

PRODUCT INFORMATION

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

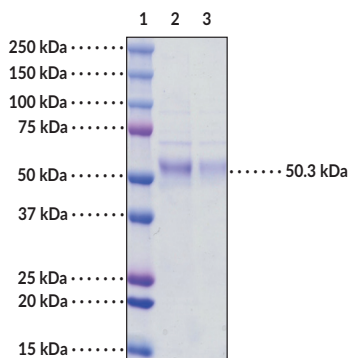
Item No. 30590

Overview and Properties

| | |
|-------------------------------|---|
| Synonyms: | SARS-CoV-2 Spike Receptor Binding Domain, S1 RBD, Severe Acute Respiratory Syndrome Coronavirus 2 Spike Glycoprotein Receptor Binding Domain, Spike S1 RBD |
| Source: | Active recombinant C-terminal rabbit IgG1 Fc-tagged SARS-CoV-2 surface glycoprotein receptor binding domain expressed in HEK293 cells |
| Amino Acids: | 319-541 |
| Uniprot No.: | PODTC2 |
| Molecular Weight: | 50.3 kDa |
| Storage: | -80°C (as supplied) |
| Stability: | ≥6 months |
| Purity: | ≥85% |
| Supplied in: | PBS, pH 7.4, with 5% mannitol, 5% trehalose, 0.01% Tween 20, and 10% glycerol |
| Protein Concentration: | <i>batch specific</i> mg/ml |
| Bioactivity: | SARS-CoV-2 Spike Glycoprotein Receptor Binding Domain was captured on a Protein G Chip S series and tested for binding with gradient concentrations of ACE2 (12.5, 25, 50, 100, and 200 nM) in 10 mM HEPES pH 7.4, 150 mM NaCl, 3 mM EDTA, 0.05% surfactant P20 at 25°C. The K_D value was calculated using the 1:1 (Langmuir) binding model. |

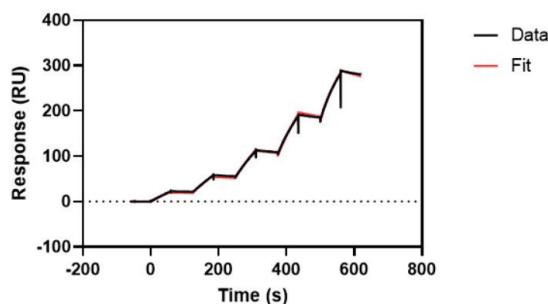
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: Marker band
Lane 2: SARS-CoV-2 spike glycoprotein (4 µg)
Lane 3: SARS-CoV-2 spike glycoprotein (2 µg)

SDS-PAGE Analysis of SARS-CoV-2 Spike Glycoprotein Receptor Binding Domain (rabbit IgG1 Fc-tagged).



SARS-CoV-2 Spike Glycoprotein Receptor Binding Domain Specifically Binds ACE2
SARS-CoV-2 Spike Glycoprotein Receptor Binding Domain was captured on a Protein G Chip S series and SPR analysis was used to determine ACE2 (human, recombinant; Item No. 30587) binding affinity on a Biacore T200, using single cycle kinetics with five concentrations of ACE2.

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, 03/18/2021

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, also known as the spike glycoprotein, is a viral structural protein encoded by the S gene in SARS-CoV-2 RNA.¹ 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 the COVID-19 respiratory illness characterized by fever, cough, and shortness of breath that can progress to pneumonia and potentially death in high-risk populations.⁸ Cayman's SARS-CoV-2 Spike Glycoprotein Receptor Binding Domain (rabbit IgG1 Fc-tagged) protein can be used for ELISA, surface plasmon resonance (SPR), and Western blot (WB) applications.

References

1. Kandeel, M., Ibrahim, A., Fayed, 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. Ahmed, S.F., Quadeer, A.A., and McKay, M.R. Preliminary identification of potential vaccine targets for the COVID-19 coronavirus (SARS-CoV-2) based on SARS-CoV immunological studies. *Viruses* **12(3)**, E254 (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).
8. 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).