

PRODUCT INFORMATION

SARS-CoV-2 Spike Glycoprotein Receptor Binding Domain (human IgG1 Fc-tagged)

Item No. 30429

Overview and Properties

Synonym: SARS-CoV-2 Surface Receptor Binding Domain, Severe Acute Respiratory Syndrome Coronavirus 2 Spike Glycoprotein Receptor Binding Domain

Amino Acids: 319-541

Molecular Weight: 52.3 kDa

Storage: -80°C (as supplied)

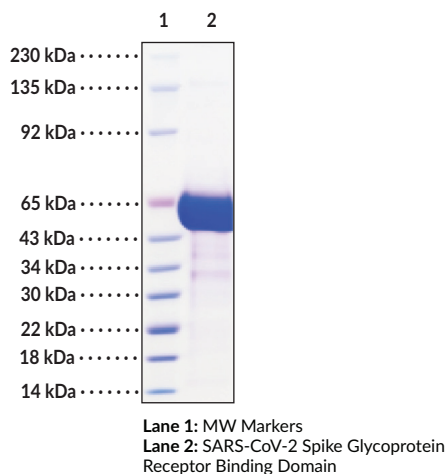
Stability: ≥1 year

Purity: ≥90% estimated by SDS-PAGE

Supplied in: 1 mg/ml in a solution of PBS, pH 7.2

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



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, 03/09/2021

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) surface glycoprotein receptor binding domain is a fragment of the SARS-CoV-2 surface glycoprotein, also known as the spike glycoprotein, which is a viral structural protein encoded by the S gene in SARS-CoV-2 RNA.¹ It contains amino acids 319-541 of the full-length SARS-CoV-2 spike glycoprotein sequence. SARS-CoV-2 is a member of the *Betacoronavirus* genus of viruses and has 88% sequence identity with two bat-derived SARS-like CoVs.² The SARS-CoV-2 genome contains approximately 30 kilobases that encode four structural proteins: spike, envelope, membrane, and nucleocapsid.^{1,3} The spike protein of SARS-CoV-2 and the related viruses SARS-CoV and Middle East respiratory syndrome coronavirus (MERS-CoV) is a transmembrane glycoprotein that assembles into homotrimers on the virus surface and is comprised 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.⁴⁻⁶ The 193-amino acid RBD of the SARS-CoV spike protein is a target for neutralizing antibodies.^{5,7} The SARS-CoV-2 RBD, which spans amino acid residues 329 to 521, is 73% identical to that of SARS-CoV 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.^{4,6,7} SARS-CoV-2 is the causative agent of COVID-19, a primarily respiratory illness characterized by fever, cough, and shortness of breath that can lead to life-threatening complications.⁸⁻¹⁰ Cayman's SARS-CoV-2 Surface Glycoprotein Receptor Binding Domain protein can be used as an antigen or for Western blot, ELISA, protein-protein interaction studies, and other *in vitro* binding and *in vivo* functional assays

References

1. Kandeel, M., Ibrahim, A., Fayed, M., et al. *J. Med. Virol.* (2020).
2. Lu, R., Zhao, X., Li, J., et al. *Lancet* **395**(10224), 565-574 (2020).
3. Ahmed, S.F., Quadeer, A.A., and McKay, M.R. *Viruses* **12**(3), E254 (2020).
4. Liu, Z., Xiao, X., Wei, X., et al. *J. Med. Virol.* (2020).
5. He, Y., Zhou, Y., Liu, S., et al. *Biochem. Biophys. Res. Commun.* **324**(2), 773-781 (2004).
6. Walls, A.C., Park, Y.-J., Tortorici, M.A., et al. *Cell* **180**, 1-12 (2020).
7. Tian, X., Li, C., Huang, A., et al. *Emerg. Microbes Infect.* **9**(1), 382-385 (2020).
8. Meo, S.A., Alhowikan, A.M., Al-Khlaiwi, T., et al. *Eur. Rev. Med. Pharmacol. Sci.* **24**(4), 2012-2019 (2020).
9. Klok, F.A., Kruij, M.J.H.A., van der Meer, N.J.M., et al. *Thromb. Res.* **S0049-3848**(20), 30120-1 (2020).
10. Yang, F., Shi, S., Zhu, J., et al. *J. Med. Virol.* (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