



## Certificate of Analysis

### Anti-Human Interferon Beta, Sheep Serum (PAb)

**Catalog No:** 31401-1

**Lot No:** 7311R

**Size:**  $\geq 2 \times 10^4$  Neutralizing Units/vial

**Description:** Sheep Polyclonal Antibody against Human Interferon Beta

**Volume:** 0.100 ml

**Activity:**  $2.7 \times 10^5$  NU/ml

**Concentration:** 0.53 mg/ml (estimate)

**Buffer:** Phosphate buffered saline (PBS)

**Antigen:** Recombinant human interferon beta 1a (expressed in mammalian cells)

**Assay Used to Measure Bioactivity:** One neutralization unit is the amount of antiserum required to neutralize one unit of human interferon beta (Hu-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 Hu-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 Hu-IFN- $\beta$ .

**Tested Applications:** Neutralization

**Suggested Applications:** ELISA; Western blot; immunoprecipitation; immunohistochemistry

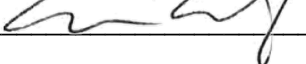
*Please note that these applications are presented for suggested use only and have not been fully evaluated by PBL.*

**Shipping Conditions:** Dry Ice

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

**Storage Conditions/Comments:** After receipt, the product may be stored at  $-20^{\circ}\text{C}$  for short-term use ( $\leq 6$  months). For long-term storage, we recommend storing the product at  $-70^{\circ}\text{C}$  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: November 16, 2023

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.