



## Certificate of Analysis

### Rat Interferon Beta

**Catalog No:** 13400-1

**Lot No:**

**Expiration:**

**Size:**  $1 \times 10^5$  units/vial

**Description:** Recombinant Rat Interferon Beta (Rat IFN- $\beta$ ). Active on rat and mouse cells.

**Contents:**  $x 10$  units/vial

**Concentration:**  $x 10$  units/ml after reconstitution with ml sterile distilled water

**Specific Activity:**  $x 10$  units/mg

**Buffer:** prior to lyophilization: phosphate buffered saline (PBS); 125mM trehalose

**Molecular Weight:** 19,680 Da

**Purity:** > 95%

**Source:** cDNA expressed in CHO cells

**Accession #:** D87919

**Assay Used to Measure Bioactivity:** 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%. As there is no international Rat IFN- $\beta$  standard, units are determined with respect to the international reference standard for mouse interferon alpha/beta (mouse 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]. Units of activity measured on rat epithelial-like (RATEC) cells with VSV.

**Shipping Conditions:** Wet Ice

**Physical State of Product During Shipping:** Lyophilized

**Storage Conditions/Comments:** Upon receipt, this product should be kept at -20°C while lyophilized, and at -70°C or below after reconstitution in sterile distilled water for retention of full activity. After reconstitution, the contents of the tube should be apportioned in separate tubes so that freezing and thawing is kept to a minimum. After thawing, this product is stable for at least four weeks at 4°C. Further dilute only 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](http://www.pbl assaysci.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.

Rev. 03