

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX PLUS Human IL-27 Kit

Kit Lot Number: K0082315

Expiration Date: 31 Mar 2025

Kit Catalog Numbers: K151WX-Series

Precision and Accuracy Specifications:

| Spot | Name | Calibrator | | TH17 Panel 1 (human) Control 1 | | TH17 Panel 1 (human) Control 2 | | TH17 Panel 1 (human) Control 3 | |
|------|-------|------------|-------|--------------------------------|-------|--------------------------------|-------|--------------------------------|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | IL-27 | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|-------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | IL-27 | 38.7 | 80% - 120% | < 20% | 13000 | 80% - 120% | < 20% |

Stock Concentration of Calibrator:


| Spot | Name | Conc. | Units |
|------|-------|-------|-------|
| 1 | IL-27 | 17700 | pg/mL |

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|---------------------------------------|------------|---------------------|-----------------|
| Human IL-27 Plate | Z0047878 | 2-8°C | 30 Nov 2025 |
| Anti-hu IL-27 Antibody | D0081712 | 2-8°C | 31 Mar 2025 |
| TH17 Panel 1 (human) Calibrator Blend | A0080324 | 2-8°C | 31 Mar 2027 |
| Diluent 3 | M0030120 | ≤ -10°C | 31 Jul 2025 |
| Diluent 43 | M0430050 | ≤ -10°C | 31 Mar 2025 |

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

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

| | Name | Function | Signature | Date |
|-----------------|---------------|----------|--|-------------|
| Review/Approval | Karen Hamilla | Quality |  | 21 JUL 2023 |