



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.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