

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IFN- γ Kit

Kit Lot Number: K0082340

Expiration Date: 31 Aug 2025

Kit Catalog Numbers: K151QO-Series, K156QO-Series

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|----------------------------|-----------------|
| Human IFN- γ Plate | Z0047897 | 2-8°C | 31 Jan 2026 |
| Anti-hu IFN- γ Antibody | D0081973 | 2-8°C | 30 Jun 2026 |
| Proinflammatory Panel 1 (human) Calibrator Blend | A0080368 | 2-8°C | 30 Jun 2028 |
| Diluent 2 | M0020126 | $\leq -10^{\circ}\text{C}$ | 31 Aug 2025 |
| Diluent 3 | M0030123 | $\leq -10^{\circ}\text{C}$ | 31 Jan 2026 |

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

Precision and Accuracy Specifications:

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


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

Stock Concentration of Calibrator:

| Spot | Name | Conc. | Units |
|------|---------------|-------|-------|
| 1 | IFN- γ | 1450 | pg/mL |

Dilution to Top of Curve = 1X

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

| | Name | Function | Signature | Date |
|-----------------|---------------|----------|--|-------------|
| Review/Approval | Karen Hamilla | Quality |  | 27 Sep 2023 |