



Certificate of Analysis

Mouse Mab to Human IFN Omega

Catalog No: 75398-9

Lot No: 7158

Size: 0.1 mg/vial

Expiration Date: June 30, 2022

Description: Monoclonal Antibody against Human Interferon Omega

Volume: 0.100 ml

Concentration: 1 mg/ml

Buffer: Phosphate Buffered Saline, pH 7.2-7.4

Antigen: human interferon omega

Isotype: IgG1

Suggested Applications: 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. 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: January 3, 2020

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.