



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX and V-PLEX Plus Human IFN- $\gamma$  Kit**Lot Number:** K0082972**Expiration Date:** 31 JUL 2028**Catalog Number:** K151QO-Series, K156QO-Series**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

| Description                                         | Lot Number | Storage Temperature        | Expiration Date |
|-----------------------------------------------------|------------|----------------------------|-----------------|
| Anti-hu IFN- $\gamma$ Antibody                      | D0082568   | 2-8°C                      | 30 JUN 2029     |
| Human IFN- $\gamma$ 96-Well Small Spot SECTOR Plate | Z0048277   | 2-8°C                      | 30 NOV 2028     |
| Proinflammatory Panel 1 (human) Calibrator Blend    | A0080461   | 2-8°C                      | 31 MAR 2031     |
| Diluent 2                                           | M0020138   | $\leq -10^{\circ}\text{C}$ | 30 SEP 2028     |
| Diluent 3                                           | M0030148   | $\leq -10^{\circ}\text{C}$ | 31 JUL 2028     |

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- $\gamma$ | 1600  | 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  |  | 07 AUG 2026 |