

PRODUCT INFORMATION

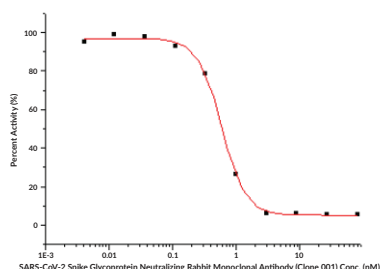
SARS-CoV-2 Spike Glycoprotein Neutralizing Rabbit Monoclonal Antibody (Clone 001)

Item No. 31803

Overview and Properties

Contents:	This vial contains 100 µg of protein A-affinity purified monoclonal antibody.
Synonyms:	SARS-CoV-2 Spike RBD, SARS-CoV-2 Spike Receptor Binding Domain, SARS-CoV-2 Surface Glycoprotein RBD, SARS-CoV-2 Surface Glycoprotein Receptor Binding Domain, Severe Acute Respiratory Syndrome Coronavirus 2 Spike Glycoprotein Receptor Binding Domain, Severe Acute Respiratory Syndrome Coronavirus 2 Surface Glycoprotein Receptor Binding Domain
Immunogen:	Recombinant SARS-CoV-2 spike glycoprotein RBD (C-terminal mouse IgG1 Fc-tagged)
Cross Reactivity:	See page 2
Species Reactivity:	See page 2
Form:	Liquid
Storage:	-80°C (as supplied)
Stability:	≥1 year
Storage Buffer:	0.2 µm filtered solution in histidine and arginine with 120 mM sodium chloride, 0.02% Tween 80, and pH 6.0
Clone:	001
Host:	Rabbit
Isotype:	IgG
Applications:	ELISA, Flow cytometry (FC), and Microneutralization (MN) assays; the recommended starting concentration is 0.1-0.2 µg/ml for ELISA and 10-40 µg/ml for FC. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

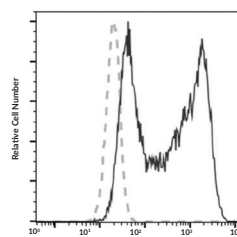
Images



Log dilutions of SARS-CoV-2 Spike Glycoprotein Neutralizing Rabbit Monoclonal Antibody (Clone 001) were detected by SARS-CoV-2 (2019-nCoV) Inhibitor Screening ELISA Kit. The IC_{50} value is typically 0.5-5 nM for the wild-type (WT) protein.

WT		Delta	
Conc. (µg/mL)	Inhibition%	Conc. (µg/mL)	Inhibition%
100	99.99%	100	100.00%
20	99.96%	20	100.00%
4	99.91%	4	99.99%
0.8	71.98%	0.8	97.57%
0.16	36.62%	0.16	50.76%
0.032	43.77%	0.032	17.36%
0.0064	24.65%	0.0064	2.53%

The neutralization activity is measured by microneutralization (MN) assay *in vitro*. The virus MN test was performed on 293T-ACE2 cells infected with SARS-CoV-2 Spike Pseudovirus, SARS-CoV-2 Spike Omicron (B.1.1.529) Pseudovirus, or SARS-CoV-2 Delta (B.1.617.2) variant Spike Pseudovirus, under treatment of log dilutions of neutralizing antibody. The infection was neutralized by increasing concentrations of SARS-CoV-2 Spike Glycoprotein Neutralizing Rabbit Monoclonal Antibody (Clone 001). Rate of inhibition was determined by comparing the Relative Light Unit (RLU) of luciferase reporter in different antibody concentrations. The IC_{50} values are 0.29 and 0.15 µg/ml for WT and Delta (B.1.617.2) variant spike pseudovirus, respectively. No neutralization activity to SARS-CoV-2 Spike Omicron (B.1.1.529) Pseudovirus was observed.



Flow cytometric analysis of SARS-CoV-2 spike glycoprotein overexpressed in HEK293 cells. Cells were stained with purified SARS-CoV-2 Spike Glycoprotein Neutralizing Rabbit Monoclonal Antibody (Clone 001), followed by 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.

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, 08/06/2024

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



Reactivity

Cross Reactivity: (+) SARS-CoV-2 (BA.2.75) spike glycoprotein S1+S2 trimer,
SARS-CoV-2 Delta (B.1.617.2) spike glycoprotein S1+S2 trimer,
SARS-CoV-2 Delta (B.1.617.2) spike glycoprotein S1 subunit,
SARS-CoV-2 spike glycoprotein S1 subunit;

