

PRODUCT INSERT
**F(ab')₂ AFFINITY-ISOLATED ANTIBODIES
SPECIFIC TO MOUSE IMMUNOGLOBULINS
FLUOROCHROME CONJUGATES**

Product Code	Description	Form	Volume	Antibody*	Excitation (nm)	Emission (nm)
M35006	Goat F(ab') ₂ Anti-Mouse IgG (H+L)	TC‡	0.5 ml	100 µg	488	670
M35018	Goat F(ab') ₂ Anti-Mouse IgG (H+L)	PE-Cy [®] 5.5§	0.5 ml	100 µg	488	694
M35011	Goat F(ab') ₂ Anti-Mouse IgG (H+L)	Cy [®] 5	1.0 ml	200 µg	600-650	670
M35201	Goat F(ab') ₂ Anti-Mouse IgG+IgM (H+L)	FITC	1.0 ml	1000 µg	488	525
M31601	Goat F(ab') ₂ Anti-Mouse IgM (µ)	FITC	1.0 ml	700 µg	488	575
M31604	Goat F(ab') ₂ Anti-Mouse IgM (µ)	R-PE	0.5 ml	100 µg	488	575

Lot No.: See label

Expiration: See label

Buffer: Phosphate buffered saline (PBS)

Preparation: Antibodies were affinity purified by column chromatography and adsorbed to remove cross reactivity to human immunoglobulins. Specificity and cross reactivities were verified by ELISA. Purified antibodies were digested with pepsin. Any aggregates, undigested goat immunoglobulins and Fc fragments were removed by column chromatography. The resulting antibody was conjugated with the indicated fluorochrome to provide the optimal fluorochrome to protein ratio for flow cytometry, cell and tissue staining use. Any free dye was removed by column chromatography.

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 when using 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 QUALITY CONTROL

Every lot is tested by flow cytometry using human peripheral blood leukocytes (PBL) prestained with an unlabeled mouse monoclonal antibody directed to a human leukocyte cell surface marker. Based on this testing it is recommended that between 0.25 and 1.0 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.




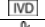




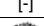
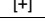


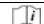
* Antibody value assigned is based on the Optical Density at 280 nm.

‡ TC, TRI-COLOR[®], PE-Cy[®]5

§ 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-Cy[®]7, 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.

Explanation of symbols

Symbol	Description	Symbol	Description
	Catalogue Number		Batch code
	Research Use Only		In vitro diagnostic medical device
	Use by		Temperature limitation
	Manufacturer		European Community authorised representative
	Without, does not contain		With, contains
	Protect from light		Consult accompanying documents
	Directs the user to consult instructions for use (IFU), accompanying the product.		

For research use only. CAUTION: Not for human or animal therapeutic or diagnostic use.

www.invitrogen.com

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