PRODUCT INFORMATION



SARS-CoV-2 (human) Neutralizing Recombinant Antibody

Item No. 32526

Overview and Properties

Contents:	This vial contains 100 μg of protein A-purified recombinant monoclonal antibody.
Synonym:	Severe Acute Respiratory Syndrome Coronavirus 2 (human) Neutralizing
Preparation:	This antibody was produced as a recombinant protein expressed in CHO cells using engineered constructs based on antibodies from COVID-19 patients
Species Reactivity:	(+) Human
Form:	Solid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 0.1% BSA when reconstituted with 100 μ l deionized water
Host:	CHO cells
Applications:	ELISA, Immunofluorescence (IF), and Western blot (WB); the recommended starting dilution is 1:1,000 for ELISA, IF, and WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



SARS-CoV-2 (human) Neutralizing Recombinant Antibody ELISA. Rabbit Fc-tagged receptor-binding domain (RBD) subunit 1 (S1 RBD) protein (100 ng/ml) and mouse anti-rabbit Fc-HRP conjugate (200 ng/ml) were incubated with serial dilutions of the SARS-CoV-2 (human) Neutralizing Recombinant Antibody in ACE2-expressing HEK293T/17 cells. Inhibition of S1 RBD binding on the cells by the SARS-CoV-2 (human) Neutralizing Recombinant Antibody was revealed by a reduction in absorbance at 450 nm after the addition of TMB substrate (IC₅₀ = 91.78 ng/ml).

ACE2 Transfected + SARS-CoV-2 (human) Neutralizing Recombinant Antibody ACF2 Transfected



Contro

Inhibition of S1 RBD-ACE2 interaction. Twenty-four hours after seeding HEK293T/17 cells on reverse transfection complexes containing ACE2 DNA (middle and right) or control DNA (left), the cells were incubated with a mixture of rabbit Fc-tagged S1 RBD and DyLight550-conjugated anti-rabbit Fc antibody (left and middle) or with the addition of SARS-CoV-2 (human) Neutralizing Recombinant Antibody (right) for 1 hour at room temperature. Cell nuclei were counterstained with DAPI in blue. The presence of SARS-CoV-2 (human) Neutralizing Recombinant Antibody significantly inhibited the interaction of S1 RBD with ACE2 on the cell surface and abolished the DyLight550-S1 RBD immunostaining in red.

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

SAFFTY 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, 12/12/2023

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) 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 surface glycoprotein, also known as the spike 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 (human) Neutralizing Recombinant Antibody disrupts the S1 RBD-ACE2 interaction and can be used for ELISA, immunofluorescence (IF), and Western blot (WB) applications. The antibody recognizes the SARS-CoV-2 spike glycoprotein at approximately 65 kDa.

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).
- 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).
- 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