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 »

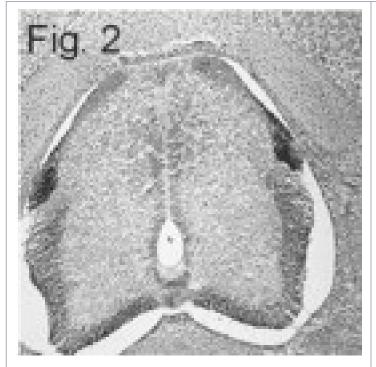


Fig. 1

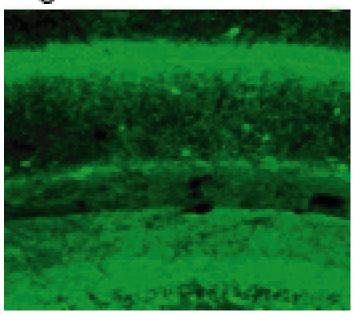


Figure 2: Immunohistochemistry of CB₁ in transverse spinal cord sections of E6 chick embryo using PAb to Cannabinoid Receptor 1 (Prod. No. <u>ALX-210-316</u>).""",Cannabinoid receptor 1 polyclonal antibody Immunohistochemistry chick,Cannabinoid receptor 1 polyclonal antibody Immunohistochemistry chick,yes 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. <u>ALX-210-327</u>).

"""""

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

Detects a band of ~60kDa by Western blot representing **Application Notes**

CB₄; 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.

Rabbit Host

Recombinant rat CB1 (cannabinoid receptor 1) (aa 1-77). **Immunogen**

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

Phone: 800.942.0430 Last modified: May 29, 2024 om

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 eu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 fr@enzolifesciences.com

Phone: +49 7621 5500 526 de@enzolifesciences.com

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