

Certificate of Analysis

Biotinylated Mab to Human IFN Beta

Catalog No: 710992-9

Lot No: 7834

Size: 0.2 mg/vial

Description: Monoclonal Antibody against Human Interferon Beta, biotin labeled

Volume: 0.400 ml

Concentration: 0.50 mg/ml

Buffer: Aqueous, protein-containing diluent and stabilizer

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

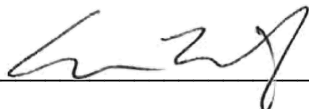
Shipping Conditions: Dry ice

Physical State of Product During Shipping: Frozen

Storage Conditions/Comments: After receipt, the product may be stored at -20°C for short-term use (≤ 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.

Authorization

Released by: _____



Date: December 18, 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.00