

Certificate of Analysis

FITC Conjugated Anti-Mouse Interferon Alpha, Clone RMMA-1 (MAb)

Catalog No: 22100-3

Lot No:

Expiration: 2.5 years from date of manufacturing of new stock [input actual date]

Size: 25 µg/vial

Description: Fluorescein Isothiocyanate (FITC) conjugated Rat Monoclonal Antibody against Mouse Interferon Alpha

Clone: RMMA-1

Volume: 0.5 ml

Antibody Concentration: 0.05 mg/ml

F/P Ratio:

Buffer: Phosphate buffered saline (PBS) containing 1% BSA and 0.05% Kathon, pH 7.2

Antigen: Mouse Interferon Alpha

Isotype: IgG₁

Tested Applications: Flow Cytometry: µg/ml of antibody is sufficient for labeling of 1×10^6 cells in 100 µl total volume of B16 mouse melanoma cell line inducing interferon alpha. It is recommended that each investigator determine appropriate dilutions for other cell lines.

Immunohistochemistry: It is recommended that each investigator determine appropriate dilutions for individual use.

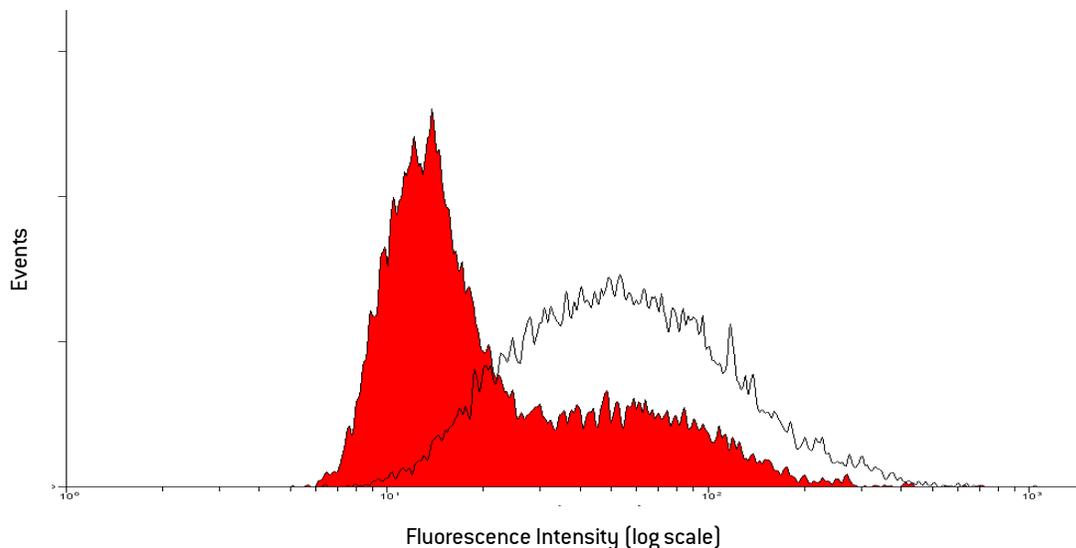


Figure 1: Log Fluorescence Intensity . FITC Mouse IFN- . Expression of endogenous Mouse Interferon Alpha in B16 cell lines. Flow Cytometry analysis on non-induced cells (closed histogram) and induced cells (open histogram).

Shipping Conditions: Wet Ice

Physical State of Product During Shipping: Liquid

Storage Conditions/Comments: Photo-sensitive material; protect from prolonged exposure to light. Store at 2-8°C. **DO NOT FREEZE.** For more information on protein handling, visit our Resource Library at www.pbl assaysci.com.



pbl assay science

131 Ethel Road West, Suite 6 | Piscataway, NJ 08854 USA | T:+1.732.777.9123 | F:+1.732.777.9141 | E: info@pbl assaysci.com | W: pbl assaysci.com

Authorization

Released by: _____

Date:

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL Assay Science. For research use only. Not for diagnostic or clinical use in, or administration to, humans. Not for resale in original or any modified form, including inclusion in a kit, for any purpose. Not for use in the preparation of any commercial product.

Rev. 04

A decorative graphic at the bottom of the page consisting of a wavy, multi-colored shape that transitions from light green on the left to bright blue on the right.