

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus NHP IL-7 Kit

**Kit Catalog Numbers:** K156RC-Series

**Kit Lot Number:** K0081124

**Expiration Date:** 30 Apr 2020

### Acceptance Data/ Percent Recovery Specifications:

| Spot | Name     | Calibrator |       | Cytokine Panel 1 (human) Control 1 |       | Cytokine Panel 1 (human) Control 2 |       | Cytokine Panel 1 (human) Control 3 |       |
|------|----------|------------|-------|------------------------------------|-------|------------------------------------|-------|------------------------------------|-------|
|      |          | Recovery   | CV    | Recovery                           | CV    | Recovery                           | CV    | Recovery                           | CV    |
| 1    | NHP IL-7 | 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-7 | 0.546 | 80% - 120% | < 20% | 563   | 80% - 120% | < 20% |

### Stock Concentration of Material Tested :

| Spot | Name     | Item                               | Conc | Units |
|------|----------|------------------------------------|------|-------|
| 1    | NHP IL-7 | Calibrator                         | 844  | pg/mL |
|      |          | Cytokine Panel 1 (human) Control 1 | 346  | pg/mL |
|      |          | Cytokine Panel 1 (human) Control 2 | 58   | pg/mL |
|      |          | Cytokine Panel 1 (human) Control 3 | 7.79 | 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-7 - Spot 1 |            |     |
|-------------------|------------|-----|
| Conc. pg/mL       | ECL Signal | %CV |
| 844               | 230,000    | 1.2 |
| 211               | 57,900     | 2.1 |
| 52.8              | 13,700     | 1.5 |
| 13.2              | 3,620      | 2.1 |
| 3.30              | 901        | 2.0 |
| 0.824             | 285        | 2.3 |
| 0.206             | 116        | 8.4 |
| 0.00              | 71         | 7.1 |

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

| Description                               | Lot Number       | Storage Temperature | Expiration Date |
|---|------------------|---------------------|-----------------|
| Human IL-7 Plate                          | Z0046945         | 2-8°C               | 31 Mar 2021     |
| Anti-hu IL-7 Antibody                     | D0081055         | 2-8°C               | 30 Apr 2020     |
| Cytokine Panel 1 (human) Calibrator Blend | A0080121         | 2-8°C               | 28 Feb 2022     |
| Diluent 43                                | M0430028         | ≤ -10°C             | 31 Jan 2021     |
| Diluent 3                                 | M0030085         | ≤ -10°C             | 28 Feb 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 (20X), 100mL                      | Not Kit Specific |
| Cytokine Panel 1 (human) Control 1            | Not Kit Specific |
| Cytokine Panel 1 (human) Control 2            | Not Kit Specific |
| Cytokine 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 |  | 28 NOV 2018 |