

Certificate of Analysis

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

Catalog No: 21886-1

Lot No: 7681

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.

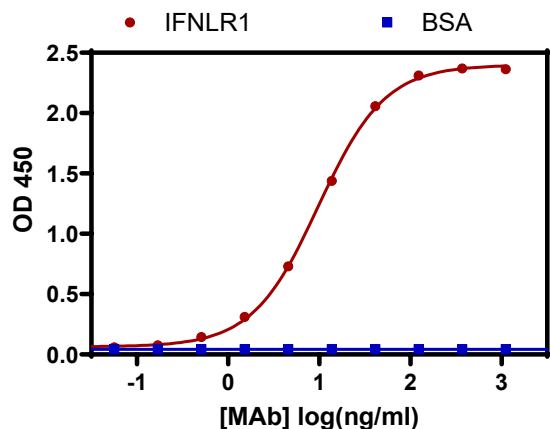


Figure 1. Representative binding curve of antibody to mammalian expressed human interferon lambda receptor (circle) and to 0.1% BSA/PBS (square) in a Direct Binding ELISA. Medium-binding polystyrene plates were coated with either 1 µg/ml human interferon lambda receptor 1 or 0.1%BSA/PBS. Thereafter, 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 H₂SO₄/HCl solution.

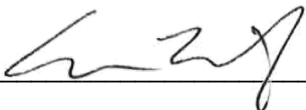


Shipping Conditions: Dry Ice

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.

Authorization

Released by: _____


Date: April 1, 2024

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