
MMD2 Polyclonal Antibody

(Catalog #A66126)

Background

perinuclear region of cytoplasm, protein kinase activity, positive regulation of neuron differentiation, positive regulation of protein kinase activity, positive regulation of Ras protein signal transduction, protein phosphorylation, regulation of protein localization

Description

MMD2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q8IY49

Purification

Protein G purified

Immunogen

Recombinant Human Monocyte to macrophage differentiation factor 2 protein (1-38AA)

Storage

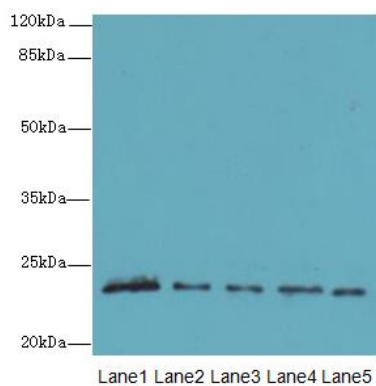
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Monocyte to macrophage differentiation factor 2, Progestin and adipoQ receptor family member 10, Progestin and adipoQ receptor family member X, MMD2, PAQR10

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:2000, IHC:1:20-1:200, IF:1:50-1:200



Western blot

All Lanes: MMD2 Polyclonal Antibody at 3ug/ml

Lane 1: Mouse brain tissue

Lane 2: Mouse kidney tissue

Lane 3: U251 whole cell lysate

Lane 4: 293T whole cell lysate

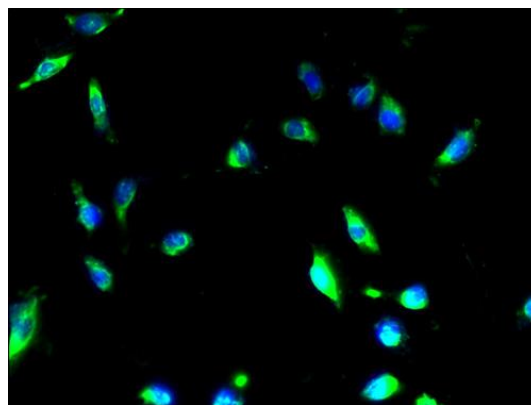
Lane 5: Mouse liver tissue

Secondary

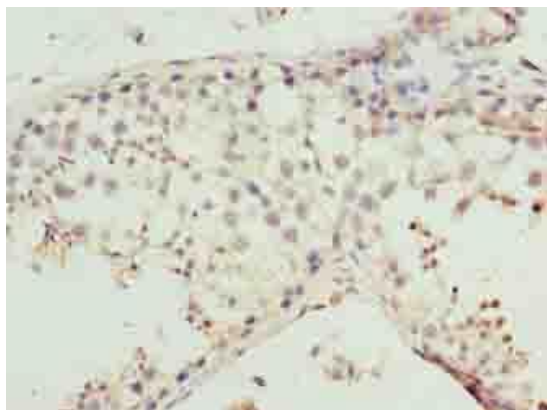
Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 31 kDa

Observed band size: 23 kDa



Immunofluorescent analysis of U251 cells using MMD2 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunohistochemistry of paraffin-embedded human testis tissue using MMD2 Polyclonal Antibody at dilution of 1:100