

## RAB25 Polyclonal Antibody

(Catalog # A63258)

### Background

Involved in the regulation of cell survival. Promotes invasive migration of cells in which it functions to localize and maintain integrin alpha-V/beta-1 at the tips of extending pseudopodia (PubMed:17925226). Involved in the regulation of epithelial morphogenesis through the control of CLDN4 expression and localization at tight junctions. May selectively regulate the apical recycling pathway. Together with MYO5B regulates transcytosis.

### Description

RAB25 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Human, Mouse

### Isotype

IgG

### Uniprot ID

P57735

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Ras-related protein Rab-25 protein (1-210AA)

### Storage

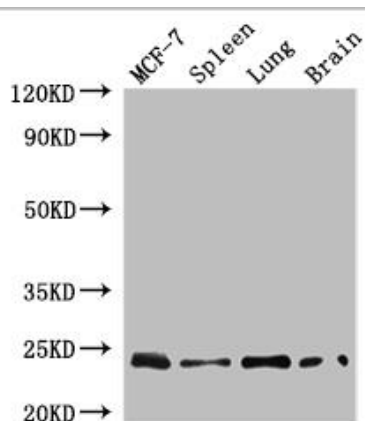
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Ras-related protein Rab-25, CATX-8, RAB25

### Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200, IF:1:50-1:200



### Western Blot

Positive WB detected in: MCF-7 whole cell lysate, Mouse spleen tissue, Mouse lung tissue, Mouse brain tissue

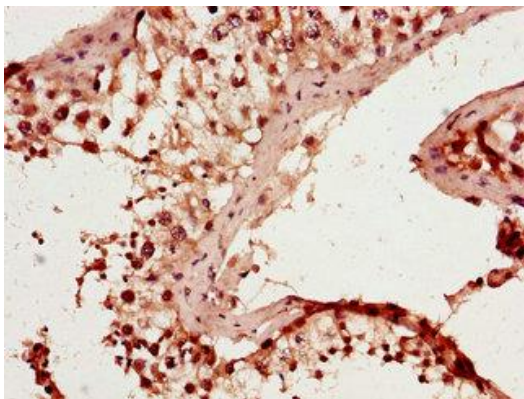
All lanes: RAB25 Polyclonal Antibody at 4ug/ml

Secondary

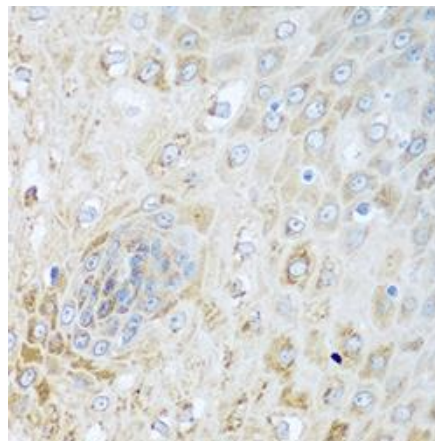
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 24 kDa

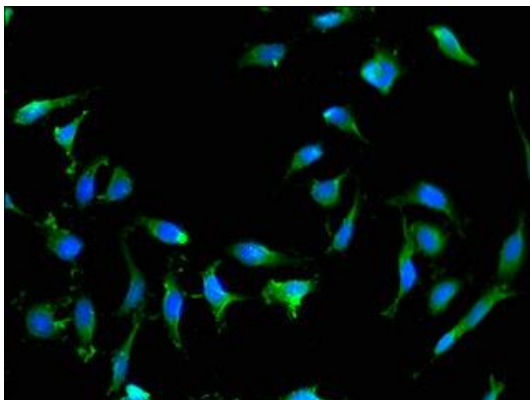
Observed band size: 24 kDa



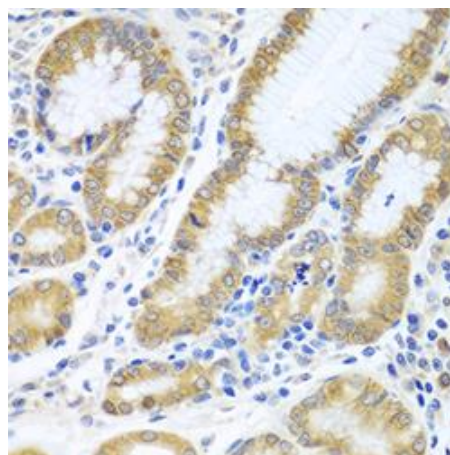
Immunohistochemistry of paraffin-embedded human testis tissue at dilution of 1:100



Immunohistochemistry of paraffin-embedded human esophagus using SIRT3 Polyclonal Antibody at dilution of 1:100 (40x lens).



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



Immunohistochemistry of paraffin-embedded human stomach using SIRT3 Polyclonal Antibody at dilution of 1:100 (40x lens).