
SECTM1 Polyclonal Antibody, FITC Conjugated

(Catalog #A60852)

Background

May be involved in thymocyte signaling.

Description

SECTM1 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

Q8WVN6

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Secreted and transmembrane protein 1 protein (29-145AA)

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

Secreted and transmembrane protein 1, SECTM1, K12, Protein K-12

Application

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