

## Certificate of Analysis

### Mouse Monoclonal Antibody against Human Interferon Alpha/Beta Receptor 1, Clone MMHAR-3 (MAb)

**Catalog No:** 21370-1

**Lot No:**

**Size:** 50 µg/vial

**Description:** Mouse Anti-Human Interferon Alpha/Beta R1 (IFNAR1), non-neutralizing

**Clone:** MMHAR-3

**Concentration:** 0.5 mg/ml

**Buffer:** Phosphate-buffered saline (PBS)

**Endotoxin:** < 1 EU/µg

**Antigen:** U266 Myeloma Cells

**Isotype:** Mouse IgG<sub>1</sub> kappa

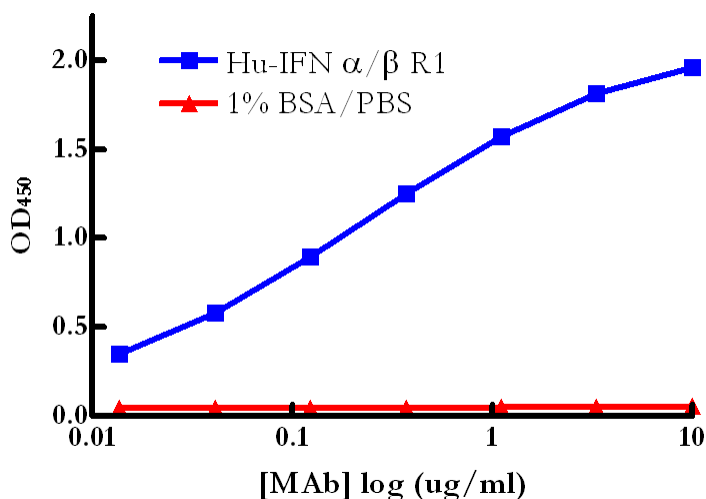
**Purity:** > 95%

**Purification Method:** Protein G affinity chromatography

**Specificity:** Detects human interferon alpha/beta R1

**Tested Applications:** *Optimal dilutions should be determined by each laboratory for each application.*

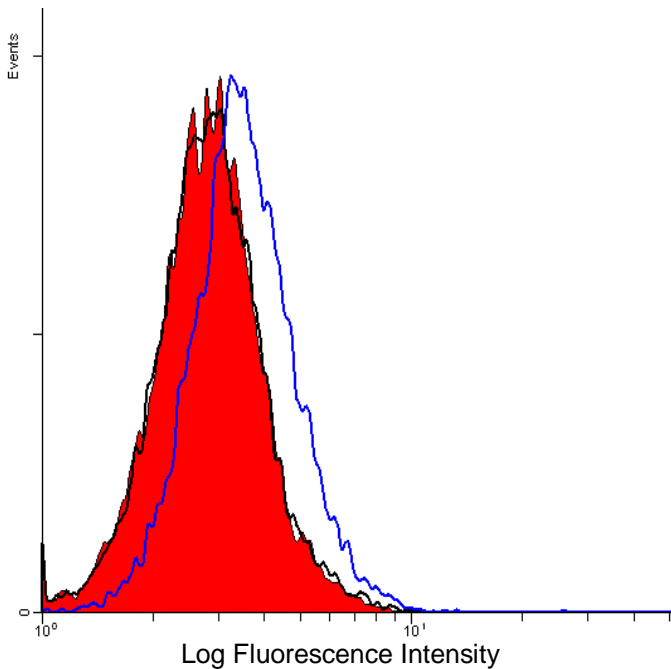
#### Direct Binding ELISA



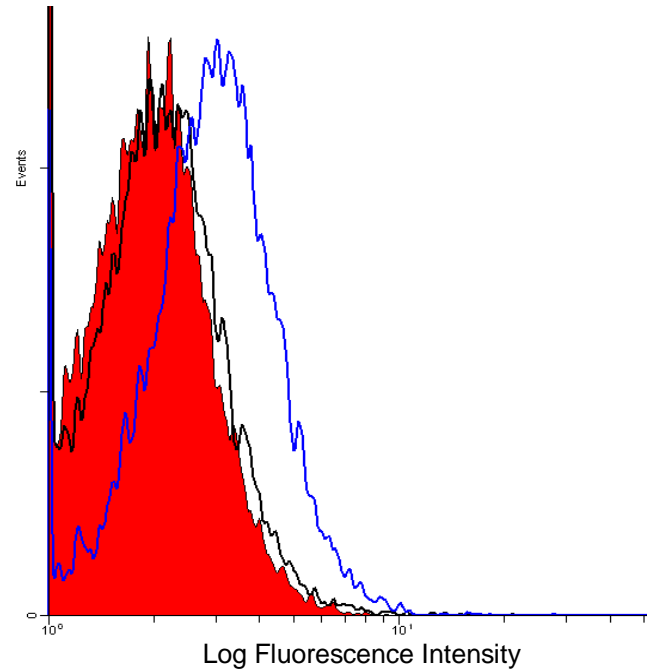
**Figure 1.** Representative binding curve of antibody to mammalian expressed human interferon alpha/beta R1 protein and to 1% BSA/PBS in a Direct Binding ELISA. High-binding polystyrene plates were coated with either 1 µg/ml of human interferon alpha/beta R1 protein or 1% BSA/PBS. Thereafter, titrations of the product were added to the wells. Goat anti-mouse IgG conjugated to HRP was used as the detection antibody. Colorimetric detection was performed using 3,3',5,5'-Tetramethylbenzidine (TMB) substrate. The HRP-TMB reaction was stopped using a diluted H<sub>2</sub>SO<sub>4</sub>/HCL solution.

