

## Certificate of Analysis

### Mouse Monoclonal Antibody Against Human IFN Beta

**Catalog No:** 710906-9

**Lot No:** 7818

**Size:** 1 mg/vial

**Description:** Mouse Anti-Human IFN-Beta

**Volume:** 0.650 ml

**Concentration:** 1.54 mg/ml

**Buffer:** 20 mM Tris pH 9, 100 mM NaCl, 10% Glycerol

**Endotoxin:** < 1 EU/ $\mu$ g

**Isotype:** Mouse IgG<sub>1</sub>, kappa

**Suggested Applications:** Western blot; immunoprecipitation; immunohistochemistry; ELISA

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 ( $\leq$  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. Re-freezing 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](http://www.pblassaysci.com).

### Authorization

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

Date: October 8, 2025

**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. 01