

PRODUCT INSERT
MONOCLONAL ANTIBODIES TO RAT B-CELLS

| Product Code | Form | Volume | Antibody* | Excitation (nm) | Peak Emission (nm) | Fluorochrome/ Antibody Molar Ratio | Matching Isotype Controls |
|--------------|----------|--------|-----------|-----------------|--------------------|--|-----------------------------------|
| MR5400 | Pur. Ab. | 1.0 ml | 200 µg | N/A | N/A | N/A | Mouse IgG-2a Purified Code MG2a00 |
| MR5401 | FITC | 1.0 ml | 100 µg | 488 | 520 | 3 to 5 | Mouse IgG-2a FITC Code MG2a01 |
| MR5404 | R-PE | 0.5 ml | 50 µg | 488 | 575 | 1 | Mouse IgG-2a R-PE Code MG2a04 |

Description: Mouse monoclonal antibody to rat B cells.
Clone RLN-9D3 Mouse IgG-2a

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

Antigen Specificity: By flow cytometry, the cellular reactivities are specific to rat B cells.

Use In Flow Cytometry:

These antibodies are provided in liquid form for application in single, dual or three color flow cytometry. 1 µg is the maximum amount of reagent required for 1×10^6 cells. Lesser amounts of antibody may be sufficient and it is recommended that the customer determine the optimum amount of antibody for each application.

Flow Cytometry Testing:

Each lot is tested in flow cytometry. Histograms are available upon request.

Preservatives:

This reagent contains *sodium azide*. Sodium azide yields hydrazoic acid under acid conditions, which is an extremely toxic compound. Azide compounds should be flushed with running water after being discarded. This procedure is recommended to avoid deposits in plumbing where explosive conditions can develop.

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

Storage & Handling:

FITC, Purified: Fluorochrome store in the dark. +2° to +8°C for short term (to 3 month) and -20°C or colder for long term. Aliquot and avoid repeated freezing and thawing. Mix thoroughly after thawing before use.

R-PE: Store in the dark. +2° to +8°C
...**DO NOT FREEZE**

For fluorochrome conjugates, light exposure should be avoided. Use dim light during handling, incubation with cells and prior to analysis.

If the reagent is being diluted, it is recommended that only the quantity to be used within one week be diluted.

References:

1. Takami, T., Oi, C-F., Kimura, H., Kawata, K., Yamada, T., Ojima, S., Koizumi, M., Uede, T. and Kikuchi, K. 1988 Analysis of rat B lymphocyte by using monoclonal antibodies Trans. Soc. Pathol. Jpn. 77:178.

* 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: L11300

(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.