

## ACAD8 Polyclonal Antibody

(Catalog # A68138)

### Background

Has very high activity toward isobutyryl-CoA. Is an isobutyryl-CoA dehydrogenase that functions in valine catabolism. Plays a role in transcriptional coactivation within the ARC complex.

### Description

ACAD8 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

Q9UKU7

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Isobutyryl-CoA dehydrogenase, mitochondrial protein (201-415AA)

### Storage

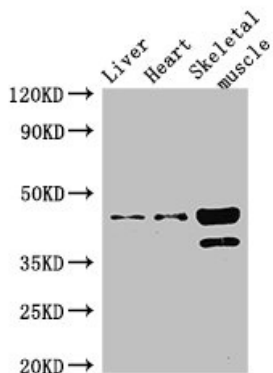
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Isobutyryl-CoA dehydrogenase, mitochondrial, Activator-recruited cofactor 42 kDa component, ARC42, Acyl-CoA dehydrogenase family member 8, ACAD-8, ACAD8, ARC42, IBD

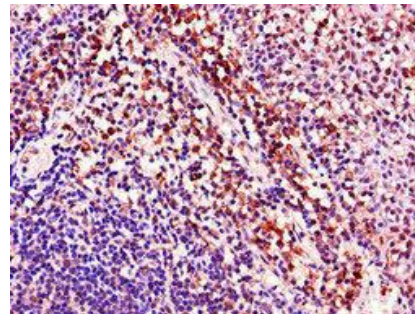
### Application

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



#### ◀ Western Blot

Positive WB detected in: Mouse liver tissue, Mouse heart tissue, Mouse skeletal muscle tissue  
All lanes: ACAD8 Polyclonal Antibody at 3ug/ml  
Secondary  
Goat polyclonal to rabbit IgG at 1/50000 dilution  
Predicted band size: 46, 32, 39 kDa  
Observed band size: 46, 39 kDa



◀ Immunohistochemistry of paraffin-embedded human tonsil tissue using ACAD8 Polyclonal Antibody at dilution of 1:100