### Flow Cytometry

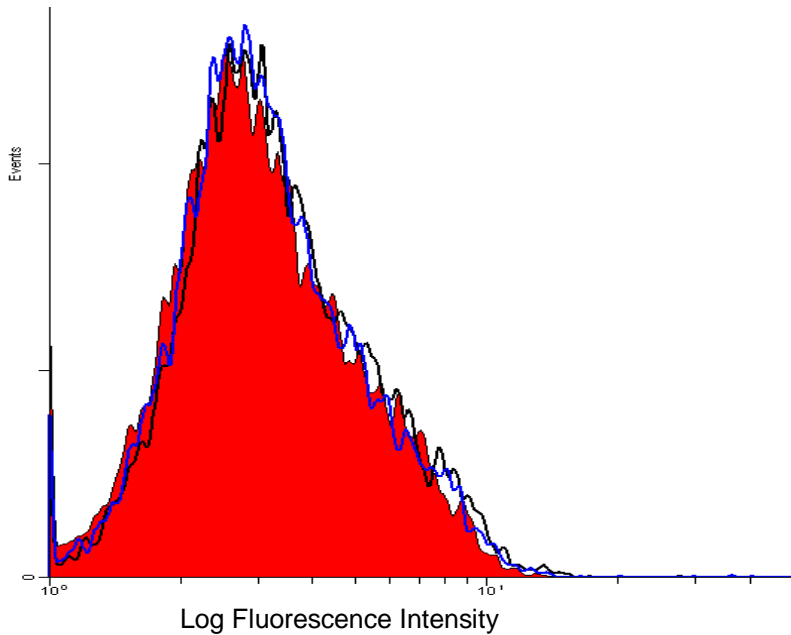
1  $\mu$ g/ml of antibody is sufficient for binding  $1 \times 10^6$  cells in 100  $\mu$ l total volume. The binding of the unlabeled antibody may be visualized by adding 10  $\mu$ l of a 0.5  $\mu$ g/ml stock solution of a secondary developing reagent such as goat anti-mouse IgG conjugated to fluorochrome.



**Figure 2.** Flow analysis on U937 (human lung lymphoblast) cells: Negative control (closed histogram; no primary or secondary antibodies), Isotype Control (black open histogram; Mouse IgG1 Kappa primary with FITC-conjugated secondary) and Anti-Human IFN- $\gamma$  / R1 (blue open histogram; #21370-1 primary with FITC-conjugated secondary)



**Figure 3.** Flow analysis on Daudi (human blood lymphoblast) cells: Negative control (closed histogram; no primary or secondary antibodies), Isotype Control (black open histogram; Mouse IgG1 Kappa primary with FITC-conjugated secondary) and Anti-Human IFN- $\gamma$  / R1 (blue open histogram; #21370-1 primary with FITC-conjugated secondary)



**Figure 4.** Flow analysis on RAW (mouse macrophage) cells: Negative control (closed histogram; no primary or secondary antibodies), Isotype Control (black open histogram; Mouse IgG1 Kappa primary with FITC-conjugated secondary) and Anti-Human IFN- $\gamma$  R1 (blue open histogram; #21370-1 primary with FITC-conjugated secondary)

**Shipping Conditions:** Dry Ice

**Physical State of Product During Shipping:** Frozen

**Storage Conditions/Comments:** After receipt, this product may be stored at  $-20^{\circ}\text{C}$  for short-term use (6 months). For long-term storage, we recommend storing the product at  $-70^{\circ}\text{C}$  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. Re-freezing 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](http://www.pbl assaysci.com).

## Authorization

Released by: \_\_\_\_\_

Date:

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. 01