

# CERTIFICATE OF ANALYSIS

Page 1 of 2

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

**Kit Catalog Numbers:** K151RC-Series

**Kit Lot Number:** K0081019

**Expiration Date:** 31 Jul 2019

## 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    | hu 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    | hu IL-7 | 0.546 | 80% - 120% | < 20% | 563   | 80% - 120% | < 20% |

## Stock Concentration of Material Tested :

| Spot | Name    | Item                               | Conc | Units |
|------|---------|------------------------------------|------|-------|
| 1    | hu IL-7 | Calibrator                         | 881  | pg/mL |
|      |         | Cytokine Panel 1 (human) Control 1 | 348  | pg/mL |
|      |         | Cytokine Panel 1 (human) Control 2 | 56.4 | pg/mL |
|      |         | Cytokine Panel 1 (human) Control 3 | 8.49 | pg/mL |

Controls tested may not be the same lot as shipped.

## Standard Curve Data:

| hu IL-7 - Spot 1 |               |     |
|------------------|---------------|-----|
| Conc.<br>pg/mL   | ECL<br>Signal | %CV |
| 881              | 219,000       | 4.0 |
| 220              | 53,700        | 2.2 |
| 55.1             | 12,800        | 2.2 |
| 13.8             | 3,230         | 1.1 |
| 3.44             | 892           | 2.0 |
| 0.860            | 280           | 1.4 |
| 0.215            | 127           | 3.8 |
| 0.00             | 80            | 4.2 |

## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Components:

| Description                               | Lot Number       | Storage Temperature | Expiration Date |
|-------------------------------------------|------------------|---------------------|-----------------|
| Human IL-7 Plate                          | Z0046841         | 2-8°C               | 30 Nov 2020     |
| Anti-hu IL-7 Antibody                     | D0080888         | 2-8°C               | 31 Jul 2019     |
| Cytokine Panel 1 (human) Calibrator Blend | A0080144         | 2-8°C               | 31 Aug 2022     |
| Diluent 43                                | M0430026         | ≤ -10°C             | 31 Aug 2020     |
| Diluent 3                                 | M0030082         | ≤ -10°C             | 31 Oct 2020     |
| MSD Read Buffer T                         | Not Kit Specific | Not Kit Specific    | N/A             |

| Additional Components Included in V-PLEX Plus |                  |
|-----------------------------------------------|------------------|
| Wash Buffer (20X), 100 mL                     | Not Kit Specific |
| Plate Seals                                   | 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        |
|-----------------|--------------|-------------------|--------------------------------------------------------------------------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance |  | 14 AUG 2018 |