

PRODUCT INSERT

GOAT ANTI-RABBIT IMMUNOGLOBULINS
FLUOROCHROME CONJUGATES

Product Code	Description	Form	Volume	Antibody*	Excitation (nm)	Peak Emission (nm)
L42001	Goat Anti-Rabbit IgG (H+L), Hu Ads	FITC	2.0 ml	2.0 mg	488	525
L42010	Goat Anti-Rabbit IgG (H+L), Hu Ads	Cy3	1.0 ml	200 µg	488	575
L42018	Goat Anti-Rabbit IgG (H+L), Hu Ads	PE-Cy5.5 [†]	0.5 ml	50 µg	488	694

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

Preparation: Antibodies were purified by affinity chromatography and adsorbed to remove cross reactivity to human immunoglobulins. Specificity and cross reactivities were verified by ELISA. The resulting antibody was conjugated with the indicated fluorochrome to provide the optimal fluorochrome to protein ratio for flow cytometry. Any free dye was removed by column chromatography.

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 & HANDLING

Store reagents at 2-8°C. 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 QUALITY CONTROL

Every lot is tested by flow cytometry using either human peripheral blood leukocytes (PBL) prestained with a rabbit polyclonal antibody directed to a human leukocyte cell surface marker, or rabbit lymphocytes. Based on this testing, it is recommended that between 0.1 and 0.25 µg of antibody be used per 1x 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.

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

[†] 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 reactions and while awaiting analysis, e.g. cover with aluminum foil.

Cy is a registered trademark of GE/Amersham Biosciences.

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 L11013

(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.