

## Certificate of Analysis

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

**Catalog No:** 22100-3

**Lot No:** 7398

**Expiration:** January 29, 2024

**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:** 5.02

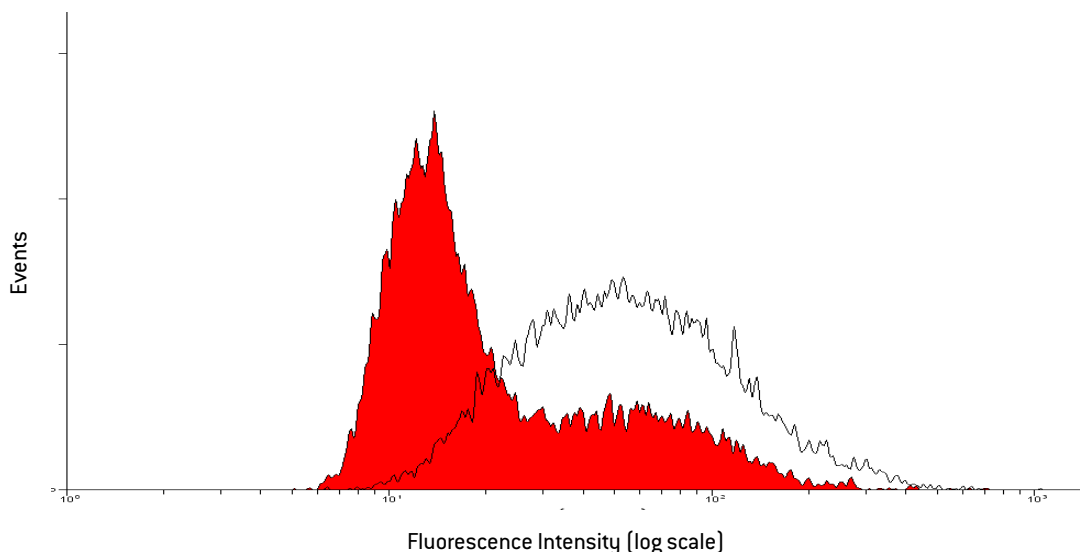
**Buffer:** Phosphate buffered saline (PBS) containing 1% BSA and 0.05% Kathon, pH 7.2

**Antigen:** Mouse Interferon Alpha

**Isotype:** IgG<sub>1</sub>

**Tested Applications:** Flow Cytometry: 0.5 µg/ml of antibody is sufficient for labeling of  $1 \times 10^6$  cells in 100 µl total volume of B16 $\alpha$  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- $\alpha$ . Expression of endogenous Mouse Interferon Alpha in B16 $\alpha$  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](http://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: August 31, 2021

**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. 05

A decorative footer at the bottom of the page consisting of a wavy, multi-colored gradient bar transitioning from light green on the left to bright blue on the right.