



Certificate of Analysis

R-PHYCOERYTHRIN (PE)-CONJUGATED MOUSE MONOCLONAL ANTIBODY AGAINST HUMAN INTERFERON ALPHA/BETA RECEPTOR 1

Catalog No: 21370-3

Lot No: 6417R

Expiration: May 8, 2017

Size: 50 tests/vial

Description: Mouse Anti-Human Interferon Alpha/Beta Receptor 1, labeled with PE

Clone: MMHAR-3

Volume: 1 ml

Buffer: phosphate-buffered saline (PBS) containing 1.5% BSA and 0.01% Kathon as a preservative, pH 7.4

Antigen: Human Interferon Alpha/Beta Receptor 1

Isotype: IgG₁

Tested Application: Flow Cytometry: 20 µl of antibody is sufficient for labeling of 1x10⁶ cells in 100 µl total volume

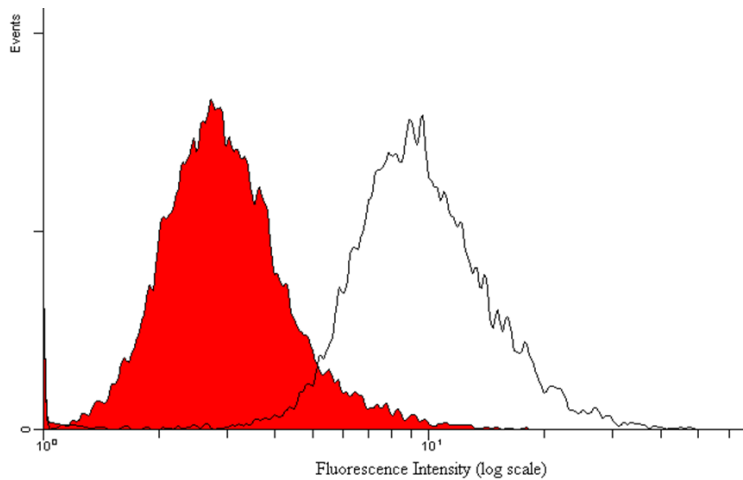


Figure 1: Flow cytometry analysis of KG-1 cells showed expression of Human Interferon Alpha/Beta Receptor 1. Isotype control (shaded histogram) and human interferon alpha/beta receptor (un-shaded histogram)

Shipping Conditions: Wet Ice

Physical State of Product During Shipping: Liquid

Storage Conditions/Comments: Store at 2-8°C, DO NOT FREEZE. Photo-sensitive material protect from prolonged exposure to light. Approximate concentration: 50ug/ml. For more information on protein handling, visit our Resource Library at www.pblassaysci.com.

Authorization

Released by: _____

Date: **January 10, 2016**

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.