
PTAFR Polyclonal Antibody

(Catalog #A63250)

Background

Receptor for platelet activating factor, a chemotactic phospholipid mediator that possesses potent inflammatory, smooth-muscle contractile and hypotensive activity. Seems to mediate its action via a G protein that activates a phosphatidylinositol-calcium second messenger system.

Description

PTAFR Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P25105

Purification

Protein G purified

Immunogen

Recombinant Human Platelet-activating factor receptor protein (297-342AA)

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

Platelet-activating factor receptor, PAF-R, PAFr, PTAFR, PAFR

Application

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



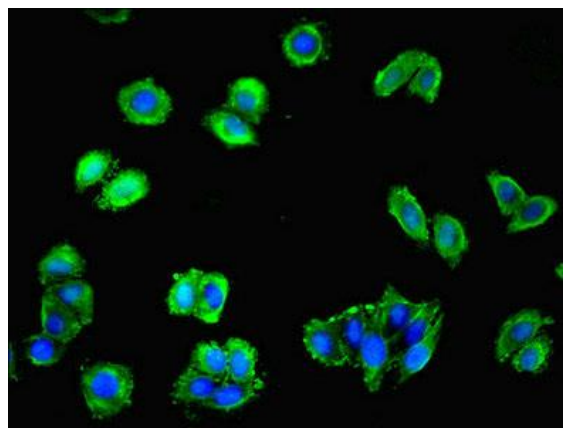
Western blot

All lanes: PTAFR Polyclonal Antibody at 4ug/ml + Mouse lung tissue

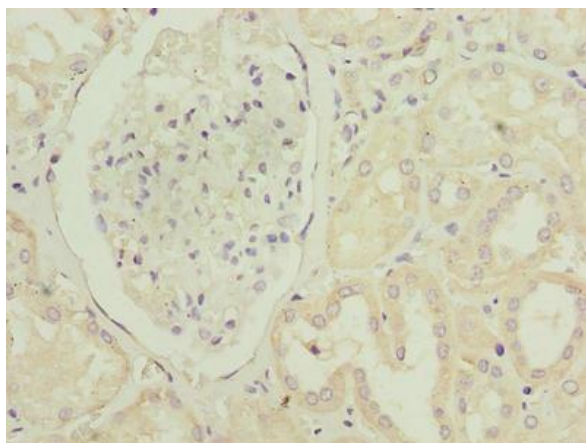
Secondary: Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 40 kDa

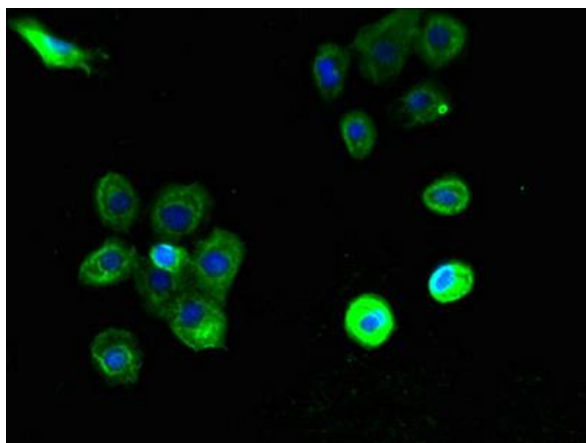
Observed band size: 40 kDa



Immunofluorescent analysis of HepG2 cells using PTAFR 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 kidney tissue using PTAFR Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of A431 cells using PTAFR Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)