# MRP1 (human) monoclonal antibody (MRPm6)

This antibody is covered by our Worry-Free Guarantee.

Citations: 25

View Online »

# **Ordering Information**

Order Online »

ALX-801-013-C250	250μg
ALX-801-013-C500	500µg

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 1, ABCC1

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

**Application Notes** Flow Cytometry: cell permeabilization required

Immunocytochemistry: cytospin preparations

Immunohistochemistry: Suitable for frozen sections, weak

in paraffin sections.

MRPm6 Clone

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

glycoprotein or human MRP2 (cMOAT) to MRP5.

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

BSA and 0.1% sodium azide.

Host Mouse

**Immunogen** Recombinant human MRP1 (multidrug resistance-

associated protein 1) (aa 1294-1531).

Isotype lgG1

**Recommendation Dilutions/Conditions** Suggested starting dilution is 1:20Suggested

dilutions/conditions may not be available for all

applications. Optimal conditions must be determined

individually for each application.

Human **Species Reactivity** 

**Specificity** 

Recognizes the internal epitope (P<sup>1511</sup>SDLLQQRGL<sup>1520</sup>) of human MRP1.

**UniProt ID** 

P33527

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-