
BCKDHA Polyclonal Antibody

(Catalog # A73493)

Background

The branched-chain alpha-keto acid (BCAA) dehydrogenase (BCKD) complex is an inner mitochondrial enzyme complex that catalyzes the second major step in the catabolism of the branched-chain amino acids leucine, isoleucine, and valine. The BCKD complex consists of three catalytic components: a heterotetrameric (alpha2-beta2) branched-chain alpha-keto acid decarboxylase (E1), a dihydrolipoyl transacylase (E2), and a dihydrolipoamide dehydrogenase (E3). This gene encodes the alpha subunit of the decarboxylase (E1) component. Mutations in this gene result in maple syrup urine disease, type IA. Multiple transcript variants encoding different isoforms have been found for this gene.

Description

BCKDHA Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P12694

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 276-445 of human BCKDHA (NP_000700.1).

Storage

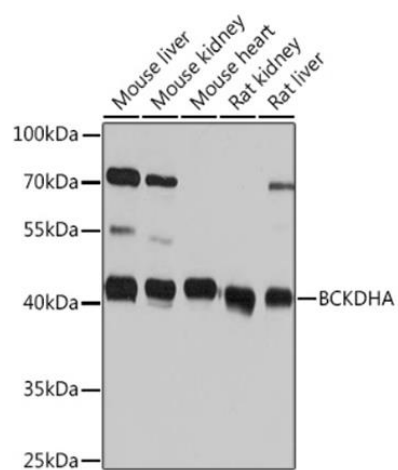
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

BCKDHA; BCKDE1A; MSU; MSUD1; OVD1A

Application

WB; Recommended dilution: WB 1:500 - 1:2000



Western blot analysis of extracts of various cell lines, using the BCKDHA Polyclonal Antibody at 1:1000 dilution.

Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

Blocking buffer: 3% nonfat dry milk in TBST.

Exposure time: 10s.