

# PRODUCT INFORMATION



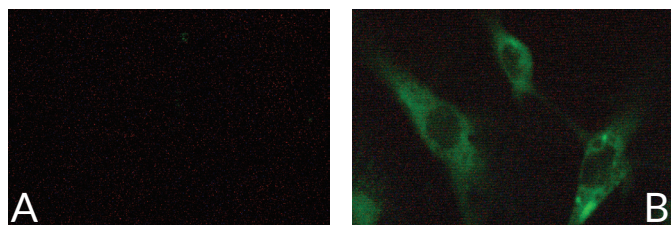
## GPR17 (C-Term) Polyclonal Antibody (Immunofluorescence)

Item No. 17087

### Overview and Properties

<b>Contents:</b>	This vial contains a 500 µl of peptide affinity-purified polyclonal antibody.
<b>Synonyms:</b>	G Protein-Coupled Receptor 17, P2Y-Like Receptor, Uracil Nucleotide/Cysteinyl Leukotriene Receptor
<b>Immunogen:</b>	Peptide from the C-terminal region of human GRP17
<b>Species Reactivity:</b>	(+) Human
<b>Uniprot No.:</b>	Q13304
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Storage Buffer:</b>	PBS, pH 7.2, with 50% glycerol, and 0.02% sodium azide
<b>Host:</b>	Rabbit
<b>Applications:</b>	Immunofluorescence (IF); the recommended starting dilution is 1:20. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



**Immunofluorescent staining of HUVEC cells.** HUVEC cells were fixed with 4% paraformaldehyde, permeabilized, and then blocked with 1% fetal bovine serum. Cells were probed with the indicated antibodies, washed between steps, and images were captured using a Leica DMIL inverted fluorescence microscope (40X objective).

**Panel A:** Control with Cayman's Goat Anti-Rabbit IgG FITC (secondary antibody; Item No. 10006588) alone (1:200)

**Panel B:** GPR17 (C-Term) Polyclonal Antibody (Immunofluorescence) (1:20)

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/17/2023

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

GPR17 is a G protein-coupled receptor that has been identified as a dualistic receptor recognizing signals from two unrelated chemical families: nucleotides and cysteinyl leukotrienes (CysLTs).<sup>1</sup> The deorphanization of GPR17 supports the suggested crosstalk between nucleotides and CysLTs during inflammation and injury. mRNA transcripts encoding this transmembrane receptor are most strongly expressed in the brain, kidney, and heart. Upon ischemic injury, GPR17 is upregulated in these tissues and its inhibition has been shown to decrease ischemic damage. This finding suggests GPR17 as a pharmacological target of ischemic injury.<sup>1</sup> Cayman's GPR17 (C-Term) Polyclonal Antibody (Immunofluorescence) is a peptide affinity-purified polyclonal antibody that recognizes the C-terminal region of human GPR17. This protein exists in two isoforms, differing by 28 amino acids at the receptor N-terminus. The longer form of the protein consists of 367 amino acids with a calculated molecular weight of 41 kDa. The human and rat proteins share 89% amino acid identity.<sup>1</sup>

## Reference

---

1. Ciana, P., Fumagalli, M., Trincavelli, M.L., *et al.* The orphan receptor GPR17 identified as a new dual uracil nucleotides/cysteinyl-leukotrienes receptor. *EMBO J.* 25, 4615-4627 (2006).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM