



Certificate of Analysis

Anti-Human Interferon Lambda Receptor 1, Clone MMHLR-1 (MAb)

Catalog No: 21885-1

Lot No:

Size: 50 µg/vial

Description: Mouse Monoclonal Antibody against Human Interferon Lambda Receptor Chain 1, neutralizing

Clone: MMHLR-1

Volume: 0.1 ml

Concentration: 0.5 mg/ml

Activity (EC₅₀): 10-60 ng/ml (measured against IFN-lambda 1)

Buffer: Phosphate-buffered saline (PBS) containing 0.1% bovine serum albumin (BSA)

Endotoxin: < 1 EU/µg

Antigen: Human Interferon Lambda Receptor 1 extracellular domains

Isotype: Mouse IgG₁, Kappa

Purity: > 95%

Purification Method: A combination of affinity, ion exchange and size exclusion chromatography

Specificity: Binds to extracellular domains of human interferon lambda receptor. Blocks biological action of Type III interferons.

Assay Used to Measure Bioactivity: Activity is reported as the concentration of antibody required to neutralize 100 pg of interferon lambda 1 to a 50% endpoint. Interferon was titrated with the use of a modified cytopathic effect inhibition assay [Rubinstein, S., Familletti, P.C., and Pestka, S. (1981) "Convenient Assay for Interferons," *J. Virol.* 37, 755-758]. In this antiviral assay, approximately 500 pg/ml of interferon lambda 1 is the quantity necessary to reduce the cytopathic effect on A549 cells challenged with EMCV to 50%.

Tested Applications: Neutralization, Flow Cytometry

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

References: Donnelly, R.P., *et al.* (2003) *Nat Immunol* 4: 69-77.

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. Thaw product vial by incubation in cold tap water until just thawed – 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.