

HMGCN Polyclonal Antibody

(Catalog # A52932)

Background

Transmembrane glycoprotein that is the rate-limiting enzyme in cholesterol biosynthesis as well as in the biosynthesis of nonsterol isoprenoids that are essential for normal cell function including ubiquinone and geranylgeranyl proteins.

Description

HMGCN 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

P04035

Purification

Protein G purified

Immunogen

Recombinant Human 3-hydroxy-3-methylglutaryl-coenzyme A reductase protein (588-887AA)

Storage

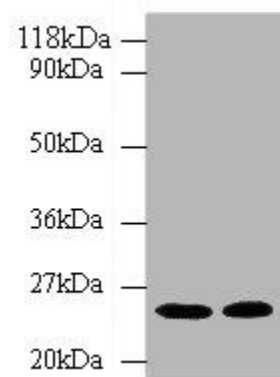
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

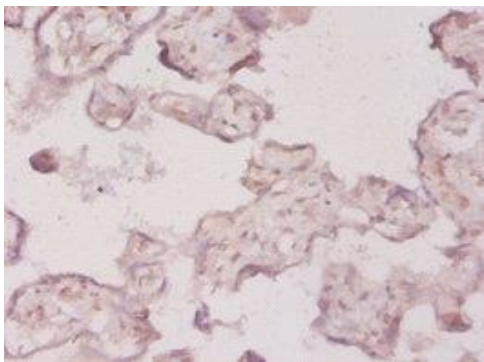
HMGCN

Application

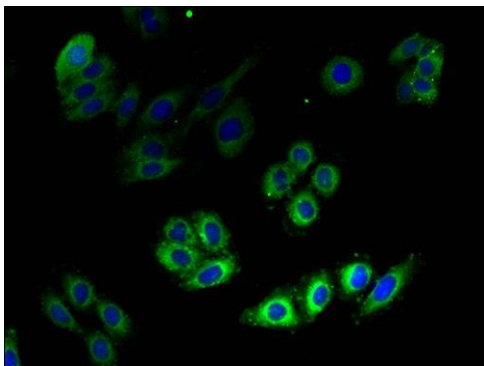
ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:2000, IHC:1:200-1:500, IF:1:50-1:200



Western blot
All lanes: HMGCN Polyclonal Antibody at 2ug/ml
Lane 1: Hela whole cell lysate
Lane 2: 293T whole cell lysate
Secondary
Goat polyclonal to rabbit at 1/10000 dilution
Predicted band size: 98, 93, 100 kDa
Observed band size: 25 kDa



Immunohistochemistry of paraffin-embedded human placenta at dilution of 1:20



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