

MRP5 (human) monoclonal antibody (M5I-1)

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

Citations: 2

[View Online »](#)

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 5, ABCC5 |
| Application | Flow Cytometry, ICC, IHC (FS), WB |
| Application Notes | Flow Cytometry: weak, cell permeabilization required Immunocytochemistry: cytospin preparations |
| Clone | M5I-1 |
| Crossreactivity | Does not cross-react with human P-glycoprotein, MDR3 P-glycoprotein, MRP1, MRP2 (cMOAT) or MRP3. |
| Formulation | Liquid. In serum free culture supernatant containing 0.7% BSA and 0.1% sodium azide. |
| Host | Rat |
| Immunogen | Recombinant human MRP5 (multidrug resistance-associated protein 5) (aa 82-168). |
| Isotype | IgG2a |

Recommendation Dilutions/Conditions

Immunocytochemistry: Use 1:20-1:50 dilution on acetone fixed cytospin preparations. Immunohistochemistry (frozen sections, weak): Use 1:20 dilution on acetone fixed frozen sections followed by incubation with rabbit anti-rat Ig and a monoclonal rat APAAP complex. Alternatively, after incubation with the primary antibody and washing, slides can be incubated with biotinylated rabbit anti-rat Ig and streptavidin conjugated to horseradish peroxidase. Western Blot: Use 1:20-1:50 dilution and anti-rat-HRP. Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity

Human

Specificity

Recognizes an internal epitope of human MRP5.

UniProt ID

O15440

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)