

Fcmmr Polyclonal Antibody

(Catalog # A56202)

Background

May play a role in the immune system processes. Protects cells from FAS-, TNF alpha- and FADD-induced apoptosis without increasing expression of the inhibitors of apoptosis BCL2 and BCLXL. Seems to activate an inhibitory pathway that prevents CASP8 activation following FAS stimulation, rather than blocking apoptotic signals downstream. May inhibit FAS-induced apoptosis by preventing CASP8 processing through CFLAR up-regulation.

Description

Fcmmr Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse, Rat

Isotype

IgG

Uniprot ID

A1KXC4

Purification

>95%, Protein G purified

Immunogen

Recombinant Mouse Fas apoptotic inhibitory molecule 3 protein (18-262AA)

Storage

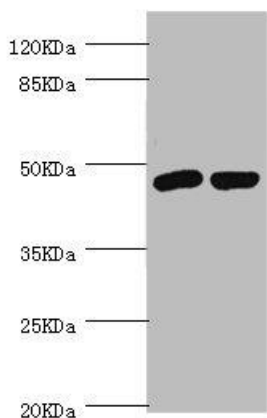
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Regulator of Fas-induced apoptosis Toso, Faim3, Toso

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western blot

All lanes: Fcmmr Polyclonal Antibody at 2ug/ml

Lane 1: rat spleen tissue

Lane 2: rat kidney tissue

Secondary

Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 47kDa

Observed band size: 47kDa