

# PRODUCT INFORMATION



## SARS-CoV/SARS-CoV-2 Spike Glycoprotein RBD Chimeric Monoclonal Antibody (Clone D004)

Item No. 31991

### Overview and Properties

|                            |   |
|----------------------------|---|
| <b>Contents:</b>           | This vial contains 50 or 100 µl of protein A-affinity purified monoclonal antibody.   |
| <b>Synonyms:</b>           | SARS-CoV/SARS-CoV-2 Spike RBD, SARS-CoV/SARS-CoV-2 Spike Receptor Binding Domain, SARS-CoV/SARS-CoV-2 Surface Glycoprotein RBD, SARS-CoV/SARS-CoV-2 Surface Glycoprotein Receptor Binding Domain, Severe Acute Respiratory Syndrome Coronavirus/Severe Acute Respiratory Syndrome Coronavirus 2 Spike Glycoprotein Receptor Binding Domain  |
| <b>Immunogen:</b>          | Recombinant C-terminal His-tagged SARS-CoV spike glycoprotein RBD   |
| <b>Cross Reactivity:</b>   | (+) SARS-CoV-2 Delta (B.1.617.2) spike glycoprotein S1+S2 trimer, SARS-CoV-2 Delta (B.1.617.2) spike glycoprotein S1+S2, SARS-CoV-2 Delta (B.1.617.2) spike glycoprotein S1 subunit, SARS-CoV-2 (BA.2.75) spike glycoprotein S1+S2 trimer, SARS-CoV-2 Omicron (BA.1.1) spike glycoprotein S1+S2 trimer, SARS-CoV-2 Omicron (B.1.1.529) S1+S2 trimer, SARS-CoV-2 Omicron (B.1.1.529) spike glycoprotein S1 subunit, SARS-CoV-2 Omicron (BA.2) spike glycoprotein S1 subunit, SARS-CoV-2 Omicron (BA.2) spike glycoprotein S1+S2 trimer, SARS-CoV-2 Omicron (BA.2.75.2) 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 spike glycoprotein S1 subunit, SARS-CoV spike glycoprotein S1 subunit; (-) SARS-CoV-2 Omicron (BA.2) spike glycoprotein S1 subunit NTD, SARS-CoV-2 Delta (B.1.617.2) spike glycoprotein S1 subunit NTD, 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 |
| <b>Species Reactivity:</b> | (+) SARS-CoV, SARS-CoV-2, SARS-CoV-2 Delta (B.1.617.2), SARS-CoV-2 Omicron (B.1.1.529), SARS-CoV-2 Omicron (BA.1.1), SARS-CoV-2 Omicron (BA.2), SARS-CoV-2 Omicron (BA.2.12.1), SARS-CoV-2 (BA.2.3.20), SARS-CoV-2 (BA.2.75), 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); other species not tested   |
| <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>              | D004  |
| <b>Host:</b>               | Chimeric monoclonal antibody combining the constant domains of human IgG1κ with variable regions from a mouse immunized with purified recombinant SARS-CoV spike glycoprotein RBD   |
| <b>Isotype:</b>            | Human IgG1κ   |
| <b>Applications:</b>       | ELISA, Flow cytometry (FC), Immunocytochemistry (ICC), Immunofluorescence (IF); the recommended starting dilution is 1:5,000-1:10,000 for ELISA, 1:25-1:100 for FC, and 1:20-1:100 for IF. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.   |

**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.

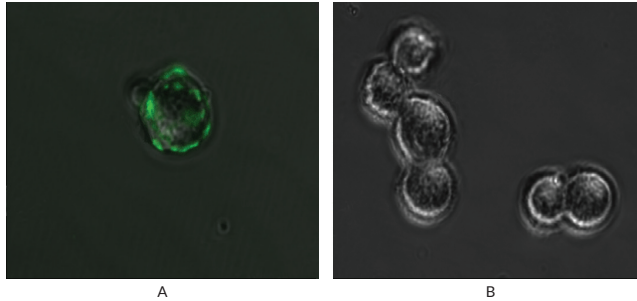
Copyright Cayman Chemical Company, 1/11/2022

