

# MRP1 (human) monoclonal antibody (QCRL-1)

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

Citations: 16

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-801-010-L001	1ml
------------------	-----

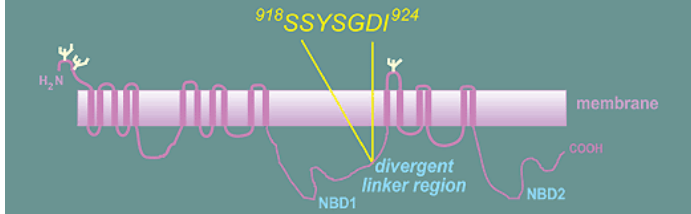
Manuals, SDS & CofA

[View Online »](#)

## Human MRP1-Specific MAb QCRL-1

Hipfner *et al.*, Cancer Res. 56: 3307 (1996)

- generated against H69AR cell membranes
- IgG<sub>1</sub>; MRP1 detection by blotting, cytochemistry, precipitation, flow cytometry and histochemistry of formalin-fixed tissues
- cytoplasmic heptapeptide epitope **NOT** found in any other ABC transporter including mouse mrp1



**Figure:** Human MRP1-specific MAb QCRL-1 (Prod.No. [ALX-801-010](#)). """,MRP1 (human) monoclonal antibody (QCRL-1) image,MRP1 (human) monoclonal antibody (QCRL-1) image,yes

ALX-801-013,801-

013.gif,[http://static.enzolifesciences.com/fileadmin/files/image/801-](http://static.enzolifesciences.com/fileadmin/files/image/801-013.gif)

013.gif,**Figure:** Epitope Mapping of MAbs MRPr1

## Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
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 (PS), IP, WB
Application Notes	Flow Cytometry: cell permeabilization required Immunocytochemistry: cytospin preparations
Clone	QCRL-1
Crossreactivity	Does not cross-react with human P-glycoprotein, MDR3 P-glycoprotein or human MRP2 (cMOAT) to MRP5, or mouse MRP gene products.
Formulation	Liquid. 0.22µm-filtered solution in PBS containing 0.7% BSA and 0.1% sodium azide.
Host	Mouse
Immunogen	Membranes prepared from the human small cell lung cancer cell line H69AR.
Isotype	IgG1
Recommendation Dilutions/Conditions	Flow Cytometry (1:20)Immunocytochemistry (1:20-1:50)Western Blot (0.02-0.2µg/ml)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 linear internal epitope (S <sup>918</sup> SYSGDI <sup>924</sup> ) of human MRP1 (multidrug resistance-associated protein 1), overexpressed in various human non-P-glycoprotein MDR tumor cell lines.

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)