

Certificate of Analysis

Anti-Human Interferon Alpha, Clone MMHA-8 (MAb)

Catalog No: 21110-1

Lot No: 7742

Size: 50 µg/vial

Description: Mouse Monoclonal Antibody against Human Interferon Alpha, neutralizing

Clone: MMHA-8

Volume: 0.1 ml

Concentration: 0.5 mg/ml

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

Endotoxin: < 1 EU/µg

Antigen: Human Leukocyte Interferon

Isotype: Mouse IgG_{2b} kappa

Purity: > 95%

Purification Method: A combination of ion exchange, hydrophobic interaction and size exclusion chromatography

Specificity: Neutralizes human interferon alpha; binds to human interferon alpha with high affinity ($K_A \approx 10^9 \text{ M}^{-1}$)

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

Tested Applications: ELISA, Neutralization

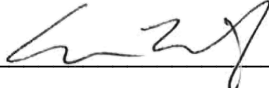
Optimal dilutions should be determined by each laboratory for each application.

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. When thawing, the contents of the tube should be apportioned in separate tubes so that freezing and thawing is kept to a minimum. Re-freezing 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: February 17, 2025

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