

MRP1 monoclonal antibody (MRPr1)

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

Citations: 51

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-801-0070	125µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

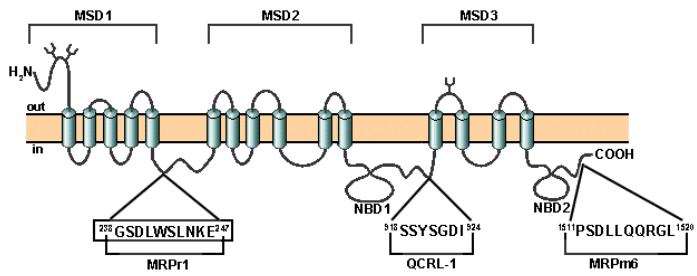


Figure: Epitope Mapping of MAbs MRPr1, QCRL-1 & MRPr6. Topological model of MRP depicting the approx. locations of the MRPr1, QCRL-1 and MRPr6 epitopes. The model shown is only one of several possible configurations. MSD=membrane spanning domain; NBD=nucleotide binding domain. *From R. Hipfner, et al.; Br. J. Cancer* **78**,1134 (1998).

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), IHC (PS), WB
Application Notes	Flow Cytometry: cell permeabilization required Immunocytochemistry: cytospin preparations
Clone	MRPr1
Crossreactivity	Does not cross-react with human P-glycoprotein, MDR3 P-glycoprotein or human MRP2 (cMOAT) to MRP 5.
Formulation	Liquid. In serum free culture supernatant containing 0.7% BSA and 0.1% sodium azide.
Host	Rat
Immunogen	Recombinant human MRP1 (multidrug resistance-associated protein 1) (aa 192-360).
Isotype	IgG2a
Recommendation Dilutions/Conditions	Suggested starting dilution is 1:20Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse
Specificity	Recognizes the internal epitope (G ²³⁸ SDLWSLNKE ²⁴⁷) of MRP1, 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](#)

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)