

TECHNICAL INSTRUCTIONS

Description	Clone	Species	Ig-isotype	Catalogue No.
Monoclonal antibody to human P-glycoprotein (P-gp) multidrug resistance related	JSB-1	mouse	IgG1	SB- JSB-1-MAB

Specificity: JSB-1 was selected after immunization with multidrug resistant Chinese hamster ovary cells (ChrC5) cells. JSB-1 reacts with a conserved cytoplasmic epitope of the plasma membrane-associated 170 kD P-gp, member of the superfamily of transmembrane transporters. JSB-1 detects P-gp overexpression in human tumour cells of all different histogenetic derivations.

Applications: Immunoblotting, Immunocytochemistry, Immunohistochemistry and Flow cytometry

Immunoblotting: Western Blot (Fig. 1)

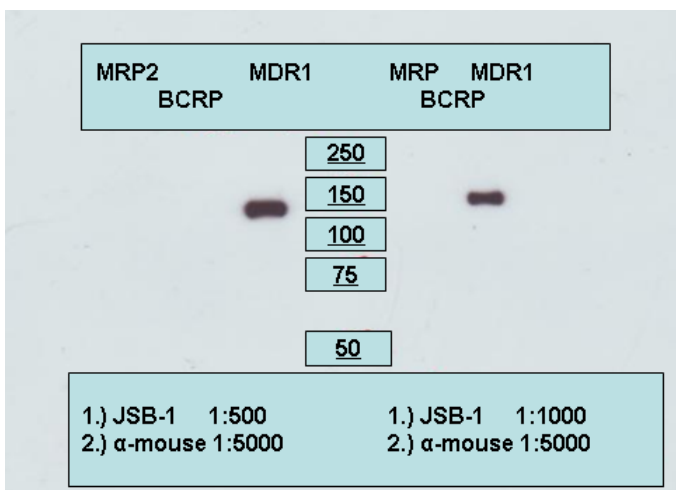


Figure 1. Western Blot detection of P-glycoprotein overexpression using JSB-1 on SOLVO membrane preparation

Immunocytochemistry: cytopsin preparations (Fig. 2)



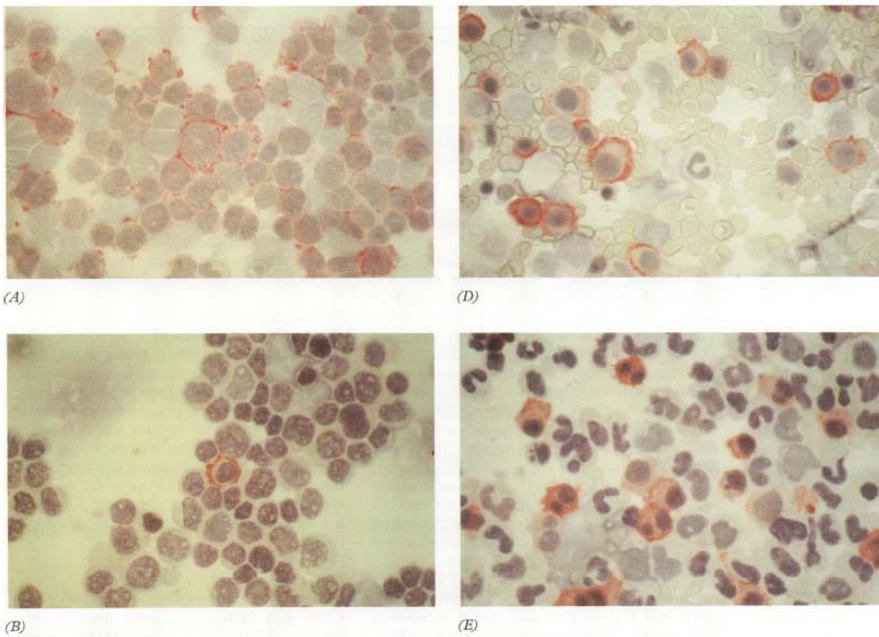


Figure 2.

Immunocytochemical (APAAP) detection of P-glycoprotein overexpression using JSB-1: (A) CEMIVLB8; (B) acute myeloid leukaemia, single positive cell; (C) acute non-lymphoblastic leukaemia-presence of P-gp intracytoplasmic localisation (arrow) (D-E)

Immunohistochemistry: the antibody can be diluted 1:20 for immuno-histochemical staining of frozen sections, air-dried or acetone fixed cells, paraffin embedded tissue sections

Flow cytometry: 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: Scheper RJ et al., Int. J. Cancer 42: 389, 1988;
Dalton WS et al., J. Clin. Oncol. 7: 415, 1989;
Tiirikainen M et al, Ann. Haematol. 65, 124, 1992;
Grogan T et al, Blood 81, 490, 1993;
Itsubo M et al, Cancer 73: 298, 1994;
Toth K et al, Am. J. Pathol 144: 227, 1994;
Gala et al, J. Clin. Pathol. 47: 619, 1994;
Sinicrope et al, Cancer 74: 2908, 1994;
Balaini et al., New Engl. J. Med. 333:1380, 1995;
Chevallard et al, Cancer 77:292, 1996;
Moran et al, Eur. J. Cancer 33: 652, 1997.



Correspondence to:

Name: **Tünde Nagy, Ph.D.**
Production Manager

Mail: **SOLVO Biotechnology**
H-6726 Szeged
Közep 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.

