

PRODUCT INSERT

RAT IgG2b ISOTYPE CONTROLS

Product Code	Form	Description	Volume	Amount*	Tests	Excitation (nm)	Peak Emission (nm)
R2b00	Purified	Rat IgG2b	0.5 ml	50 µg	50 min.	N/A	N/A
R2b15	Biotin	Rat IgG2b	0.5 ml	50 µg	50 min.	N/A	N/A
R2b26	Alexa Fluor® 405	Rat IgG2b	0.5 ml	50 µg	50 min.	405	421
R2b28	Pacific Blue™	Rat IgG2b	0.5 ml	50 µg	50 min.	405	455
R2b30	Pacific Orange™	Rat IgG2b	0.5 ml	50 µg	50 min.	405	551
R2b20	Alexa Fluor® 488	Rat IgG2b	0.5 ml	50 µg	50 min.	488	519
R2b01	FITC	Rat IgG2b	0.5 ml	50 µg	50 min.	488	525
R2b04	R-PE	Rat IgG2b	0.5 ml	50 µg	50 min.	488	575
R2b17	PE-TR†	Rat IgG2b	0.5 ml	50 µg	50 min.	488	615
R2b22	PE-Alexa Fluor® 610	Rat IgG2b	0.5 ml	50 µg	50 min.	488	628
R2b06	TC‡	Rat IgG2b	0.5 ml	50 µg	50 min.	488	670
R2b18	PE-Cy5.5	Rat IgG2b	0.5 ml	50 µg	50 min.	488	694
R2b24	PE-Alexa Fluor® 700	Rat IgG2b	0.5 ml	50 µg	50 min.	488	723
R2b05	APC	Rat IgG2b	0.5 ml	50 µg	50 min.	600-650	660
R2b21	Alexa Fluor® 647	Rat IgG2b	0.5 ml	50 µg	50 min.	600-650	668
R2b19	APC-Cy5.5	Rat IgG2b	0.5 ml	50 µg	50 min.	600-650	694
R2b29	Alexa Fluor® 700	Rat IgG2b	0.5 ml	50 µg	50 min.	630-702	723
R2b27	APC-Alexa Fluor® 750	Rat IgG2b	0.5 ml	50 µg	50 min.	600-650	775

PRODUCT DESCRIPTION

Rat IgG2b isotype controls

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.

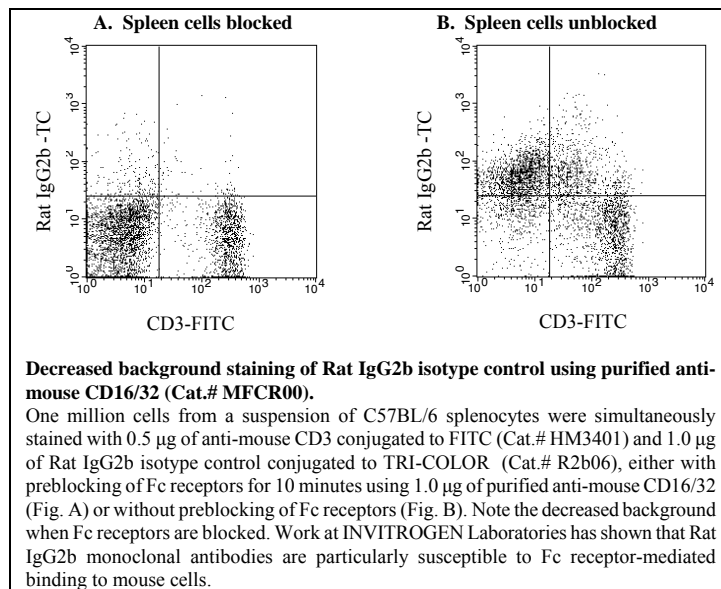
**Stabilizer:** For conjugated products only, a highly purified grade of BSA has been added as a stabilizing agent.

STORAGE & HANDLING

Store reagents at 2-8°C. Light exposure should be avoided with fluorochrome-conjugated reagents. 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 QUALITY CONTROL

Every lot is tested by flow cytometry using freshly harvested mouse splenocytes and/or human peripheral blood leukocytes (PBL). When using isotype controls as negative controls, it is recommended that the amount of isotype control and testing antibody used be equivalent.



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

† TR, Texas Red®

‡ TC, TRI-COLOR®, PE-Cy5

The efficiency of energy transfer in tandem dyes can be significantly decreased by exposure to visible light. We recommend that longer wavelength fluorochrome conjugates, e.g. PE-Cy7, PE-Alexa Fluor® 700, be protected from light during staining and while awaiting analysis, e.g. cover with aluminum foil.

**FIX & PERM®** and **COMBI-IC** reagents are produced for Invitrogen Laboratories by An Der Grub Bio Research GmbH, Austria.

The Texas Red®, Alexa Fluor® and Pacific Blue® dye conjugates in this product are sold under license from Molecular Probes, Inc., for research use only or as analyte specific reagents, except for use in combination with microarrays or high content screening, and are covered by pending and issued patents.

Cy™ is a trademark of GE/Amersham Biosciences.

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

[www.invitrogen.com](http://www.invitrogen.com)

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

PI: L11515

(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](http://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.