

PRODUCT INFORMATION



CB₂ Receptor Polyclonal FITC Antibody

Item No. 10010712

Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified antibody.
Synonyms:	Cannabinoid Receptor 2, CNR ₂
Immunogen:	Synthetic peptide from the N-terminal region of human CB ₂ receptor
Species Reactivity:	(+) Human; other species not tested
Uniprot No.:	P34972
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Applications:	Flow cytometry (FC) and immunofluorescence (IF); the recommended starting dilution for FC and IF is 1:40. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Description

The central cannabinoid (CB₁) and peripheral cannabinoid (CB₂) receptors are G protein-coupled receptors (GPCRs) 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-terminal and the first transmembrane domain of the protein of the human CB₂ receptor.¹ Conserved amino acids between the CB₁ and CB₂ receptors in this region are minimal.² Human and mouse CB₂ receptors exhibit 82% homology at the amino acid level over the complete protein sequence.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).

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, 10/25/2019

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