MRP2 (human) monoclonal antibody (M2I-4)

This antibody is covered by our Worry-Free Guarantee.

Citations: 9

View Online »

Ordering Information

Order Online »

ALX-801-015-C250

250µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Multidrug resistance-associated protein 2, cMOAT, ABCC2

Application ICC, IHC (FS), WB

Application Notes Immunohistochemistry: Suitable for frozen sections, weak

in paraffin sections.

Clone M2I-4

Crossreactivity Does not cross-react with human P-glycoprotein, MDR3 P-

glycoprotein, MRP1, MRP3 or MRP5. Does not cross-react

with mouse, rat or guinea pig MRP2.

Formulation Liquid. In serum free culture supernatant containing 0.7%

BSA and 0.1% sodium azide.

Host Mouse

Immunogen Recombinant human MRP2 (multidrug resistance-

associated protein 2; cMOAT) (aa 215-310).

lsotype lgG1

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 rabbit anti-mouse IgG and a monoclonal mouse APAAP complex. Or, after incubation with the primary antibody and washing, slides can be incubated with biotinylated rabbit anti-mouse IgG and streptavidin conjugated to horseradish peroxidase.Western Blot: Use 1:20-1:50 dilution and anti-mouse-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 MRP2.

UniProt ID

Q92887

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024