



## Certificate of Analysis

### Anti-Human Interferon Gamma Receptor Chain 2, Clone MMHGR-2 (MAb)

**Catalog No:** 21585-1

**Lot No:** 7432

**Size:** 250 µg/vial

**Description:** Mouse Monoclonal Antibody against Human Interferon Gamma Receptor Chain 2, non-neutralizing

**Clone:** MMHGR-2

**Volume:** 0.129 ml

**Concentration:** 1.94 mg/ml

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

**Endotoxin:** < 1 EU/µg

**Antigen:** Human Interferon Gamma Receptor Chain 2

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

**Purity:** > 95%

**Purification Method:** A combination of ion exchange, hydrophobic interaction and size exclusion chromatography

**Specificity:** Binds to but does not neutralize human interferon gamma receptor

**Tested Applications:** ELISA (direct binding)

*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. 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: February 8, 2022

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.