



Certificate of Analysis

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

Catalog No: 22100-1

Lot No: 7477

Size: 250 µg/vial

Description: Rat Monoclonal Antibody against Mouse Interferon Alpha, neutralizing

Clone: RMMA-1

Volume: 0.1 ml

Concentration: 2.5 mg/ml

Buffer: 0.3M Sodium Bicarbonate; 0.2M Sodium Chloride, pH 8.0

Endotoxin: < 1 EU/µg

Antigen: Mouse Interferon Alpha

Isotype: Rat IgG₁

Purity: > 95%

Purification Method: A combination of affinity, ion exchange, hydrophobic interaction and size exclusion chromatography

Specificity: Neutralizes mouse interferon alpha; binds to mouse interferon alpha

Assay Used to Measure Bioactivity: One neutralization unit is the amount of antiserum required to neutralize one unit of mouse interferon alpha (Mu-IFN-α) to a 50% endpoint. Interferon was titrated with the use of the cytopathic effect inhibition assay using L929 cells as described [Rubinstein, S., Familletti, P.C., and Pestka, S. (1981) "Convenient Assay for Interferons," *J. Virol.* 37, 755-758] with the exception that EMCV was used as the challenge virus. In this antiviral assay for interferon, about 5 unit/ml of interferon is the quantity necessary to produce a cytopathic effect of 50%. The units are determined with respect to the international reference standard for Mu-IFN-α provided by the National Institutes of Health (Ga02-901-511).

Tested Applications: ELISA, Neutralization

Optimal dilutions should be determined by each laboratory for each application.

Suggested Applications: Immunoprecipitation

Please note that these applications are presented for suggested use only and have not been fully evaluated by PBL.

Selected References:

- (1) Asselin-Paturel *et al.* (2005). Type I interferon dependence of plasmacytoid dendritic cell activation and migration. *J. Exp. Med.* 201:1157. Used this antibody for FACS analysis by intracellular staining of dendritic cells for IFN-Alpha production.
- (2) Mellor *et al.* (2005). CpG oligonucleotides induce splenic CD19+ dendritic cells to acquire potent indoleamine 2,3-dioxygenase-dependent T cell regulatory functions via IFN Type 1 signaling. *J. Immunol.* 175:5601. Used this antibody to block IFN-alpha activity in the stimulation of IDO expression of dendritic cells.

Shipping Conditions: Dry Ice

Physical State of Product During Shipping: Frozen

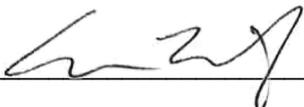
Storage Conditions/Comments: After receipt, the product may be stored at -20°C for short-term use (≤ 6 months). For long-term storage, we recommend storing the product at -70°C or below for retention of full activity. When thawing, the contents of the tube should be apportioned in separate tubes so that freezing and thawing is kept to a minimum. Re-freezing should be done on dry ice or in a dry ice/alcohol bath. Further dilution of the product should be in buffers containing protein such as 0.1% bovine serum albumin (BSA). 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: February 1, 2022

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. 06

