



Certificate of Analysis

Biotinylated Mouse Mab to Human IFN Alpha

Catalog No: 710745-9

Lot No: 7521

Size: 0.1 mg/vial

Description: Biotin Conjugated Mouse Monoclonal Antibody against Human Interferon Alpha

Volume: 0.105 ml

Concentration: 0.956 mg/ml

Buffer: Stabilzyme Select

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. For more information on protein handling, visit our Resource Library at www.pblassaysci.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.

Rev. 00