

## Certificate of Analysis

### Rat Mab to Mouse Interferon Beta, Clone RMMB-1 (MAb)

**Catalog No:** 22400-9

**Lot No:** 7173

**Size:** 0.2 mg/vial

**Description:** Rat Monoclonal Antibody against Mouse Interferon Beta

**Clone:** RMMB-1

**Volume:** 0.050 ml

**Concentration:** 4.04 mg/ml

**Buffer:** 0.3 Sodium Bicarbonate, 0.2 M Sodium Chloride

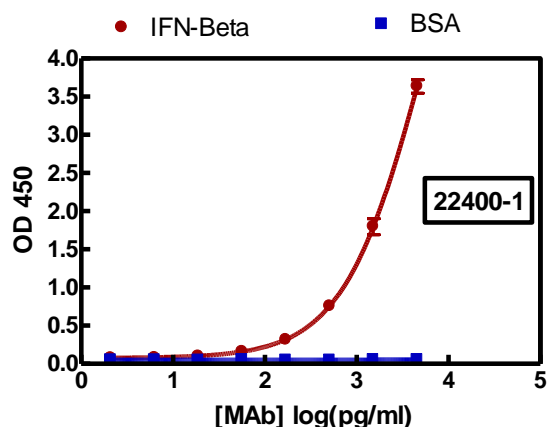
**Antigen:** Mouse Interferon Beta

**Isotype:** IgG<sub>2a,K</sub>

**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- $\beta$ ) 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- $\beta$  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- $\beta$ .

**Tested Applications:** Direct Binding ELISA



**Figure 1.** Representative binding curve of antibody to recombinant Mu-IFN- $\beta$  (circle) and to 1% BSA/PBS (square) in a Direct Binding ELISA. High-binding polystyrene plates were coated with either 1  $\mu$ g/ml Mu-IFN- $\beta$  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, 5'-Tetramethylbenzidine (TMB) substrate. The HRP-TMB reaction was stopped using a diluted H<sub>2</sub>SO<sub>4</sub>/HCl solution.



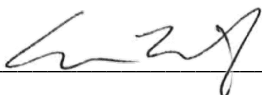
**Note:** PBL has not tested the use of this product in western blot, flow cytometry, immunoprecipitation, or immunohistochemistry.

**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 ( $\leq$  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. 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](http://www.pblassaysci.com).

**Authorization**

Released by: \_\_\_\_\_ 

Date: January 17, 2020

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.