

MRP2 monoclonal antibody (M2III-5)

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

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-801-0370	250µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Multidrug resistance-associated protein 2, cMOAT, ABCC2

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

Application Notes Immunocytochemistry: cytospin preparations
Immunohistochemistry: Does not recognize mouse MRP2 in IHC.

Clone M2III-5

Crossreactivity Does not cross-react with human P-glycoprotein, MDR3 P-glycoprotein, MRP1, MRP3 or MRP5.

Formulation Liquid. In serum free culture supernatant containing 0.7% BSA and 0.1% sodium azide.

Host Mouse

Immunogen Recombinant C-terminal rat MRP2 (multidrug resistance-associated protein 2; cMOAT) (aa 1339-1541).

Isotype IgG2b

Recommendation Dilutions/Conditions

Immunocytochemistry: Use 1:20-1:50 dilution on acetone fixed cytospin preparations. Immunohistochemistry (frozen sections, paraffin sections): Use 1:20 dilution on acetone fixed frozen sections followed by incubation with rabbit anti-mouse IgG and a monoclonal mouse APAAP complex or on paraffin sections after pretreatment with 0.01 M citric acid, pH 6.0 in distilled water at 100°C for 5 min 3 times. After washing, slides can be incubated with biotinylated rabbit anti-mouse IgG and streptavidin conjugated to horseradish peroxidase. Western Blot: Use 1:20-1:50 dilution and anti-mouse-HRP. Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity

Human, Mouse, Rat

Specificity

Recognizes an internal epitope of MRP2.

UniProt ID

Q63120

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)