



## Certificate of Analysis

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

**Catalog No:** 32120-1

**Lot No:** 7337

**Size:** 500 µg/vial

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

**Volume:** 0.100 ml

**Concentration:** 5 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](http://www.pbl assaysci.com).

### Authorization

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

Date: January 20, 2021

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.