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.

lgG2a

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/mlSuggested 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