# LAMP2 monoclonal antibody (GL2A7)

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: 3

**View Online »** 

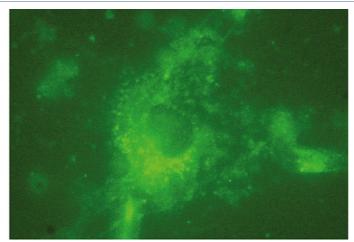
# **Ordering Information**

**Order Online** »

ADI-VAM-EN003-D	50µl
ADI-VAM-EN003-F	200µl

Manuals, SDS & CofA

**View Online** »



Mouse astrocytes stained with VAM-EN003. Cells were fixed using 4% paraformaldehyde and permeabilized using 0.01% saponin.

## **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 2

**Application** ICC, IP, WB

Clone GL2A7

Formulation Liquid. Rat Ascites.

GenBank ID J05287

**Host** Rat

**Immunogen** Mouse liver lysosomal membrane preparation.

**Isotype** IgG

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 Mouse, Rabbit

UniProt ID P17047

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com