CD39L1 (mouse) polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ALX-215-044-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 Ectonucleoside triphosphate diphosphohydrolase 2,

NTPDase 2, ENTPD2

Application Flow Cytometry, ICC

Formulation Liquid. Antiserum diluted with PBS containing 0.1% BSA

and 0.02% sodium azide.

Host Rabbit

Immunogen Vector containing the cDNA of mouse CD39L1.

Positive Control HEK cells transfected with an expression plasmid

for mouse CD39L1 can be used as positive control.

Recommendation Dilutions/Conditions 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.

Species Reactivity Mouse

UniProt ID O55026

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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com