



## Certificate of Analysis

### Anti-Human Interferon Alpha/Beta Receptor 1, Clone MMHAR-1 (MAb)

**Catalog No:** 21375-1

**Lot No:** 7277

**Expiration:** July 31, 2021

**Size:** 500 µg/vial

**Description:** Mouse Monoclonal Antibody against Human Interferon Alpha/Beta R1

**Clone:** MMHAR-1

**Concentration:** 0.5 mg/ml; after reconstitution with 1 ml sterile PBS

**Buffer:** Lyophilized from 125µL of a 0.2 µm filtered solution in phosphate-buffered saline (PBS) containing 5% trehalose

**Endotoxin:** < 1 EU/µg

**Antigen:** S21-derived Hu-IFN-α/β R1 extracellular domain

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

**Purity:** > 95%

**Purification Method:** Protein G affinity chromatography

**Specificity:** Detects human interferon alpha/beta R1; in direct ELISAs, this antibody shows no cross reactivity with Hu-IFN-γ R1, Hu-IFN-γ R2, Mu-IFN-α/β R2, Hu-IL-10 Rβ, Hu-IL-10 Rα or Hu-IL-20 Rβ

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

**Western Blot:** 1 - 2 µg/ml with the appropriate secondary reagents to detect human IFN-α/β R1.

**Flow Cytometry:** Dilute antibody to 25 µg/ml and add 10 µl of this solution to 1 – 2.5 x 10<sup>5</sup> cells in a total reaction volume not exceeding 200 µl. The binding of the unlabeled antibody may be visualized by adding 10 µl of a 25 µg/ml stock solution of a secondary developing reagent such as goat anti-mouse IgG conjugated to fluorochrome.

**Direct ELISA:** 0.5 - 1.0 µg/ml with the appropriate secondary reagents to detect human IFN-α/β R1.

**Shipping Conditions:** Wet Ice

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

**Storage Conditions/Comments:** Upon receipt, this product should be kept at -20 to -70°C for retention of full activity. After reconstitution, the contents of the tube should be apportioned in separate tubes so that freezing and thawing is kept to a minimum. This product may be kept at 2 to 8°C for up to 1 month and -20 to -70°C for up to 6 months after reconstitution. For more information on protein handling, visit our Resource Library at [www.pbl assaysci.com](http://www.pbl assaysci.com).

### Authorization

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

Date: August 6, 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.