

Cannabinoid receptor 1 polyclonal antibody

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

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

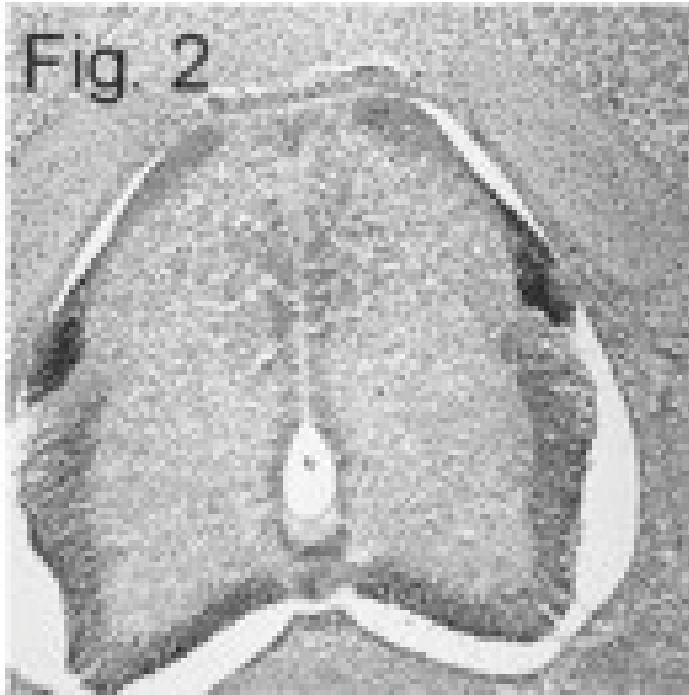


Figure 2: Immunohistochemistry of CB₁ in transverse spinal cord sections of E6 chick embryo using PAb to Cannabinoid Receptor 1 (Prod. No. [ALX-210-316](#)). Cannabinoid receptor 1 polyclonal antibody Immunohistochemistry chick, Cannabinoid receptor 1 polyclonal antibody Immunohistochemistry chick, yes 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](#)).

“”””””

Fig. 1

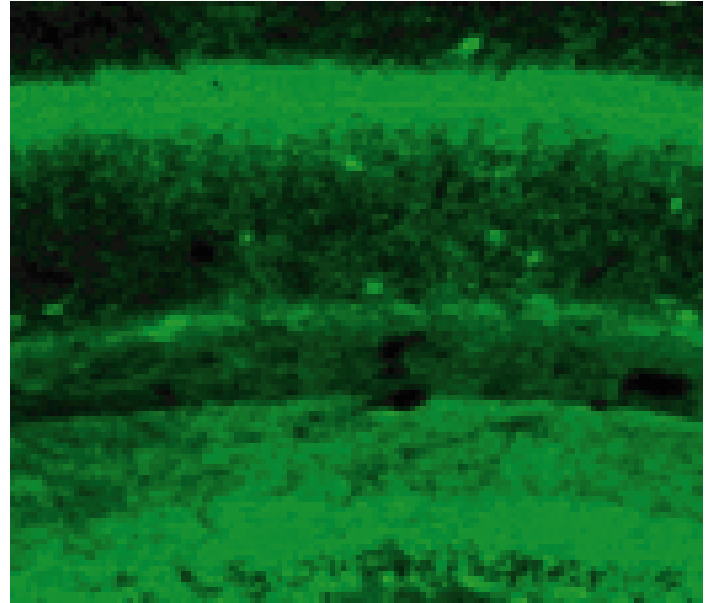


Figure 1: Immunolocalization of CB₁ in rat hippocampus using PAb to Cannabinoid Receptor 1 (Prod. No. ALX-210-316).

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 CB₁; 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 [Worry-Free Guarantee](#)



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

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

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

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

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

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

Last modified: May 29, 2024