



Certificate of Analysis

Anti-Mouse Interferon Alpha, Chicken IgY, Purified (PAb)

Catalog No: 32120-1

Lot No:

Size: 500 µg/vial

Description: Chicken Polyclonal Antibody against Mouse Interferon Alpha, non neutralizing

Volume: ml

Concentration: mg/ml

Buffer: Phosphate buffered saline (PBS) containing 0.1% BSA

Antigen: Recombinant mouse interferon alpha

Isotype: Chicken IgY

Specificity: Binds to but does not neutralize mouse IFN alpha

Tested Applications: Western Blot (1:500), ELISA (direct binding)

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. 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