

PRODUCT INFORMATION



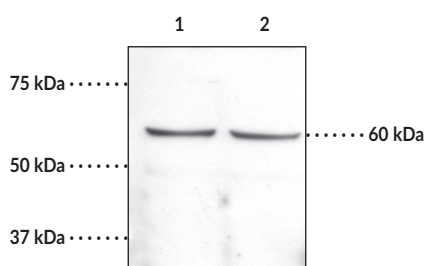
CB₁ Receptor (C-Term) Polyclonal Antibody

Item No. 10006590

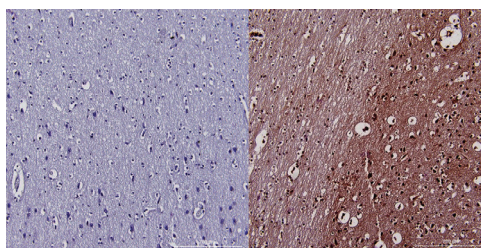
Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified IgG
Synonyms:	Cannabinoid Receptor 1, CNR1
Immunogen:	Synthetic peptide from the C-terminal region of human protein CB ₁ receptor
Species Reactivity:	(+) Human, mouse, rat
Uniprot No.:	P21554
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Applications:	Immunofluorescence (IF), Immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution for IF and IHC is 1:100 and 1:200 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

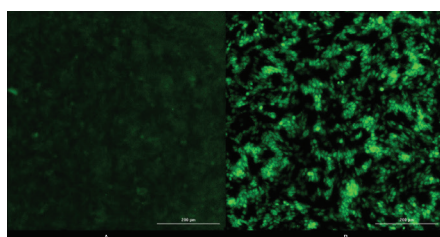
Images



Lane 1: RAW cell pellet (25 µg)
Lane 2: RAW cell supernatant (50 µg)



Immunohistochemical staining of formalin-fixed and paraffin-embedded human brain tissue after heat-induced antigen retrieval in citrate buffer, pH 6.0, using CB₁ Receptor (C-Term) Polyclonal Antibody (Item No. 10006590) at a dilution of 1:100 (left panel, secondary alone)



Immunofluorescent staining of SH-SY5Y cells. SH-SY5Y cells were fixed with 3.7% PFA and blocked with 1% FBS in PBS. Cells were probed with an anti-rabbit FITC secondary antibody alone (A) or CB₁ Receptor (C-Term) Polyclonal Antibody (Item No. 10006590) (B) at a dilution of 1:100, which was followed by Goat Anti-Rabbit IgG FITC (Item No. 10006588)

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, 12/11/2023

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

The CB₁ receptor is a G protein-coupled receptor that binds the active component of cannabis, Δ⁹-tetrahydrocannabinol. This antibody has been raised against the C-terminal (amino acids 461-472) intracellular region of the human CB₁ receptor.^{1,2} Human and rat CB₁ receptors exhibit 97.3% homology at the amino acid level over the complete protein, and 100% homology within the peptide sequence used to make this antibody.^{3,4} This peptide exhibits no homology with the CB₂ receptor. Based on the amino acid sequence, the CB₁ receptor has a molecular weight of approximately 52,800.⁴ The CB₁ receptor and the splice variant CB_{1a} are localized mainly in the brain, whereas the CB₂ receptor is localized predominantly in peripheral tissues, including the spleen and hemopoietic cells.³⁻⁶

References

1. Howlett, A.C., Song, C., Berglund, B.A., *et al.* Characterization of CB₁ cannabinoid receptors using receptor peptide fragments and site-directed antibodies. *Mol. Pharmacol.* **53(3)**, 504-510 (1998).
2. McIntosh, H.H., Song, C., and Howlett, A.C. CB₁ cannabinoid receptor: Cellular regulation and distribution in N18TG2 neuroblastoma cells. *Mol. Brain Res.* **53(1-2)**, 163-173 (1998).
3. Gérard, C.M., Mollereau, C., Vassart, G., *et al.* Molecular cloning of a human cannabinoid receptor which is also expressed in testis. *Biochem. J.* **279(Pt 1)**, 129-134 (1991).
4. Matsuda, L.A., Lolait, S.J., Brownstein, M.J., *et al.* Structure of a cannabinoid receptor and functional expression of the cloned cDNA. *Nature* **346(6284)**, 561-564 (1990).
5. Shire, D., Carillon, C., Kaghad, M., *et al.* An amino-terminal variant of the central cannabinoid receptor resulting from alternative splicing. *J. Biol. Chem.* **270(8)**, 3726-3731 (1995).
6. Shire, D., Calandra, B., Rinaldi-Carmona, M., *et al.* Molecular cloning, expression and function of the murine CB₂ peripheral cannabinoid receptor. *Biochim. Biophys. Acta* **1307(2)**, 132-136 (1996).

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