
ACSBG2 Polyclonal Antibody

(Catalog # A57978)

Background

Mediates activation of long-chain fatty acids for both synthesis of cellular lipids, and degradation via beta-oxidation. Able to activate long-chain fatty acids. Also able to activate very long-chain fatty acids; however, the relevance of such activity is unclear in vivo. Has increased ability to activate oleic and linoleic acid. May play a role in spermatogenesis.

Description

ACSBG2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q5FVE4

Purification

Protein G purified

Immunogen

Recombinant Human Long-chain-fatty-acid--CoA ligase ACSBG2 protein (301-600AA)

Storage

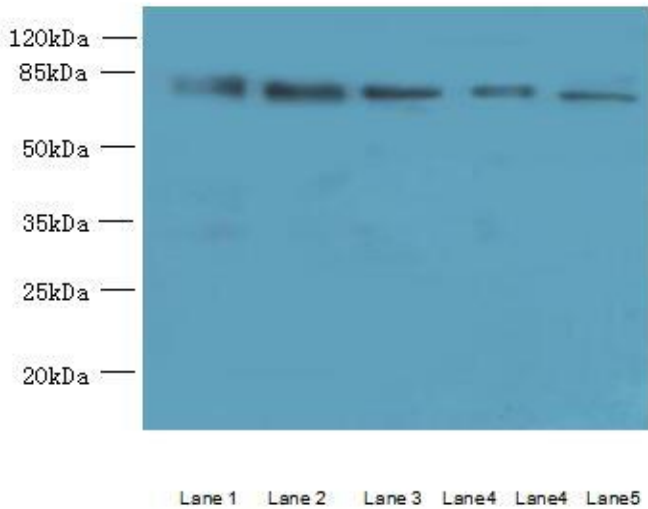
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Long-chain-fatty-acid--CoA ligase ACSBG2 (EC:6.2.1.3), ACSBG2, BGR, UNQ2443/PRO5005, Acyl-CoA synthetase bubblegum family member 2, Bubblegum-related protein, PRTD-NY3

Application

ELISA, WB, IHC; Recommended dilution: WB:1:500-1:2000, IHC:1:20-1:200



Western blot

All Lanes: ACSBG2 Polyclonal Antibody at 5ug/ml

Lane 1: Hela whole cell lysate

Lane 2: K562 whole cell lysate

Lane 3: A549 whole cell lysate

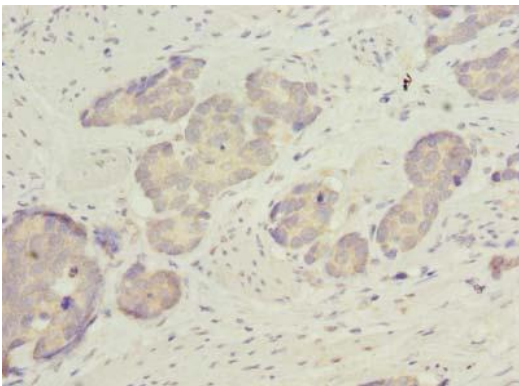
Lane 4: Human high value serum
Lane 5: A431 whole cell lysate

Secondary

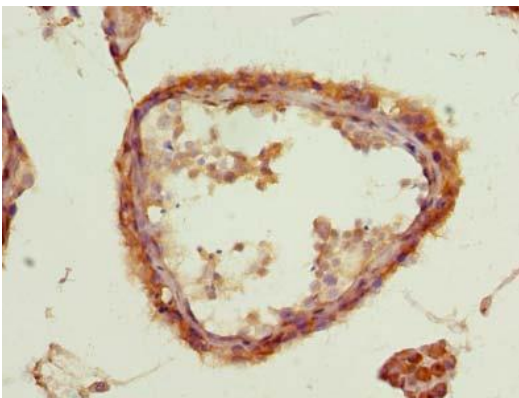
Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 74 kDa

Observed band size: 74 kDa



Immunohistochemistry of paraffin-embedded human gastric cancer using ACSBG2 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human testis tissue using ACSBG2 Polyclonal Antibody at dilution of 1:100