



Certificate of Analysis

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

Catalog No: 21370-1

Lot No: 7557

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 IgG₁ kappa

Purity: > 95%

Purification Method: Protein G affinity chromatography

Specificity: Detects human interferon alpha/beta R1

Tested Application: Direct Binding ELISA.

Optimal dilutions should be determined by each laboratory for each application.

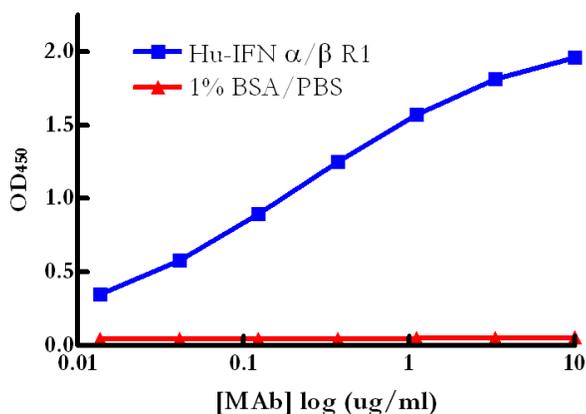


Figure 1. Representative binding curve of antibody to mammalian expressed human interferon alpha/beta R1 protein and to 1% BSA/PBS in a Direct Binding ELISA. High-binding polystyrene plates were coated with either 1 µg/ml of human interferon alpha/beta R1 protein or 1% BSA/PBS. Thereafter, titrations 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'-Tetramethyl-benzidine (TMB) substrate. The HRP-TMB reaction was stopped using a diluted H₂SO₄/HCL solution.

Shipping Conditions: Dry Ice

Physical State of Product During Shipping: Frozen

Storage Conditions/Comments: After receipt, this product may be stored at -20°C for short-term use (≤ 6 months). For long-term storage, we recommend storing the product at -70°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.pblassaysci.com.

Authorization

Released by: _____

Date: October 19, 2022

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.