

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Mouse IFN- $\gamma$  Kit  
**Kit Lot Number:** K0082313  
**Expiration Date:** 30 Sep 2025  
**Kit Catalog Numbers:** K152QO-Series

**Precision and Accuracy Specifications:**

| Spot | Name          | Calibrator |       | Proinflammatory Panel 1 (mouse) Control 1 |       | Proinflammatory Panel 1 (mouse) Control 2 |       | Proinflammatory Panel 1 (mouse) Control 3 |       |
|------|---------------|------------|-------|---|-------|---|-------|---|-------|
|      |               | Recovery   | CV    | Recovery                                  | CV    | Recovery                                  | CV    | Recovery                                  | CV    |
| 1    | IFN- $\gamma$ | 80% - 120% | < 20% | 90% - 140%                                | < 25% | 90% - 140%                                | < 25% | 90% - 140%                                | < 25% |

| Spot | Name          | LLOQ  |            |       | ULOQ  |            |       |
|------|---------------|-------|------------|-------|-------|------------|-------|
|      |               | Value | Recovery   | CV    | Value | Recovery   | CV    |
| 1    | IFN- $\gamma$ | 0.390 | 75% - 125% | < 25% | 570   | 75% - 125% | < 25% |

**Stock Concentration of Calibrator:**


| Spot | Name          | Conc. | Units |
|------|---------------|-------|-------|
| 1    | IFN- $\gamma$ | 1030  | pg/mL |

**Lot-Specific Components:**

| Description                                      | Lot Number | Storage Temperature | Expiration Date |
|--|------------|---------------------|-----------------|
| Mouse IFN- $\gamma$ Plate                        | Z0047877   | 2-8°C               | 30 Nov 2025     |
| Anti-ms IFN- $\gamma$ Antibody                   | D0081787   | 2-8°C               | 30 Sep 2025     |
| Proinflammatory Panel 1 (mouse) Calibrator Blend | A0080351   | 2-8°C               | 31 Dec 2027     |
| Diluent 41                                       | M0410052   | $\leq$ -10°C        | 30 Sep 2025     |
| Diluent 45                                       | M0450052   | $\leq$ -10°C        | 30 Sep 2025     |

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

|                 | Name          | Function | Signature  | Date        |
|-----------------|---------------|----------|--|-------------|
| Review/Approval | Karen Hamilla | Quality  |  | 22 AUG 2023 |