

MRP4 monoclonal antibody (M4I-10)

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

Citations: 17

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-801-038-C075	75µg
ALX-801-038-C150	150µg

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 4, ABCC4

Application ICC, IHC (FS), WB

Application Notes Immunocytochemistry: cytospin preparations

Clone M4I-10

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

Host Rat

Immunogen Recombinant human MRP4 (multidrug resistance-associated protein 4) (aa 372-431).

Isotype IgG2a

Recommendation Dilutions/Conditions Immunocytochemistry: Use 1:20-1:50 dilution on acetone fixed cytospin preparations.Immunohistochemistry (frozen sections): Use 1:20 dilution on acetone fixed frozen sections followed by incubation with biotin-labelled rabbit anti-rat IgG and HRP-labelled streptavidin.Western Blot: Use 1:20-1:50 dilution and HRP-anti-rat-IgG.Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse

UniProt ID O15439

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)