

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-17A/F Kit
Kit Catalog Numbers: K152WN-Series
Kit Lot Number: K0081234
Expiration Date: 30 Jun 2021

Acceptance Data/ Percent Recovery Specifications:

| Spot | Name | Calibrator | | Cytokine Panel 1 (mouse) Control 1 | | Cytokine Panel 1 (mouse) Control 2 | | Cytokine Panel 1 (mouse) Control 3 | |
|------|----------|------------|-------|------------------------------------|-------|------------------------------------|-------|------------------------------------|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | IL-17A/F | 80% - 120% | < 20% | 75% - 125% | < 20% | 75% - 125% | < 20% | 75% - 125% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|----------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | IL-17A/F | 1.39 | 75% - 125% | < 20% | 1620 | 75% - 125% | < 20% |

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|----------|------------------------------------|------|-------|
| 1 | IL-17A/F | Calibrator | 2370 | pg/mL |
| | | Cytokine Panel 1 (mouse) Control 1 | 2220 | pg/mL |
| | | Cytokine Panel 1 (mouse) Control 2 | 212 | pg/mL |
| | | Cytokine Panel 1 (mouse) Control 3 | 18.6 | pg/mL |

Controls tested may not be the same lot as shipped.

Standard Curve Data:

| IL-17A/F - Spot 1 | | |
|-------------------|---------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 2370 | 401,000 | 3.8 |
| 593 | 90,800 | 4.5 |
| 148 | 22,700 | 3.0 |
| 37.0 | 5,450 | 4.1 |
| 9.26 | 1,470 | 4.1 |
| 2.31 | 423 | 5.1 |
| 0.579 | 167 | 5.4 |
| 0.00 | 87 | 6.5 |

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|---|------------------|---------------------|-----------------|
| Mouse IL-17A/F Plate | Z0047074 | 2-8°C | 30 Sep 2021 |
| Anti-ms IL-17A/F Antibody | D0081115 | 2-8°C | 31 Dec 2021 |
| Cytokine Panel 1 (mouse) Calibrator blend | A0080156 | 2-8°C | 31 Mar 2023 |
| Diluent 41 | M0410027 | ≤ -10°C | 31 Aug 2021 |
| Diluent 45 | M0450026 | ≤ -10°C | 30 Jun 2021 |
| 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 (mouse) Control 1 | Not Kit Specific |
| Cytokine Panel 1 (mouse) Control 2 | Not Kit Specific |
| Cytokine Panel 1 (mouse) 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 |  | 13 JUN 2019 |