

## ZC3H12A Polyclonal Antibody

(Catalog # A61850)

### Background

Has RNase activity and selectively degrades specific target mRNA species. Modulates the immune response and inflammation by regulating the decay of specific mRNA molecules. Recognizes the 3'-untranslated region (UTR) of the mRNA for IL6, CALCR and IL12B. Required for normal decay of IL6 mRNA. Triggers apoptosis and promotes angiogenesis in response to the binding of CCL2 to CCR2. Regulates expression of CDH12 and CHD19.

### Description

ZC3H12A 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

Q5D1E8

### Purification

Protein G purified

### Immunogen

Recombinant Human Endoribonuclease ZC3H12A protein (1-260AA)

### Storage

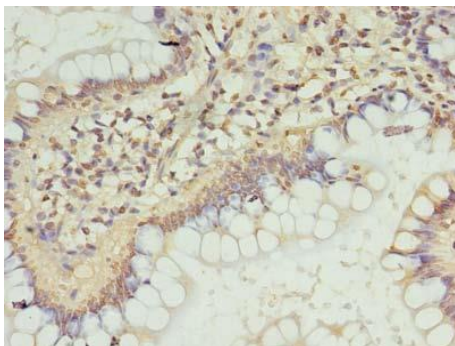
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Ribonuclease ZC3H12A (EC:3.1.-.-), ZC3H12A, MCP-induced protein 1, Zinc finger CCCH domain-containing protein 12A

### Application

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



Immunohistochemistry of paraffin-embedded human small intestine tissue using ZC3H12A Polyclonal Antibody at dilution of 1:100