
LAMP2 Polyclonal Antibody, Biotin Conjugated

(Catalog A50333)

Background

Implicated in tumor cell metastasis. May function in protection of the lysosomal membrane from autodigestion, maintenance of the acidic environment of the lysosome, adhesion when expressed on the cell surface (plasma membrane), and inter- and intracellular signal transduction. Protects cells from the toxic effects of methylating mutagens.

Description

LAMP2 Polyclonal Antibody, Biotin Conjugated. Biotin. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

P13473

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Lysosome-associated membrane glycoprotein 2 protein (54-268AA)

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

LAMP-2, Lysosome-associated membrane protein 2, CD107 antigen-like family member B, LAMP2

Application

ELISA; Recommended dilution: ELISA 1:500-1:1000
