

Anti-Human Interferon α/β Receptor 1 (IFNAR1) Antibodies

Non-conjugated and PE-conjugated

Catalog No. 21370-1 and 21370-3

Mouse Monoclonal Antibody against Human IFNAR1 (21370-1) is a non-neutralizing unconjugated monoclonal antibody which provides scientists with a new tool for Direct Binding ELISAs and flow cytometry (FACS) studies. This antibody is reported to work in immunoprecipitation and immunohistochemistry applications.

PE Conjugated Mouse Monoclonal Antibody against Human IFNAR1 (21370-3) is a conjugated monoclonal antibody highly recommended for use in flow cytometry (FACS) studies.

21370-3 Recommended for use in Flow Cytometry

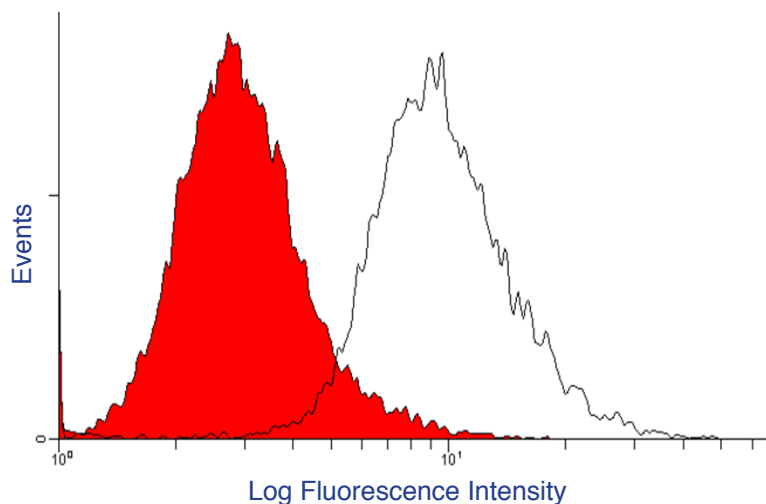


Figure 1. Flow cytometry analysis of KG-1 cells showed expression of Human Interferon Alpha/Beta Receptor 1. Isotype control (shaded histogram) and human IFN Alpha/Beta Receptor (unshaded histogram).

PBL Assay Science (formerly PBL InterferonSource)
131 Ethel Road West, Suite 6 | Piscataway, NJ 08854

Phone: (732) 777-9123

Fax: (732) 777-9141

E-mail: info@pblassaysci.com

Web: www.pblassaysci.com

21370-1 Recommended for use in Direct Binding ELISA

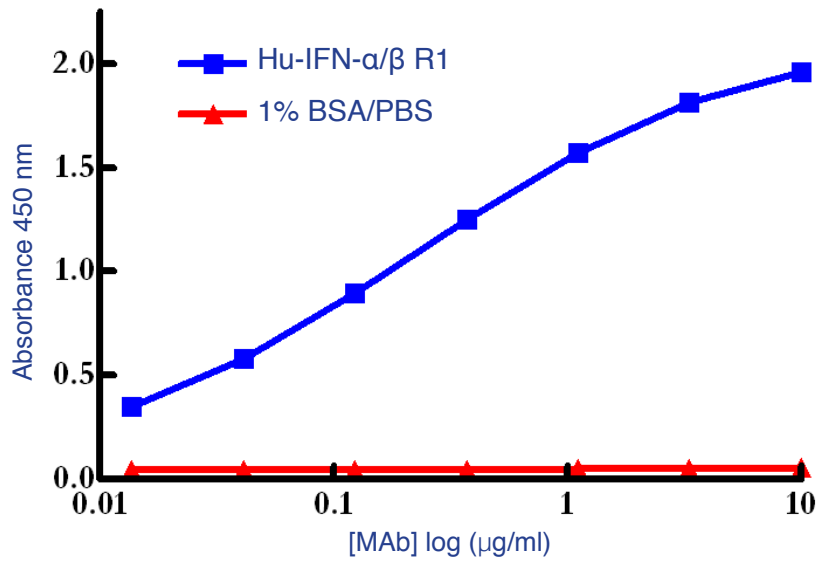


Figure 2. Representative binding curve of antibody to mammalian expressed human interferon alpha/beta R1 protein and to 1% BSA/PBS in a Direct Binding ELISA.

Product Information

Catalog Number 21370-1

21370-3

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

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

Size 50 µg

50 tests



For more information on these products, including a general cell surface staining procedure, visit our website.