

## Certificate of Analysis

### Anti-Human Interferon Lambda Receptor 1 (MAb), Clone HLR14

**Catalog No:** 21886-1

**Lot No:**

**Size:** 50 µg/vial

**Description:** Mouse Monoclonal Antibody against Human Interferon Lambda Receptor 1

**Clone:** HLR14

**Volume:** 0.1 ml

**Concentration:** 0.5 mg/ml

**Buffer:** 0.3 M Sodium Bicarbonate, 0.2 M Sodium Chloride, pH 8.0

**Endotoxin:** < 1 EU/µg

**Antigen:** Human Interferon Lambda Receptor 1

**Isotype:** Mouse IgG1

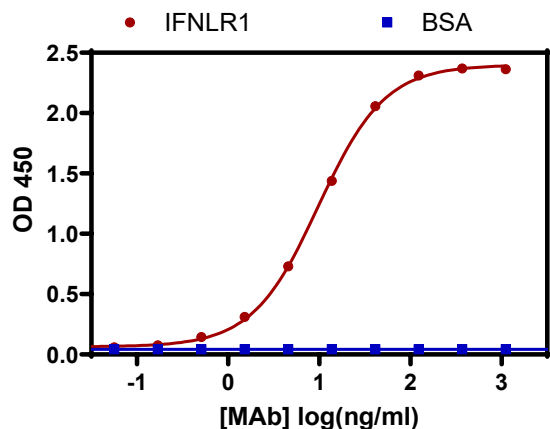
**Purity:** > 95%

**Purification Method:** Protein G

**Specificity:** Binds to human interferon lambda receptor 1 with high affinity

**Tested Applications:** Direct Binding ELISA, Flow Cytometry

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



**Figure 1.** Representative binding curve of antibody to mammalian expressed human interferon lambda receptor (circle) and to 0.1% BSA/PBS (square) in a Direct Binding ELISA. Medium-binding polystyrene plates were coated with either 1 µg/ml human interferon lambda receptor 1 or 0.1%BSA/PBS. Thereafter, titrations of the product were added to the wells. Donkey anti-mouse IgG conjugated to HRP was used as the detection antibody. Colorimetric detection was performed using 3, 3', 5, 5'-Tetramethylbenzidine (TMB) substrate. The HRP-TMB reaction was stopped using a diluted H<sub>2</sub>SO<sub>4</sub>/HCl solution.



**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 ( $\leq$  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](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