

Nucleocapsid-Fc

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

Catalog code: fc-sars2-n

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

For research use only, not for diagnostic or therapeutic use

Version 2019-NJ

PRODUCT INFORMATION

Contents:

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

Protein construction:

Full-length Nucleocapsid [M1-A419] with a human IgG1 Fc tag in C-terminus

Accession sequence: PODTC9

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

Tag: C-terminal human IgG1 Fc

Total protein size: 669 a.a. (secreted form)

Molecular weight: ~79 kDa (SDS-PAGE)

Purification: Protein A 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.
- Nucleocapsid-Fc has been functionally validated by ELISA using an Anti-SARS Nucleocapsid antibody.
- 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 Nucleocapsid (N) is an important structural protein that plays important roles in the viral life cycle including replication, transcription, and genome packaging¹. The SARS-CoV-2 N features two important NTD and CTD functional domains in N-terminal and C-terminal, respectively¹⁻⁶. NTD interacts with both the RNA genome and M proteins to form virion particles. The N protein interaction with the RNA forms the virus ribonucleoprotein core which is packed as a helical “beads-on-a-string” conformation. CTD allows RNA synthesis through binding of the replication-transcription complexes (RTCs), oligomerization of multiple N proteins through its dimerization domain, and genome incorporation into the new virion. N is a major immunogen of SARS-CoV-2. Indeed, elevated Anti-SARS-CoV-2 N IgG and IgM antibody titers have been reported in COVID-19 patients’ sera⁷⁻⁹. These observations make SARS-CoV-2 N an attractive tool for early diagnosis⁷⁻⁹ and a potential therapeutic drug-target³.

PRODUCT DESCRIPTION

Nucleocapsid-Fc is a soluble protein generated by fusing the full-length SARS-CoV-2 nucleocapsid [M1-A419] to a C-terminal human IgG1 Fc tag with a TEV (Tobacco Etch Virus) sequence linker. This fusion protein has a molecular weight of ~79 kDa on a SDS PAGE gel. Nucleocapsid-Fc has been generated by recombinant DNA technology, produced in HEK293 cells, and purified by protein G affinity chromatography.

APPLICATIONS

- **Vaccination studies:** using combinations of Nucleocapsid protein antigens and adjuvants.
- **Antibody screening:** finding anti-Nucleocapsid antibodies in COVID-19 patients’ sera.
- **Inhibitor screening:** finding small molecules able to block the SARS-CoV-2 nucleocapsid interaction with replication-transcription complexes (RTCs).

METHODS

Nucleocapsid-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

PROTEIN SEQUENCE

MEIKVLFALICIAVAEAKPTELEMSDNGPQNQRNA
PRITFGGSPSDSTGSNQNNGERSGARSKQRRPQGLP
NNTASWFTALTQHGKEDLKFRGGQGVPIINTNSSP
DDQIGYYRRATRRIIRGGDGKMKDLSRPRWYFYLLG
TGPEAGLPYGANKDGIIVVATEGALNTPKDHIGT
RNPANNAAIVLQLPQGTTLPKGFYAEGSRGGSQA
SSRSSSRSRNSSRNS TPGSSRGTSPARMAGNGGD
AALALLLLDRLNQLLESKMSGKQQQQGQTVTKKS
AAEASKKPRQKRATKAYNVTQAFGRRGPEQTQG
NFGDQELIRQGTDYKHWPQIAQFAPSASAFFGMS
RIGMEVTPSGTWLTYTGAIKLDDKDPNFKDQVILL
NKHIDAYKTFFPTEPKKDKKKKADETTALPQRQK
KQQTVTLLPAADLDDFSKQLQQSMSSADSTQART
ENLYFQGSSEPKSSDKTHTCPPCPAPEAEGGPSV
FLFPPKPKDQLMISRTPEVTCVVVDVSHEDPEVKF
NWyVDGVEVHNAKTKPREEQYNSTYRVVSVLTVL
HQDWLNGKEYKCKVSNKALPASIIEKTISKAKGQP
REPQVYTLPPSREEMTKNQVSLTCLVKGFYPSDIA
VEWESNGQPENNYKTTPPVLDSDGSFFLYSKLTVD
KSRWQQGNVFCSSVLEALHNHYTQKSLSLSPGK

Green: signal sequence

Purple: stabilizing amino acid sequence

Blue: Nucleocapsid sequence

Black: TEV cleavage sequence

Red: human IgG1 Fc sequence

REFERENCES

1. Mu, J. *et al.*, 2020. SARS-CoV-2-encoded nucleocapsid protein acts as a viral suppressor of RNA interference in cells. *Sci China Life Sci* 63, 1-4. 2. Chang C. *et al.*, 2006. Modular organization of SARS coronavirus nucleocapsid protein. *J. Biom. Sci.* 13:59-72. 3. Krokhn O. *et al.*, 2003. Mass spectrometric characterization of proteins from the SARS virus. *Mol. & Cell. Prot.* 2:346-356. 4. Cubuk, J. *et al.*, 2020. The SARS-CoV-2 nucleocapsid protein is dynamic, disordered, and phase separates with RNA. *bioRxiv*. doi:10.1101/2020.06.17.158121. 5. Kang, S. *et al.*, 2020. Crystal structure of SARS-CoV-2 nucleocapsid protein RNA binding domain reveals potential unique drug targeting sites. *Acta Pharm Sin B*. doi:10.1016/j.apsb.2020.04.009. 6. Khan, M.T. *et al.*, 2020. SARS-CoV-2 nucleocapsid and Nsp3 binding: an in silico study. *Arch Microbiol*. doi: 10.1007/s00203-020-01998-6. 7. Liu, W. *et al.*, 2020. Evaluation of Nucleocapsid and Spike Protein-Based Enzyme-Linked Immunosorbent Assays for Detecting Antibodies against SARS-CoV-2. *J Clin Microbiol* 58. 8. Guo L. *et al.*, 2020. Profiling Early Humoral Response to Diagnose Novel Coronavirus Disease (COVID-19). *Clinical Infectious Diseases*. 71(15) :778-785. 9. To K. K-W. *et al.*, 2020. Temporal profiles of viral load in posterior oropharyngeal saliva samples and serum antibody responses during infection by SARS-CoV-2: an observational cohort study. *The Lancet Infectious Diseases*. 20(5):565-574.

RELATED PRODUCTS

Product	Catalog Code
Nucleocapsid-His	his-sars2-n
Spike-S1-Fc	fc-sars2-s1
Spike-S1-His	his-sars2-s1
Spike-RBD-Fc	fc-sars2-srbd
Spike-RBD-His	his-sars2-srbd

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