

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus NHP Eotaxin-3 Kit
Kit Catalog Numbers: K156NU-series
Kit Lot Number: K0080728
Expiration Date: 30 Jun 2019

Acceptance Data/ Percent Recovery Specifications:

| Spot | Name | Calibrator | | Chemokine Panel 1 (human) Control 1 | | Chemokine Panel 1 (human) Control 2 | | Chemokine Panel 1 (human) Control 3 | |
|------|---------------|------------|-------|-------------------------------------|-------|-------------------------------------|-------|-------------------------------------|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | NHP Eotaxin-3 | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|---------------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | NHP Eotaxin-3 | 10.2 | 80% - 120% | < 20% | 3750 | 80% - 120% | < 20% |

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|---------------|-------------------------------------|------|-------|
| 1 | NHP Eotaxin-3 | Calibrator | 7120 | pg/mL |
| | | Chemokine Panel 1 (human) Control 1 | 4670 | pg/mL |
| | | Chemokine Panel 1 (human) Control 2 | 794 | pg/mL |
| | | Chemokine Panel 1 (human) Control 3 | 127 | pg/mL |

Controls tested may not be the same lot as shipped.

The calibrator and controls for this assay contain a human recombinant cytokine calibrator sequence that is highly homologous to the non-human primate cytokine. This assay has been verified to cross-react with homologs of rhesus and cynomolgus monkeys. See Product Insert for experimental data.

CERTIFICATE OF ANALYSIS

Standard Curve Data:

| NHP Eotaxin-3 - Spot 1 | | |
|------------------------|------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 7120 | 69,000 | 3.1 |
| 1780 | 21,500 | 3.0 |
| 445 | 5,850 | 3.1 |
| 111 | 1,530 | 1.8 |
| 27.8 | 429 | 1.3 |
| 6.95 | 149 | 3.3 |
| 1.74 | 78 | 1.5 |
| 0.00 | 57 | 5.4 |


Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------------|---------------------|-----------------|
| Human Eotaxin-3 Plate | Z0046404 | 2-8°C | 31 Jul 2019 |
| Anti-hu Eotaxin-3 Antibody | D0080862 | 2-8°C | 30 Jun 2019 |
| Chemokine Panel 1 (human) Calibrator Blend | A0080122 | 2-8°C | 28 Feb 2022 |
| Diluent 43 | M0430020 | ≤ -10°C | 30 Jun 2019 |
| Diluent 3 | M0030074 | ≤ -10°C | 30 Jun 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), 100mL | Not Kit Specific |
| Chemokine Panel 1 (human) Control 1 | Not Kit Specific |
| Chemokine Panel 1 (human) Control 2 | Not Kit Specific |
| Chemokine 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 |  | 27 MAR 2017 |