



Certificate of Analysis

Human Interferon Gamma

Catalog No: 11500-1

Lot No:

Size: 50 µg/vial

Description: Recombinant Human Interferon Gamma (Hu IFN-)

Volume: ml

Concentration: mg/ml

Specific Activity: x 10 units/mg

Buffer: phosphate buffered saline containing 0.1 % bovine serum albumin (BSA)

Endotoxin: < 1 EU/ g

Molecular Weight: 16.8 kDa

Purity: > 95%

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

Source: human T lymphocyte Interferon cDNA expressed in *E. coli*

Synonyms: None

Accession #: V00543

Human Gene: IFNG

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%. The units are determined with respect to the international reference standard for human interferon gamma (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]. Unit of activity measured on A549 cells with Encephalomyocarditis Virus (EMCV)].

Shipping Conditions: Dry ice

Physical State of Product during Shipping: Frozen

Storage Conditions/Comments: After receipt, this product should be kept 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.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.