

## FPR2 Polyclonal Antibody

(Catalog # A59154)

### Background

Low affinity receptor for N-formyl-methionyl peptides, which are powerful neutrophils chemotactic factors. Binding of FMLP to the receptor causes activation of neutrophils. This response is mediated via a G-protein that activates a phosphatidylinositol-calcium second messenger system. The activation of LXA4R could result in an anti-inflammatory outcome counteracting the actions of proinflammatory signals such as LTB4 (leukotriene B4).

### Description

FPR2 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

P25090

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human N-formyl peptide receptor 2 protein (307-351AA)

### Storage

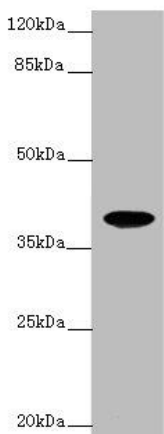
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

N-formyl peptide receptor 2, FPR2, FPRH1, FPRL1, LXA4R, FMLP-related receptor I, FMLP-R-I, Formyl peptide receptor-like 1, HM63, Lipoxin A4 receptor, LXA4 receptor, RFP

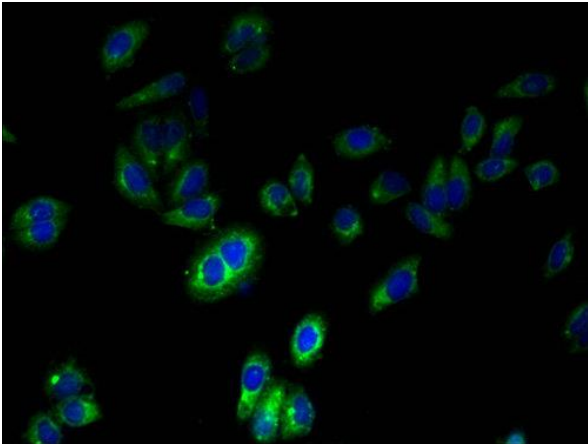
### Application

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



### Western blot

All lanes: FPR2 Polyclonal Antibody at 1ug/ml+ K562 whole cell lysate  
Goat polyclonal to rabbit at 1/10000 dilution  
Predicted band size: 39 kDa  
Observed band size: 39 kDa



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