



Certificate of Analysis

Anti-Rat Interferon Beta, Rabbit IgG, Protein A purified (PAb)

Catalog No: 33400-1

Lot No:

Expiration:

Size: 500 µg/vial

Description: Rabbit Polyclonal Antibody against Rat Interferon Beta, Protein A purified, neutralizing

Concentration: 1 mg/ml after reconstitution with 0.5 ml sterile distilled water

Reconstitution: Dissolve contents of the vial by injection of 0.5 ml sterile distilled water

Buffer: 0.5 ml Phosphate-buffered saline (PBS) + 125 mM trehalose prior to lyophilization

Purification Method: Antibodies were sequentially purified by ammonium sulphate precipitation and Protein A-affinity chromatography

Antigen: Recombinant rat interferon beta derived from Chinese hamster ovary (CHO) cells

Activity: > 5 x 10³ neutralizing units/mg

Specificity: Neutralizes both natural and recombinant rat interferon beta. Cross-reactivity with human interferon beta has not been tested.

Assay Used to Measure Bioactivity: One neutralization unit is the amount of antiserum required to neutralize one unit of recombinant rat interferon beta (rat IFN- β) to a 50% endpoint (one unit corresponds to approximately 15 pg of purified rat IFN- β). 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%. As there is no international rat IFN- β standard, the units are determined with respect to the international reference standard for mouse interferon beta (mouse IFN- β) 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 rat IFN- β .

Suggested Applications: Western blot; ELISA; Neutralization

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

Reference:

1. Ruuls, S.R. *et al.* (1996) *J. Immunol.* 157: 5721;
2. Zou, L.P. *et al.* 1999. *Neuroscience Res.* 56: 123

Shipping Conditions: Wet Ice

Physical State of Product During Shipping: Lyophilized

Storage Conditions/Comments: Upon receipt, this product is stable at 4°C until the expiration date listed above. After reconstitution, the contents can be safely stored at 4°C for one month or at -20°C for one year 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. Add 0.02% sodium azide to prevent bacterial growth. For more information on protein handling, visit our Resource Library at www.pblassaysci.com.



pbl assay science

131 Ethel Road West, Suite 6 | Piscataway, NJ 08854 USA | T:+1.732.777.9123 | F:+1.732.777.9141 | E: info@pbl assaysci.com | W: 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.

Rev. 03

A decorative graphic at the bottom of the page consisting of a wavy, multi-colored shape that transitions from light green on the left to bright blue on the right.