

Purinergic receptor P2X7 (mouse) monoclonal antibody (HANO43)

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

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-802-027-L001	1ml
------------------	-----

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 P2X purinoceptor 7, P2RX7

Application Flow Cytometry, ICC, IP

Clone HANO43

Formulation Liquid. Hybridoma supernatant containing 0.02% sodium azide.

Host Rat

Immunogen Vector containing the cDNA of mouse purinergic receptor P2X7.

Isotype IgG2b

Positive Control HEK cells transfected with an expression plasmid for mouse purinergic receptor P2X7 can be used as positive control.

Recommendation Dilutions/Conditions Flow Cytometry (1:1-1:10)Immunocytochemistry (1:1-1:10)Immunoprecipitation (1:1-1:10)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Mouse

UniProt ID Q9Z1M0

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



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-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@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)