

LAMP1 monoclonal antibody (Ly1C6)

Lysosome-associated membrane proteins (LAMP1 and LAMP2) are type I integral membrane protein that are transported from the trans-Golgi network to endosomes and then lysosomes. Following activation of platelets, T-cells, neutrophils, or endothelium, LAMP1 and LAMP2 are expressed on the cell surface. Cell surface LAMP1 and LAMP2 have been shown to promote adhesion of human peripheral blood mononuclear cells (PBMC) to vascular endothelium, and loss of LAMP2 expression is associated with impaired autophagy observed in Danons disease.

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

Citations: 40

[View Online »](#)

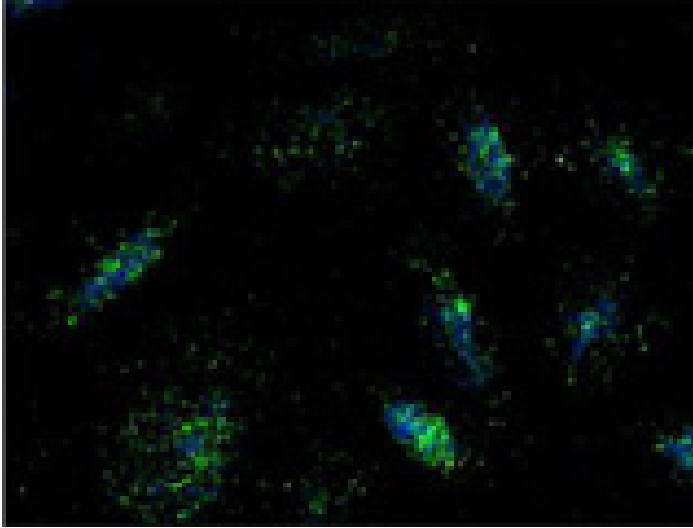
Ordering Information

[Order Online »](#)

ADI-VAM-EN001-D	50µg
ADI-VAM-EN001-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Immunocytochemistry analysis of CHO K1 cells incubated with LAMP1, mAb (Ly1C6) at 1:100 and goat anti-mouse Alexa 488 secondary antibody (Image provided by Dr. Guy Groblewski, University of Wisconsin-Madison).

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Lysosomal-associated membrane protein 1
-------------------------	---

Application	Electron microscopy, ICC, IF, IP, WB
--------------------	--------------------------------------

Application Notes	Detects a band of ~120kDa by Western blot.
--------------------------	--

Clone	LY1C6
-------	-------

Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
--------------------	--

GenBank ID M34959

Host _____ Mouse _____

Immunogen	Rat liver lysosomal membrane preparation.
------------------	---

Isotype IgG1

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions	Immunocytochemistry (1:100)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
--	--

Source	Purified from ascites.
---------------	------------------------

Species Reactivity	Hamster, Human, Rat
---------------------------	---------------------

UniProt ID P14562

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