

Coronavirus (SARS-CoV NP) antigen (full length), (recombinant) (His-tag)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Formulation	Lyophilized. Sterile filtered from 0.5 mg/ml in 0.05M Acetate buffer pH 4.0.
MW	47.8 kDa
Purity	≥ 95%
Purity Detail	Purified.
Reconstitution	Add 0.2 ml of 0.1M Acetate buffer pH4.0 and let the lyophilized pellet dissolve completely. For conversion into higher pH value, we recommend intensive dilution by relevant buffer to a concentration of 10µg/ml. In higher concentrations the solubility of this antigen is limited.
Source	Produced in <i>E. coli</i> . Recombinant protein containing 422 amino acid residues of the SARS-CoV Nucleocapsid protein and 15 additional amino acid residues.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com