

# MRP2 monoclonal antibody (M2III-6)

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

Citations: 20

[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)