

LAMP3 (human) monoclonal antibody (104G4) (Alexa Fluor[®] 546)

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

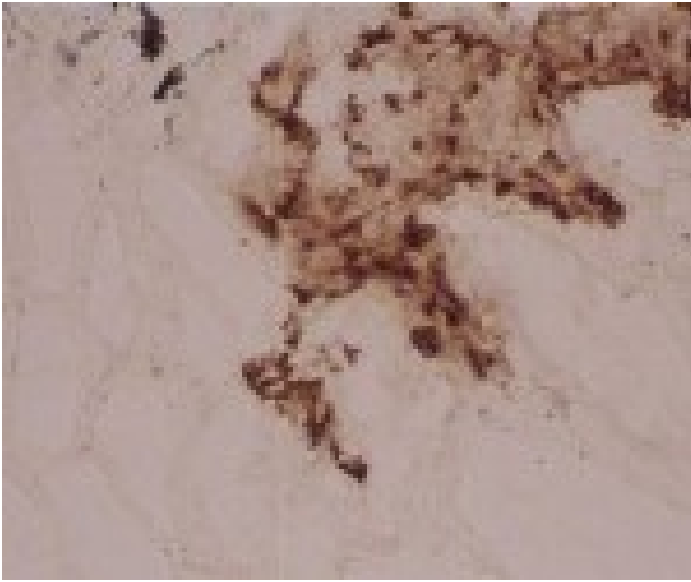
Ordering Information

[Order Online »](#)

ENZ-ABS185-0100	100µg
-----------------	-------

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 IF, 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

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](#)

.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
Last modified: May 29, 2024
usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@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