

# PRODUCT INFORMATION



## SARS-CoV-2 Spike Glycoprotein Receptor Binding Domain K417N, E484K, N501Y variant (rabbit IgG1 Fc-tagged)

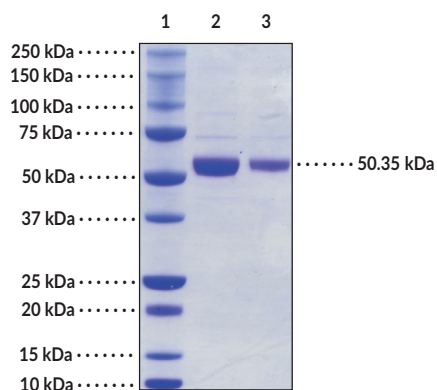
Item No. 33868

### Overview and Properties

<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, South African Variant, Spike S1 RBD, WHO Label: Beta
<b>Source:</b>	Active recombinant C-terminal rabbit IgG1 Fc-tagged SARS-CoV-2 Spike Glycoprotein Receptor Binding Domain K417N, E484K, N501Y variant expressed in HEK293 cells
<b>Amino Acids:</b>	21-243, 319-541 of P0DTC2
<b>Uniprot No.:</b>	P0DTC2
<b>Molecular Weight:</b>	50.35 kDa
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Purity:</b>	<i>batch specific</i> (≥90% estimated by SDS-PAGE)
<b>Supplied in:</b>	PBS, pH 7.4, with 5% mannitol, 5% trehalose, 0.01% polysorbate 20, and 10% glycerol
<b>Protein Concentration:</b>	<i>batch specific</i> mg/ml
<b>Bioactivity:</b>	See figure for details

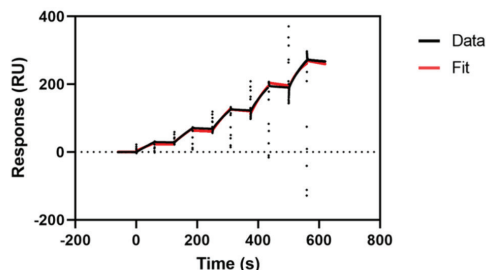
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Images



Lane 1: MW Markers  
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 K417N, E484K, N501Y variant (rabbit IgG1 Fc-tagged).



SARS-CoV-2 Spike Glycoprotein Receptor Binding Domain K417N, E484K, N501Y variant (rabbit IgG1 Fc-tagged) was captured on a Protein G Chip 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, 09/17/2021

**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) 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> In this way, ACE2 acts as the functional receptor for SARS-CoV-2. The SARS-CoV-2 variant of concern (VOC) B.1.351, originally identified in South Africa, contains three substitutions located within the spike glycoprotein RBD.<sup>11,12</sup> The lysine-to-asparagine substitution at position 417 (K417N) reduces SARS-CoV-2 affinity for human ACE2, whereas the glutamate-to-lysine substitution at position 484 (E484K) and the asparagine-to-tyrosine substitution at position 501 (N501Y) increase SARS-CoV-2 affinity for human ACE2.<sup>11</sup> Collectively, the K417N, E484K, and N501Y substitutions are associated with enhanced viral infectivity and resistance to antibody-mediated neutralization.<sup>12</sup> Cayman's SARS-CoV-2 Spike Glycoprotein Receptor Binding Domain K417N, E484K, N501Y variant (rabbit IgG1 Fc-tagged) can be used for ELISA, surface plasmon resonance (SPR), and Western blot 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. 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).
11. Boehm, E., Kronig, I., Neher, R.A., *et al.* Novel SARS-CoV-2 variants: The pandemics within the pandemic. *Clin. Microbiol. Infect.* (2021).
12. Kuzmina, A., Khalaila, Y., Voloshin, O., *et al.* SARS-CoV-2 spike variants exhibit differential infectivity and neutralization resistance to convalescent or post-vaccination sera. *Cell Host Microbe* **29(4)**, 522-528e2 (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