**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

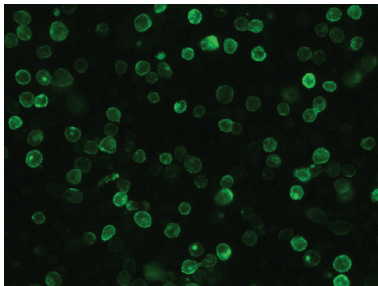
# PRODUCT INFORMATION



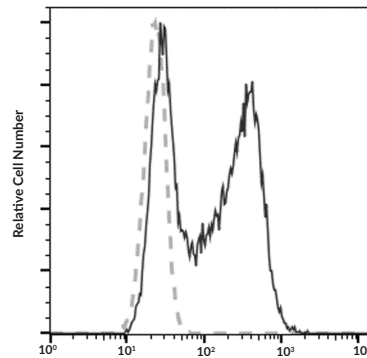
## Images



Immunofluorescence labeling of the SARS-CoV-2 spike glycoprotein in ACE2-overexpressed 293T cells infected (A) or noninfected (B) by 2019-nCoV-spike pseudovirus (PSV). Cells were blocked with 10% serum and incubated with SARS-CoV/SARS-CoV-2 Spike Glycoprotein RBD Chimeric Monoclonal Antibody (Clone D004) at a dilution of 1:60 at 37°C for 1 hour. Cells were then labeled with a FITC-conjugated goat anti-human IgG secondary antibody (green).



Immunofluorescent analysis of HEK293 cells overexpressing the spike glycoprotein. Cells were labeled with purified SARS-CoV/SARS-CoV-2 Spike Glycoprotein RBD Chimeric Monoclonal Antibody (Clone D004) followed by a FITC-conjugated secondary antibody.



SARS-CoV/SARS-CoV-2 Spike Glycoprotein RBD Chimeric Monoclonal Antibody (Clone D004) with negative control

Flow cytometric analysis of the SARS-CoV-2 spike glycoprotein overexpressed in HEK293 cells. Cells were labeled with purified SARS-CoV/SARS-CoV-2 Spike Glycoprotein RBD Chimeric Monoclonal Antibody (Clone D004) followed by a FITC-conjugated secondary antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.

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) spike glycoprotein, also known as the surface glycoprotein, is a viral structural protein encoded by the S gene in SARS-CoV RNA that contains the receptor binding domain (RBD).<sup>1</sup> SARS-CoV is a member of the *Betacoronavirus* genus of viruses and has an approximately 79% sequence identity with SARS-CoV-2, the causative agent of COVID-19.<sup>2,3</sup> SARS-CoV spike glycoprotein is a transmembrane glycoprotein that assembles into homotrimers on the virus surface and is composed of an N-terminal S1 subunit, which contains the receptor binding domain (RBD), and a C-terminal S2 subunit, which facilitates fusion between viral and host cell membranes.<sup>4-6</sup> The 193-amino acid RBD of the SARS-CoV spike protein is a target for neutralizing antibodies.<sup>5,7</sup> The SARS-CoV RBD, which spans amino acid residues 318 to 510, is 73% identical to that of SARS-CoV-2 and can bind to human angiotensin-converting enzyme 2 (ACE2), which is the host cell surface receptor for both SARS-CoV and SARS-CoV-2.<sup>4-7</sup> SARS-CoV is the causative agent of SARS, a primarily respiratory illness characterized by fever, cough, shortness of breath, and an approximately 10% fatality rate.<sup>3</sup> Cayman's SARS-CoV/SARS-CoV-2 Spike Glycoprotein RBD Chimeric Monoclonal Antibody (Clone D004) is composed of human IgG1k constant domains and variable regions from a mouse immunized with purified recombinant SARS-CoV spike glycoprotein RBD. It can be used for ELISA, flow cytometry (FC), immunocytochemistry (ICC), and immunofluorescence (IF) applications.

## 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. 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).
5. He, Y., Zhou, Y., Liu, S., *et al.* Receptor-binding domain of SARS-CoV spike protein induces highly potent neutralizing antibodies: Implication for developing subunit vaccine. *Biochem. Biophys. Res. Commun.* **324(2)**, 773-781 (2004).
6. 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).
7. Tian, X., Li, C., Huang, A., *et al.* Potent binding of 2019 novel coronavirus spike protein by a SARS coronavirus-specific human monoclonal antibody. *Emerg. Microbes Infect.* **9(1)**, 382-385 (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