

PRODUCT INFORMATION



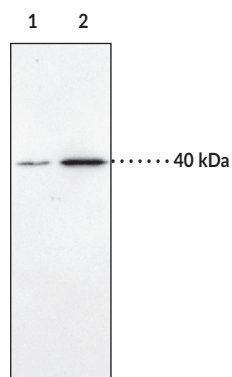
CB₂ Receptor Polyclonal Antibody

Item No. 101550

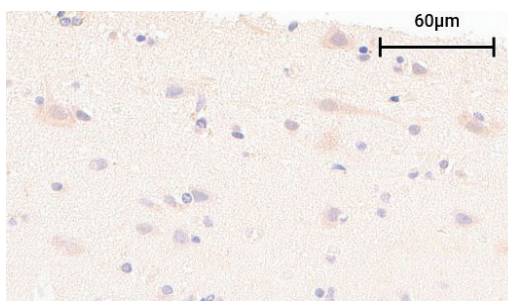
Overview and Properties

Contents: This vial contains 500 µl of peptide-affinity purified polyclonal antibody.
Synonyms: Cannabinoid Receptor 2, CNR₂
Immunogen: Synthetic peptide from the N-terminal region of human CB₂ receptor
Cross Reactivity: (+) CB₂ receptor
Species Reactivity: (+) Human and mouse; other species not tested
Uniprot No.: P34972
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: Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:50 for IHC and 1:200 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: Jurkat (human T cell leukemia) lysate (50 µg)
Lane 2: Jurkat (human T cell leukemia) lysate (100 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 CB₂ Receptor Polyclonal Antibody (Item No. 101550) 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, 11/01/2023

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 CB₂ receptor is localized predominantly in peripheral tissues, including the spleen and hemopoietic cells.¹ The CB₁ and CB₂ receptors are G-protein coupled receptors that bind the active component of cannabis, Δ⁹-tetrahydrocannabinol, as well as anandamide which is an endogenous CB receptor ligand. This antibody has been raised against a sequence between the N-terminus and the first transmembrane domain of the protein of the human CB₂ receptor.¹ It can be used for western blotting and immunohistochemistry applications. Conserved amino acids between the CB₁ and CB₂ receptors in this region are minimal.² Human and murine CB₂ receptors exhibit 82% homology at the amino acid level over the complete protein.² The CB₂ receptor is localized predominantly in peripheral tissues, including the spleen and hemopoietic cells.¹

References

1. Munro, S., Thomas, K.L., and Abu-Shaar, M. Molecular characterization of a peripheral receptor for cannabinoids. *Nature* **365(6441)**, 61-65 (1993).
2. 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