



Certificate of Analysis

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

Catalog No: 21375-1

Lot No: 7300

Expiration: October 8, 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 IgG₁

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 10⁵ 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.

Authorization

Released by: _____

Date: October 9, 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.