

# PRODUCT INFORMATION



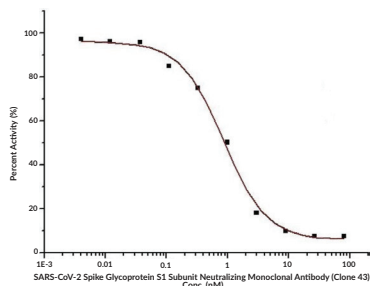
## SARS-CoV-2 Spike Glycoprotein S1 Subunit Neutralizing Monoclonal Antibody (Clone 43)

Item No. 31999

### Overview and Properties

|                            |   |
|----------------------------|---|
| <b>Contents:</b>           | This vial contains 100 µg of protein A-affinity purified monoclonal antibody.   |
| <b>Synonyms:</b>           | SARS-CoV-2 Spike S1 Subunit, SARS-CoV-2 Surface Glycoprotein S1 Subunit, Severe Acute Respiratory Syndrome Coronavirus 2 Spike Glycoprotein S1 Subunit, Severe Acute Respiratory Syndrome Coronavirus 2 Surface Glycoprotein S1 Subunit   |
| <b>Immunogen:</b>          | Recombinant SARS-CoV-2 spike glycoprotein S1 subunit (C-terminal mouse IgG1 Fc-tagged)  |
| <b>Cross Reactivity:</b>   | See page 2  |
| <b>Species Reactivity:</b> | See page 2  |
| <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>              | 43  |
| <b>Host:</b>               | Mouse   |
| <b>Isotype:</b>            | IgG1  |
| <b>Applications:</b>       | ELISA, Flow cytometry (FC), Immunohistochemistry (IHC; paraffin), and Microneutralization (MN) assays; the recommended starting concentration is 0.5-1 µg/ml for ELISA, 10-40 µg/ml for FC, and 10-50 µg/ml for IHC; paraffin. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically. |

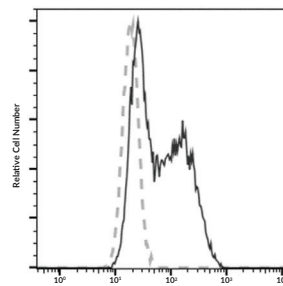
### Images



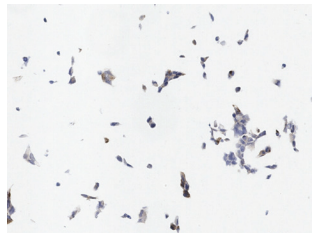
Serial dilutions of SARS-CoV-2 Spike Glycoprotein S1 Subunit Neutralizing Monoclonal Antibody (Clone 43) was detected by SARS-CoV-2 (2019-nCoV) Inhibitor Screening ELISA Kit. The IC<sub>50</sub> value is typically 0.5-5 nM for the wild-type protein.

| Conc. (µg/mL) | Inhibition% |
|---------------|-------------|
| 100           | 96.39%      |
| 10            | 94.57%      |
| 1             | 42.24%      |
| 0.1           | -5.53%      |

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 under treatment of log dilutions of neutralizing antibody. The infection was neutralized by increasing concentrations of SARS-CoV-2 Spike Glycoprotein S1 Subunit Neutralizing Monoclonal Antibody (Clone 43). Rate of inhibition was determined by comparing the Relative Light Unit (RLU) of luciferase reporter in different antibody concentrations. The IC<sub>50</sub> value is typically 1.41 µg/ml for the wild-type protein.



Flow cytometric analysis of SARS-CoV-2 spike overexpressed in HEK293 cells. Cells were stained with purified SARS-CoV-2 Spike Glycoprotein S1 Subunit Neutralizing Monoclonal Antibody (Clone 43), 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.



