

# Cannabinoid receptor 1 polyclonal antibody

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

Citations: 1

[View Online »](#)

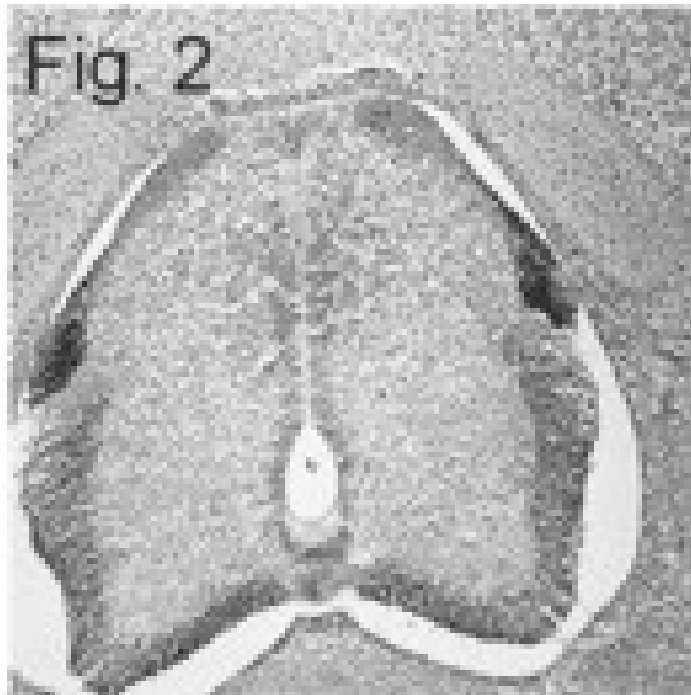
## Ordering Information

[Order Online »](#)

ALX-210-316-R100	100µl
------------------	-------

Manuals, SDS & CofA

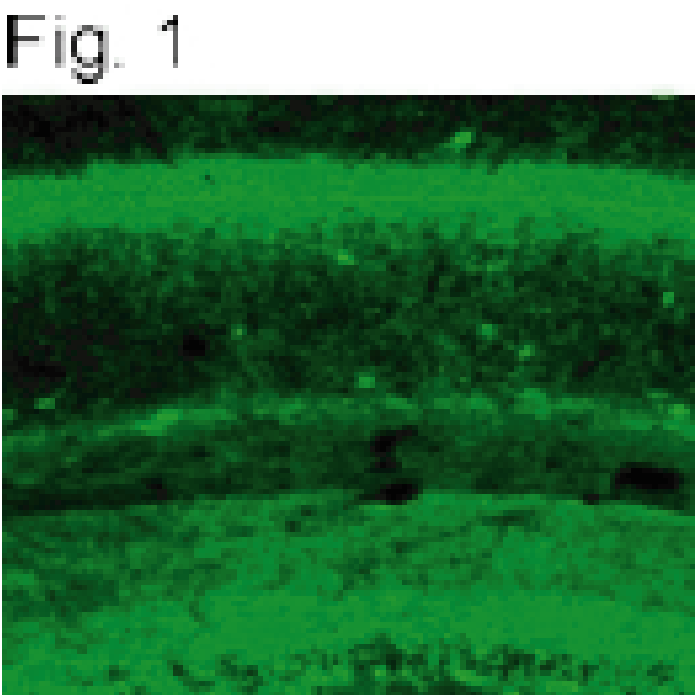
[View Online »](#)



**Figure 2:** Immunohistochemistry of CB<sub>1</sub> in transverse spinal cord sections of E6 chick embryo using PAb to Cannabinoid Receptor 1 (Prod. No. [ALX-210-316](#)).  
 Cannabinoid receptor 1 polyclonal antibody  
 Immunohistochemistry chick, Cannabinoid receptor 1 polyclonal antibody Immunohistochemistry chick, yes  
 ALX-210-327, 210-327.gif, <http://static.enzolifesciences.com/fileadmin/files/image/210-327.gif>,

**Figure:** Western blot analysis of XIAP expression in Jurkat cell lysate using PAb to XIAP (Prod. No. [ALX-210-327](#)).

“””””””””



**Figure 1:** Immunolocalization of CB<sub>1</sub> in rat hippocampus using PAb to Cannabinoid Receptor 1 (Prod. No. ALX-210-316).

## Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	CB1, CNR1, CBR
Application	Flow Cytometry, ICC, IHC (PS), WB
Application Notes	Detects a band of ~60kDa by Western blot representing CB <sub>1</sub> ; also detects less intense bands of ~23kDa, ~72kDa and ~180kDa from rat brain homogenate.
Formulation	Liquid. In PBS containing 50% glycerol, 1mg/ml BSA and 0.05% sodium azide.
Host	Rabbit
Immunogen	Recombinant rat CB1 (cannabinoid receptor 1) (aa 1-77).
Purity Detail	Affinity purified IgG.
Recommendation Dilutions/Conditions	Immunocytochemistry (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Amphibian, Chicken, Fish, Human, Monkey, Mouse, Rat
UniProt ID	P20272
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



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-eu@enzolifesciences.com)

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