(-) SARS-CoV-2 Omicron (BA.1.1) spike glycoprotein S1+S2 trimer,
SARS-CoV-2 Omicron (B.1.1.529) spike glycoprotein S1 subunit,
SARS-CoV-2 Omicron (B.1.1.529) S1+S2 trimer,
SARS-CoV-2 Omicron (BA.2) spike glycoprotein S1+S2 trimer,
SARS-CoV-2 Omicron (BA.2) spike glycoprotein S1 subunit,
SARS-CoV-2 Omicron (BA.2) spike glycoprotein S1 subunit NTD,
SARS-CoV-2 Omicron (BA.2.75.2) spike glycoprotein S1+S2 trimer,
SARS-CoV-2 XD (BA.1 x AY.4) spike glycoprotein S1+S2 trimer,
SARS-CoV-2 (BA.4.6) spike glycoprotein S1+S2 trimer,
SARS-CoV-2 Omicron (BF.7) spike glycoprotein S1+S2 trimer,
SARS-CoV-2 Omicron (BQ.1.1) spike glycoprotein S1+S2 trimer,
SARS-CoV-2 Delta (B.1.617.2) spike glycoprotein S1 subunit NTD,
SARS-CoV spike glycoprotein S1 subunit,
MERS-CoV spike glycoprotein S1 subunit,
HCoV-HKU1 (isolate N1) spike glycoprotein S1 subunit,
HCoV-HKU1 (isolate N5) spike glycoprotein S1 subunit,
HCoV-NL63 spike glycoprotein S1 subunit,
HCoV-229E spike glycoprotein S1 subunit,
HCoV-OC43 spike glycoprotein S1+S2 ECD

Species Reactivity: (+) SARS-CoV-2,
SARS-CoV-2 Delta (B.1.617.2),
SARS-CoV-2 (BA.2.75),
SARS-CoV-2 (BA.2.3.20),
SARS-CoV-2 Omicron (BA.1.1);

(-) SARS-CoV-2 Omicron (BA.1.1.529),
SARS-CoV-2 Omicron (BA.2),
SARS-CoV-2 Omicron (BA.2.12.1),
SARS-CoV-2 Omicron (BA.2.75.2),
SARS-CoV-2 Omicron (BA.4),
SARS-CoV-2 Omicron (BA.4.6/BF.7),
SARS-CoV-2 Omicron (BA.5),
SARS-CoV-2 Omicron (BQ.1.1),
SARS-CoV-2 Omicron (XBB)

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) is an enveloped positive-stranded RNA virus, a member of the *Betacoronavirus* genus, and the causative agent of COVID-19.¹⁻⁵ The SARS-CoV-2 spike glycoprotein, also known as the surface glycoprotein, is located on the outer envelope of the virion.¹ It is composed of an S1 and S2 subunit divided by a furin S-cleavage site not found in other SARS-CoVs.^{6,7} The S1 subunit contains the receptor-binding domain (RBD), which binds to the carboxypeptidase angiotensin-converting enzyme 2 (ACE2), and the S1 and S2 subunits are cleaved by the protease TMPRSS2 to facilitate viral fusion with the host cell membrane.⁸⁻¹⁰ In this way, ACE2 acts as the functional receptor for SARS-CoV-2. SARS-CoV-2 infection can result in the production of neutralizing antibodies, which bind to the SARS-CoV-2 spike RBD preventing further viral entry and infection, starting approximately 4-10 days after symptom onset.^{11,12} Cayman's SARS-CoV-2 Spike Glycoprotein Neutralizing Rabbit Monoclonal Antibody (Clone 001) disrupts the spike glycoprotein S1 subunit-ACE2 interaction and can be used for ELISA, and flow cytometry (FC) applications, as well as microneutralization (MN) assays.

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., 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).
5. Yang, F., Shi, S., Zhu, J., *et al.* Analysis of 92 deceased patients with COVID-19. *J. Med. Virol.* **92**(11), 2511-2515 (2020).
6. Liu, Z., Xiao, X., Wei, X., *et al.* Composition and divergence of coronavirus spike proteins and host ACE2 receptors predict potential intermediate hosts of SARS-CoV-2. *J. Med. Virol.* **92**(6), 595-601 (2020).
7. Walls, A.C., Park, Y.-J., Tortorici, M.A., *et al.* Structure, function, and antigenicity of the SARS-CoV-2 spike glycoprotein. *Cell* **181**(2), 281-292 (2020).
8. Hoffmann, M., Kleine-Weber, H., Schroeder, S., *et al.* SARS-CoV-2 cell entry depends on ACE2 and TMPRSS2 and is blocked by a clinically proven protease inhibitor. *Cell* **181**(2), 271-280 (2020).
9. Yan, R., Zhang, Y., Li, Y., *et al.* Structural basis for the recognition of the SARS-CoV-2 by full-length human ACE2. *Science* **267**(6485), 1444-1448 (2020).
10. Wrapp, D., Wang, N., Corbett, K.S., *et al.* Cryo-EM structure of the 2019-nCoV spike in the prefusion conformation. *Science* **367**(6483), 1260-1263 (2020).
11. Wang, A., Zhang, L., Sang, L., *et al.* Kinetics of viral load and antibody response in relation to COVID-19 severity. *J. Clin. Invest.* **130**(10), 5235-5244 (2020).
12. Xiang, F., Wang, X., He, X., *et al.* Antibody detection and dynamic characteristics in patients with coronavirus disease 2019. *Clin. Infect. Dis.* **71**(8), 1930-1934 (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