

PRODUCT INSERT

RAT anti-MOUSE CD45

Product Code	Form	Volume	Antibody*	Excitation (nm)	Peak Emission (nm)	Matching Isotype Controls	
RM6400	Purified	1.0 ml	200 µg	N/A	N/A	Rat IgG2b Purified	Code R2b00
RM6404	PE	0.5 ml	50 µg	488	575	Rat IgG2b R-PE	Code R2b04
RM6404-3	PE	3.0 ml	300 µg				

PRODUCT DESCRIPTION

Rat monoclonal antibody to mouse CD45

Clone: YW62.3

Isotype: Rat IgG2b

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

STORAGE AND HANDLING

Store reagents at 2-8°C. Light exposure should be avoided with fluorochrome-conjugated antibodies. 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: The YW62.3 monoclonal antibody (mAb) reacts with CD45 antigen (LCA, Ly-5), a transmembrane tyrosine phosphatase required for activation of T cells through B and T cell antigen receptor signaling. CD45 antigen is found on all leukocytes and cells of hematopoietic origin, but not erythrocytes. The YW62.3 mAb recognizes all isoforms of CD45 antigen and can be used in immunostaining for flow cytometry.

PRODUCT QUALITY CONTROL

Every lot is tested by flow cytometry using freshly harvested mouse splenocytes. From this testing it is recommended that between 0.10 and 0.25 µg of antibody be used per 1 x 10⁶ 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 each application.

* The amount of antibody is determined by measuring the optical density using a spectrophotometer. The antibody titer is verified by immunofluorescent staining and flow cytometric analysis.

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: L11366

(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.