

PRODUCT INSERT
CYTO-IC MONOCLONAL ANTIBODIES
SPECIFIC TO MURINE CYTOKINES

Product Code	Form	Volume	Antibody*	Excitation (nm)	Peak Emission (nm)	Matching Isotype Controls	
RM9060	Purified	1.0 ml	200 µg	N/A	N/A	Rat IgG1 Purified	Code R100
RM9061	FITC	1.0 ml	100 µg	488	520	Rat IgG1 FITC	Code R101
RM9061-3	FITC	3.0 ml	300 µg				
RM9064	R-PE	0.5 ml	50 µg	488	575	Rat IgG1 R-PE	Code R104
RM9064-3	R-PE	3.0 ml	300 µg				
RM90624	PE-Alexa Fluor® 700	1.0 ml	100 µg	488	723	Rat IgG1 PE-Alexa Fluor® 700	Code R124

PRODUCT DESCRIPTION

Rat monoclonal antibody to Mouse IL-5

Clone TRFK5

Isotype: Rat IgG1

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, highly purified grade of BSA has been added as a stabilizing agent.

STORAGE & HANDLING

Store reagents at 2-8°C. Light exposure should be avoided with fluorochrome conjugated reagents. It is recommended that investigators dilute only the quantity of antibody to be used within one week.

PRODUCT CHARACTERISATION

Antigen Specificity: Mouse IL-5

PRODUCT QUALITY CONTROL

Every lot of fluorochrome conjugated antibody is tested by flow cytometry using activated mouse splenocytes. Based on this testing it is recommended that a maximum of 1.0 µg of antibody be used per 1×10^6 cells in a 100 µl staining volume. Because conditions may vary, it is recommended that each investigator determine the optimal amount of antibody to be used for their application.

REFERENCES:

Schumacher, J. H., A. O'Garra, et al. 1988 The characterization of four monoclonal antibodies specific for mouse IL-5 and development of mouse and human IL-5 enzyme-linked immunosorbent. *J. Immunol.* 141(5): 1576-81.

* Antibody value assigned is based on the extinction coefficient of IgG: Optical Density at 280 nm of 1.4 equals 1 mg antibody.

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

PI: L11400

(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.