

## CYP2U1 Polyclonal Antibody

(Catalog # A65794)

### Background

Catalyzes the hydroxylation of arachidonic acid, docosahexaenoic acid and other long chain fatty acids. May modulate the arachidonic acid signaling pathway and play a role in other fatty acid signaling processes.

### Description

CYP2U1 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

Q7Z449

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Cytochrome P450 2U1 protein (363-496AA)

### Storage

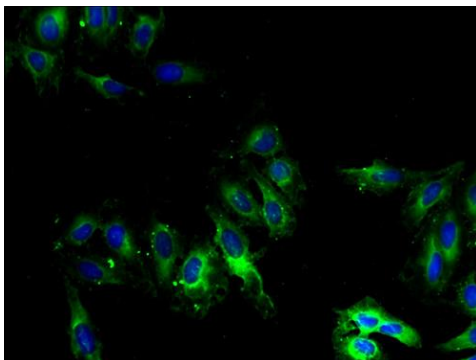
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Cytochrome P450 2U1, CYP2U1

### Application

ELISA, IF; Recommended dilution: IF:1:50-1:200



Immunofluorescent analysis of HeLa cells using CYP2U1 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)