

---

## HACD2 Polyclonal Antibody

(Catalog # A65662)

---

### Background

Catalyzes the third of the four reactions of the long-chain fatty acids elongation cycle. This endoplasmic reticulum-bound enzymatic process, allows the addition of two carbons to the chain of long- and very long-chain fatty acids/VLCFAs per cycle. This enzyme catalyzes the dehydration of the 3-hydroxyacyl-CoA intermediate into trans-2,3-enoyl-CoA, within each cycle of fatty acid elongation. Thereby, it participates to the production of VLCFAs of different chain lengths that are involved in multiple biological processes as precursors of membrane lipids and lipid mediators.

### Description

HACD2 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

Q6Y1H2

### Purification

Protein G purified

### Immunogen

Recombinant Human Very-long-chain (3R)-3-hydroxyacyl-CoA dehydratase 2 protein (7-41AA)

### Storage

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

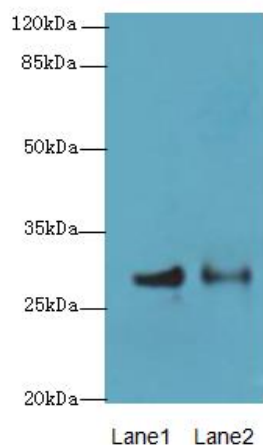
### Alternative Names

Very-long-chain (3R)-3-hydroxyacyl-CoA dehydratase 2, 3-hydroxyacyl-CoA dehydratase 21, HACD21

### Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:2000, IHC:1:20-1:200, IF:1:50-1:200

---



#### Western blot

All Lanes: HACD2 Polyclonal Antibody at 4ug/ml

Lane 1: MCF7 whole cell lysate

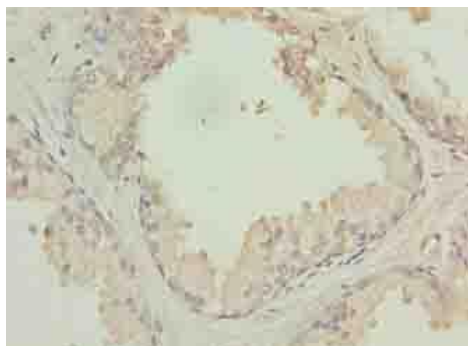
Lane 2: LO2 whole cell lysate

Secondary

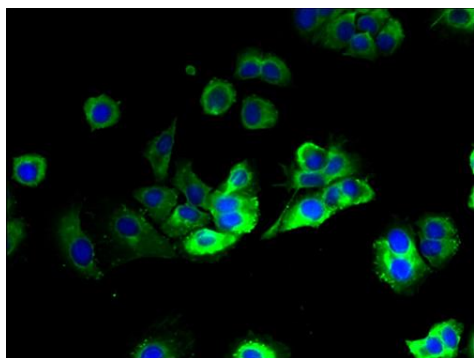
Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 28 kDa

Observed band size: 28 kDa



Immunohistochemistry of paraffin-embedded human prostate cancer using HACD2 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of MCF-7 cells using HACD2 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)