



promising results

GTCTCTTACC CCGATGTTCA ACCAAAAG ACTTACTACC TTTATTTTAT GTTTACTTTT TATAGATTTG CTTTTTATC
TCCTTACTGC CCGTCAACAA ACACTAAA AAAAAAGTCA AATACTACTA CATCAAAACG CATATTCCCT AGAAAAAA

Cynomolgus IFN Alpha 2 (ILE 16) (mammalian)

Product Number: 16105-1

Lot Number: 5852

Expiration Date: August 13, 2014

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

Description: Recombinant Cynomolgus IFN Alpha 2 (ILE 16), mammalian

Volume: 0.100 ml

Activity: 2.71×10^6 units/ml

Specific Activity: 1.69×10^8 units/mg

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

Endotoxin: < 1 EU/ μ 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 ~ 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- 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]. 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

Special Conditions/Comments: After receipt, this product should be kept 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 the PBL website at www.interferonsource.com

Authorization

Released by: _____

Date: September 10, 2013

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL InterferonSource. 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.00

131 Ethel Rd West, Suite 6, Piscataway, NJ 08854 • T: +1-732-777-9123 • F: +1-732-777-9141
E: info@interferonsource.com • W: www.interferonsource.com