

# CD39L1 (human) polyclonal antibody

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

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

|                  |       |
|------------------|-------|
| ALX-215-045-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

|  |  |
|--|--|
| <b>Alternative Name</b>                    | Ectonucleoside triphosphate diphosphohydrolase 2, NTPDase 2, ENTPD2  |
| <b>Application</b>                         | Flow Cytometry, ICC  |
| <b>Formulation</b>                         | Liquid. Antiserum diluted with PBS containing 0.1% BSA and 0.02% sodium azide.   |
| <b>Host</b>                                | Rabbit   |
| <b>Immunogen</b>                           | Vector containing the cDNA of human CD39L1.  |
| <b>Positive Control</b>                    | HEK cells transfected with an expression plasmid for human CD39L1 can be used as positive control.   |
| <b>Recommendation Dilutions/Conditions</b> | Flow Cytometry (1:100-1:400)Immunocytochemistry (1:100-1:400)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application. |
| <b>Species Reactivity</b>                  | Human  |
| <b>UniProt ID</b>                          | Q9Y5L3   |
| <b>Worry-free Guarantee</b>                | This antibody is covered by our <a href="#">Worry-Free Guarantee</a>   |

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

France  
Phone: +33 472 440 655  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)