

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-
be@enzolifesciences.com](mailto:info-be@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)