

# MRP1 (human) monoclonal antibody (MRPm6)

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

Citations: 25

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-801-013-C250	250µg
ALX-801-013-C500	500µg

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 1, ABCC1

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

Application Notes Flow Cytometry: cell permeabilization required  
Immunocytochemistry: cytospin preparations  
Immunohistochemistry: Suitable for frozen sections, weak in paraffin sections.

Clone MRPm6

Crossreactivity Does not cross-react with human P-glycoprotein, MDR3 P-glycoprotein or human MRP2 (cMOAT) to MRP5.

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

Host Mouse

Immunogen Recombinant human MRP1 (multidrug resistance-associated protein 1) (aa 1294-1531).

Isotype IgG1

Recommendation Suggested starting dilution is 1:20Suggested dilutions/conditions may not be available  
Dilutions/Conditions for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human

Specificity Recognizes the internal epitope (P<sup>1511</sup>SDLLQQRGL<sup>1520</sup>) of human MRP1.

UniProt ID P33527

Worry-free Guarantee

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



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)