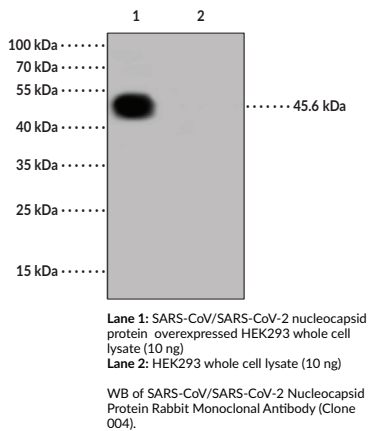
**SAFETY DATA**  
 This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your 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, 01/10/2023

**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



Flow cytometric analysis of SARS-CoV-2 nucleocapsid protein overexpressed in HEK293 Cells. Cells were stained with purified SARS-CoV/SARS-CoV-2 Nucleocapsid Protein Rabbit Monoclonal Antibody (Clone 004), then a FITC-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.

## Reactivity

**Cross Reactivity:** (+) Nucleocapsid protein,  
SARS-CoV-2 Delta (B.1.617) nucleocapsid protein (D377Y),  
SARS-CoV-2 Delta (B.1.617) nucleocapsid protein (R230M, D377Y),  
SARS-CoV-2 Delta (B.1.617.3) nucleocapsid protein (P67S, R203M, D377Y),  
SARS-CoV-2 Alpha (B.1.1.7) nucleocapsid protein (R203K, G204R),  
SARS-CoV-2 Alpha (B.1.1.7) nucleocapsid protein (I292T),  
SARS-CoV-2 Alpha (B.1.1.7) nucleocapsid protein (D3L, R203K, G204R, S235F),  
SARS-CoV-2 Alpha (B.1.1.7) nucleocapsid protein (D3L, S235F),  
SARS-CoV-2 Alpha (B.1.1.7/B.1.237) nucleocapsid protein (S194L),  
SARS-CoV-2 Alpha/Beta (B.1.1.7/B.1.351/A.2.2) nucleocapsid protein (P13L),  
SARS-CoV-2 Beta (B.1.351/B.1.351.2/B.1.351.3/B.1.427/B.1.429) nucleocapsid protein (T205I),  
SARS-CoV-2 Gamma (P.1/P.1.1/P.1.2) nucleocapsid protein (P80R),  
SARS-CoV-2 Eta (B.1.525) nucleocapsid protein (A12G, T205I)

**Species Reactivity:** (+) SARS-CoV,  
SARS-CoV-2,  
SARS-CoV-2 Delta (B.1.617),  
SARS-CoV-2 Delta (B.1.617.2),  
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);

(-) MERS-CoV,  
HCoV-229E,  
HCoV-NL63,  
HCoV-HKU1 (isolate N5),  
HCoV-OC43

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 (SARS-CoV) and SARS-CoV-2 nucleocapsid proteins are encoded by the N gene in SARS-CoV and SARS-CoV-2 RNA.<sup>1,2</sup> SARS-CoV and SARS-CoV-2 are members of the *Betacoronavirus* genus of viruses that have approximately 79% sequence identity and share 27 T cell epitopes in common.<sup>3-5</sup> The SARS-CoV-2 nucleocapsid protein has greater than 90% similarity to the SARS-CoV nucleocapsid protein and contains two unique B cell epitopes and two T cell epitopes that are structurally stable, non-allergenic, and induce production of IFN- $\gamma$ .<sup>2,5</sup> SARS-CoV and SARS-CoV-2 nucleocapsid proteins package the viral RNA into a helical ribonucleoprotein complex (RNP), which is a template for viral replication, and are integral for viral self-assembly and involved in regulation of the host cell cycle.<sup>2,6</sup> SARS-CoV and SARS-CoV-2 are the causative agents of SARS and COVID-19, respectively, both of which are primarily respiratory illnesses characterized by fever, cough, and shortness of breath that can lead to life-threatening complications.<sup>7-9</sup> Cayman's SARS-CoV/SARS-CoV-2 Nucleocapsid Protein Rabbit Monoclonal Antibody (Clone 004) can be used for ELISA, flow cytometry (FC), immunocytochemistry (ICC), immunofluorescence (IF), immunohistochemistry-paraffin (IHC-P), and Western blot (WB) applications. This recombinant antibody recognizes wild-type and various mutant nucleocapsid proteins at 45.6 kDa from SARS-CoV, SARS-CoV-2, and several SARS-CoV-2 Alpha, Beta, Gamma, Delta, Eta, and Omicron subvariants.

## 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. Kwarteng, A., Asiedu, E., Sakyi, S.A., *et al.* Targeting the SARS-CoV2 nucleocapsid protein for potential therapeutics using immuno-informatics and structure-based drug discovery techniques. *Biomed. Pharmacother.* **132**, 110914 (2020).
3. 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).
4. 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).
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. 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).
7. 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).
8. 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