

## ALOX5 Polyclonal Antibody

(Catalog #A58062)

### Background

Catalyzes the first step in leukotriene biosynthesis, and thereby plays a role in inflammatory processes.

### Description

ALOX5 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

0.03% Proclin 300. 50% Glycerol, 0.01M PBS, pH 7.4.

### Specificity

Human

### Isotype

IgG

### Uniprot ID

P09917

### Purification

Protein G purified

### Immunogen

Recombinant Human Arachidonate 5-lipoxygenase protein (2-149AA)

### Storage

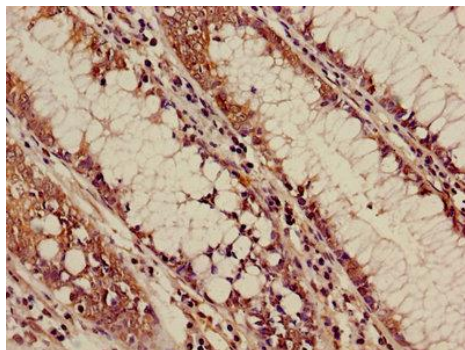
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

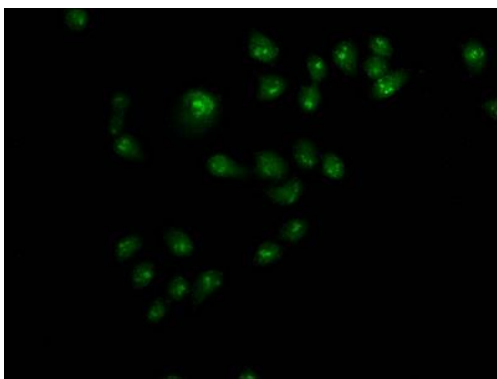
Arachidonate 5-lipoxygenase (EC:1.13.11.34), 5-LO, 5-lipoxygenase, ALOX5, LOG5

### Application

ELISA, IHC, IF; Recommended dilution: IHC 1:20 - 1:200, IF 1:50 - 1:200



Immunohistochemistry of paraffin-embedded human colon cancer using ALOX5 Antibody at dilution of 1:100



Immunofluorescence staining of HeLa cells with ALOX5 Antibody at 1:110, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L).