

# PRODUCT INFORMATION



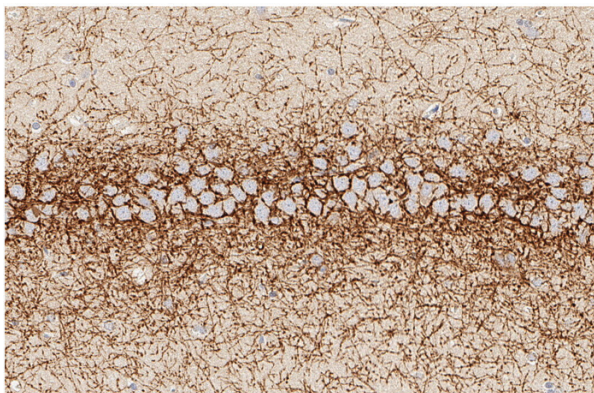
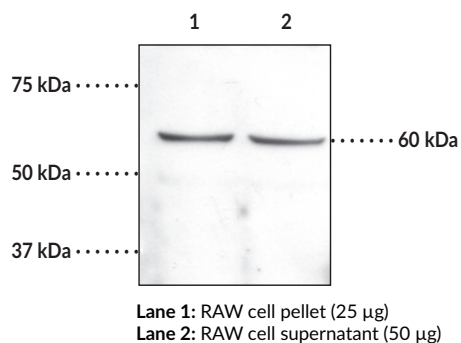
## CB<sub>1</sub> Receptor (C-Term) Polyclonal Antibody

Item No. 10006590

### Overview and Properties

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

### Images



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) rat brain tissue after heat induced antigen retrieval in pH 6.0 citrate buffer. After incubation with CB<sub>1</sub> receptor (C-term) polyclonal antibody (Item No. 10006590), at a 1:200 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, 09/17/2018

**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 CB<sub>1</sub> receptor is a G protein-coupled receptor that binds the active component of cannabis,  $\Delta^9$ -tetrahydrocannabinol. This antibody has been raised against the C-terminal (amino acids 461-472) intracellular region of the human CB<sub>1</sub> receptor.<sup>1,2</sup> Human and rat CB<sub>1</sub> 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.<sup>3,4</sup> This peptide exhibits no homology with the CB<sub>2</sub> receptor. Based on the amino acid sequence, the CB<sub>1</sub> receptor has a molecular weight of approximately 52,800.<sup>4</sup> The CB<sub>1</sub> receptor and the splice variant CB<sub>1a</sub> are localized mainly in the brain, whereas the CB<sub>2</sub> receptor is localized predominantly in peripheral tissues, including the spleen and hemopoietic cells.<sup>3-6</sup>

## References

---

1. Howlett, A.C., Song, C., Berglund, B.A., *et al.* Characterization of CB<sub>1</sub> cannabinoid receptors using receptor peptide fragments and site-directed antibodies. *Mol. Pharmacol.* **53**, 504-510 (1998).
2. McIntosh, H.H., Song, C., and Howlett, A.C. CB<sub>1</sub> cannabinoid receptor: Cellular regulation and distribution in N18TG2 neuroblastoma cells. *Mol. Brain Res.* **53**, 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**, 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**, 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**, 3726-3731 (1995).
6. Shire, D., Calandra, B., Rinaldi-Carmona, M., *et al.* Molecular cloning, expression and function of the murine CB<sub>2</sub> peripheral cannabinoid receptor. *Biochim. Biophys. Acta* **1307**, 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