

Purinergic receptor P2X7 (mouse) monoclonal antibody (HANO43)

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

Citations: 9

[View Online »](#)

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

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-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)