

LAMP2 polyclonal antibody

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](#).

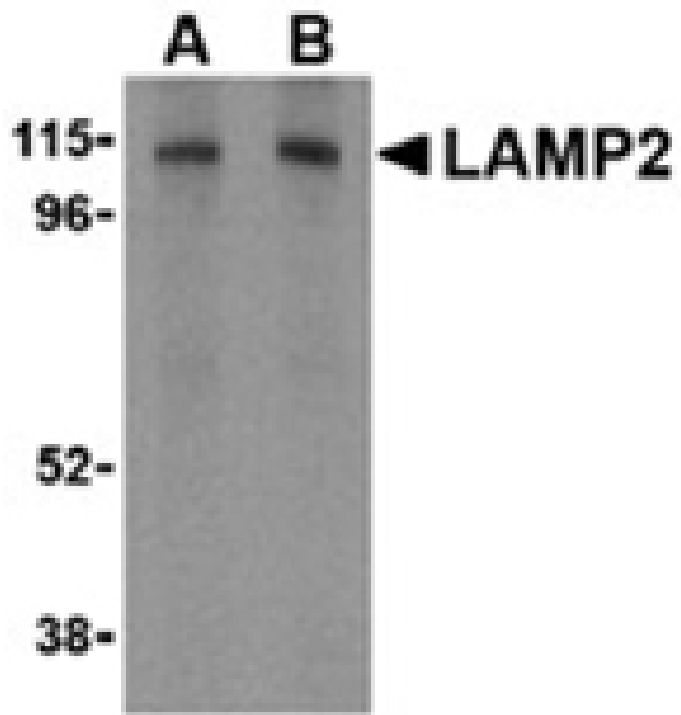
Ordering Information

[Order Online »](#)

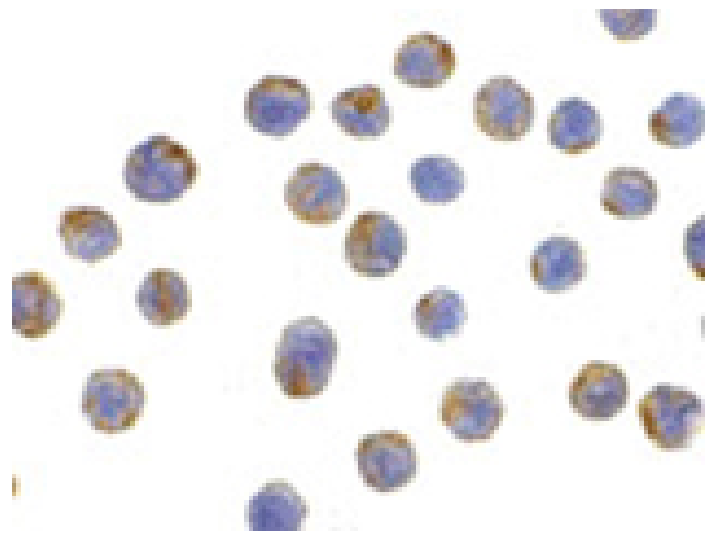
ADI-905-723-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of LAMP-2 in HepG2 cell lysate with LAMP-2 antibody at (A) 1 and (B) 2 μ g/ml.



Immunocytochemistry analysis of LAMP-2 in HepG2 cells with LAMP-2 antibody at 10 μ g/ml.

Handling & Storage

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, WB

Formulation Liquid. In PBS containing 0.02% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the C-terminus of human LAMP2.

Purity Detail Peptide affinity purified.

Species Reactivity Human, Mouse

UniProt ID P13473

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)