

---

## UBIAD1 Polyclonal Antibody

(Catalog # A73155)

---

### Background

This gene encodes a protein thought to be involved in cholesterol and phospholipid metabolism. Mutations in this gene are associated with Schnyder crystalline corneal dystrophy.

### Description

UBIAD1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

### Specificity

Human, Mouse, Rat

### Isotype

IgG

### Uniprot ID

Q9Y5Z9

### Purification

Affinity Purified

### Immunogen

A synthetic peptide corresponding to a sequence within amino acids 1-100 of human UBIAD1 (NP\_037451.1).

### Storage

Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

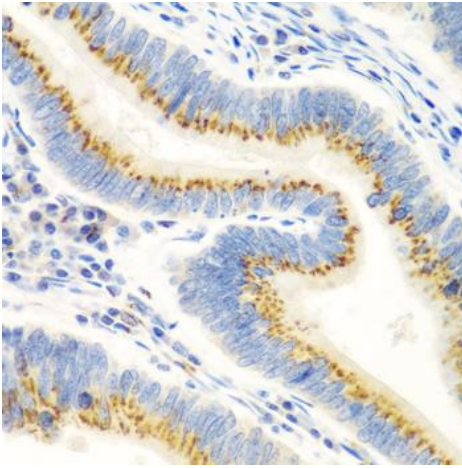
### Alternative Names

UBIAD1; SCCD; TERE1; UbiA prenyltransferase domain containing 1

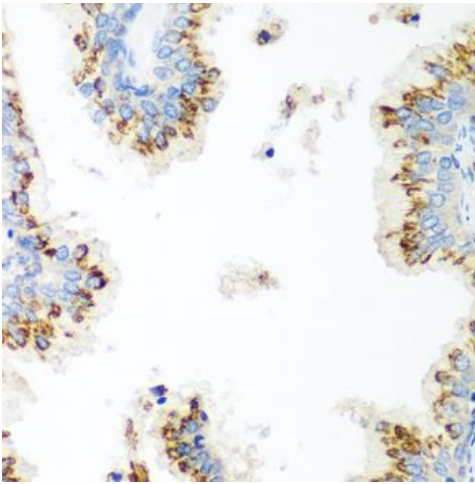
### Application

WB, IHC; Recommended dilution: WB 1:500 - 1:2000, IHC 1:100 - 1:200

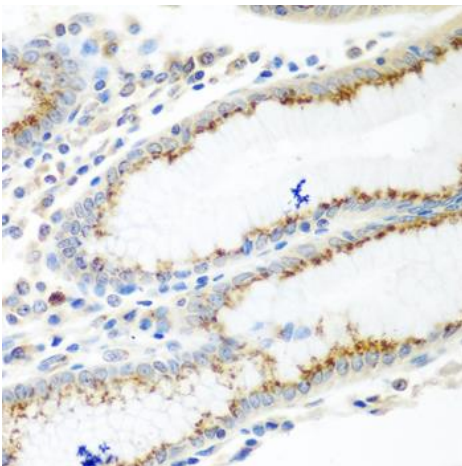
---



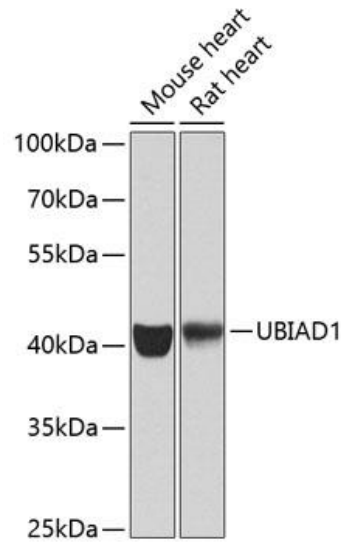
Immunohistochemistry of paraffin-embedded human colon carcinoma using UBIAD1 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded human prostate using UBIAD1 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded human stomach using UBIAD1 Polyclonal Antibody at dilution of 1:100 (40x lens).



Western blot analysis of extracts of various cell lines, using UBIAD1 Polyclonal Antibody at 1:1000 dilution. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug per lane. Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 60s.