



## Certificate of Analysis

### Cynomolgus Interferon Alpha 2 (Ile 16), mammalian

**Catalog No:** 16105-1

**Lot No:** 7708

**Expiration:** July 2, 2026

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

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

**Volume:** 0.100 ml

**Activity:**  $1.89 \times 10^6$  units/ml

**Specific Activity:**  $1.18 \times 10^8$  units/mg

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

**Endotoxin:**  $< 1$  EU/ $\mu$ 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  $EC_{50}$  for interferon in this assay is  $\sim 5$  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- $\alpha$ 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 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  $-20^{\circ}\text{C}$  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. 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.pblassaysci.com](http://www.pblassaysci.com).

### Authorization

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

Date: August 7, 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.