

# PRODUCT INFORMATION

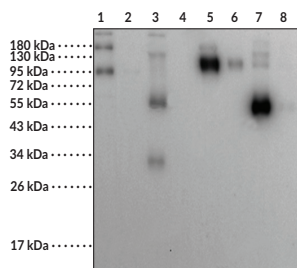
## SARS-CoV-2 Spike Glycoprotein Receptor Binding Domain Polyclonal Antibody

Item No. 31997

### Overview and Properties

<b>Contents:</b>	This vial contains 50 or 100 µl of protein A-affinity purified polyclonal antibody.
<b>Synonyms:</b>	SARS-CoV-2 Spike RBD, SARS-CoV-2 Spike Receptor Binding Domain, Severe Acute Respiratory Syndrome Coronavirus 2 Spike Glycoprotein Receptor Binding Domain, Spike S1 RBD
<b>Immunogen:</b>	Recombinant SARS-CoV-2 spike glycoprotein RBD
<b>Cross Reactivity:</b>	(+) SARS-CoV-2 spike glycoprotein RBD (mouse Fc-tagged), SARS-CoV spike glycoprotein RBD (His-tagged), SARS-CoV-2 spike glycoprotein S1 subunit (His-tagged), SARS-CoV spike glycoprotein S1 subunit (His-tagged)
<b>Species Reactivity:</b>	(+) SARS-CoV-2, SARS-CoV; 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>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Applications:</b>	ELISA and Western blot (WB); the recommended starting dilution is 1:5,000-1:10,000 for ELISA and 1:1,000-1:5,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically

### Image



Lane 1: SARS-CoV Spike Glycoprotein S1 Subunit (His Tag) (30 ng)  
 Lane 2: SARS-CoV Spike Glycoprotein S1 Subunit (His Tag) (5 ng)  
 Lane 3: SARS-CoV Spike Glycoprotein RBD (His Tag) (30 ng)  
 Lane 4: SARS-CoV Spike Glycoprotein RBD (His Tag) (5 ng)  
 Lane 5: SARS-CoV-2 Spike Glycoprotein S1 Subunit (His Tag) (30 ng)  
 Lane 6: SARS-CoV-2 Spike Glycoprotein S1 Subunit (His Tag) (5 ng)  
 Lane 7: SARS-CoV-2 Spike Glycoprotein RBD (mFc Tag) (30 ng)  
 Lane 8: SARS-CoV-2 Spike Glycoprotein RBD (mFc Tag) (5 ng)

WB of SARS-CoV-2 Spike Glycoprotein Receptor Binding Domain Polyclonal Antibody at 1:2,000 dilution.

**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, 03/09/2023

**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.<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> Cayman's SARS-CoV-2 Spike Glycoprotein Receptor Binding Domain Polyclonal Antibody can be used for ELISA and Western blot (WB) applications. The antibody recognizes the spike glycoprotein RBD from SARS-CoV-2 and SARS-CoV.

## 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).