

Certificate of Analysis

Anti-Mouse Interferon Beta, Clone RMMB-1 (MAb)

Catalog No: 22400-1

Lot No:

Size: 250 µg/vial

Description: Rat Monoclonal Antibody against Mouse Interferon Beta

Clone: RMMB-1

Volume: ml

Concentration: mg/ml

Buffer: Phosphate buffered saline (PBS) containing 0.1% bovine serum albumin (BSA)

Antigen: Mouse Interferon Beta

Isotype: IgG_{2a,κ}

Bioactivity: Weakly neutralizes mouse interferon beta, but not recommended for neutralization assays (for this application products 32400-1 or 32401-1 are recommended); binds to mouse interferon beta

Assay Used to Measure Bioactivity: One neutralization unit is the amount of antiserum required to neutralize one unit of mouse interferon beta (Mu-IFN-) to a 50% endpoint. Interferon was titrated with the use of the cytopathic effect inhibition assay as described [Rubinstein, S., Familletti, P.C., and Pestka, S. (1981) "Convenient Assay for Interferons," *J. Virol.* 37, 755-758; Familletti, P.C., Rubinstein, S., and Pestka, S. (1981) "A Convenient and Rapid Cytopathic Effect Inhibition Assay for Interferon," in *Methods in Enzymology*, Vol. 78 (S. Pestka, ed.), Academic Press, New York, 387-394]. In this antiviral assay for interferon about 1 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- provided by the National Institutes of Health [see Pestka, S. (1986) "Interferon Standards and General Abbreviations," in *Methods in Enzymology* (S. Pestka, ed.), Academic Press, New York 119, 14-23]. This material is prepared specifically for effective neutralization of Mu-IFN- .

Tested Applications: Direct Binding ELISA

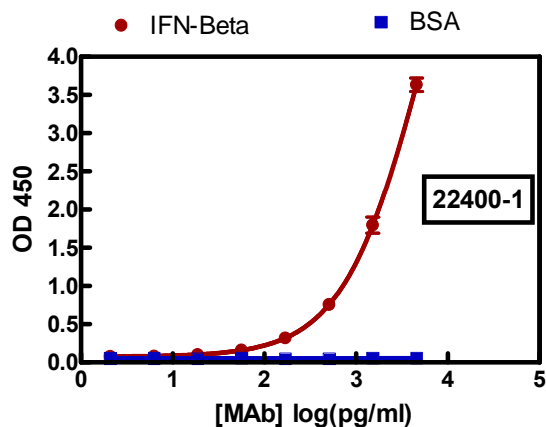


Figure 1. Representative binding curve of antibody to recombinant Mu-IFN- (circle) and to 1% BSA/PBS (square) in a Direct Binding ELISA. High-binding polystyrene plates were coated with either 1 µg/ml Mu-IFN- or 1% BSA/PBS. Thereafter, titrations of the product were added to the wells. Donkey anti-rat IgG conjugated to HRP was used as the detection antibody. Colorimetric detection was performed using 3, 3', 5'-Tetramethylbenzidine (TMB) substrate. The HRP-TMB reaction was stopped using a diluted H₂SO₄/HCl solution.



Note: PBL has not tested the use of this product in western blot, flow cytometry, immunoprecipitation, or immunohistochemistry.

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 (m6 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. Refreezing 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.pblassaysci.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.