

# 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 5×10<sup>5</sup> 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

<b>Specificity</b>	Recognizes an external epitope of human P-glycoprotein. Also recognizes P-glycoprotein of human cortical adenomas.
<b>UniProt ID</b>	P08183
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



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)