

# MRP1 (human) monoclonal antibody (MRPm6)

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

Citations: 25

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-801-0130	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: cyospin 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 Dilutions/Conditions** Suggested starting dilution is 1:20. Suggested dilutions/conditions may not be available 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](#)

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)