

Certificate of Analysis

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

Catalog No: 21376-3

Lot No: 7276

Expiration: July 31, 2021

Size: 100 tests/vial

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

Clone: MMHAR-1

Volume: 1 ml

Antibody Concentration: 0.05 µg/ml

Antigen: Human interferon alpha/beta receptor 1

Isotype: Mouse IgG₁

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

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

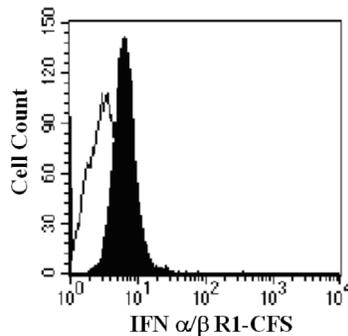


Figure 1. U937 cells were stained with anti-human IFN- α/β R1-CFS (filled histogram). Isotype control (open histogram).

Product Information: Type I interferons (IFN- α , IFN- β , IFN- ω) bind to the type I IFN receptor, also called IFN- α/β receptor. This receptor is composed of 2 chains, IFN- α/β R1 and R2.

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: 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.
Rev. 01