



Certificate of Analysis

Mouse Interferon Alpha All-Subtype Neutralizing Viral Protein

Catalog No: 12185-1

Lot No:

Size: 10 µg/vial

Description: Mouse Interferon Alpha All-Subtype Neutralizing Viral Protein

Volume: 0.1 ml

Concentration: 0.1 mg/ml

Molecular Weight: 50kDa

Activity (EC₅₀): 0.3 – 11.5 ng/ml (varies by subtype)

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

Endotoxin: < 1 EU/µg

Purity: > 95%

Purification Method: A combination of ion exchange and size exclusion chromatography

Specificity: Neutralizes all mouse interferon alpha subtypes

Cross-Reactivity:

- Neutralizes all human interferon alpha subtypes, human interferon beta, human interferon omega, rat interferon alpha 1 and mouse limitin
- Does not neutralize mouse interferon beta, human interferon gamma, human interferon lambda 1, human interferon lambda 2 or human interferon lambda 3 when tested at ≥ 100X the typical neutralizing dose

Assay Used to Measure Bioactivity: Activity is reported as the concentration of protein required to neutralize one unit of mouse interferon alpha 4 (Mu-IFN-α4) to a 50% endpoint. Interferon was titrated with the use of a modified cytopathic effect inhibition assay [Rubinstein, S., Familletti, P.C., and Pestka, S. (1981) "Convenient Assay for Interferons," *J. Virol.* 37, 755-758]. In this antiviral assay, approximately 1 unit/ml of interferon is the quantity necessary to reduce the cytopathic effect on L929 cells challenged with EMCV to 50%. The units are determined with respect to the international reference standard for Mu-IFN-α provided by the National Institutes of Health (Ga02-901-511).

Tested Applications: Neutralization, Western Blot

Optimal dilutions should be determined by each laboratory for each application.

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 (≤ 6 months). For long-term storage, we recommend storing the product 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:

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.