

Cannabinoid receptor 2 polyclonal antibody

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

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-3150	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 Notes By 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



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-
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)