

Certificate of Analysis

Mouse Interferon Beta, mammalian

Catalog No: 12405-1

Lot No: 7757R

Expiration: July 15, 2027

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

Description: Recombinant Mouse Interferon Beta, mammalian

Volume: 0.100 ml

Activity: 3.30×10^6 units/ml

Specific Activity: 6.60×10^8 units/mg

Buffer: 20 mM Sodium Phosphate, pH 6.4; 300 mM NaCl; 5% Glycerol; 0.1% bovine serum albumin (BSA)

Endotoxin: < 1 EU/ μ g

Molecular Weight: 19.6 kDa

Purity: $> 95\%$

Purification Method: Purified from mammalian tissue culture media by a combination of ion exchange, hydrophobic interaction and size exclusion chromatography.

Source: Gene was obtained from mouse DNA expressed in mammalian cells

Synonyms: Mouse Fibroblast Interferon

Accession #: K00020

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

Mouse (L929/EMCV) – performed as described [Rubinstein, *et al.* (1981) *J. Virol.* 37(2):755]. The EC_{50} for interferon in this assay is ~ 2.5 U/ml. Lot Activity was derived from multiple determinations in the above assay.

The units are determined by use of a Mu-IFN- β laboratory standard calibrated to the international reference standard for mouse interferon beta 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 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 our Resource Library at www.pbl assaysci.com.

Authorization

Released by: _____

Date: July 21, 2025

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. 05