



Certificate of Analysis

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

Catalog No: 33100-1

Lot No: 7646

Expiration: January 31, 2028

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: > 10⁴ 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- α) in the antiviral bioassay. As there is no international rat IFN- α standard, the units are determined with respect to a laboratory standard preparation of recombinant rat IFN- α (one unit corresponds to approximately 5 pg of purified rat IFN- α). This material is prepared specifically for effective neutralization of rat IFN- α .

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.pblassaysci.com.

Authorization

Released by: _____

Date: August 4, 2023

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.