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

# **Certificate of Analysis**

## Anti-Mouse Interferon Beta, Rabbit Serum (PAb)

**pbl** assay science

Catalog No: 32400-1 Lot No: Size:  $\ge 2 \times 10^4$  Neutralizing Units/vial

Description: Rabbit polyclonal antibody against Mouse Interferon BetaVolume:mlActivity using *E. coli* interferon:x 10NU/mlActivity using mammalian interferon:x 10NU/mlConcentration:mg/ml (estimate)Buffer:Neat serum diluted in phosphate buffered saline (PBS)

Antigen: Recombinant mouse interferon beta (mammalian expressed)

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 [Rubinstein, S., Familletti, P.C., and Pestka, S. (1981) *J. Virol.* 37, 755-758] using L929 cells and EMCV as the challenge virus. In this antiviral assay for interferon, about 2.5 unit/ml of interferon is the quantity necessary to produce a cytopathic effect of 50%. The units are determined by use of a laboratory standard calibrated with respect to the international reference standard for Mu-IFN- $\beta$  provided by the National Institutes of Health [Gb02-902-511]. This material is prepared specifically for effective neutralization of Mu-IFN- $\beta$ .

## Tested Applications: Neutralization

**Suggested Applications:** ELISA; Western blot; immunoprecipitation; immunohistochemistry *Please note that these applications are presented for suggested use only and have not been fully evaluated by PBL.* 

## Shipping Conditions: Dry ice

#### Physical State of Product During Shipping: Frozen

**Storage Conditions/Comments:** After receipt, this product should be kept at -70°C or below for retention of full activity. Thaw product vial by incubation in cold tap water until just thawed – 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 <u>www.pblassaysci.com</u>.

#### Selected References:

- (1) Asselin-Paturel *et al.* (*J Exp. Med*, 2005) used 32400 in combination with Rabbit anti-mouse IFN-alpha (32100) to explore the role of Type I IFN in the migration and activation of dendritic cells.
- (2) Kamath *et al.* (*J. Immunol.* 2005) used a similar cocktail to demonstrate that dendritic cell derived Type I IFN stimulates bystander T-cells.
- (3) Seimon *et al.* (*PNAS*, 2006) used 32400 to examine the role of IFN-Beta in protecting LPS treated macrophages from apoptosis. Zheng *et al.* (*J. Biol. Chem*, 2006) used 32400 to examine the role of autocrine IFN-Beta in RANKL stimulated iNOS expression in macrophages.

### 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