



Certificate of Analysis

Anti-Human Interferon Beta, Rabbit IgG, protein A purified (PAb)

Catalog No: 31405-1

Lot No:

Size: 2×10^4 Neutralizing Units/vial

Description: Rabbit Polyclonal Antibody against Human Interferon Beta, protein A purified

Volume: ml

Activity: x 10 NU/ml

Concentration: mg/ml

Buffer: 0.3 M Sodium Bicarbonate; 0.2 M Sodium Chloride

Antigen: Recombinant human interferon beta (expressed in *E. coli*)

Assay Used to Measure Bioactivity: One neutralization unit is the amount of antiserum required to neutralize one unit of human interferon beta 1a (Hu-IFN- β 1a) to a 50% endpoint. Interferon was titrated with the use of the cytopathic effect inhibition assay [Rubinstein, S., Familletti, P.C., and Pestka, S. (1981) *J. Virol.* 37:75] using A549 cells and EMCV was used as the challenge virus [Budd *et al.* (1985) *Canc. Chem. Pharm.* 12:39]. 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 by use of a laboratory standard calibrated with respect to the international reference standard for Hu-IFN- β provided by the National Institutes of Health [Gb23-902-531].

Tested Applications: Neutralization

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

Suggested Applications: ELISA; Western blot; 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:

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. 01