

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus NHP IL-1 $\beta$  Kit  
**Kit Catalog Numbers:** K156QP-Series  
**Kit Lot Number:** K0080942  
**Expiration Date:** 31 Jul 2019

### 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    | NHP IL-1 $\beta$ | 80% - 120% | < 20% | 80% - 120%                                | < 20% | 80% - 120%                                | < 20% | 80% - 120%                                | < 20% |

| Spot | Name             | LLOQ  |            |       | ULOQ  |            |       |
|------|------------------|-------|------------|-------|-------|------------|-------|
|      |                  | Value | Recovery   | CV    | Value | Recovery   | CV    |
| 1    | NHP IL-1 $\beta$ | 2.14  | 80% - 120% | < 20% | 375   | 80% - 120% | < 20% |

### Stock Concentration of Material Tested :

| Spot | Name             | Item                                      | Conc | Units |
|------|------------------|-------------------------------------------|------|-------|
| 1    | NHP IL-1 $\beta$ | Calibrator                                | 566  | pg/mL |
|      |                  | Proinflammatory Panel 1 (human) Control 1 | 215  | pg/mL |
|      |                  | Proinflammatory Panel 1 (human) Control 2 | 52   | pg/mL |
|      |                  | Proinflammatory Panel 1 (human) Control 3 | 11   | 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 IL-1 $\beta$ - Spot 1 |            |      |
|---------------------------|------------|------|
| Conc. pg/mL               | ECL Signal | %CV  |
| 566                       | 1,130,000  | 1.9  |
| 142                       | 304,000    | 2.0  |
| 35.4                      | 78,200     | 1.9  |
| 8.84                      | 20,000     | 2.0  |
| 2.21                      | 5,250      | 2.0  |
| 0.553                     | 1,480      | 1.6  |
| 0.138                     | 547        | 2.2  |
| 0.00                      | 238        | 13.0 |

**CERTIFICATE OF ANALYSIS**

**Components:**

| Description                                      | Lot Number       | Storage Temperature | Expiration Date |
|--------------------------------------------------|------------------|---------------------|-----------------|
| Human IL-1 $\beta$ Plate                         | Z0046726         | 2-8°C               | 30 Jun 2020     |
| Anti-hu IL-1 $\beta$ Antibody                    | D0080915         | 2-8°C               | 31 Jul 2019     |
| Proinflammatory Panel 1 (human) Calibrator Blend | A0080147         | 2-8°C               | 31 Aug 2022     |
| Diluent 2                                        | M0020096         | $\leq$ -10°C        | 30 Nov 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), 100 mL                 | Not Kit Specific |
| Proinflammatory Panel 1 (human) Control 1 | Not Kit Specific |
| Proinflammatory Panel 1 (human) Control 2 | Not Kit Specific |
| Proinflammatory 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 | <i>Dawn Stollar</i> | 06 APR 2018 |