

PRODUCT INSERT
CYTO-IC MONOCLONAL ANTIBODIES
SPECIFIC TO HUMAN IL-2

Product Code	Form	Volume	Antibody*	Tests	Excitation (nm)	Peak Emission (nm)	Matching Isotype Controls
RHCIL200	Purified	1.0 ml	200 µg		N/A	N/A	Rat IgG2a Purified Code R2a00
RHCIL220	Alexa Fluor ^{®†} 488	0.5 ml		100 min.	488	519	Rat IgG2a Alexa Fluor 488 Code R2a20
RHCIL201	FITC	1.0 ml		100 min.	488	525	Rat IgG2a FITC Code R2a01
RHCIL201-3	FITC	3.0 ml		300 min.			
RHCIL204	R-PE	0.5 ml		50 min.	488	575	Rat IgG2a R-PE Code R2a04
RHCIL204-3	R-PE	3.0 ml		300 min.			
RHCIL205	APC	0.5 ml		100 min.	600-650	660	Rat IgG2a APC Code R2a05

PRODUCT DESCRIPTION

Rat monoclonal antibody to human IL-2

Clone: MQ1-17H12

Isotype: Rat IgG2a

Lot No.: See label **Expiration:** See label

Buffer: Phosphate buffered saline (PBS)

Preservative: 0.1% *sodium azide*. Sodium azide is an extremely toxic and dangerous compound particularly when combined with acids or metals. Solutions containing sodium azide should be disposed of properly.

Stabilizer: For conjugated products only, a highly purified grade of BSA has been added as a stabilizing agent.

STORAGE & HANDLING

Store reagents at 2-8°C. For fluorochrome-conjugated antibodies only, light exposure should be avoided. Use dim light during handling, incubation with cells and prior to analysis. It is recommended that cells be analyzed within 18 hours of staining. If the reagent is being diluted, it is recommended that only the quantity to be used within one week be diluted.

PRODUCT CHARACTERIZATION

Antigen Specificity: According to the literature this antibody recognizes human interleukin-2 (IL-2)^{1,2}. This antibody has been demonstrated to inhibit the biological function of IL-2.

Staining Protocol: This antibody has been optimized for use with Invitrogen's FIX & PERM[®] fixation and permeabilization kits (catalog numbers GAS-003 and GAS-004). The staining protocol is given in the product insert for these kits.

PRODUCT QUALITY CONTROL

Each lot is tested by flow cytometry using activated human PBMCs (HiCK-1 cells, Becton Dickinson Biosciences). Based on this testing it is recommended that 5 µl of antibody be used per 1 x 10⁶ cells in a 100 µl staining volume. Because conditions may vary, it is suggested that each investigator determine the optimal amount of antibody to be used for each application.

REFERENCES:

- Andersson, J., J. Abrams, L. Bjork, K. Funai, M. Litton, K. Agren, and U. Andersson. 1994. *Immunology* 83: 16.
- Fernandez, V., J. Andersson, U. Andersson, and M. Troye-Blomberg. 1994. *Eur. J. Immunol.* 24: 1808.
- Prussin, C. and D. Metcalfe. 1995. *J. Immunol. Meth.* 188: 117.
- Abrams J. S., M.-G. Roncarolo, H. Yssel, U. Andersson, G. J. Gleich, and J. Silver. 1992. *Immunol. Rev.* 127: 5.
- Gaines Das R. E., and S. Poole. 1993. *J. Immunol. Meth.* 160: 147.
- McKenzie A. N. J. and G. Zurawski. 1994. Measurement of IL-13. In *Current Protocols in Immunology*. J. E. Coligan, A. M. Kruisbeek, D. H. Margulies, E. M. Shevach, and W. Strober, eds. John Wiley & Sons. New York, NY, p. 6.18.1.

* Antibody value assigned is based on the Optical Density at 280 nm.

† The Alexa Fluor dye conjugate in this product is sold under license from Molecular Probes, Inc., and is covered by pending and issued patents.

FOR RESEARCH USE ONLY. . . NOT FOR THERAPEUTIC OR IN VITRO DIAGNOSTIC USE

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PF: L13016

(Rev 12/08) DCC-08-1818

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.