

ACADM Polyclonal Antibody

(Catalog # A62038)

Background

Acyl-CoA dehydrogenase specific for acyl chain lengths of 4 to 16 that catalyzes the initial step of fatty acid beta-oxidation. Utilizes the electron transfer flavoprotein (ETF) as an electron acceptor to transfer electrons to the main mitochondrial respiratory chain via ETF-ubiquinone oxidoreductase (ETF dehydrogenase).

Description

ACADM Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P11310

Purification

Protein G purified

Immunogen

Recombinant Human Medium-chain specific acyl-CoA dehydrogenase, mitochondrial protein (26-421AA)

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

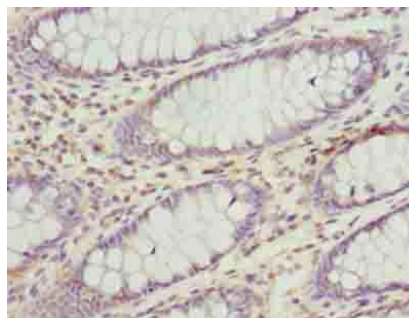
Medium-chain specific acyl-CoA dehydrogenase, mitochondrial, MCAD, ACADM

Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human kidney tissue at dilution of 1:100



Immunohistochemistry of paraffin-embedded human colon cancer at dilution of 1:100