

PRODUCT INFORMATION



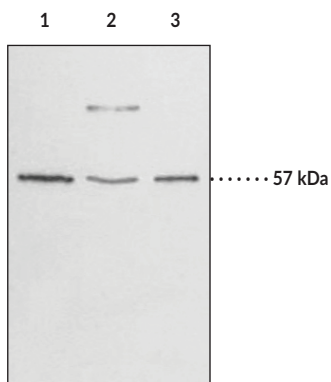
GPR17 (C-Term) Polyclonal Antibody

Item No. 10136

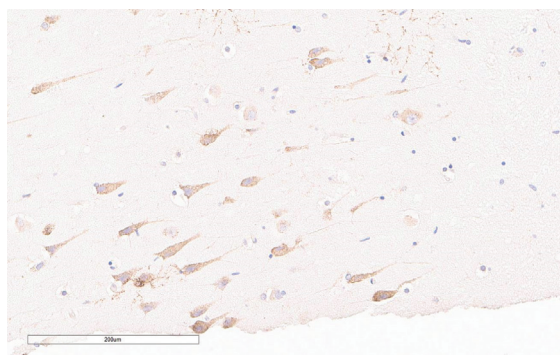
Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified antibody.
Synonym:	G Protein-Coupled Receptor 17
Immunogen:	A synthetic peptide from the C-terminal region of human GPR17
Species Reactivity:	(+) Human, mouse, and rat; other species not tested
Uniprot No.:	Q13304
Form:	Liquid
Storage:	-20°C (as supplied)
Storage Buffer:	PBS, pH 7.2, with 50% glycerol, and 0.02% sodium azide
Stability:	≥1 year
Host:	Rabbit
Applications:	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, 08/13/2019

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.¹ 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.¹ 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.¹ 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**, 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