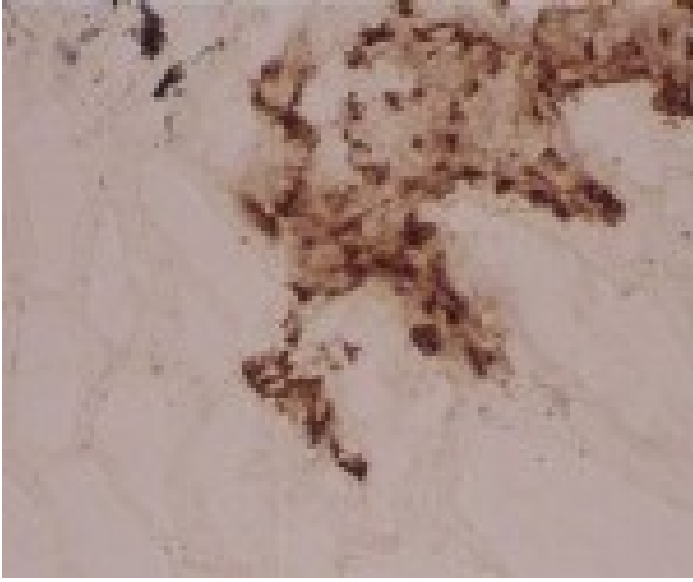


LAMP3 (human) monoclonal antibody (104G4) (biotin conjugate)

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

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human lung frozen section with LAMP3 (human) mAb, clone 104G4.

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	CD208, Lysosome-associated membrane glycoprotein 3, DC LAMP, Protein TSC403
Application	Flow Cytometry, IHC (FS)
Clone	104G4
Formulation	Liquid. In PBS containing 50% glycerol.
Host	Mouse
Immunogen	Human DC-LAMP alkaline phosphatase fusion protein.
Isotype	IgG1
Purity Detail	Purified using QMA Hyper D ion exchange chromatography
Recommendation Dilutions/Conditions	Flow Cytometry (intracellular)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Sheep
Specificity	Recognizes an epitope in the intracytoplasmic domain of DC-LAMP.
UniProt ID	Q9UQV4
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](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)