



## Certificate of Analysis

### Mouse Interferon Alpha 11

**Catalog No:** 12125-9

**Lot No:** 7197

**Size:** 0.5 mg/vial

**Description:** Recombinant Mouse Interferon Alpha 11 (Mu-IFN- $\alpha$ 11), carrier free

**Volume:** 0.439 ml

**Activity:**  $9.69 \times 10^6$  units/ml

**Specific Activity:**  $8.5 \times 10^6$  units/mg

**Buffer:** 50 mM HEPES pH 6.5, 250 mM NaCl, 6% Glycerol

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

**Molecular Weight:** 19.1 kDa

**Purity:** > 95%

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

**Source:** Murine Leukocyte Interferon cDNA expressed in *E. coli*

**Synonyms:** Mu-IFN- $\alpha$ 11

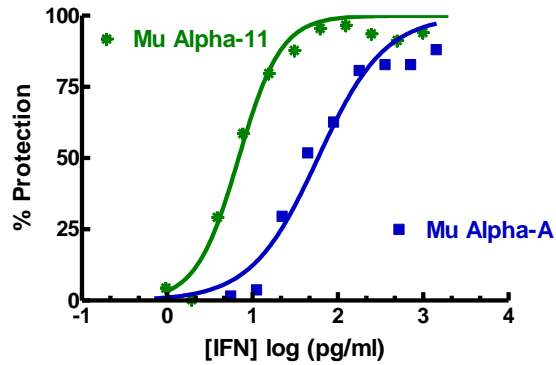
**Gene:** IFNA11

**Accession #:** AY225954

**Assay Used to Measure Bioactivity:** Interferon was titrated with the use of the cytopathic effect inhibition assay as described [Familletti, *et al.* (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] with the exception that EMCV rather than VSV was used as the challenge. The activity was determined relative to a lab standard of Mu-IFN- $\alpha$ A which was calibrated to the NIH Murine IFN- $\alpha$  standard (Ga02-901-511). Mouse (L929/EMCV) in this assay the EC<sub>50</sub> for IFN is ~5 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].

**Product Information:** Most mammalian species have multiple IFN- $\alpha$  subtypes. Although the reasons for these multiple subtypes are not fully known, there are clear cell type and temporal differences in their expression. A study established a nomenclature for the murine IFN- $\alpha$  subtypes (van Pesch, *et al.* 2004) and determined relative activities of the subtypes with protein quantification by phosphorimaging of metabolically-labeled protein. In this study, Mu-IFN- $\alpha$ A was defined to have average antiviral activity in order to compare the potencies of the other subtypes.

**Comparison of Mu Alpha-A and Mu Alpha-11  
Antiviral Activity**



**Figure 1:** The activity of Mu Alpha-A and Mu Alpha-11 was compared in an L929/EMCV CPE assay. The EC<sub>50</sub> for Mu Alpha A in this experiment was 58 pg/ml while the EC<sub>50</sub> for Mu Alpha-11 was 7 pg/ml. This relationship was confirmed for several batches of Mu Alpha-11.

*Results are representative and may vary depending upon experimental conditions.*

**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.pbl assaysci.com](http://www.pbl assaysci.com).

**Authorization**

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

Date: March 3, 2020

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.