



## Certificate of Analysis

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

**Catalog No:** 33100-1

**Lot No:** 6230

**Expiration:** June 30, 2017

**Size:** 500 µg/vial

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

**Concentration:** 1.0 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

**Antigen:** Recombinant rat interferon alpha derived from baculovirus

**Isotype:** Rabbit Ig

**Purification Method:** Ammonium sulfate precipitation and protein-A affinity chromatography

**Activity:** > 10<sup>4</sup> neutralizing units/mg

**Specificity:** Neutralizes both natural and recombinant rat interferon alpha

**Assay Used to Measure Bioactivity:** One neutralization unit is the amount of antiserum required to neutralize one unit of recombinant rat interferon alpha (rat IFN- $\alpha$ ) in the antiviral bioassay. As there is no international rat IFN- $\alpha$  standard, the units are determined with respect to a laboratory standard preparation of recombinant rat IFN- $\alpha$  (one unit corresponds to approximately 5 pg of purified rat IFN- $\alpha$ ). This material is prepared specifically for effective neutralization of rat IFN- $\alpha$ .

**Suggested Applications:** Western blot; *in vitro* neutralization

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

#### Reference:

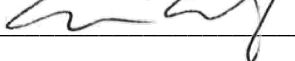
1. Van der Meide, P.H. *et al.* (1986) *Methods Enzymol.* 119: 441.

**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.pblsaysci.com](http://www.pblsaysci.com).

#### Authorization

Released by: \_\_\_\_\_ 

Date: March 12, 2015

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.