

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-Plex Plus NHP MIP-1 $\alpha$  Kit  
**Kit Catalog Numbers:** K156NQ-Series  
**Kit Lot Number:** K0080889  
**Expiration Date:** 31 Jul 2019

### Acceptance Data/ Percent Recovery Specifications:

| Spot | Name               | Calibrator |       | Chemokine Panel 1 (human) Control 1 |       | Chemokine Panel 1 (human) Control 2 |       | Chemokine Panel 1 (human) Control 3 |       |
|------|--------------------|------------|-------|-------------------------------------|-------|-------------------------------------|-------|-------------------------------------|-------|
|      |                    | Recovery   | CV    | Recovery                            | CV    | Recovery                            | CV    | Recovery                            | CV    |
| 1    | NHP MIP-1 $\alpha$ | 80% - 120% | < 20% | 80% - 120%                          | < 20% | 80% - 120%                          | < 20% | 80% - 120%                          | < 20% |

| Spot | Name               | LLOQ  |            |       | ULOQ  |            |       |
|------|--------------------|-------|------------|-------|-------|------------|-------|
|      |                    | Value | Recovery   | CV    | Value | Recovery   | CV    |
| 1    | NHP MIP-1 $\alpha$ | 13.8  | 80% - 120% | < 20% | 743   | 80% - 120% | < 20% |

### Stock Concentration of Material Tested :

| Spot | Name               | Item                                | Conc | Units |
|------|--------------------|-------------------------------------|------|-------|
| 1    | NHP MIP-1 $\alpha$ | Calibrator                          | 1080 | pg/mL |
|      |                    | Chemokine Panel 1 (human) Control 1 | 824  | pg/mL |
|      |                    | Chemokine Panel 1 (human) Control 2 | 371  | pg/mL |
|      |                    | Chemokine Panel 1 (human) Control 3 | 152  | pg/mL |

Controls tested may not be the same lot as shipped.

The calibrator and controls for this assay contain a human recombinant cytokine calibrator sequence that is highly homologous to the non-human primate cytokine. This assay has been verified to cross-react with homologs of rhesus and cynomolgus monkeys. See Product Insert for experimental data.

### Standard Curve Data:

| NHP MIP-1 $\alpha$ - Spot 1 |               |     |
|-----------------------------|---------------|-----|
| Conc.<br>pg/mL              | ECL<br>Signal | %CV |
| 1080                        | 238,000       | 2.4 |
| 270                         | 26,800        | 4.4 |
| 67.5                        | 2,580         | 4.9 |
| 16.9                        | 273           | 3.3 |
| 4.22                        | 92            | 3.4 |
| 1.05                        | 73            | 3.4 |
| 0.264                       | 70            | 5.5 |
| 0.00                        | 69            | 7.6 |

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

| Description                                | Lot Number       | Storage Temperature | Expiration Date |
|--|------------------|---------------------|-----------------|
| Human MIP-1 $\alpha$ Plate                 | Z0046645         | 2-8°C               | 28 Feb 2020     |
| Anti-hu MIP-1 $\alpha$ Antibody            | D0080902         | 2-8°C               | 31 Jul 2019     |
| Chemokine Panel 1 (human) Calibrator Blend | A0080122         | 2-8°C               | 28 Feb 2022     |
| Diluent 43                                 | M0430021         | $\leq$ -10°C        | 30 Sep 2019     |
| Diluent 3                                  | M0030077         | $\leq$ -10°C        | 30 Nov 2019     |
| 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 (20X), 100mL                      | Not Kit Specific |
| Chemokine Panel 1 (human) Control 1           | Not Kit Specific |
| Chemokine Panel 1 (human) Control 2           | Not Kit Specific |
| Chemokine Panel 1 (human) Control 3           | Not Kit Specific |

### Statement:

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

|                        | Name         | Function          | Signature  | Date        |
|------------------------|--------------|-------------------|--|-------------|
| <b>Review/Approval</b> | Dawn Stollar | Quality Assurance |  | 18 OCT 2017 |