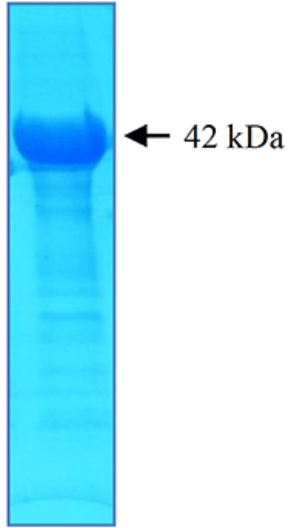


Coronavirus (hCoV-229E NP) antigen, (recombinant) (His-tag)

Manuals, SDS & CofA

[View Online »](#)



Coomassie-stained SDS-PAGE showing purified recombinant HCoV-229E nucleoprotein.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Formulation Liquid. In PBS with 3M Urea.

Purity ≥ 95%

Purity Detail Purified.

Source Produced in *E. coli*. Recombinant protein (359 aa).

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