

P-glycoprotein (human) monoclonal antibody (MRK16)

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

Citations: 25

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-801-008-C150	150µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	ABCB1, CD243, P-glycoprotein 1
Application	Flow Cytometry, FUNC, IHC (FS), IP
Application Notes	Functional Application: Inhibits P-glycoprotein function. Not recommended for Western blot (epitope recognized by this antibody may be lost in the presence of protein denaturing agents such as SDS).
Clone	MRK16
Crossreactivity	Does not cross-react with mouse P-glycoprotein.
Formulation	Liquid. In PBS containing 1% BSA and 0.1% sodium azide.
Host	Mouse
Immunogen	Adriamycin-resistant human myelogenous leukemia K-562/ADM cells.
Isotype	IgG2a
Purity Detail	Protein A affinity purified.

Recommendation Dilutions/Conditions

Flow Cytometry: For 5×10⁵ pelleted cells use 50µl of 10-100µg/ml. Immunohistochemistry (frozen sections): 0.1-10µg/ml. Use mild fixation treatment such as 4% paraformaldehyde. Do not use acetone, paraffin, methanol or ethanol as a fixative. Iodine labelling can also be performed on MRK16. Immunoprecipitation: 25-50µg/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 an external epitope of human P-glycoprotein. Also recognizes P-glycoprotein of human cortical adenomas.

UniProt ID

P08183

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com