

---

## GCDH Polyclonal Antibody

(Catalog # A66898)

---

### Background

Catalyzes the oxidative decarboxylation of glutaryl-CoA to crotonyl-CoA and CO<sub>2</sub> in the degradative pathway of L-lysine, L-hydroxylysine, and L-tryptophan metabolism. It uses electron transfer flavoprotein as its electron acceptor. Isoform Short is inactive.

### Description

GCDH 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

Q92947

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Glutaryl-CoA dehydrogenase, mitochondrial protein (45-300AA)

### Storage

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

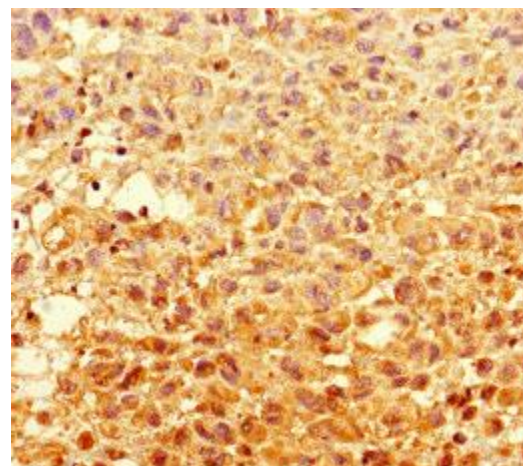
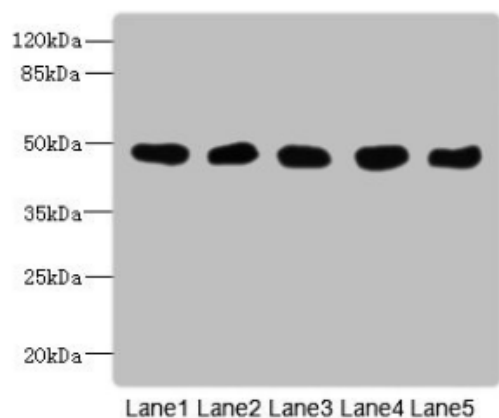
### Alternative Names

Glutaryl-CoA dehydrogenase, mitochondrial, GCD, GCDH

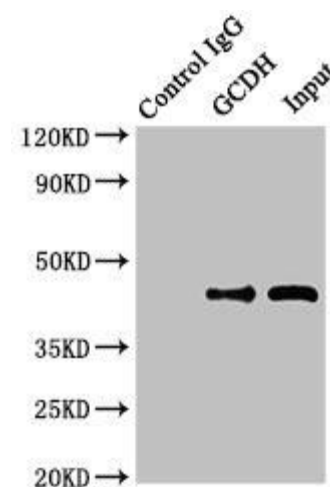
### Application

ELISA, WB, IHC, IP; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200, IP:1:200-1:2000

---



Immunohistochemistry of paraffin-embedded human melanoma cancer using GCDH Polyclonal Antibody at dilution of 1:100



Immunoprecipitating GCDH in HeLa whole cell lysate  
 Lane 1: Rabbit monoclonal IgG (1 ug) in HeLa whole cell lysate. For western blotting, a HRP-conjugated Protein G antibody was used as the secondary antibody (1/2000)  
 Lane 2: GCDH Polyclonal Antibody (8 ug) + HeLa whole cell lysate (500 ug)  
 Lane 3: HeLa whole cell lysate (10 ug)

#### Western Blot

All lanes: GCDH Polyclonal Antibody at 8ug/ml

Lane 1: Mouse kidney tissue

Lane 2: Mouse liver tissue

Lane 3: HeLa whole cell lysate

Lane 4: MCF7 whole cell lysate

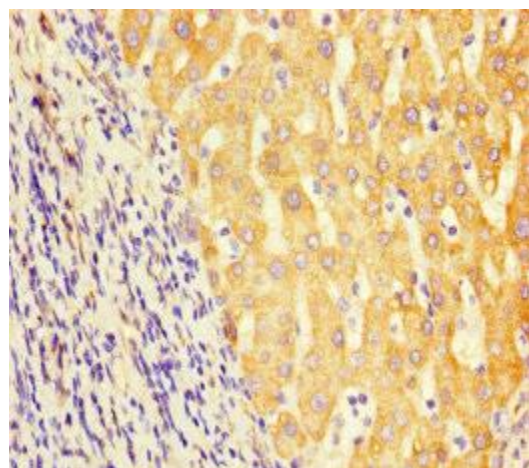
Lane 5: LO2 whole cell lysate

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 49, 48 kDa

Observed band size: 48 kDa



Immunohistochemistry of paraffin-embedded human liver cancer using GCDH Polyclonal Antibody at dilution of 1:100