



Certificate of Analysis

Porcine Interferon Alpha

Catalog No: 17105-1

Lot No:

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

Description: Recombinant Porcine Interferon Alpha

Volume: 0.1 ml

Activity: $\times 10$ units/ml (MDBK/VSV)

Specific Activity: $\times 10$ units/mg (MDBK/VSV)

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

Endotoxin: < 1 EU/ g

Molecular Weight: 19.1 kDa

Purity: $\sim 95\%$

Purification Method: Purified from mammalian tissue culture media by a combination of ion exchange and size exclusion chromatography

Source: cDNA obtained from porcine leukocyte mRNA expressed in mammalian cells

Porcine Gene: IFNA1

Synonyms: Pig IFN-

Accession #: NM_214393.1

Reference: Lefevre, *et al.* (1990) *J. Gen. Virol.* 71:1057-1063.

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 EC₅₀ for interferon in this assay is ~ 5 U/ml. Lot Activity was derived from multiple determinations in the above assay.

The units are determined by use of a Human IFN Alpha A (Hu IFN- A [2a]) laboratory standard calibrated to the international reference standard for Human Interferon Alpha A (Hu IFN- A [2a]) provided by the National Institutes of Health [Meager, *et al.* (2001) *J. Immunol Methods.* 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°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) or Tissue Culture Media with serum. Dilution for freezing of material below 2×10^5 units/ml is not recommended. One freeze-thaw cycle is equivalent to thawing an aliquot prepared from the material received. For more information on protein handling, visit our Resource Library at www.pblassaysci.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.

Rev. 01