

Certificate of Analysis

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

Catalog No: 21375-3

Lot No:

Expiration:

Size: 100 tests/vial

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

Clone: MMHAR-1

Volume: ml

Antibody Concentration: µg/ml

Antigen: Human interferon alpha/beta receptor 1

Isotype: Mouse IgG₁

Tested Applications: Flow Cytometry: use material as is; no preparation needed.

Recommended Concentration for Use: 10 µl/1x10⁶ cells

It is recommended that each investigator determine appropriate dilutions for individual use.

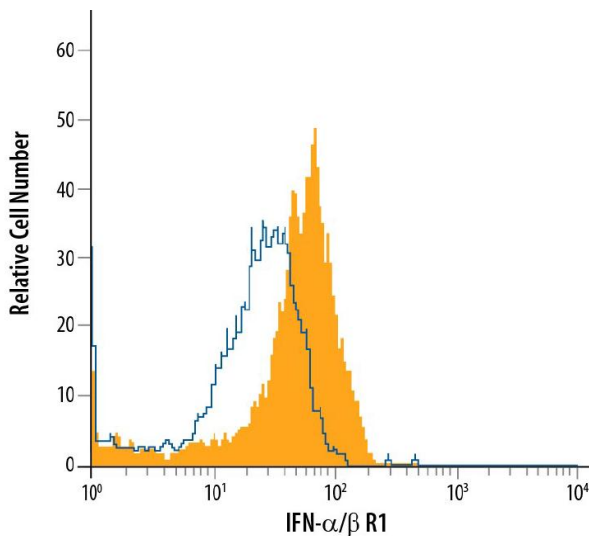


Figure 1. Human peripheral monocytes were stained with anti-human IFN- / R1-PE (filled histogram). Isotype control (open histogram).

Product Information: IFN- / R1, also known as IFNAR1, belongs to the class II cytokine receptor family of proteins. Class II cytokine receptors form heterodimeric receptor complexes that mediate class II cytokine signals. Subunits of the different receptor complexes are shared and serve multiple functions. IFN- / R1, in association with IFN- / R2, is required for propagating antiviral signal transduction triggered by IFN- and IFN-. IFN- / R1 interacts very weakly or not at all with type 1 interferons and does not stably interact with IFN- / R2. Ligands associate with IFN- / R2, and this complex subsequently forms a stable ternary assembly with IFN- / R1. IFN- / R1 also associates with IFN- R2 even in the absence of IFN- stimulation. Human- / R1 contains a nuclear localization signal in its extracellular domain that is required for receptor translocation to the nucleus following interaction with ligand.

Note: Contains sodium azide as a preservative. Sodium azide may react with lead and copper plumbing to form explosive metal azides. Flush with large volumes of water during disposal.



Shipping Conditions: Wet Ice

Physical State of Product During Shipping: Liquid

Storage Conditions/Comments: Photo-sensitive material . protect from prolonged exposure to light. **Store at 2-8°C. DO NOT FREEZE.** For more information on protein handling, visit our Resource Library at 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. 03

A decorative graphic at the bottom of the page consisting of a wavy, horizontal band with a color gradient from light green on the left to bright blue on the right.