

Spike-S1-His

Soluble SARS-CoV-2 Spike S1 protein fused to a poly-histidine tag

Catalog code: his-sars2-s1

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

For research use only, not for diagnostic or therapeutic use

Version 20I23-NJ

PRODUCT INFORMATION

Contents:

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

Protein construction:

Codon-optimized spike glycoprotein S1 domain [V16-R685] with a C-terminal poly-histidine tag

Accession sequence: YP_009724390 (native sequence)

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

Tag: C-terminal poly-histidine (6 x His)

Total protein size: 681 a.a. (secreted form)

Molecular weight: ~123 kDa

Purification: Ni²⁺ 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-His 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)¹⁻³. In its resting conformation, S1 exerts a physical constraint on the Spike fusion subunit³. 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^{4,5}.

PRODUCT DESCRIPTION

Spike-S1-His is a soluble protein generated by fusing the SARS-CoV-2 Spike S1 domain [V16-R685] to a C-terminal poly-histidine (6 x Histidine) tag with a tri-amino acid linker. This fusion protein has a molecular weight of ~123 kDa on an SDS PAGE gel. Spike-S1-His has been generated by recombinant DNA technology, produced in HEK293 cells, and purified by Ni²⁺ 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-His 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

PROTEIN SEQUENCE

MEIKVLFALICIAVAEAL **E**VNLTTRTQLPPAYTNSF
TRGVVYYPDKVFRSSVLHSTQDLFLPFFSNVTWFHA
IHVSGTNGTKRFDNPVLPFNDGVYFASTEKSNIIIR
GWIFGTTLD SKTQSL LIVNNATNVVIK VCEFQFCN
DPFLGVYYHKNNKSWMESEFRVYSSANNCTFEYV
SQPF LMDLEGKQGNFKNLREFVFKNIDGYFKIYSK
HTPINLV RDLPQGFSALEPLVDLP IGINITRFQTL L
ALHRSYLT PGDSSSGW TAGAAAYVGYLQPRTFLL
KYNENGTITDAVDCALDPLSETKCTLSFTVEKGI
YQTSNFRVQPTESIVRFPNITNLCPFGGEVFNATRFA
SVYAWNRKRISNCVADYSVLYNSASFSTFKCYGVS
PTKLN DL CFTNVYADSFVIRGDEV RQIAPGQTGKI
ADYNYKLPDDFTGCVIAWNSNNLDSKVGGNYNYL
YRLF RKSNLKP FERDISTEIQAGSTPCNGVEGFN
CYFPLQSYGFQPTNGVGYQPYRVVLSFELLHAPA
TVCGP KKSTNLVKNKCVNFNFNGLTGTGVLTESN
KKFLPFQQFGRDIADTTDAVRDPQTLEILDITPCS
FGGVS VITPGTNTSNQVAVLYQDVNCTEVPVAIHA
DQLTPTWRVYSTGSNVFQTRAGCLIGAEHVNNSY
ECDIPIGAGICASYQTQ TNSPRRAR GSG **H H H H H H**

Green: signal sequence

Purple: stabilizing amino acid sequence

Blue: Spike S1 sequence

Black: tri-amino acid short linker

Red: 6 x Histidine

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-Fc	fc-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