

## ZDHHC23 Polyclonal Antibody

(Catalog No. A61866)

### Background

Palmitoyltransferase that mediates palmitoylation of KCNMA1, regulating localization of KCNMA1 to the plasma membrane. May be involved in NOS1 regulation and targeting to the synaptic membrane.

### Description

ZDHHC23 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Human, Mouse

### Isotype

IgG

### Uniprot ID

Q8IYP9

### Purification

Caprylic Acid Ammonium Sulfate Precipitation Purified

### Immunogen

Recombinant human Palmitoyltransferase ZDHHC23 protein

### Storage

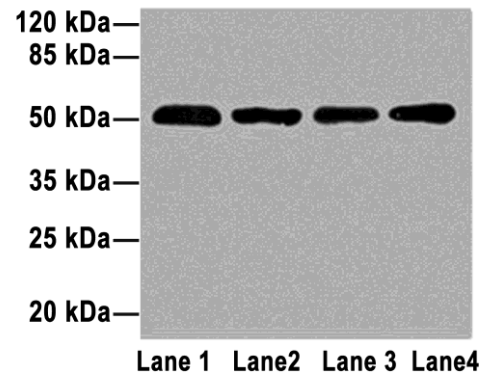
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Application

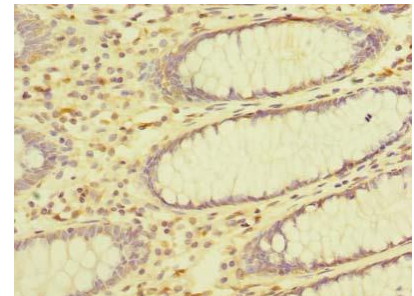
ELISA, WB, IHC, IF

### Ordering Information

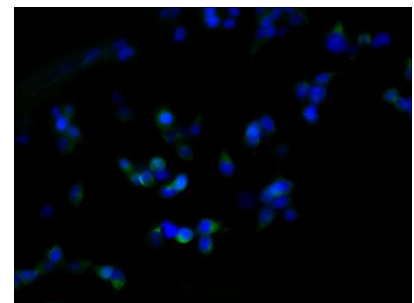
Products	Size	Cat. No.
ZDHHC23 Polyclonal Antibody	50 µg	A61866-050
	100 µg	A61866-100



- Western blot - All Lanes: ZDHHC23 Polyclonal Antibody at 8µg/ml  
Lane 1: 293T whole cell lysate  
Lane 2: Mouse brain tissue  
Lane 3: A549 whole cell lysate  
Lane 4: Human placenta tissue  
Secondary  
Goat polyclonal to Rabbit IgG at 1/10000 dilution  
Predicted band size: 46 kDa  
Observed band size: 46 kDa



- Immunohistochemistry of paraffin-embedded human colon cancer using ZDHHC23 Polyclonal Antibody at dilution of 1:100.



- Immunofluorescent analysis of 293T cells using ZDHHC23 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L).

*This product is for research purposes only. Not intended for use in diagnostic procedures.*