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

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

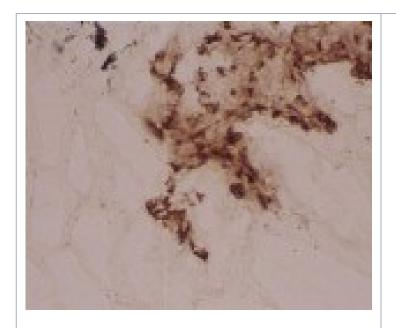
Order Online »

ENZ-ABS210-0100

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** Flow Cytometry, IHC (FS)

**Clone** 104G4

**Formulation** Liquid. In PBS containing 50% glycerol.

**Host** Mouse

Immunogen Human DC-LAMP alkaline phosphatase fusion protein.

lgG1

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



info-

eu@enzolifesciences.com