Immunohistochemical analysis of SARS-CoV-2 spike overexpressed in HEK293 cells. Cells were stained with purified SARS-CoV-2 Spike Glycoprotein S1 Subunit Neutralizing Monoclonal Antibody (Clone 43), followed by a HRP-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, 01/05/2023

**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



## Reactivity

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

(-) SARS-CoV-2 Omicron (BA.1.1) spike glycoprotein S1+S2 trimer, SARS-CoV-2 Omicron (BA.1.1) spike glycoprotein RBD, SARS-CoV-2 Omicron (B.1.1.529) S1+S2 trimer, SARS-CoV-2 Omicron (B.1.1.529) spike glycoprotein RBD, SARS-CoV-2 Omicron (BA.2) spike glycoprotein S1+S2 trimer, SARS-CoV-2 Omicron (BA.2) spike glycoprotein RBD, SARS-CoV-2 Omicron (BA.2) spike glycoprotein S1 subunit-NTD, SARS-CoV-2 Omicron (BA.2.12.1) spike glycoprotein RBD, SARS-CoV-2 Omicron (BA.4) spike glycoprotein RBD, SARS-CoV-2 Omicron (BA.5) spike glycoprotein RBD, SARS-CoV-2 XD (BA.1 x AY.4) spike glycoprotein S1+S2 trimer, SARS-CoV-2 Delta (B.1.617.2) spike glycoprotein S1 subunit-NTD, SARS-CoV-2 spike glycoprotein S1 subunit-NTD (C-terminal human IgG1 Fc- and AVI-tagged), SARS-CoV spike glycoprotein S1 subunit (mouse Fc-tagged), SARS-CoV spike glycoprotein RBD (His-tagged), 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 Delta (B.1.617.2) spike glycoprotein S1+S2 trimer, SARS-CoV-2 Delta (B.1.617.2) spike glycoprotein RBD, SARS-CoV-2 spike glycoprotein RBD;

(-) SARS-CoV-2 Omicron (BA.1.1) spike glycoprotein S1+S2 trimer, SARS-CoV-2 Omicron (BA.1.1) spike glycoprotein RBD, SARS-CoV-2 Omicron (B.1.1.529) S1+S2 trimer, SARS-CoV-2 Omicron (B.1.1.529) spike glycoprotein RBD, SARS-CoV-2 Omicron (BA.2) spike glycoprotein S1+S2 trimer, SARS-CoV-2 Omicron (BA.2) spike glycoprotein RBD, SARS-CoV-2 Omicron (BA.2) spike glycoprotein S1 subunit-NTD, SARS-CoV-2 Omicron (BA.2.12.1) spike glycoprotein RBD, SARS-CoV-2 Omicron (BA.4) spike glycoprotein RBD, SARS-CoV-2 Omicron (BA.5) spike glycoprotein RBD, SARS-CoV-2 XD (BA.1 x AY.4) spike glycoprotein S1+S2 trimer, SARS-CoV-2 Delta (B.1.617.2) spike glycoprotein S1 subunit-NTD, SARS-CoV-2 spike glycoprotein S1 subunit-NTD (C-terminal human IgG1 Fc- and AVI-tagged), SARS-CoV spike glycoprotein S1 subunit (mouse Fc-tagged), SARS-CoV spike glycoprotein RBD (His-tagged), 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

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.<sup>1-5</sup> The SARS-CoV-2 spike glycoprotein, also known as the surface glycoprotein, is located on the outer envelope of the virion.<sup>1</sup> It is composed of an S1 and S2 subunit divided by a furin S-cleavage site not found in other SARS-CoVs.<sup>6,7</sup> 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.<sup>8-10</sup> In this way, ACE2 acts as the functional receptor for SARS-CoV-2. The SARS-CoV-2 spike glycoprotein S1 subunit induces inflammatory gene expression in the frontal cortex, hippocampus, and hypothalamus, as well as activates toll-like receptor 2 (TLR2) and TLR4 signaling and increases social avoidance in the juvenile social exploration test in rats.<sup>11</sup> 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.<sup>12,13</sup> Cayman's SARS-CoV-2 Spike Glycoprotein S1 Subunit Neutralizing Monoclonal Antibody (Clone 43) disrupts the spike glycoprotein S1 subunit-ACE2 interaction and can be used for ELISA, flow cytometry (FC), and immunohistochemistry (IHC; paraffin) 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. Frank, M.G., Ngyuen, K.H., Ball, J.B., *et al.* SARS-CoV-2 spike S1 subunit induces neuroinflammatory, microglial and behavioral sickness responses: Evidence of PAMP-like properties. *Brain Behav. Immun.* **100**, 267-277 (2022).
12. 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).
13. 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