

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human IFN- $\gamma$  Kit  
**Kit Catalog Numbers:** K151QO-Series  
**Kit Lot Number:** K0081193  
**Expiration Date:** 31 Mar 2021

### Acceptance Data/ Percent Recovery 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    | hu IFN- $\gamma$ | 80% - 120% | < 20% | 80% - 120%                                | < 20% | 80% - 120%                                | < 20% | 80% - 120%                                | < 20% |

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

### Stock Concentration of Material Tested :

| Spot | Name             | Item                                      | Conc | Units |
|------|------------------|-------------------------------------------|------|-------|
| 1    | hu IFN- $\gamma$ | Calibrator                                | 1230 | pg/mL |
|      |                  | Proinflammatory Panel 1 (human) Control 1 | 544  | pg/mL |
|      |                  | Proinflammatory Panel 1 (human) Control 2 | 146  | pg/mL |
|      |                  | Proinflammatory Panel 1 (human) Control 3 | 37.8 | pg/mL |

Controls tested may not be the same lot as shipped.

### Standard Curve Data:

| hu IFN- $\gamma$ - Spot 1 |               |     |
|---------------------------|---------------|-----|
| Conc.<br>pg/mL            | ECL<br>Signal | %CV |
| 1230                      | 67,500        | 2.9 |
| 308                       | 16,700        | 2.4 |
| 76.9                      | 4,180         | 2.4 |
| 19.2                      | 1,070         | 2.3 |
| 4.80                      | 316           | 2.2 |
| 1.20                      | 126           | 3.4 |
| 0.300                     | 81            | 5.1 |
| 0.00                      | 65            | 8.4 |

# CERTIFICATE OF ANALYSIS

Page 2 of 2

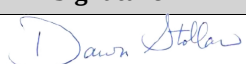
## Components:

| Description                                      | Lot Number       | Storage Temperature | Expiration Date |
|--------------------------------------------------|------------------|---------------------|-----------------|
| Human IFN- $\gamma$ Plate                        | Z0047031         | 2-8°C               | 31 Jul 2021     |
| Anti-hu IFN- $\gamma$ Antibody                   | D0081180         | 2-8°C               | 31 Mar 2022     |
| Proinflammatory Panel 1 (human) Calibrator Blend | A0080187         | 2-8°C               | 31 Dec 2023     |
| Diluent 2                                        | M0020103         | $\leq$ -10°C        | 31 Mar 2021     |
| Diluent 3                                        | M0030087         | $\leq$ -10°C        | 30 Jun 2021     |
| MSD Read Buffer T                                | Not Kit Specific | Not Kit Specific    | N/A             |

| Additional Components Included in V-PLEX Plus |                  |
|-----------------------------------------------|------------------|
| Plate Seals                                   | Not Kit Specific |
| Wash Buffer                                   | Not Kit Specific |
| Proinflammatory Panel 1 (human) Control 1     | A00C0640         |
| Proinflammatory Panel 1 (human) Control 2     | A00C0641         |
| Proinflammatory Panel 1 (human) Control 3     | A00C0642         |

## Statement:

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

|                 | Name         | Function          | Signature                                                                            | Date        |
|-----------------|--------------|-------------------|--------------------------------------------------------------------------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance |  | 25 MAR 2019 |