

Certificate of Analysis

FITC Conjugated Anti-Human Interferon Beta, Clone MMHB-3 (MAb)

Catalog No: 21400-3

Lot No: 7225

Expiration: April 24, 2021

Size: ≥ 20 µg/vial

Description: Fluorescein Isothiocyanate (FITC) Conjugated Mouse Monoclonal Antibody against Human Interferon Beta

Clone: MMHB-3

Volume: 0.5 ml

Antibody Concentration: 0.05 mg/ml

F/P Ratio: 9.21

Buffer: Phosphate buffered saline (PBS) containing 1% BSA and 0.05% Kathon as a preservative, pH 7.2

Antigen: Human Interferon Beta

Isotype: Mouse IgG1 κ

Tested Applications: Flow Cytometry: 0.25 µg/ml of antibody is sufficient for labeling of 1×10^6 cells in 100 µl total volume of A549-human lung epithelial cells. It is recommended that each investigator determine appropriate dilutions for other cell lines. Immunocytochemistry: It is recommended that each investigator determine appropriate dilutions for individual use.

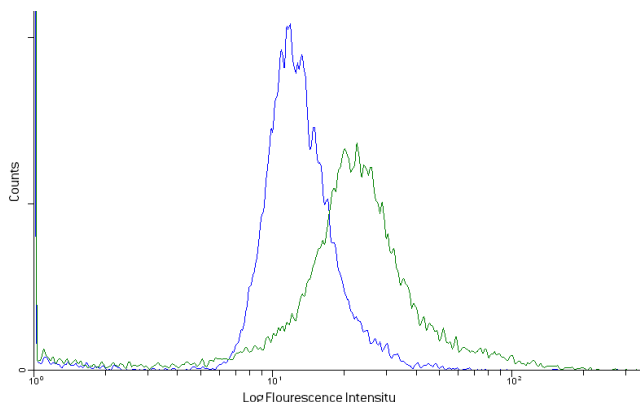


Figure 1. Log Fluorescence Intensity-FITC Human Interferon Beta. Expression of endogenous Human Interferon Beta in A549 human lung epithelial cells infected with Encephalomyocarditis virus (EMCV). Flow cytometry analysis on non-induced cells (blue) and induced cells (green).

Shipping Conditions: Wet Ice

Physical State of Product during Shipping: Liquid

Storage Conditions/Comments: Photo-sensitive material protect from prolonged exposure to light. Store at 2-8°C, DO NOT FREEZE. For more information on protein handling, visit our Resource Library at www.pbl assaysci.com.

Authorization

Released by: _____

Date: May 21, 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.