

---

## TMEM55A Polyclonal Antibody

(Catalog # A61430)

---

### Background

Catalyzes the hydrolysis of the 4-position phosphate of phosphatidylinositol 4,5-bisphosphate. Does not hydrolyze phosphatidylinositol 3,4,5-trisphosphate, phosphatidylinositol 3,4-bisphosphate, inositol 3,5-bisphosphate, inositol 3,4-bisphosphate, phosphatidylinositol 5-monophosphate, phosphatidylinositol 4-monophosphate and phosphatidylinositol 3-monophosphate.

### Description

TMEM55A 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

Q8N4L2

### Purification

Protein G purified

### Immunogen

Recombinant Human Type 2 phosphatidylinositol 4,5-bisphosphate 4-phosphatase protein (1-191AA)

### Storage

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

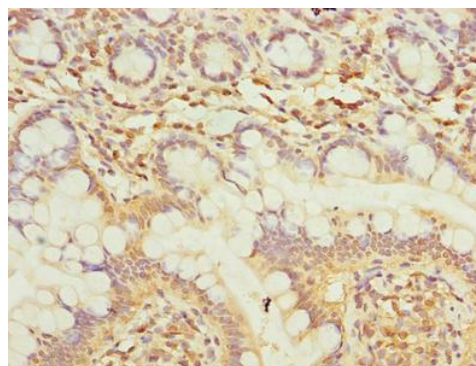
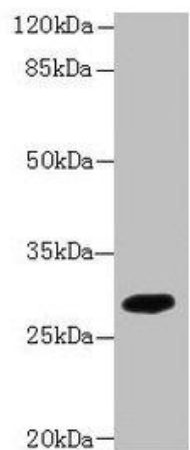
### Alternative Names

Type 2 phosphatidylinositol 4, 5-bisphosphate 4-phosphatase (EC:3.1.3.78), Type 2 PtdIns-4, 5-P2 4-Ptase, TMEM55A, PtdIns-4, 5-P2 4-Ptase II, Transmembrane protein 55A, PIP4P2

### Application

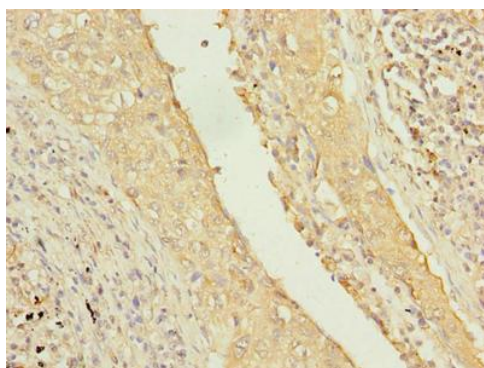
ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:2000, IHC:1:20-1:200, IF:1:50-1:200

---

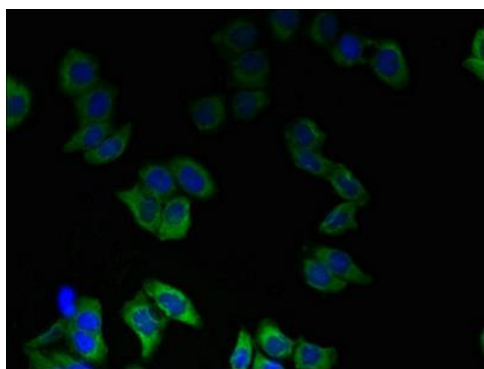


Immunohistochemistry of paraffin-embedded human small intestine tissue using TMEM55A Polyclonal Antibody at dilution of 1:100

Western blot  
All lanes: TMEM55A Polyclonal Antibody at 5ug/ml + Mouse liver tissue  
Secondary  
Goat polyclonal to rabbit IgG at 1/10000 dilution  
Predicted band size: 29 kDa  
Observed band size: 29 kDa



Immunohistochemistry of paraffin-embedded human lung cancer using TMEM55A Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of LO2 cells using TMEM55A Polyclonal Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)