

Validation data for Nucleocapsid-Fc

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

For research use only

Version 20I22-NJ

Nucleocapsid-Fc is a soluble SARS-CoV-2 fusion protein generated by fusing the full-length 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 (Figure 1). The recognition of the Nucleocapsid-His protein by an Anti-SARS-CoV-Nucleocapsid antibody has been verified by ELISA (Figure 2).

Nucleocapsid-Fc purity analysis by SDS-PAGE

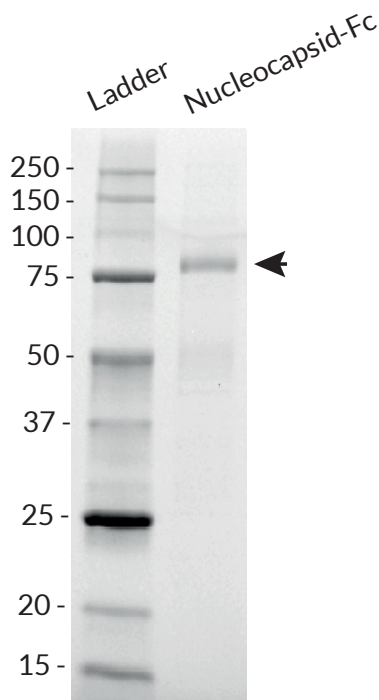


Figure 1: SDS-PAGE analysis of the SARS-CoV-2 Nucleocapsid-Fc protein. 2 µg of the fusion protein was loaded onto a 12% Mini-PROTEAN® TGX Stain-Free™ Precast Gel (Bio-Rad). Detection was performed as per manufacturer's instructions.

Recognition of Nucleocapsid-Fc by an Anti-SARS-CoV-Nucleocapsid mouse IgG1

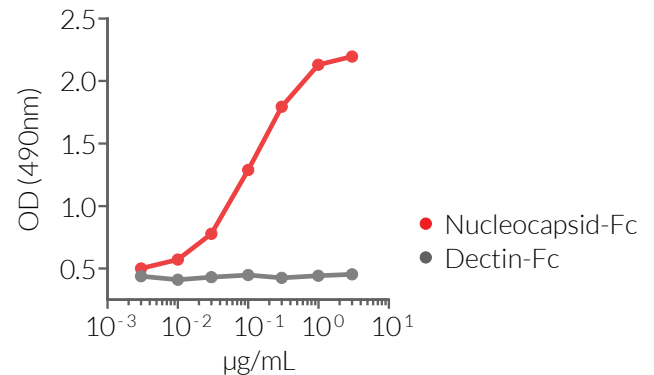


Figure 2: ELISA detection of Nucleocapsid-Fc fusion protein with an Anti-SARS-CoV-Nucleocapsid mAb. Anti-SARS-CoV-Nucleocapsid antibody (2 µg/ml) was coated onto ELISA plates overnight. Following this, a 3-fold serial dilution of Nucleocapsid-Fc (red curve) or control protein (Dectin-Fc; grey curve) were added and incubated for 1 hour. Binding was detected using a HRP-labelled anti-His antibody (1/1000 dilution) and the HRP substrate OPD (o-phenylenediamine dihydrochloride). Absorbance was read at 490 nm.

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 3-622-34-80
E-mail: info@invivogen.com