

Certificate of Analysis

Human Interferon Alpha A (Alpha 2a)

Catalog No: 11100-9

Lot No: 7855

Size: 0.065 mg/vial

Description: Recombinant Human Interferon Alpha A (Alpha 2a) (Hu-IFN- α -2a)

Volume: 0.050 ml

Concentration: 1.29 mg/ml

Activity: 4.11×10^8 units/ml (MDBK/VSV)

Specific Activity: 3.18×10^8 units/mg (MDBK/VSV)

Buffer: 20 mM Tris pH 8.0; 200 mM NaCl; 6% Glycerol

Endotoxin: < 1 EU/ μ g

Molecular Weight: 19.2 kDa

Purity: > 95%

Purification Method: A combination of ion exchange, hydrophobic interaction and size exclusion chromatography.

Source: cDNA obtained from human leukocyte mRNA expressed in *E. coli*

Synonyms: None

Accession #: V00549

Assay Used to Measure Bioactivity: Interferon was titrated with the use of the cytopathic effect inhibition assay as described:

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- 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

Storage Conditions/Comments: After receipt, the product may be stored at -20°C for short-term use (\leq 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. For more information on protein handling, visit our Resource Library at www.pblassaysci.com.

Authorization

Released by: _____ 

Date: March 3, 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. 03