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: 39

View Online »

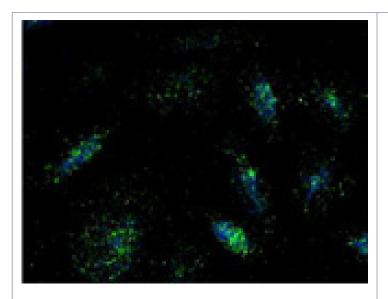
Ordering Information

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 & Storage

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 NotesDetects 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.

lsotype 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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com