

---

## ACKR1 Polyclonal Antibody

(Catalog # A50945)

---

### Background

Atypical chemokine receptor that controls chemokine levels and localization via high-affinity chemokine binding that is uncoupled from classic ligand-driven signal transduction cascades, resulting instead in chemokine sequestration, degradation, or transcytosis. Also known as interceptor (internalizing receptor) or chemokine-scavenging receptor or chemokine decoy receptor. Has a promiscuous chemokine-binding profile, interacting with inflammatory chemokines of both the CXC and the CC subfamilies but not with homeostatic chemokines. Acts as a receptor for chemokines including CCL2, CCL5, CCL7, CCL11, CCL13, CCL14, CCL17, CXCL5, CXCL6, IL8/CXCL8, CXCL11, GRO, RANTES, MCP-1, TARC and also for the malaria parasites *P. vivax* and *P. knowlesi*. May regulate chemokine bioavailability and, consequently, leukocyte recruitment through two distinct mechanisms: when expressed in endothelial cells, it sustains the abluminal to luminal transcytosis of tissue-derived chemokines and their subsequent presentation to circulating leukocytes; when expressed in erythrocytes, serves as blood reservoir of cognate chemokines but also as a chemokine sink, buffering potential surges in plasma chemokine levels.

### Description

ACKR1 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

Q16570

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Atypical chemokine receptor 1 protein (1-63AA)

### Storage

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

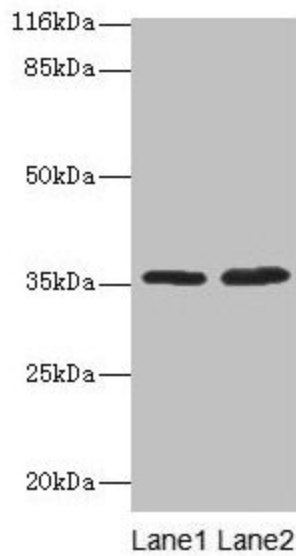
### Alternative Names

Atypical chemokine receptor 1, Duffy antigen/chemokine receptor, Glycoprotein D, Plasmodium vivax receptor, ACKR1, DARC, FY, GPD

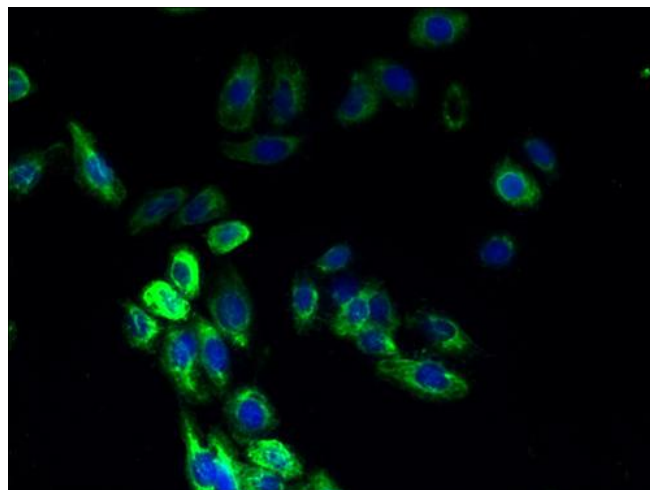
### Application

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

---



Western blot  
All lanes: ACKR1 Polyclonal Antibody at 2ug/ml  
Lane 1: mouse kidney tissue  
Lane 2 mouse spleen tissue  
Secondary  
Goat polyclonal to rabbit at 1/10000 dilution  
Predicted band size: 36kDa  
Observed band size: 36kDa



Immunofluorescent analysis of HepG2 cells using ACKR1 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)