



Certificate of Analysis

Anti-Mouse Interferon Gamma, Rabbit Serum (PAb)

Catalog No: 32500-1

Lot No:

Size: 2×10^4 Neutralizing Units/vial

Description: Rabbit polyclonal antibody against Mouse Interferon Gamma

Volume: ml

Activity: x10 NU / ml

Concentration: mg/ml (estimate)

Buffer: Neat serum

Antigen: Mouse interferon gamma

Assay Used to Measure Bioactivity: One neutralization unit is the amount of antiserum required to neutralize one unit of mouse interferon gamma (mouse IFN- γ) to a 50% endpoint. Interferon was titrated with the use of the cytopathic effect inhibition assay as described [Rubinstein, S., Familletti, P.C., and Pestka, S. (1981) "Convenient Assay for Interferons," *J. Virol.* 37, 755-758; Familletti, P.C., Rubinstein, S., and Pestka, S. (1981) "A Convenient and Rapid Cytopathic Effect Inhibition Assay for Interferon," in *Methods in Enzymology*, Vol. 78 (S. Pestka, ed.), Academic Press, New York, 387-394]. In this antiviral assay for interferon about 1 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 mouse IFN- γ provided by the National Institutes of Health [see Pestka, S. (1986) "Interferon Standards and General Abbreviations," in *Methods in Enzymology* (S. Pestka, ed.), Academic Press, New York 119, 14-23]. This material is prepared specifically for effective neutralization of mouse IFN- γ .

Tested Applications: Neutralization

Suggested Applications: Western blot (1:500); immunoprecipitation; immunohistochemistry; ELISA

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, this 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.pblassaysci.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.

Rev. 02