



promising results

GTCTCTTACC CGGATGTTCA ACCAAAAG ACTTACTACC TTTATTTTAT GTTTACTTTT TATAGATTGT CTTTTTATC
TCCTACTGCG CGTGCAACAA ACACTAAA AAAAAAGTGA AATACTACTA CATCAAAACG CATATTCCCT AGAAAAAAA

R-PHYCOERYTHRIN (PE)-CONJUGATED MOUSE MONOCLONAL ANTIBODY AGAINST HUMAN INTERFERON ALPHA/BETA RECEPTOR CHAIN 2

Product Number: 21385-3

Lot Number:

Expiration: 2.5 years from date of manufacturing of new stock (input actual date)

Size: 50 tests

Description: Mouse Anti-Human Interferon Alpha/Beta Receptor Chain 2, labeled with PE

Clone: MMHAR-2

Contents: 50 tests

Volume: 1 ml

Buffer: Phosphate buffered saline (PBS) containing 1.5% bovine serum albumin (BSA) and 0.1% Kathon as a preservative, pH 7.4

Antigen: Human Interferon Alpha/Beta Receptor Chain 2

Isotype: IgG_{2a}

Special Condition/comments: Photo-sensitive material protect from prolonged exposure to light. Store at 2-8 °C, DO NOT FREEZE.

Tested Applications: Flow Cytometry: 20uL is sufficient for labeling of 1×10^6 cells in 100uL total volume.

Figure1: Flow cytometry analysis of KG-1 cells showed expression of Human Interferon Alpha/Beta Receptor Chain 2. Isotype control (shaded histogram) and human interferon alpha receptor (un-shaded histogram).

Shipping Conditions: Wet ice

Physical State of Product during Shipping: Liquid

Authorization

Released by: _____

Date: October 1, 2013

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL InterferonSource. 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.03

131 Ethel Rd West, Suite 6, Piscataway, NJ 08854 • T: +1-732-777-9123 • F: +1-732-777-9141
E: info@interferonsource.com • W: www.interferonsource.com