

## GPR17 Polyclonal Antibody

(Catalog # A59290)

### Background

Dual specificity receptor for uracil nucleotides and cysteinyl leukotrienes (CysLTs). Signals through G(i) and inhibition of adenylyl cyclase. May mediate brain damage by nucleotides and CysLTs following ischemia.

### Description

GPR17 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

Q13304

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Uracil nucleotide/cysteinyl leukotriene receptor protein (1-36AA)

### Storage

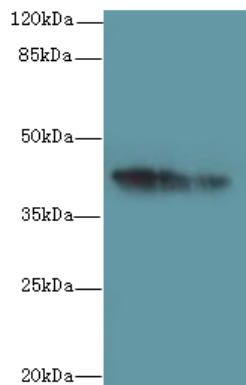
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Uracil nucleotide/cysteinyl leukotriene receptor, UDP/CysLT receptor, GPR17, G-protein coupled receptor 17, P2Y-like receptor, R12

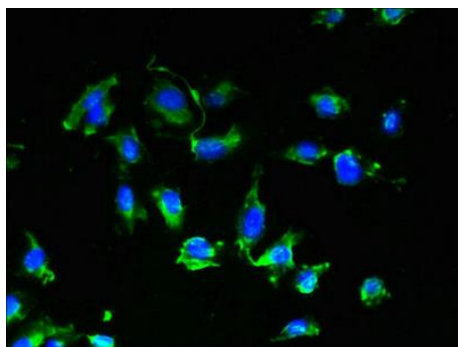
### Application

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



#### Western blot

All Lanes: GPR17 Polyclonal Antibody at 10ug/ml  
Lane 1: Mouse brain tissue  
Lane 2: Human high value serum  
Secondary  
Goat polyclonal to Rabbit IgG at 1/10000 dilution  
Predicted band size: 41 kDa  
Observed band size: 41 kDa



Immunofluorescent analysis of U251 cells using GPR17 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)