

**PRODUCT INSERT**

**CYTO-IC MONOCLONAL ANTIBODIES  
RAT anti-MOUSE TNF- $\alpha$**

Product Code	Form	Amount	Excitation (nm)	Peak Emission (nm)	Matching Isotype Controls
RM90128	Pacific Blue <sup>™</sup>	1.0 ml	405	455	
RM9011	FITC	1.0 ml	488	520	Rat IgG1 FITC Code R101

To see a complete list of our antibody offerings, please visit [www.lifetechnologies.com/antibodies](http://www.lifetechnologies.com/antibodies)

**PRODUCT DESCRIPTION**

Rat monoclonal antibody to Mouse TNF- $\alpha$

**Clone** MP6-XT22

**Isotype:** Rat IgG1

**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, 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. It is recommended that investigators dilute only the quantity of antibody to be used within one week.








**PRODUCT USE**

Every lot is tested by flow cytometry using activated mouse splenocytes. Because conditions may vary, it is recommended that each investigator determine the optimal amount of antibody to be used for each application.

**REFERENCES:**

1. Abrams, J.S., M.G. Roncaolo, I. Yssel, U. Anderson, G.J. Gleich, and J. Silver. 1992. Strategies of anti-cytokine monoclonal antibody development. *Immunol. Rev.* 127: 5-24.
2. Litton, M.J., B. Sander, E. Murphy, A.O. O'Garra, and J.S. Abrams. 1994. Early expression of cytokines in lymph nodes after treatment *in vivo* with *Staphylococcus* enterotoxin B. *J. Immunol. Meth.* 175: 47-58.

**Explanation of symbols**

Symbol	Description	Symbol	Description
	Manufacturer		Catalog number
	Use by		Temperature limitation
	Consult instructions for use		Caution, consult accompanying documents
	Batch code		

**For Research Use Only. Not for use in diagnostic procedures.**

© 2013 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners. For support visit [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support) or email [techsupport@lifetech.com](mailto:techsupport@lifetech.com)

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact [outlicensing@lifetech.com](mailto:outlicensing@lifetech.com) or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.