



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX and V-PLEX Plus Human IL-17A Kit  
**Lot Number:** K0082991  
**Expiration Date:** 29 FEB 2028  
**Catalog Number:** K151RF-Series, K156RF-Series

**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

| Description                                  | Lot Number | Storage Temperature | Expiration Date |
|----------------------------------------------|------------|---------------------|-----------------|
| Anti-hu IL-17A Antibody                      | D0082350   | 2-8°C               | 29 FEB 2028     |
| Cytokine Panel 1 (human) Calibrator Blend    | A0080454   | 2-8°C               | 31 DEC 2030     |
| Human IL-17A 96-Well Small Spot SECTOR Plate | Z0048226   | 2-8°C               | 31 JUL 2028     |
| Diluent 3                                    | M0030148   | ≤ -10°C             | 31 JUL 2028     |
| Diluent 43                                   | M0430060   | ≤ -10°C             | 30 JUN 2028     |

Additional components may be provided to execute the assay. Additional components are not lot-specific.

**Precision, Accuracy, and Sensitivity Specifications:**

|      |        | Calibrator |       | Cytokine Panel 1 (human)<br>Control 1 |       | Cytokine Panel 1 (human)<br>Control 2 |       | Cytokine Panel 1 (human)<br>Control 3 |       |
|------|--------|------------|-------|---------------------------------------|-------|---------------------------------------|-------|---------------------------------------|-------|
| Spot | Name   | Recovery   | CV    | Recovery                              | CV    | Recovery                              | CV    | Recovery                              | CV    |
| 1    | IL-17A | 80% - 120% | < 20% | 90% - 130%                            | < 20% | 90% - 130%                            | < 20% | 90% - 130%                            | < 20% |

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

**Stock Concentration of Calibrator:**

| Spot | Name   | Conc. | Units |
|------|--------|-------|-------|
| 1    | IL-17A | 6310  | pg/mL |

Dilution to top of Curve: 1X

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

|                 | Name          | Function | Signature                                                                            | Date        |
|-----------------|---------------|----------|--------------------------------------------------------------------------------------|-------------|
| Review/Approval | Karen Hamilla | Quality  |  | 12 AUG 2026 |