

SECTM1 Polyclonal Antibody

(Catalog # A60850)

Background

May be involved in thymocyte signaling.

Description

SECTM1 Polyclonal Antibody. Unconjugated. 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

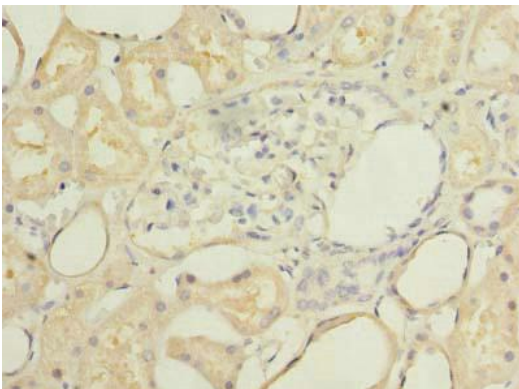
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

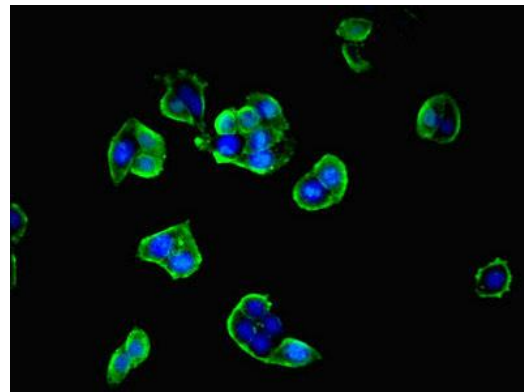
Secreted and transmembrane protein 1, SECTM1, K12, Protein K-12

Application

ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human kidney tissue using SECTM1 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of PC3 cells using SECTM1 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)