

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-17A Kit

Kit Catalog Numbers: K151RF-Series; K156RF-Series

Kit Lot Number: K0081434

Expiration Date: 30 Jun 2022

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 | IL-17A | 80% - 120% | < 20% | 90% - 130% | < 20% | 90% - 130% | < 20% | 90% - 130% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|--------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | IL-17A | 3.19 | 80% - 120% | < 20% | 3650 | 80% - 120% | < 20% |

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|--------|------------------------------------|------|-------|
| 1 | IL-17A | Calibrator | 5770 | pg/mL |
| | | Cytokine Panel 1 (human) Control 1 | 2220 | pg/mL |
| | | Cytokine Panel 1 (human) Control 2 | 305 | pg/mL |
| | | Cytokine Panel 1 (human) Control 3 | 46.4 | pg/mL |

Controls tested may not be the same lot as shipped.

Standard Curve Data:

| IL-17A - Spot 1 | | |
|-----------------|------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 5770 | 691,000 | 3.8 |
| 1440 | 190,000 | 3.7 |
| 361 | 48,300 | 2.8 |
| 90.2 | 12,600 | 3.4 |
| 22.5 | 3,220 | 2.1 |
| 5.63 | 876 | 4.6 |
| 1.41 | 259 | 4.4 |
| 0.00 | 58 | 7.4 |

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|---|------------------|---------------------|-----------------|
| Human IL-17A Plate | Z0047274 | 2-8°C | 31 Jul 2022 |
| Anti-hu IL-17A Antibody | D0081208 | 2-8°C | 31 Mar 2023 |
| Cytokine Panel 1 (human) Calibrator Blend | A0080227 | 2-8°C | 31 Oct 2024 |
| Diluent 43 | M0430037 | ≤ -10°C | 31 Aug 2022 |
| Diluent 3 | M0030096 | ≤ -10°C | 30 Jun 2022 |
| 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 |
|------------------------|---------------|----------|--|-------------|
| Review/Approval | Karen Hamilla | Quality |  | 08 APR 2020 |