

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IFN- γ Kit
Kit Lot Number: K0081798
Expiration Date: 31 Oct 2023
Kit Catalog Numbers: K151QO-Series; K156QO-Series

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% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 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 | 1480 | pg/mL |

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|---------------------|-----------------|
| Human IFN- γ Plate | Z0047570 | 2-8°C | 31 Dec 2023 |
| Anti-hu IFN- γ Antibody | D0081686 | 2-8°C | 31 Oct 2024 |
| Proinflammatory Panel 1 (human) Calibrator Blend | A0080291 | 2-8°C | 30 Apr 2026 |
| Diluent 2 | M0020117 | \leq -10°C | 31 Oct 2023 |
| Diluent 3 | M0030108 | \leq -10°C | 30 Nov 2023 |

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 | Xiomara Talavera-Herrera | Quality |  | 07 OCT 2021 |