

# PRODUCT INFORMATION



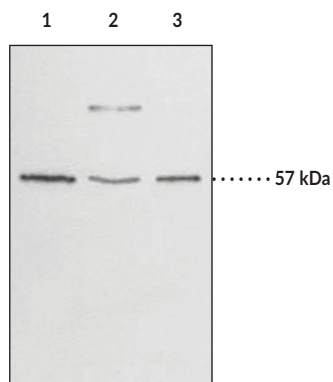
## GPR17 (C-Term) Polyclonal Antibody

Item No. 10136

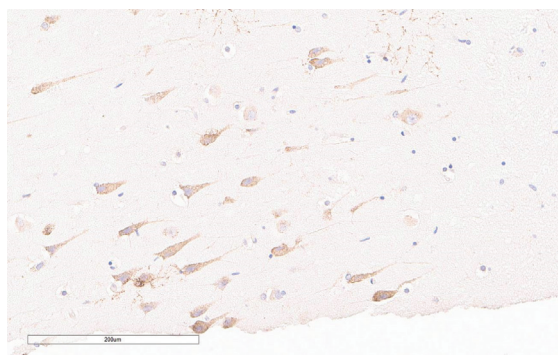
### Overview and Properties

<b>Contents:</b>	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
<b>Synonym:</b>	G Protein-Coupled Receptor 17
<b>Immunogen:</b>	A synthetic peptide from the C-terminal region of human GPR17
<b>Species Reactivity:</b>	(+) Human, mouse, and rat; other species not tested
<b>Uniprot No.:</b>	Q13304
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Storage Buffer:</b>	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
<b>Stability:</b>	≥3 years
<b>Host:</b>	Rabbit
<b>Applications:</b>	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:40 and 1:200, respectively. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



Lane 1: C6 (mouse) cell lysate (30 µg)  
Lane 2: HUVEC cell lysate (30 µg)  
Lane 3: HL60 cell lysate (25 µg)



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) human brain tissue after heat induced antigen retrieval in pH 6.0 citrate buffer. After incubation with GPR17 (C-term) Polyclonal Antibody (Item No. 10136) at a 1:40 dilution, slides were incubated with biotinylated secondary antibody, followed by alkaline phosphatase-streptavidin and chromogen (DAB).

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 via 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, 11/02/2023

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 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 peptide affinity-purified polyclonal antibody 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> Post-translational modifications may explain the observed SDS-PAGE gel-migration to 57 kDa on immunoblot.

## 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(19)**, 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