



## Certificate of Analysis

### Mouse Mab to Human IFN Alpha

**Catalog No:** 710636-9

**Lot No:** 7520

**Size:** 0.1 mg/vial

**Description:** Monoclonal Antibody against Human Interferon Alpha

**Volume:** 0.014 ml

**Concentration:** 7.13 mg/ml

**Buffer:** 0.2 M NaHCO<sub>3</sub>, 0.3 M NaCl, pH 8

**Endotoxin:** <1 EU/μg

**Isotype:** mouse IgG1 kappa

**Suggested Applications:** ELISA, Neutralization

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

**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. When thawing, 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: Michael Skawinski

Date: May 16, 2022

**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.**