



## Certificate of Analysis

### Human Interferon Alpha 4a (Alpha M1)

**Catalog No:** 11177-1

**Lot No:** 7724

**Size:**  $\geq 1 \times 10^5$  units

**Description:** Recombinant Human Interferon Alpha 4a (Alpha M1) (Hu-IFN- $\alpha$  M1)

**Volume:** 0.100 ml

**Activity:**  $3.16 \times 10^6$  units/ml (MDBK/VSV)

**Specific Activity:**  $3.32 \times 10^8$  units/mg (MDBK/VSV)

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

**Endotoxin:**  $< 1$  EU/ $\mu$ g

**Molecular Weight:** 19.4 kDa

**Purity:**  $\geq 95\%$

**Purification Method:** Purified from *E. coli* by a combination of Ion Exchange, Hydrophobic Interaction and Size Exclusion Chromatography

**Source:** cDNA obtained from human leukocyte mRNA expressed in *E. coli*

**Human Gene:** IFNA4

**Synonyms:** Hu-IFN- $\alpha$  M1

**Accession Number:** NM\_021068

**Assay Used to Measure Bioactivity:** Interferon was titrated with the use of the cytopathic effect inhibition assay as listed

*Bovine (MDBK/VSV)* – Performed as described [Rubinstein, *et al.* (1981) *J. Virol.* 37 (2):755]. The EC<sub>50</sub> for interferon in this assay is  $\sim 5$  U/ml. Lot Activity was derived from multiple determinations in the above assay.

*Human (A549/EMCV)* – Performed as described [Budd, *et al.* (1985) *Canc. Chem. Pharm.* 12:39]. The EC<sub>50</sub> for interferon in this assay is  $\sim 1$  U/ml.      **Activity:**  $2.28 \times 10^6$  units/ml      **Specific Activity:**  $2.40 \times 10^8$  units/mg

The units are determined by use of a Human IFN Alpha A (Hu-IFN- $\alpha$  A [2a]) laboratory standard calibrated to the international reference standard for Human Interferon Alpha A (Hu-IFN- $\alpha$  A [2a]) provided by the National Institutes of Health [Meager, *et al.* (2001) *J. Immunol. Meth.* 257 (1-2):17-33]. 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:** Dry Ice

**Physical State of Product During Shipping:** Frozen

**Storage Conditions/Comments:** After receipt, the product may be stored at  $-20^{\circ}\text{C}$  for short-term use ( $\leq 6$  months). For long-term storage, we recommend storing the product at  $-70^{\circ}\text{C}$  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. 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) or tissue culture media with serum. Dilution of material below  $2 \times 10^5$  units/ml for freezing is not recommended. For more information on protein handling, visit our Resource Library at [www.pbl assaysci.com](http://www.pbl assaysci.com).

### Authorization

Released by: \_\_\_\_\_

Date: October 15, 2024

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. 04