



Certificate of Analysis

Anti-Human Interferon Alpha (All Subtype), Rabbit Serum (PAb)

Catalog No: 31135-1

Lot No:

Size: $\geq 1 \times 10^5$ Neutralizing Units/vial

Description: Rabbit Polyclonal Antibody against Human Interferon Alpha

Volume: ml

Activity: x 10 NU/ml

Buffer: Neat serum diluted in phosphate buffered saline (PBS)

Antigen: Mixture of all recombinant human interferon alpha subtypes (IFNA2, IFNA8, IFNA10, IFNA1, IFNA21, IFNA5, IFNA14, IFNA17, IFNA7, IFNA6, IFNA4, IFNA16)

Specificity: Binds to and blocks biological activity of all human interferon alpha subtypes.

Assay Used to Measure Bioactivity: One neutralization unit is the amount of antiserum required to neutralize one unit of human interferon alpha 2a (Hu-IFN- α 2a) 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- α 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- α .

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°C for short-term use (≤ 6 months). For long-term storage, we recommend storing the product 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 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.

Rev. 00