
APOC3 Polyclonal Antibody

(Catalog # A55307)

Background

Component of triglyceride-rich very low-density lipoproteins (VLDL) and high density lipoproteins (HDL) in plasma. Plays a multifaceted role in triglyceride homeostasis. Intracellularly, promotes hepatic very low-density lipoprotein 1 (VLDL1) assembly and secretion; extracellularly, attenuates hydrolysis and clearance of triglyceride-rich lipoproteins (TRLs). Impairs the lipolysis of TRLs by inhibiting lipoprotein lipase and the hepatic uptake of TRLs by remnant receptors. Formed of several curved helices connected via semiflexible hinges, so that it can wrap tightly around the curved micelle surface and easily adapt to the different diameters of its natural binding partners

Description

APOC3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Rat

Isotype

IgG

Uniprot ID

P02656

Purification

Protein G purified

Immunogen

Recombinant Human Apolipoprotein C-III protein (21-99AA)

Storage

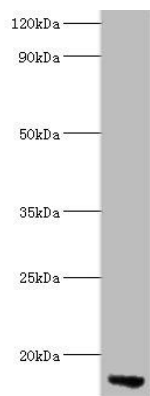
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Apolipoprotein C3 APOC3

Application

ELISA, WB, IHC; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200



Western blot

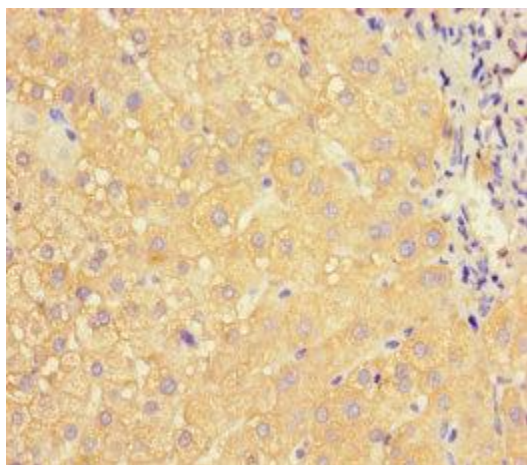
All lanes: APOC3 Polyclonal Antibody at 2ug/ml+rat liver tissue

Secondary

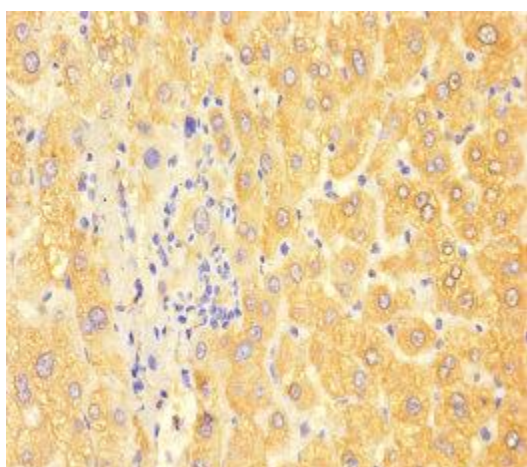
Goat polyclonal to Rabbit IgG at 1/15000 dilution

Predicted band size: 11kDa

Observed band size: 11kDa



Immunohistochemistry of paraffin-embedded human liver cancer at dilution of 1:100



Immunohistochemistry of paraffin-embedded human liver tissue at dilution of 1:100