

Certificate of Analysis

Human Interferon Alpha 1 (Alpha D)

Catalog No: 11175-9

Lot No: 7869

Size: 0.010 mg/vial

Expiration Date: April 16, 2027

Description: Recombinant Human Interferon Alpha 1 (Alpha D) (Hu-IFN- α 1; Hu-[Ala¹¹⁴] IFN- α D)

Volume: 0.100 ml

Specific Activity of Source Material: 1.14 x 10⁸ units/mg (MDBK/VSV)

Concentration: 0.100 mg/ml

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

Molecular Weight: 19.4 kDa

Purity: \geq 95%

Purification Method: Purified from *E. coli* by a combination of ion exchange, hydrophobic interaction and size exclusion chromatography

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

Human Gene: IFNA1

Synonyms: Hu-[Ala¹¹⁴] IFN- α D

Accession #: J00210

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. 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) or tissue culture media with serum. Dilution of material below 2 x 10⁵ units/ml for freezing is not recommended. One freeze-thaw cycle is equivalent to thawing an aliquot prepared from the material received. For more information on protein handling, visit our Resource Library at www.pblassaysci.com.

Authorization

Released by: _____

Date: April 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. 02