

GOT1 Polyclonal Antibody

(Catalog # A52800)

Background

Plays a key role in amino acid metabolism.

Description

GOT1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P17174

Purification

Protein G purified

Immunogen

Recombinant Human Aspartate aminotransferase, cytoplasmic protein (2-411AA)

Storage

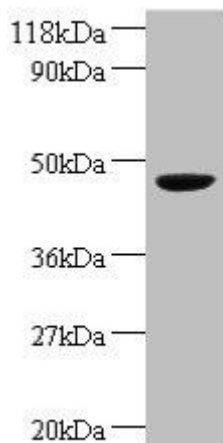
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

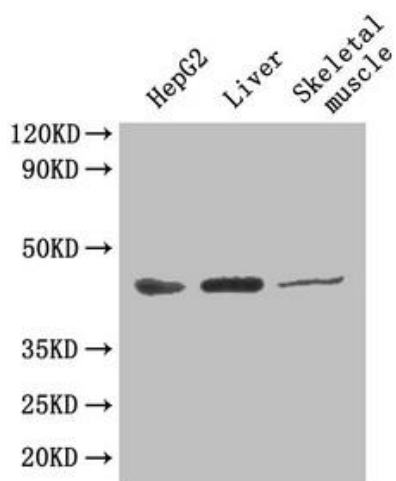
Glutamate oxaloacetate transaminase 1, Transaminase, GOT1

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:200-1:500, IF:1:50-1:200



All lanes: GOT1 Polyclonal Antibody at 2ug/ml+mouse brain tissue
Secondary
Goat polyclonal to Rabbit IgG at 1/10000 dilution
Predicted band size: 47, 45 kDa
Observed band size: 46 kDa



Western Blot

Positive WB detected in: hepG2 cell, mouse liver, mouse skeletal muscle

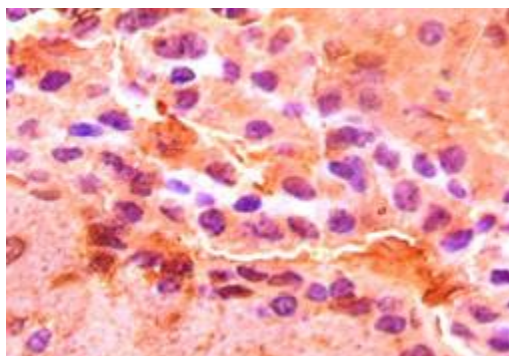
All lanes: GOT1 Polyclonal Antibody at 4.5ug/ml

Secondary

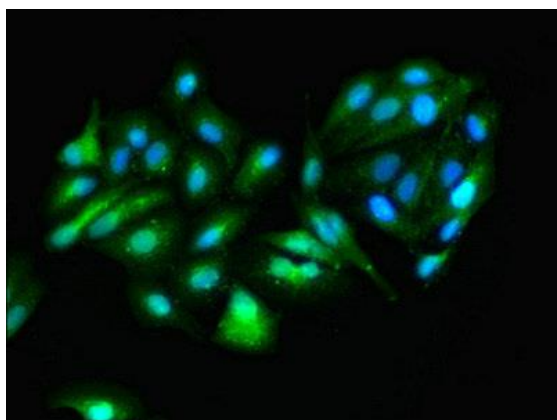
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 47, 45 kDa

Observed band size: 46 kDa



Immunohistochemistry of paraffin-embedded mouse brain using GOT1 Polyclonal Antibody at dilution of 1:50



Immunofluorescent analysis of HepG2 cells using GOT1 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)