
BCAT1 Polyclonal Antibody

(Catalog #A62146)

Background

Catalyzes the first reaction in the catabolism of the essential branched chain amino acids leucine, isoleucine, and valine.

Description

BCAT1 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

P54687

Purification

Protein G purified

Immunogen

Recombinant Human Branched-chain-amino-acid aminotransferase, cytosolic protein (4-276AA)

Storage

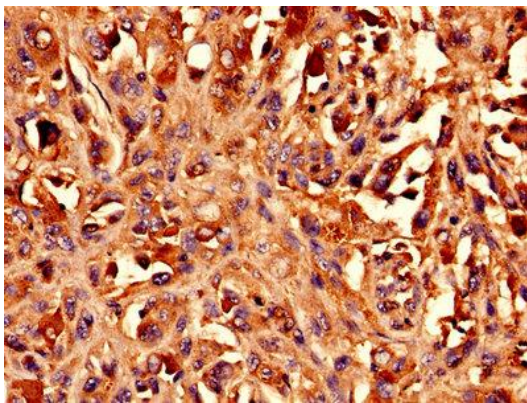
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

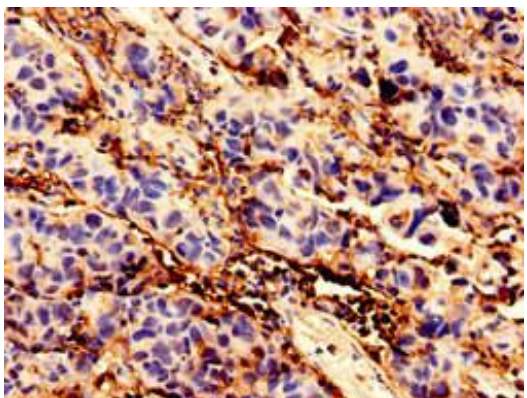
Branched-chain-amino-acid aminotransferase, cytosolic, BCAT(c), Protein ECA39, BCAT1, BCT1, ECA39

Application

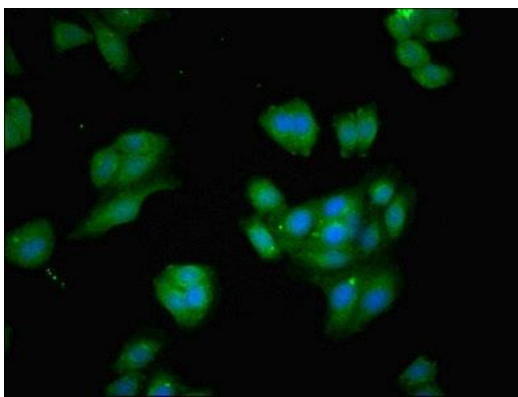
ELISA, IHC, IF; Recommended dilution: IHC 1:20 - 1:200, IF 1:50 - 1:200



Immunocytochemistry analysis of human melanoma cancer using BCAT1 Polyclonal Antibody at dilution of 1:100



Immunocytochemistry analysis of human bladder cancer using BCAT1 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HepG2 cells using BCAT1 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)