

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-27 Kit
Kit Catalog Numbers: K151WX-Series
Kit Lot Number: K0080595
Expiration Date: 31 Jan 2019

Acceptance Data/ Percent Recovery Specifications:

| Spot | Name | Calibrator | | TH17 Panel 1 (human) Control 1 | | TH17 Panel 1 (human) Control 3 | | TH17 Panel 1 (human) Control 2 | |
|------|-------|------------|-------|--------------------------------|-------|--------------------------------|-------|--------------------------------|-------|
| | | 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 Material Tested :

| Spot | Name | Item | Conc | Units |
|------|-------|--------------------------------|-------|-------|
| 1 | IL-27 | Calibrator | 17100 | pg/mL |
| | | TH17 Panel 1 (human) Control 1 | 8470 | pg/mL |
| | | TH17 Panel 1 (human) Control 2 | 1960 | pg/mL |
| | | TH17 Panel 1 (human) Control 3 | 477 | pg/mL |

Controls tested may not be the same lot as shipped.

Standard Curve Data:

| IL-27 - Spot 1 | | |
|----------------|------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 17100 | 122,000 | 2.9 |
| 4280 | 26,900 | 2.5 |
| 1070 | 6,530 | 2.4 |
| 267 | 1,630 | 1.9 |
| 66.8 | 503 | 1.9 |
| 16.7 | 193 | 2.9 |
| 4.17 | 121 | 4.3 |
| 0.00 | 92 | 6.1 |



CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|---------------------------------------|------------------|---------------------|-----------------|
| Human IL-27 Plate | Z0046180 | 2-8°C | 28 Feb 2019 |
| Human IL-27 Antibody | D0080760 | 2-8°C | 31 Jan 2019 |
| TH17 Panel 1 (human) Calibrator Blend | A0080101 | 2-8°C | 30 Sep 2021 |
| Diluent 43 | M0430018 | ≤ -10°C | 31 Jan 2019 |
| Diluent 3 | M0030073 | ≤ -10°C | 31 Mar 2019 |
| 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), 100 mL | Not Kit Specific |
| TH17 Panel 1 (human) Control 1 | Not Kit Specific |
| TH17 Panel 1 (human) Control 2 | Not Kit Specific |
| TH17 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 | | 15 FEB 2017 |