

TECHNICAL INSTRUCTIONS

Description	Clone	Species	Ig-isotype	Catalogue No.
Monoclonal antibody to the Multidrug Resistance-related Protein MRP1	MRPm5	mouse	IgG2a	SB- MRPm5-MAB

Specificity: MRPm5 reacts with an internal epitope of MRP1, a 180-195 kD transmembrane transporter protein overexpressed in various human non-P-glycoprotein MDR tumor cell lines. MRPm5 was raised against a bacterial fusion protein of MRP1, containing amino acids 986-1204 of the protein. MRPm5 does not cross-react with the human MDR1 and MDR3 gene products.

Applications: Immunoblotting, Immunohistochemistry and Flow cytometry

Immunoblotting: the antibody can be diluted 1:20.

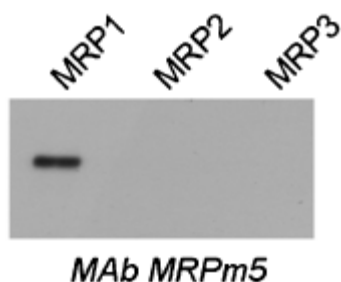


Figure 1. Immunoblots of membrane vesicles enriched for MRP1, MRP2, or MRP3. Membrane vesicles (2 lg/lane) were immunoblotted with MRP1-specific MAb MRPm5.

Immunohistochemical staining: the antibody can be diluted 1:20 for frozen sections, air-dried or acetone fixed cells, paraffin-embedded tissue sections.

Flow cytometry: the antibody can be diluted 1:20 (note: cell permeabilization required).

Note: Optimal conditions must be determined individually for each application.

Presentation: 1 ml vials (>>200 tests) containing antibody in serumfree culture supernatant, with 0.7% BSA (Roche, Mannheim, Germany) and 0.1% Sodium azide. Concentration 250 µg immunoglobulin/ml (by ELISA)

Shipping: Ambient temperature.



Storage: Maintain refrigerated at 2-8°C for up to 6 months. For long-term storage prepare small aliquots and store at -20°C.

References: Flens et al. Cancer Res. 54: 4557-4564, 1994.
Cole et al. Science 258: 1650-1654, 1992.
Nooter et al. Ann oncol 7: 75-81, 1996.
Flens et al. Am J. path. 148: 1237-1247, 1996.
Scheffer et al. Cancer Res 60: 5269, 2000

Correspondence to:

Name: Tünde Nagy, Ph.D.
Production Manager

Mail: **SOLVO Biotechnology**
H-6726 Szeged
Közép fasor 52.
Hungary

E-Mail: t.nagy@solvo.com
Phone: +36 62 424 729
Cell: +36 30 676 8640
Fax: +36 62 426 098

LIMITATIONS: This is a laboratory reagent, not to be administered to humans or animals nor used for any drug purpose.

Safety information about the cell lines and culture media used in the production of the MAb.

MAb producing cells: The hybridoma cell line was obtained by fusion of lymph node cells from an immunized mouse (Balb/c) with SP2/O mouse myeloma cells.

Culture medium: IMDM (BioWhittaker), supplemented with Nutridoma-SP (Boehringer, Indianapolis, USA), without serum or added enzymes. Antibody containing supernatant has been concentrated and filtered through a 0.22 micron filter.

NOTE: this monoclonal antibody has been produced in a clinical laboratory in which no animal viruses are being studied or cultured.

