
CLPTM1L Polyclonal Antibody, FITC Conjugated

(Catalog #A70145)

Background

Enhances cisplatin-mediated apoptosis, when overexpressed.

Description

CLPTM1L 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

Q96KA5

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Cleft lip and palate transmembrane protein 1-like protein (32-172AA)

Storage

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

Alternative Names

Cleft lip and palate transmembrane protein 1-like protein, CLPTM1-like protein, Cisplatin resistance-related protein 9, CRR9p, CLPTM1L, CRR9

Application

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