

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-8 Kit
Kit Lot Number: K0081599
Expiration Date: 31 Jul 2022
Kit Catalog Numbers: K151RA-Series, K156RA-Series

Precision and Accuracy Specifications:

| | | Calibrator | | Proinflammatory Panel 1(human) Control 1 | | Proinflammatory Panel 1(human) Control 2 | | Proinflammatory Panel 1(human) Control 3 | |
|------|------|------------|-------|--|-------|--|-------|--|-------|
| Spot | Name | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | IL-8 | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |

| | | LLOQ | | | ULOQ | | |
|------|------|-------|------------|-------|-------|------------|-------|
| Spot | Name | Value | Recovery | CV | Value | Recovery | CV |
| 1 | IL-8 | 0.591 | 80% - 120% | < 20% | 375 | 80% - 120% | < 20% |

Stock Concentration of Calibrator:


| Spot | Name | Conc. | Units |
|------|------|-------|-------|
| 1 | IL-8 | 635 | pg/mL |

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|---|------------|---------------------|-----------------|
| Human IL-8 Plate | Z0047425 | 2-8°C | 30 Apr 2023 |
| Anti-hu IL-8 Antibody | D0081255 | 2-8°C | 31 Jul 2022 |
| Proinflammatory Panel 1(human) Calibrator Blend | A0080239 | 2-8°C | 31 Oct 2024 |
| Diluent 2 | M0020114 | ≤ -10°C | 31 Mar 2023 |
| Diluent 3 | M0030104 | ≤ -10°C | 30 Jun 2023 |

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 |  | 12 MAR 2021 |