
ZDHHC6 Polyclonal Antibody

(Catalog # A61886)

Background

Palmitoylates calnexin (CALX) which is required for its association with the ribosome-translocon complex and efficient folding of glycosylated proteins.

Description

ZDHHC6 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q9H6R6

Purification

Affinity Purified

Immunogen

Recombinant Human Palmitoyltransferase ZDHHC6 protein (227-409AA)

Storage

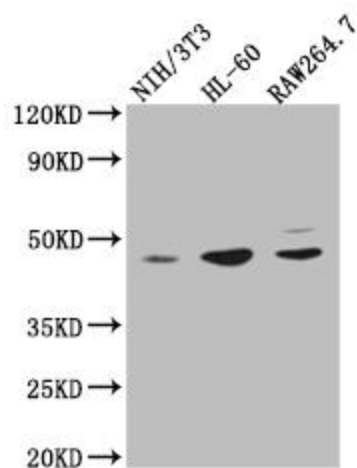
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Palmitoyltransferase ZDHHC6 (EC:2.3.1.225), ZDHHC6, ZNF376, Transmembrane protein H4, Zinc finger DHHC domain-containing protein 6, DHHC-6, Zinc finger protein 376

Application

ELISA, WB; Recommended dilution: WB:1:500-1:2000



Western Blot

Positive WB detected in: NIH/3T3 whole cell lysate, HL-60 whole cell lysate, RAW264.7 whole cell lysate

All lanes: ZDHHC6 antibody at 1:2000

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 48 kDa

Observed band size: 48 kDa.