

SARS-CoV-2 – nucleocapsid protein – 48.3 kDa (L pack)

Ordering Information

[Order Online »](#)

JBS-PR-1454-L	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name SARS-CoV-2 – coronavirus nucleocapsid protein

Formulation liquid (Supplied in Sodium Phosphate 12.5 mM, pH 8.0, Tris-HCl 20 mM, Sodium Chlorid 300 mM, Imidasol 62.5 mM, Sarcosyl 0.5 %)

MW about 48.3 kDa

Source recombinant, E. coli

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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