



## Certificate of Analysis

### Anti-Human Interleukin-28B/Interferon Lambda 3, neutralizing, clone MMHL-3 (MAb)

**Catalog No:** 21730-1

**Lot No:**

**Size:** 100 µg/vial

**Description:** Mouse Monoclonal Antibody against Human Interleukin-28B/Interferon Lambda 3, neutralizing

**Clone:** MMHL-3

**Volume:** 0.1 ml

**Concentration:** 1 mg/ml

**Activity (EC<sub>50</sub>):** 250 - 1050 ng/ml

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

**Endotoxin:** < 1 EU/ g

**Antigen:** Human IL28B/Interferon Lambda 3

**Isotype:** Mouse IgG<sub>1</sub> kappa

**Purity:** > 95%

**Purification Method:** Protein G affinity chromatography

**Specificity:** Neutralizes human IL-28B/interferon lambda 3. Less than 0.1% cross-reactivity is observed with human interferon lambda 1 and human interferon lambda 2.

**Assay Used to Measure Bioactivity:** Human interferon lambda 3 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 1.3 ng/ml of human interferon lambda 3 (Human IL-28B) 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](http://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