

Certificate of Analysis

Human Interferon Omega

Catalog No: 11395-1

Lot No: 7845

Expiration: August 31, 2028

Size: 25 µg/vial

Description: Recombinant Human Interferon Omega (Hu-IFN-ω)

Volume: 0.5 ml

Concentration: 0.05 mg/ml

Specific Activity: 1 x 10⁸ units/mg

Buffer: Phosphate buffered saline (PBS) containing 2% bovine serum albumin (BSA)

Molecular Weight: 20 kDa

Purity: > 95%

Source: Recombinant Hu-IFN-ω expressed in *E. coli*

Synonyms: IFN-ω

Accession #: X58822

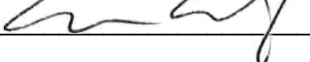
Assay Used to Measure Bioactivity: Interferon was titrated with the use of the cytopathic effect inhibition assay as described [Rubinstein, S., Familletti, P.C., and Pestka, S. (1981) "Convenient Assay for Interferons," *J. Virol.* 37, 755-758; Familletti, P.C., Rubinstein, S., and Pestka, S. (1981) "A Convenient and Rapid Cytopathic Effect Inhibition Assay for Interferon," in *Methods in Enzymology*, Vol. 78 (S. Pestka, ed.), Academic Press, New York, 387-394]. In this antiviral assay for interferon about 1 unit/ml of interferon is the quantity necessary to produce a cytopathic effect of 50%. As there is no international standard for human interferon omega, the units are determined with respect to the international reference standard for human interferon alpha A (Hu-IFN-αA) provided by the National Institutes of Health [see Pestka, S. (1986) "Interferon Standards and General Abbreviations," in *Methods in Enzymology* (S. Pestka, ed.), Academic Press, New York 119, 14-23]. Unit of activity measured on human lung carcinoma cell line A549 with encephalomyocarditis virus (EMCV).

Shipping Conditions: Dry Ice

Physical State of Product During Shipping: Frozen

Storage Conditions/Comments: After receipt, the product should be kept at -20°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 2% bovine serum albumin (BSA). For more information on protein handling, visit our Resource Library at www.pblassaysci.com.

Authorization

Released by: _____ 

Date: February 20, 2026

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. 04