

---

## UGT2A3 Polyclonal Antibody

(Catalog # A65650)

---

### Background

UDP-glucuronosyltransferases catalyze phase II biotransformation reactions in which lipophilic substrates are conjugated with glucuronic acid to increase water solubility and enhance excretion. They are of major importance in the conjugation and subsequent elimination of potentially toxic xenobiotics and endogenous compounds.

### Description

UGT2A3 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

Q6UWM9

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human UDP-glucuronosyltransferase 2A3 protein (201-491AA)

### Storage

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

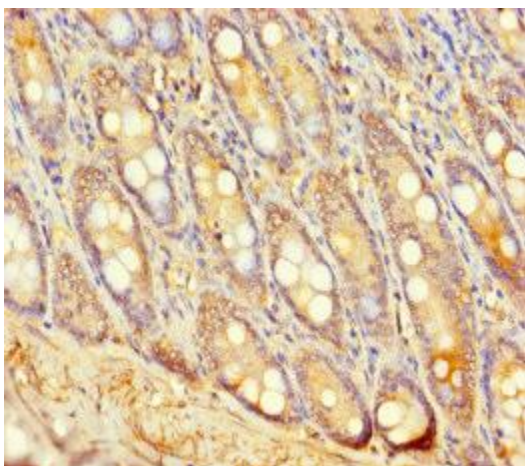
### Alternative Names

UDP-glucuronosyltransferase 2A3, UDPGT 2A3, UGT2A3, UNQ2559/PRO6239

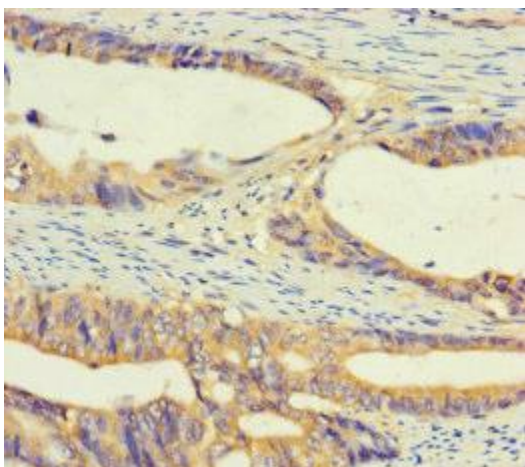
### Application

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

---



Immunohistochemistry of paraffin-embedded human small intestine tissue using UGT2A3 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human colon cancer using UGT2A3 Polyclonal Antibody at dilution of 1:100