# MRP2 monoclonal antibody (M2III-5)

This antibody is covered by our Worry-Free Guarantee.

Citations: 8

View Online »

## **Ordering Information**

Order Online »

ALX-801-037-C125	125µg
ALX-801-037-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 2, cMOAT, ABCC2

**Application** ICC, IHC (FS), IHC (PS), WB

Application Notes Immunocytochemistry: cytospin preparations

Immunohistochemistry: Does not recognize mouse MRP2

in IHC.

Clone M2III-5

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

glycoprotein, MRP1, MRP3 or MRP5.

**Formulation** Liquid. In serum free culture supernatant containing 0.7%

BSA and 0.1% sodium azide.

**Host** Mouse

Immunogen Recombinant C-terminal rat MRP2 (multidrug resistance-

associated protein 2; cMOAT) (aa 1339-1541).

**Isotype** IgG2b

#### **Recommendation Dilutions/Conditions**

Immunocytochemistry: Use 1:20-1:50 dilution on acetone fixed cytospin preparations.Immunohistochemistry (frozen sections, paraffin 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 on paraffin sections after pretreatment with 0.01 M citric acid, pH 6.0 in distilled water at 100°C for 5 min 3 times. After 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, Mouse, Rat

Specificity

Recognizes an internal epitope of MRP2.

**UniProt ID** 

Q63120

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

