

MVD Polyclonal Antibody

(Catalog # A69603)

Background

Performs the first committed step in the biosynthesis of isoprenes.

Description

MVD 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

P53602

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Diphosphomevalonate decarboxylase protein (69-221AA)

Storage

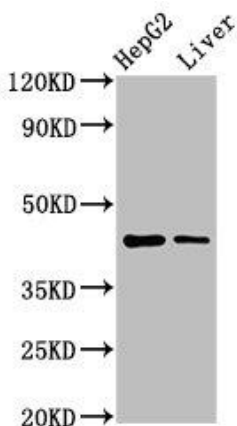
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Diphosphomevalonate decarboxylase, Mevalonate (diphospho)decarboxylase, MDDase, Mevalonate pyrophosphate decarboxylase, MVD, MPD

Application

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



Western Blot

Positive WB detected in: HepG2 whole cell lysate, Rat liver tissue

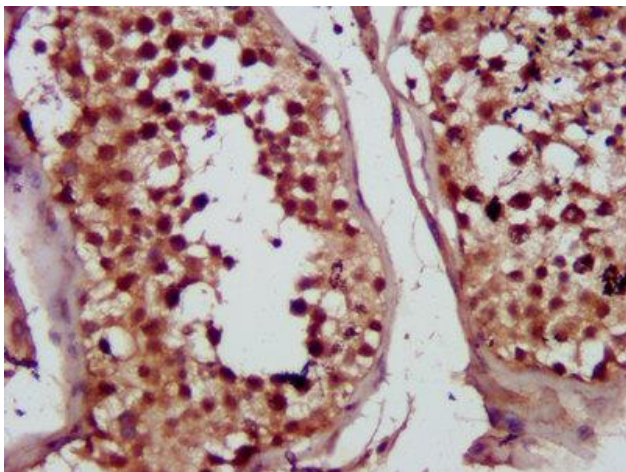
All lanes: MVD antibody at 7.4ug/ml

Secondary

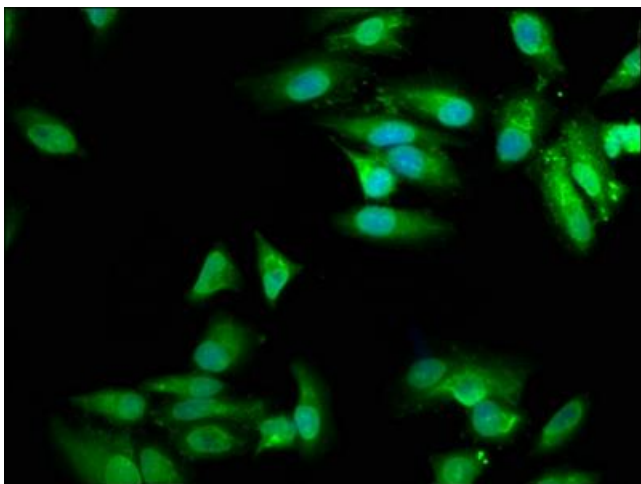
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 44 kDa

Observed band size: 44 kDa



IHC image of MVD Polyclonal Antibody diluted at 1:300 and staining in paraffin-embedded human testis tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



Immunofluorescence staining of HeLa cells with MVD Polyclonal Antibody at 1:100, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L).