
LCP1 Polyclonal Antibody, Biotin Conjugated

(Catalog #A57160)

Background

Actin-binding protein. Plays a role in the activation of T-cells in response to costimulation through TCR/CD3 and CD2 or CD28. Modulates the cell surface expression of IL2RA/CD25 and CD69.

Description

LCP1 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

P13796

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Plastin-2 protein (2-627AA)

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

L-plastin, LC64P, Lymphocyte cytosolic protein 1, LCP-1, PLS2, LCP1

Application

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