

## FAM167A Polyclonal Antibody

(Catalog #A66462)

### Description

FAM167A 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

Q96KS9

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Protein FAM167A protein (1-214AA)

### Storage

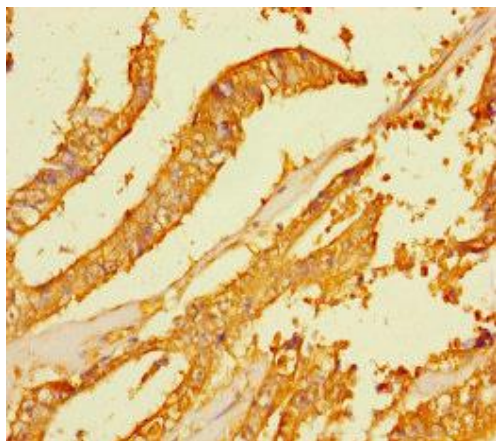
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

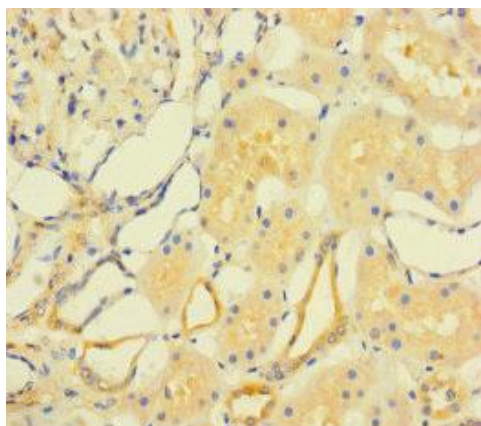
Protein FAM167A, FAM167A, C8orf13

### Application

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



Immunohistochemistry of paraffin-embedded human endometrial cancer using FAM167A Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human kidney tissue using FAM167A Polyclonal Antibody at dilution of 1:100