



Certificate of Analysis

Cynomolgus Interferon Alpha 2 (Ile 16), mammalian

Catalog No: 16105-1

Lot No:

Expiration:

Size: ~ 1 x 10⁵ units/vial

Description: Recombinant Cynomolgus Interferon Alpha 2 (Ile 16), mammalian

Volume: ml

Activity: x 10 units/ml

Specific Activity: x 10 units/mg

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

Endotoxin: < 1 EU/ g

Molecular Weight: 19.5 kDa

Purity: > 95%

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

Source: Cynomolgus (*Macaca Fascicularis*) genomic DNA expressed in mammalian cell line

Synonyms: None

Accession #: None

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). For more information on protein handling, visit our Resource Library at www.pbl assaysci.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.