

# Cannabinoid receptor

## CB<sub>2</sub> polyclonal antibody

The cannabinoid receptors CB1 and CB2 are GPCRs named for their signaling function in response to cannabinoid drugs such as Delta9-tetrahydrocannabinol (Delta9-THC). In addition to Delta9-THC and its synthetic analogs, CB1 and CB2 receptors are bound by endogenous eicosanoid ligands (e.g. anandamide), and signal via Gi/o alpha subunits to regulate adenylyl cyclase activity (generally inhibitory), ion channel activation, intracellular Ca<sup>2+</sup> levels, nitric oxide synthase activation, and immediate early gene expression. CB1 and CB2 signals in turn regulate neurological and immunological functions such as inhibition of neurotransmitter release and immune cell migration, antigen processing, and cytokine production.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

### Ordering Information

[Order Online »](#)

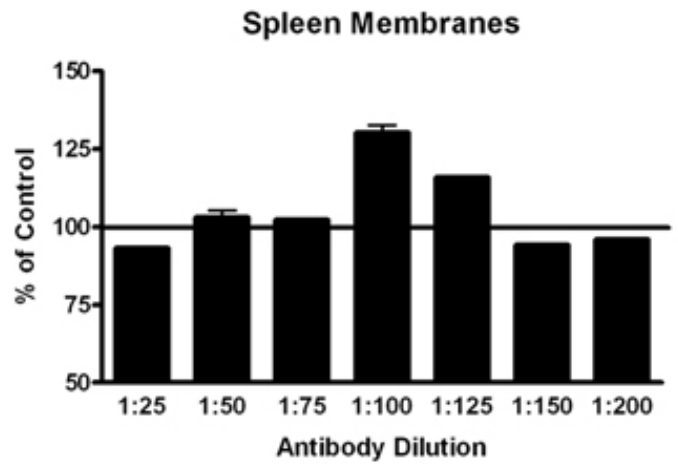
ADI-905-749-100	100µg
-----------------	-------

### Manuals, SDS & CofA

[View Online »](#)



Western blot analysis: MW marker (1) and 10µg rat brain extract (2) probed with Cannabinoid Receptor (CB2) pAb at 4µg/ml.



Membrane ELISA: Lewis rat spleen membranes (5µg/well) were treated with 1 µM concentrations of agonist ( $\Delta^9$ -THC) and probed with Cannabinoid Receptor (CB2) pAb (1:25 to 1:200 of a 0.1µg/µl stock solution) by ELISA. Data from vehicle treated cells were taken as 100%. Results are the mean  $\pm$  SEM (n=2).

## Handling & Storage

**Handling** Keep on ice at all times.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** CB2, CNR2

**Application** ELISA, IHC, WB

**Application Notes** Suitable for cell-based and membrane-based ELISA.  
Predicted MW of ~35kDa. Higher MW species (~50kDa) and intermediates also observed by Western blot which reflect post-translational modification of the receptor.

**Crossreactivity** Predicted species reactivity based on epitope sequence homology: human, mouse, bovine, horse, rhesus macaque, dog, chimpanzee.

**Formulation** Liquid. In PBS containing 50% glycerol and 0.01% sodium azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to the sequence near the N-terminus of rat cannabinoid receptor CB2.

**Purity Detail** Peptide affinity purified.

**Species Reactivity** Rat

**UniProt ID** Q9QZN9

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)