MRP5 (human) monoclonal antibody (M5I-1)

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

Ordering Information

Order Online »

ALX-801-021-C250

250µg

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 5, ABCC5

Application Flow Cytometry, ICC, IHC (FS), WB

Application Notes Flow Cytometry: weak, cell permeabilization required

Immunocytochemistry: cytospin preparations

Clone M5I-1

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

glycoprotein, MRP1, MRP2 (cMOAT) or MRP3.

Formulation Liquid. In serum free culture supernatant containing 0.7%

BSA and 0.1% sodium azide.

Host Rat

Immunogen Recombinant human MRP5 (multidrug resistance-

associated protein 5) (aa 82-168).

lsotype lgG2a

Recommendation Dilutions/Conditions

Immunocytochemistry: Use 1:20-1:50 dilution on acetone fixed cytospin preparations. Immunohistochemistry (frozen sections, weak): Use 1:20 dilution on acetone fixed frozen sections followed by incubation with rabbit anti-rat Ig and a monoclonal rat APAAP complex. Alternatively, after incubation with the primary antibody and washing, slides can be incubated with biotinylated rabbit anti-rat Ig and streptavidin conjugated to horseradish peroxidase. Western Blot: Use 1:20-1:50 dilution and anti-rat-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 MRP5.

UniProt ID

O15440

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

eu@enzolifesciences.com