

PRDM11 Polyclonal Antibody

(Catalog # A-2011)

Background

The PR domain-containing protein, PRDM11, (also known as PFM8) is an amino acid protein that produces two isoforms through alternative splicing. This gene localizes to the minimal liver tumor suppressor region within chromosome 11p11.2-p12.

Description

PRDM11 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q9NQV5

Purification

Protein G purified

Immunogen

Recombinant Human PR domain-containing protein 11 protein (63-215AA)

Storage

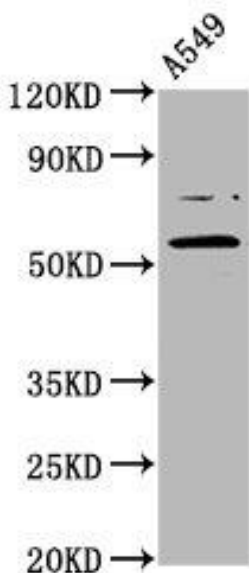
Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

Alternative Names

PFM8 antibody, PR domain containing protein 11 antibody

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:500-1:1000, IF:1:200-1:500



Western Blot

Positive WB detected in: A549 whole cell lysate

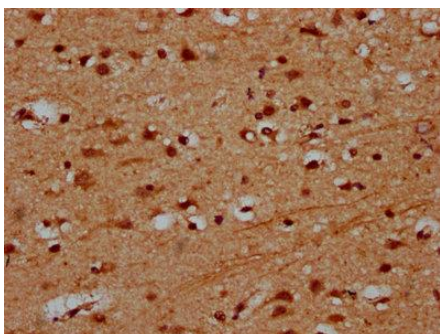
All lanes: PRDM11 antibody at 5.6µg/ml

Secondary

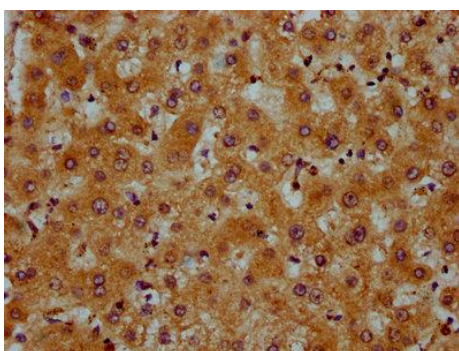
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 58, 54 kDa

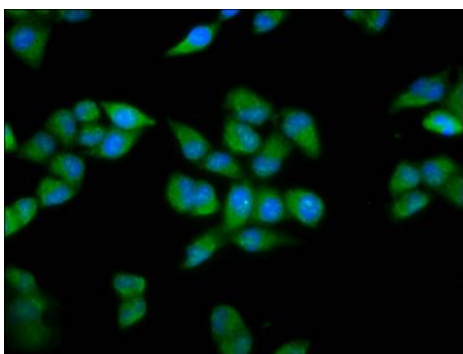
Observed band size: 58 kDa



IHC image of PRDM11 Polyclonal Antibody diluted at 1:600 and staining in paraffin-embedded human brain 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.



IHC image of PRDM11 Polyclonal Antibody diluted at 1:600 and staining in paraffin-embedded human liver 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 PRDM11 Polyclonal Antibody at 1:200, 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).