

N. gonorrhoeae native antigen

Bacterial lysate ideal for use as a broad positive control

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Sonicate/homogenize briefly before use.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Neisseria gonorrhoeae native antigen
Appearance	Cloudy liquid.
Formulation	Liquid. In PBS containing 0.09% sodium azide.
Purity	≥99%
Purity Detail	Homogenized lysate of pure bacteria.
Source	From <i>Neisseria gonorrhoeae</i> , strain ATCC 19424. Inactivated by 3 hours at 56°C.

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