

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human Eotaxin-3 Kit
Kit Lot Number: K0081553
Expiration Date: 31 Aug 2022
Kit Catalog Numbers: K151NU-Series; K156NU-Series

Precision and Accuracy 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 | Eotaxin-3 | 80% - 120% | < 20% | 75% - 125% | < 25% | 75% - 125% | < 25% | 75% - 125% | < 25% |

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

Stock Concentration of Calibrator:


| Spot | Name | Conc. | Units |
|------|-----------|-------|-------|
| 1 | Eotaxin-3 | 5405 | pg/ml |

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|---|------------|---------------------|-----------------|
| Human Eotaxin-3 plate | Z0047396 | 2-8°C | 31 Mar 2023 |
| Anti-hu Eotaxin-3 Antibody | D0081260 | 2-8°C | 31 Oct 2022 |
| Chemokine Panel 1(human) Calibrator Blend | A0080266 | 2-8°C | 31 Aug 2025 |
| Diluent 43 | M0430039 | ≤ -10°C | 30 Sep 2022 |
| Diluent 3 | M0030097 | ≤ -10°C | 31 Aug 2022 |

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 | Xiomara Talavera | Quality |  | 16 NOV 2020 |