
OXCT1 Polyclonal Antibody, FITC Conjugated

(Catalog # A54774)

Background

Key enzyme for ketone body catabolism. Transfers the CoA moiety from succinate to acetoacetate. Formation of the enzyme-CoA intermediate proceeds via an unstable anhydride species formed between the carboxylate groups of the enzyme and substrate.

Description

OXCT1 Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P55809

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Succinyl-CoA:3-ketoacid coenzyme A transferase 1, mitochondrial protein (40-520AA)

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

3-oxoacid CoA-transferase 1 Somatic-type succinyl-CoA:3-oxoacid CoA-transferase OXCT1 OXCT, SCOT

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
