

Purinergic receptor P2X4 (mouse) polyclonal antibody

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

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|------------------|-------|
| ALX-215-033-R100 | 100µl |
|------------------|-------|

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](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)