

Certificate of Analysis

Human Interferon Beta, 1a (mammalian)

Catalog No: 710785-9

Lot No: 7852

Size: 0.25 mg/vial

Description: Recombinant Human Interferon Beta 1a, mammalian (Hu IFN- β 1a)

Volume: 0.887 ml

Activity: 6.77×10^7 units/ml

Specific Activity: 2.4×10^8 units/mg

Buffer: 10mM Acetic Acid

Endotoxin: < 1 EU/ μ g

Molecular Weight: ~23 kDa

Purity: > 95%, one N-Linked glycosylation methods

Purification Method: Proprietary chromatography method

Source: cDNA obtained from human fibroblast mRNA expressed in *CHO cells*

Human Gene: IFNA2

Synonyms: Fibroblast Interferon

Accession #: V00534

Assay Used to Measure Bioactivity: Interferon was titrated with the use of the cytopathic effect inhibition assay as described [Familletti, *et al.* (1981) "A Convenience and Rapid Cytopathic Effect Inhibition Assay for Interferon," in *Methods in Enzymology*, Vol. 78 (S. Pestka, ed.), Academic Press, New York, 387-394]. Encephalomyocarditis virus (EMCV) was used to challenge A549 cells [Budd *et al.* (1985) *Canc. Chem. Pharm.* 12:39]. The activity was determined relative to a lab standard of Hu-IFN- β which was calibrated to the NIH Hu-IFN- β standard (Gb23-902-531). The EC₅₀ for IFN is ~1 U/ml. Lot Activity was derived from multiple determinations in the above assay. Please note that IFN assays vary between labs and assay systems [Meager *et al.* (2001). *J. Immunol. Meth.* 257:17. Meager and Das (2005) *J. Immunol. Meth.* 306:1]

Shipping Conditions: Wet Ice

Physical State of Product During Shipping: Liquid

Storage Conditions/Comments: After receipt, the product should be stored at 2-8°C for retention of full, activity. For more information on protein handling, visit our Resource Library at www.pbl assaysci.com.

Authorization

Released by: _____

Date: February 24, 2026

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