



Certificate of Analysis

Anti-Human Interleukin-28A/Interferon Lambda 2, neutralizing, clone MMHL-2 (MAb)

Catalog No: 21720-1

Lot No:

Size: 200 µg/vial

Description: Mouse Monoclonal Antibody against Human Interleukin-28A/Interferon Lambda 2, neutralizing

Clone: MMHL-2

Volume: 0.1 ml

Concentration: 2 mg/ml

Activity (EC₅₀): 750 - 3000 ng/ml

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

Endotoxin: < 1 EU/ g

Antigen: Human IL28A/Interferon Lambda 2

Isotype: Mouse IgG₁ kappa

Purity: > 95%

Purification Method: Protein G affinity chromatography

Specificity: Neutralizes human IL-28A/interferon lambda 2. Approximately 3% and 8% cross-reactivity is observed with human interferon lambda 1 and human interferon lambda 3, respectively.

Assay Used to Measure Bioactivity: Human interferon lambda 2 protein was titrated with the use of the cytopathic effect inhibition assay as described [Rubinstein, S., Familletti, P.C., and Pestka, S. (1981) "Convenient Assay for Interferons," *J. Virol.* 37, 755-758]. Neutralizing activity is reported as the concentration of antibody required to neutralize 50% of the protective effect of 5.7 ng/ml of human interferon lambda 2 (Human IL-28A) in a cytopathic assay of A549 cells challenged with EMCV [Budd *et al.* (1985) Phase-I Trial of Ultrapure™ Human Leukocyte Interferon in Human Malignancy, *Canc. Chem. Pharm.* 12, 39-42].

Tested Applications: Neutralization

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

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. Re-freezing 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.