

FAM96A Polyclonal Antibody

(Catalog # A59086)

Background

May play a role in chromosome segregation through establishment of sister chromatid cohesion.

Description

FAM96A Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q9H5X1

Purification

Protein G purified

Immunogen

Recombinant Human Cytosolic iron-sulfur assembly component 2A protein (1-160AA)

Storage

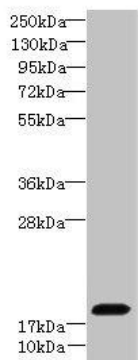
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

CIAO2A antibody; CIA2A antibody; FAM96A antibody; Cytosolic iron-sulfur assembly component 2A antibody; MIP18 family protein FAM96A antibody

Application

ELISA, WB; Recommended dilution: WB:1:500-1:2000



Western blot

All lanes: FAM96A Polyclonal Antibody at 8ug/ml + NIH/3T3 whole cell lysate

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 19, 12 kDa

Observed band size: 19 kDa