



Certificate of Analysis

Anti-Mouse Interferon Alpha, Rabbit Serum (PAb)

Catalog No: 32100-1

Lot No: 7558

Size: $\geq 2 \times 10^4$ Neutralizing Units/vial

Description: Rabbit Polyclonal Antibody against Mouse Interferon Alpha

Volume: 0.100 ml

Activity: 3.95×10^5 NU/ml

Concentration: 0.42 mg/ml (estimate)

Form: Neat serum diluted in phosphate buffered saline

Antigen: Recombinant Mouse Interferon Alpha A (expressed in *E. coli*)

Assay Used to Measure Bioactivity: One Neutralization Unit (NU) is the amount of antiserum required to neutralize one unit of mouse interferon alpha (Mu-IFN- α) to a 50% endpoint. Interferon was titrated with the use of the cytopathic effect inhibition assay using L929 cells as described [Rubinstein, S., Familletti, P.C., and Pestka, S. (1981) "Convenient Assay for Interferons," *J. Virol.* 37, 755-758] with the exception that EMCV was used as the challenge virus. In this antiviral assay for interferon, about 5 unit/ml of interferon is the quantity necessary to produce a cytopathic effect of 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). This material is prepared specifically for effective neutralization of Mu-IFN- α .

Tested Applications: Neutralization

Suggested Applications: ELISA; immunoprecipitation; immunohistochemistry

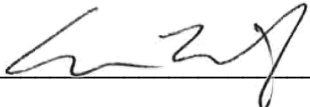
Please note that these applications are presented for suggested use only and have not been fully evaluated by PBL.

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. 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). For more information on protein handling, visit our Resource Library at www.pblassaysci.com.

Authorization

Released by: _____ 

Date: November 15, 2022

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.