

Purinergic receptor P2X4 (mouse) polyclonal antibody

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

Citations: 2

[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 4, P2RX4
Application	Flow Cytometry, ICC, IP
Formulation	Liquid. Antiserum diluted with PBS, containing 0.1% BSA and 0.02% sodium azide.
Host	Rabbit
Immunogen	Vector containing the cDNA of mouse purinergic receptor P2X4.
Positive Control	HEK cells transfected with an expression plasmid for mouse purinergic receptor P2X4 can be used as positive control.
Recommendation Dilutions/Conditions	Flow Cytometry (1:100-1:400)Immunocytochemistry (1:100-1:400)Immunoprecipitation (1:100-1:400)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Mouse
UniProt ID	Q9JJX6
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

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