

GPR34 Polyclonal Antibody

(Catalog # A68062)

Background

Orphan receptor.

Description

GPR34 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

Q9UPC5

Purification

Protein G purified

Immunogen

Recombinant Human Probable G-protein coupled receptor 34 protein (1-61AA)

Storage

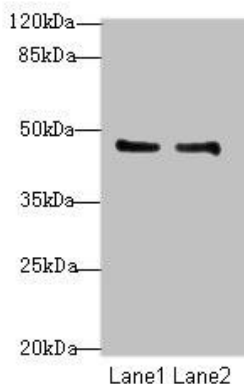
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Probable G-protein coupled receptor 34, GPR34

Application

ELISA, WB, IF; Recommended dilution: WB:1:1000-1:5000, IF:1:50-1:200



Western blot

All lanes: GPR34 antibody at 8ug/ml

Lane 1: Mouse liver tissue

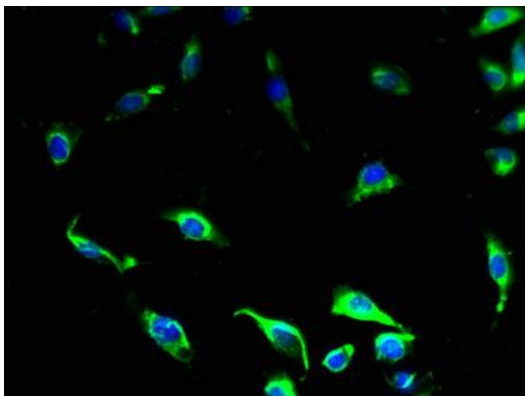
Lane 2: Mouse kidney tissue

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 44 kDa

Observed band size: 44 kDa



Immunofluorescent analysis of U251 cells using GPR34
Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated
AffiniPure Goat Anti-Rabbit IgG(H+L)