

CD39 (human) monoclonal antibody (AC2.5) (Alexa fluor[®] 647)

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Short Term Storage -20°C

Long Term Storage -80°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Ectonucleoside triphosphate diphosphohydrolase 1, NTPDase 1, ENTPD1
Application	Flow Cytometry
Clone	AC2.5
Formulation	Liquid. In PBS containing 50% glycerol.
Host	Mouse
Immunogen	EBV-transformed human B lymphoblastoid cell line.
Isotype	IgG1
Purity Detail	Purified using QMA Hyper D ion exchange chromatography
Recommendation Dilutions/Conditions	Flow Cytometry (surface)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P49961

Worry-free Guarantee

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



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-fr@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

Last modified: May 29, 2024