
FPR1 Polyclonal Antibody

(Catalog # A59150)

Background

High affinity receptor for N-formyl-methionyl peptides (fMLP), which are powerful neutrophil chemotactic factors (PubMed:2161213, PubMed:2176894, PubMed:10514456, PubMed:15153520). Binding of fMLP to the receptor stimulates intracellular calcium mobilization and superoxide anion release (PubMed:2161213, PubMed:1712023, PubMed:15153520). This response is mediated via a G-protein that activates a phosphatidylinositol-calcium second messenger system (PubMed:1712023, PubMed:10514456).

Description

FPR1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P21462

Purification

Protein G purified

Immunogen

Recombinant Human fMet-Leu-Phe receptor protein (306-350AA)

Storage

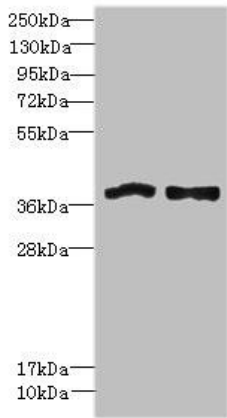
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

fMet-Leu-Phe receptor, fMLP receptor, FPR1, N-formyl peptide receptor, FPR, N-formylpeptide chemoattractant receptor

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: FPR1 antibody at 6ug/ml

Lane 1: Colo320 whole cell lysate

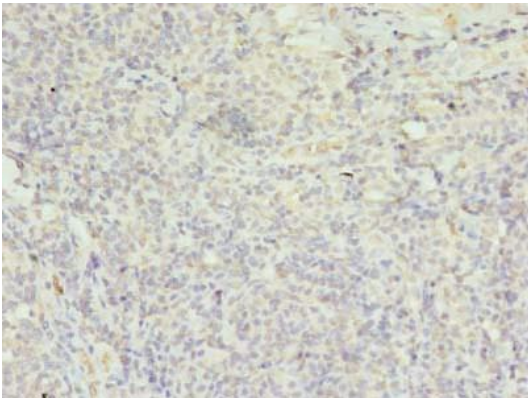
Lane 2: HT29 whole cell lysate

Secondary

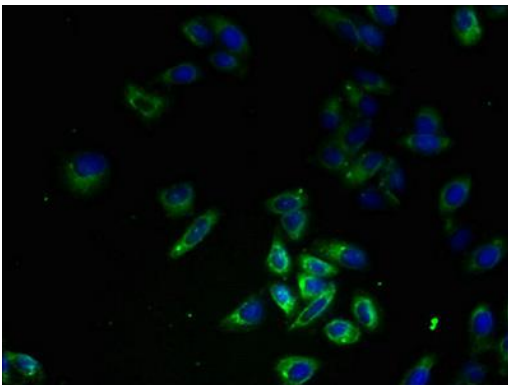
Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 39 kDa

Observed band size: 39 kDa



Immunohistochemistry of paraffin-embedded human tonsil tissue using FPR1 Antibody at dilution of 1:100



Immunofluorescent analysis of HepG2 cells using FPR1 Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)