
CALML5 Polyclonal Antibody, FITC Conjugated

(Catalog # A51315)

Background

Binds calcium. May be involved in terminal differentiation of keratinocytes.

Description

CALML5 Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q9NZT1

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Calmodulin-like protein 5 protein (2-146AA)

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

Calmodulin-like protein 5, Calmodulin-like skin protein, CALML5, CLSP

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
