

MRP2 monoclonal antibody (M2III-6)

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

Citations: 19

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-801-016A-C100	100µg
-------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
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
Clone	M2III-6
Crossreactivity	Does not cross-react with human P-glycoprotein, MDR3, MRP1, MRP3 or MRP5.
Formulation	Liquid. In PBS containing 0.1% BSA and 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant C-terminal rat MRP2 (multidrug resistance-associated protein-2; cMOAT) (aa 1339-1541).
Isotype	IgG2a

Recommendation Dilutions/Conditions

Immunocytochemistry: Use 1:20-1:50 dilution on acetone fixed cytospin preparations. Immunohistochemistry (frozen sections) (1:20): Can be use on acetone fixed frozen sections. Can be followed by incubation with rabbit anti-mouse antibody and a monoclonal mouse APAAP complex. Immunohistochemistry (paraffin sections) (1:20, weak): Can be used on formaldehyde-fixed paraffin-embedded tissues and tumors, after pretreatment with 0.01M citric acid (pH 6.0) in distilled water at 100°C for 5 min 3 times. After incubation with the primary antibody and washing, slides can be incubated with biotinylated rabbit anti-mouse antibody and streptavidin conjugated to horseradish peroxidase. Western Blot: Use 1:20-1:50 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, 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)