

---

## COL6A6 Polyclonal Antibody

(Catalog # A57485)

---

### Background

Collagen VI acts as a cell-binding protein.

### Description

COL6A6 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

A6NMZ7

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Collagen alpha-6(VI) chain protein (430-982AA)

### Storage

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

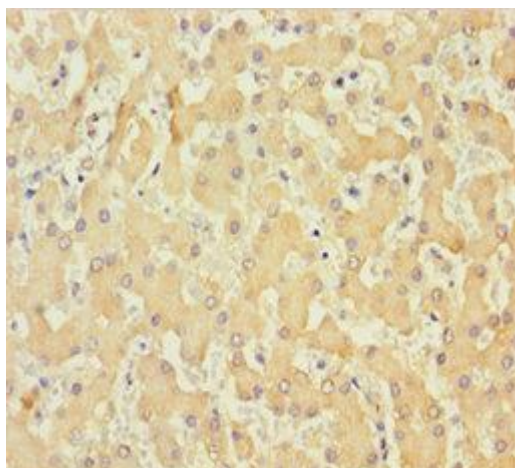
### Alternative Names

COL6A6

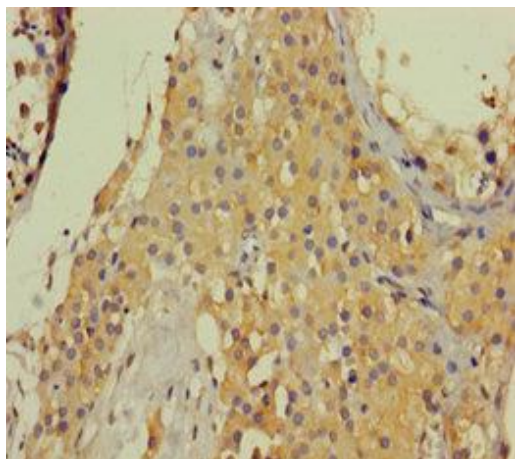
### Application

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

---



Immunohistochemistry of paraffin-embedded human liver tissue using COL6A6 Polyclonal Antibody at dilution 1:100



Immunohistochemistry of paraffin-embedded human testis tissue using COL6A6 Polyclonal Antibody at dilution 1:100