

NLRP1/NALP1 (human) (NT) polyclonal antibody (Bur 241)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-018-R050	50µl
------------------	------

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

Alternative Name Caspase recruitment domain-containing protein 7, NAC, NACHT-, LRR- and PYD-containing protein 1, DEFCAP, CARD7

Application IHC (PS), IP, WB

Crossreactivity Does not cross-react with mouse NLRP1/NALP1.

Formulation Liquid. Antiserum containing PBS and 0.05% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 161-180 (P161SSPDHESPSQESPNAPTST180) of N-terminal human NLRP1/NALP1 (NACHT-, LRR- and PYD-containing protein 1).

Species Reactivity Human

UniProt ID Q9C000

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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