

ZDHHC12 Polyclonal Antibody

(Catalog #A66298)

Background

endoplasmic reticulum, Golgi apparatus, palmitoyltransferase activity, protein-cysteine S-palmitoyltransferase activity, peptidyl-L-cysteine S-palmitoylation, protein palmitoylation, protein targeting to membrane

Description

ZDHHC12 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

Q96GR4

Purification

Affinity Purified

Immunogen

Recombinant Human Probable palmitoyltransferase ZDHHC12 protein (121-197AA)

Storage

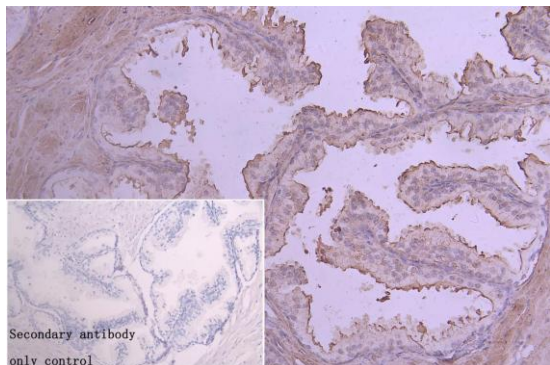
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Probable palmitoyltransferase ZDHHC12, Zinc finger DHHC domain-containing protein 12, DHHC-12, Zinc finger protein 400, ZDHHC12, ZNF400, PSEC0008

Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:100



IHC image of ZDHHC12 Polyclonal Antibody diluted at 1:20 and staining in paraffin-embedded human prostate cancer performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a Goat anti-rabbit polymer IgG labeled by HRP and visualized using 0.05% DAB. Secondary antibody only control: uses 1% BSA instead of primary antibody.