

Product Information



CB₂ Receptor Polyclonal FITC Antibody

Item No. 10010712 • Lot. No. XXXXXX

Contents:	This vial contains <i>lot specific</i> µg peptide affinity-purified IgG in <i>lot specific</i> µl PBS, pH 7.2, containing 50% glycerol, 0.5 mg/ml BSA, and 0.02% sodium azide
Synonym:	Cannabinoid Receptor 2
Antigen:	Synthetic peptide from human CB ₂ amino acids 20-33 (NPMKDYMILSGPQK); the antigen alignment with other known sequences is as follows: Human N P M K D Y M I L S G P Q K Mouse N P M K e Y M I L S s g Q q Human CB ₁ - d i e c f M v L n p s Q q Rat CB ₁ - d m e c f M I L n p s Q q
Host:	Rabbit
Cross Reactivity:	(+) Human and mouse CB ₂ receptor
Stability:	≥1 year at -20°C
Applications:	The recommended starting dilution for flow cytometry and immunofluorescence is 5 µg/ml. Cayman's CB ₂ Receptor Polyclonal Antibody (Item No. 101550) also is intended for western blot, however this product works well for that application at 2 µg/ml. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

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 thereby preventing cross-reactivity of this antibody with the CB₁ receptor.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**, 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**, 132-136 (1996).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10010712

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/19/2011

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com