



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX PLUS Human IFN- $\gamma$  Kit  
**Lot Number:** K0082454  
**Expiration Date:** 30 JUN 2026  
**Catalog Number:** K151QO-Series; K156QO-Series

**Lot-Specific Components:**

| Description                                      | Lot Number | Storage Temperature        | Expiration Date |
|--------------------------------------------------|------------|----------------------------|-----------------|
| Anti-hu IFN- $\gamma$ Antibody                   | D0082162   | 2-8°C                      | 30 APR 2027     |
| Human IFN- $\gamma$ Plate                        | Z0047968   | 2-8°C                      | 30 SEP 2026     |
| Proinflammatory Panel 1 (human) Calibrator Blend | A0080385   | 2-8°C                      | 31 DEC 2028     |
| Diluent 2                                        | M0020129   | $\leq -10^{\circ}\text{C}$ | 30 JUN 2026     |
| Diluent 3                                        | M0030127   | $\leq -10^{\circ}\text{C}$ | 30 SEP 2026     |

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

**Precision, Accuracy, and Sensitivity Specifications:**

|      |               | Calibrator |       | Proinflammatory Panel 1 (human) Control 1 |       | Proinflammatory Panel 1 (human) Control 2 |       | Proinflammatory Panel 1 (human) Control 3 |       |
|------|---------------|------------|-------|-------------------------------------------|-------|-------------------------------------------|-------|-------------------------------------------|-------|
| Spot | Name          | Recovery   | CV    | Recovery                                  | CV    | Recovery                                  | CV    | Recovery                                  | CV    |
| 1    | IFN- $\gamma$ | 80% - 120% | < 20% | 90% - 130%                                | < 20% | 90% - 130%                                | < 20% | 90% - 130%                                | < 20% |

|      |               | LLOQ  |            |       | ULOQ  |            |       |
|------|---------------|-------|------------|-------|-------|------------|-------|
| Spot | Name          | Value | Recovery   | CV    | Value | Recovery   | CV    |
| 1    | IFN- $\gamma$ | 1.76  | 80% - 120% | < 20% | 938   | 80% - 120% | < 20% |

**Stock Concentration of Calibrator:**

| Spot | Name          | Conc. | Units |
|------|---------------|-------|-------|
| 1    | IFN- $\gamma$ | 1380  | pg/mL |

Dilution to top of Curve: 1X

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

|                 | Name          | Function | Signature                                                                            | Date        |
|-----------------|---------------|----------|--------------------------------------------------------------------------------------|-------------|
| Review/Approval | Karen Hamilla | Quality  |  | 08 MAY 2024 |