



## Certificate of Analysis

### Anti-Human Interferon Alpha, Clone MMHA-14 (MAb)

**Catalog No:** 21129-1

**Lot No:**

**Size:** 50 µg/vial

**Description:** Mouse Monoclonal Antibody against Human Interferon Alpha

**Clone:** MMHA-14

**Volume:** ml

**Concentration:** mg/ml

**Buffer:** Phosphate-buffered saline (PBS) containing 0.1% bovine serum albumin (BSA)

**Antigen:** Recombinant human interferon alpha D

**Isotype:** IgG<sub>1K</sub>

**Bioactivity:** Binds to and neutralizes human interferon alpha

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

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

*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 (m6 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](http://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