

Certificate of Analysis

Anti-Mouse IFN-Alpha/Beta Receptor Chain 1 (IFNAR1), Clone MAR1-5A3 (MAb)

Catalog No: 22370-1

Lot No:

Size: 1 mg/vial

Description: MAb to Mouse Interferon Alpha/Beta Receptor 1 (MAR1-5A3), Neutralizing

Clone: MAR1-5A3

Volume: ml

Concentration: 5 mg/ml

Buffer: Phosphate Buffered Saline

Endotoxin: < 1 EU/mg

Antigen: Recombinant Mouse Interferon Alpha/Beta Receptor Chain 1

Isotype: Mouse IgG1

Purity: > 95%

Purification Method: Protein G

Specificity: Neutralizes the extracellular domain of the IFNAR1 subunit of the mouse IFN alpha/beta receptor

Antigen: Extracellular domain of mouse IFNAR-1

Bioactivity: Measured on ability to neutralize mouse IFNAR1 mediated protection of L929 cells when challenged with EMCV in the cytopathic effect assay (CPE).

Tested Applications: Neutralization

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

Suggested Applications: ELISA; Western blot; immunoprecipitation; immunohistochemistry

Please note that these applications are presented for suggested use only and have not been fully evaluated by PBL.

Shipping Conditions: Dry Ice

Physical State of Product During Shipping: Frozen

Storage Conditions/Comments: After receipt, the 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. Refreezing 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: _____

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.00