



Certificate of Analysis

Universal™ Type I Interferon

Catalog No: 11200-1

Lot No: 7324R

Size: $\geq 1 \times 10^5$ units/vial

Description: Recombinant Human Interferon Alpha hybrid protein; Universal™ Type I IFN is an alpha interferon hybrid (Human Interferon Alpha A/D [Bg/II]) constructed from recombinant Human IFN alpha A and Human IFN alpha D. Activity has been observed on many mammalian cells, including human, monkey, mouse, bovine, rat, cat, pig, rabbit, guinea pig, or hamster. This product is useful for cross-species testing with Interferon or for species where autologous interferon is not available.

Volume: 0.100 ml

Activity: 4.99×10^6 units/ml (MDBK/VSV)

Specific Activity: 3.56×10^8 units/mg (MDBK/VSV)

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

Endotoxin: < 1 EU/ μ g

Molecular Weight: 19.4 kDa

Purity: > 95%

Purification Method: A combination of hydrophobic interaction, ion exchange and gel permeation chromatography

Source: Construct described in Rehberg, E., (1982) *J. Biol. Chem.* 257: 11497 derived from *E. coli*

Synonyms: Hu-IFN- α A/D [Bg/II]

Accession #: Not applicable

Human Gene: Not applicable

Assay Used to Measure Bioactivity: Interferon was titrated in the Cytopathic Effect Inhibition (CPE) assays as listed.

Bovine (MDBK/VSV) – performed as described [Rubinstein, *et al.* (1981) *J. Virol.* 37(2):755]. The EC₅₀ for interferon in this assay is ~5 U/ml.

Human (A549/EMCV) – performed as described [Budd, *et al.* (1985) *Canc. Chem. Pharm.* 12:39]. The EC₅₀ for interferon in this assay is ~1 U/ml.

Activity: 4.15×10^6 units/ml

Specific Activity: 2.96×10^8 units/mg

Mouse (L929/EMCV) – developed and performed at PBL Biomedical Laboratories. The EC₅₀ for interferon in this assay is ~5 U/ml.

Activity: 2.31×10^5 units/ml

Specific Activity: 1.65×10^7 units/mg

The units are determined by use of a Human IFN Alpha A (Hu-IFN- α A [2a]) laboratory standard calibrated to the international reference standard for Human Interferon Alpha A (Hu-IFN- α A [2a]) provided by the National Institutes of Health [Meager, *et al.* (2001) *J. Immunol. Methods.* 257(1-2):17-33]. On L929 cells the laboratory standard has been calibrated to the reference standard for Mouse Interferon Alpha provided by the National Institutes of Health. 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: 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. 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.pblassaysci.com.

Comparison of Universal™ Type I IFN and Mouse IFN Alpha A Antiviral Activity on Mouse Cells

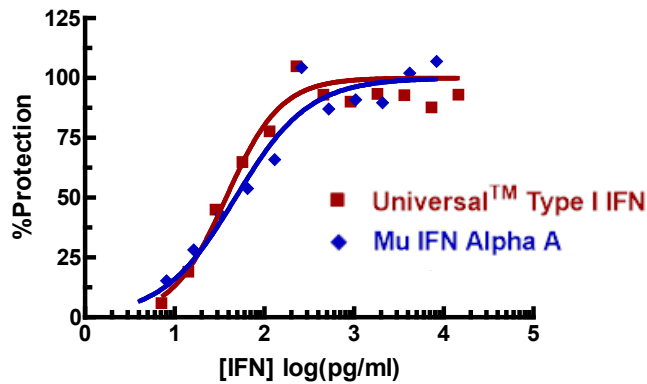
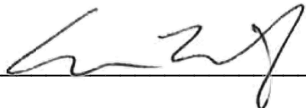


Figure 1: The activity of Mouse IFN Alpha-A and Universal™ Type-I IFN was compared in an L929/EMCV CPE assay. The EC₅₀ for Universal™ Type I IFN in this experiment was 38 pg/ml while the EC₅₀ for Mu Alpha-A was 48 pg/ml. Similar results were obtained for several batches of Mouse Alpha-A. Results are representative and may vary depending upon experimental conditions.

Authorization

Released by: _____ 

Date: May 13, 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.