Cannabinoid receptor 2 polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 3

View Online »

Ordering Information

Order Online »

ALX-210-315-R100

100µl

Manuals, SDS & CofA

View Online »

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 CB2, CNR2

Application Flow Cytometry, ICC, IHC (FS), IHC (PS), WB

Application NotesBy Western blot detects bands of ~60kDa from human

spinal cord homogenate and ~40kDa from RBL2H3 cells,

representing different glycosylated forms of CB₂.

Formulation Liquid. In PBS containing 50% glycerol, 1mg/ml BSA and

0.05% sodium azide.

Host Rabbit

Immunogen Recombinant human CB2 (cannabinoid receptor 2) (aa 1-

33).

Purity Detail Epitope-affinity purified IgG.

Recommendation Dilutions/Conditions Flow Cytometry (1:50)Immunocytochemistry

(1:100)Immunohistochemistry (paraffin sections,

1:100)Western Blot (1:250)Suggested dilutions/conditions may not be available for all applications.Optimal conditions

must be determined individually for each application.

Species Reactivity Human, Rat

UniProt ID P34972

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com