

# Spike-S1-Fc

Soluble SARS-CoV-2 Spike S1 protein fused to a human IgG1 Fc tag

Catalog code: fc-sars2-s1

<https://www.invivogen.com/sars2-spike-s1-proteins>

For research use only, not for diagnostic or therapeutic use

Version 20123-NJ

## PRODUCT INFORMATION

### Contents:

- 50 µg of lyophilized Spike-S1-Fc protein
- 1.5 ml endotoxin-free water

### Protein construction:

Codon-optimized spike glycoprotein S1 domain [V16-R685] with a C-terminal human IgG1 Fc tag

**Accession sequence:** YP\_009724390 (native sequence)

**Species:** SARS-CoV-2 (2019-nCoV); Wuhan-Hu-1 (D614) isolate

**Tag:** C-terminal human IgG1 Fc

**Total protein size:** 920 a.a. (secreted form)

**Molecular weight:** ~124 kDa

**Purification:** Protein G affinity chromatography

**Purity:** >95% (SDS-PAGE)

### Formulation:

0.2 µm filtered solution in a sodium phosphate buffer with glycine, saccharose, and stabilizing agents

### Storage:

- Product is shipped at room temperature. Store lyophilized product at -20°C. Lyophilized product is stable for at least 1 year.
- Reconstituted protein is stable for 1 month when stored at 4°C and for 1 year when aliquoted and stored at -20°C. Avoid repeated freeze-thaw cycles.

### Quality control:

- The size and purity of the protein has been confirmed by SDS-PAGE.
- Spike-S1-Fc has been functionally validated by ELISA using the Anti-SARS-CoV-Spike human IgM (clone CR3022).
- Absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and TLR4 cellular assays.

## BACKGROUND

The SARS-CoV-2 Spike S1 subunit plays a crucial role in the viral entry into the target cell. The S1 subunit features an N-term S1-NTD region and a C-term S1-CTD region. While S1-NTD is thought to mediate sugar-binding, the S1-CTD allows the virus binding to ACE2 through the receptor-binding domain (RBD)<sup>1-3</sup>. In its resting conformation, S1 exerts a physical constraint on the Spike fusion subunit<sup>3</sup>. Research is ongoing to understand the exact mechanisms that drive conformation changes in S1 allowing subsequent membrane fusion events. S1 is a candidate for subunit vaccines against SARS-CoVs<sup>4,5</sup>.

## PRODUCT DESCRIPTION

Spike-S1-Fc is a soluble SARS-CoV-2 protein generated by fusing the Spike S1 domain [V16-R685] to a C-terminal human IgG1 Fc tag with a TEV (Tobacco Etch Virus) sequence linker. This fusion protein has a molecular weight of ~124 kDa on an SDS-PAGE gel. Spike-S1-Fc has been generated by recombinant DNA technology, produced in CHO cells, and purified by protein G affinity chromatography.

## APPLICATIONS

- **Vaccination studies:** using combinations of Spike protein antigens and adjuvants.
- **Antibody screening:** finding anti-Spike antibodies that can neutralize the SARS-CoV-2 infection.
- **Inhibitor screening:** finding small molecules, or antibodies able to block the SARS-CoV-2 RBD interaction with the ACE2 receptor.
- **ACE2 cellular expression screening:** in primary isolated cells or transfected cells.

## METHODS

### Spike-S1-Fc resuspension (100 µg/ml)

*Note: Ensure you see the lyophilized pellet before resuspension.*

- Add 500 µl of endotoxin-free water to the vial and gently pipette until completely resuspended.
- Prepare aliquots and store at -20°C or 4°C.

## TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)

## PROTEIN SEQUENCE

MEIKVLFALICIAVAEAKPTELEVNLTTRTQLPPAY  
TNSFTRGVVYYPDKVFRSSVLHSTQDLFLPFFSNVT  
WFHAIHVS GTNGTKRFDNPVLPFNDGVYFASTEK  
SNIIRGWIFGTTLDLSDKTQSLIVNNATNVVIVKVEF  
QFCNDPFLGVYYHKNNKSWMESEFRVYSSANNCT  
FEYVVSQPFLMDLEGKQGNFKNLREFVFKNIDGYF  
KIYSKHTPINLVRDLPPQGFSALEPLVDLPIGINITR  
FQTLALLHRSYLT PGDSSSGWTAGAAAYVGYLQP  
RTFLLKY NENGTITDAVDCALDPLSETKCTLKSFT  
VEKGIYQTSNFRVQPTESIVRFPNITNLCPFGEVF  
NATRFASVYAWNRKRISNCVADYSVLYNSASFSTF  
KCYGVSPTKLNLDLCFTNVYADSFVIRGDEVQRQIAP  
GQTGKIADYNYKLPDDFTGCVIAWNSNNLDSKVG  
GNYNYLYRFLRKS NLKPFERDISTEIQAGSTPCN  
GVEGFNCYFPLQSYGFPQTNGVGYQPYRVVLSF  
ELLHAPATVCGPKKSTNLVKNKCVNFNFNGLTGT  
GVLTESNKKFLPFQGFGRDIADTTDAVRDPQTLEI  
LDITPCSFGGVSVITPGTNTSNQVAVLYQDVNCTE  
VPVAIHADQLTPTWRVYSTGSNVFQTRAGCLIGAE  
HVNNSYECDIPIGAGICASYQTQTN SPRRARRTEN  
LYFQGS GSEPKSSDKTHTCPPCPAPEAEGGSPVFL  
FPPKPKDQLMISRTPEVTCVVVDVSHEDPEVKFN  
WYVDGVEVHNAKTKPREEQYNSTYRVVSVLTVLH  
QDWLNGKEYKCKVSNKALPASIEKTISKAKGQPR  
EPQVYTLPPSREEMTKNQVSLTCLVKGFYPSDIAV  
EWESNGQPENNYKTTTPVLDSDGSFFLYSKLTVD  
KSRWQQGNV FSCSVLHEALHNHYTQKLSLSLSPGK

Green: signal sequence

Purple: stabilizing amino acid sequence

Blue: Spike S1 sequence

Black: TEV cleavage sequence

Red: Human IgG1 Fc sequence

## REFERENCES

1. Li F., 2016. Structure, function, and evolution of coronavirus spike proteins. *Annu. Rev. Virol.* 3:237-261. 2. Li F. *et al.*, 2005. Structure of SARS coronavirus spike receptor-binding domain complexed with receptor. *Science*. 309:1864-1868. 3. Walls A.C. *et al.*, 2020. Structure, function, and antigenicity of the SARS-CoV-2 spike glycoprotein. *Cell*. 181(2):281-292.e6. 4. Wang N. *et al.*, 2020. Subunit vaccines against emerging pathogenic human coronaviruses. *Front. Microbiol.* 11:298. DOI: 10.3389/fmicb.2020.00298. 5. Padron-Regalado E., 2020. Vaccines for SARS-CoV-2: Lessons from other coronavirus strains. *Infect. Dis. Ther.* DOI: 10.1007/s40121-020-00300-x.

## RELATED PRODUCTS

Product	Catalog Code
Anti-Spike-RBD-hIgG1	srbd-mab1
Anti-Spike-RBD-hIgM	srbd-mab5
Anti-Spike-RBD-hIgA2	srbd-mab6
hACE2-Fc	fc-hace2
Spike-S1-His	his-sars2-s1
Spike-RBD-His	his-sars2-srbd
Spike-RBD-Fc	fc-sars2-srbd
Nucleocapsid-His	his-sars2-n
Nucleocapsid-Fc	fc-sars2-n
pDUO2-hACE2-TMPRSS2a	pduo2-hace2tpsa

### TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)