

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-Plex Plus Human IL-8 Kit
Kit Lot Number: K0082228
Expiration Date: 30 Sep 2024
Kit Catalog Numbers: K151RA-Series; K156RA-Series

Precision and Accuracy Specifications:

| Spot | Name | Calibrator | | Proinflammatory Panel 1 (human) Control 1 | | Proinflammatory Panel 1 (human) Control 2 | | Proinflammatory Panel 1 (human) Control 3 | |
|------|------|------------|-------|---|-------|---|-------|---|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | IL-8 | 80% - 120% | < 20% | 90% - 130% | < 20% | 90% - 130% | < 20% | 90% - 130% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|------|-------|------------|-------|-------|------------|-------|
| | | 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 | 592 | pg/ml |

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|---------------------|-----------------|
| Human IL-8 Plate | Z0047820 | 2-8°C | 31 Jul 2025 |
| Anti-hu IL-8 Antibody | D0081885 | 2-8°C | 31 Jan 2026 |
| Proinflammatory Panel 1 (human) Calibrator Blend | A0080229 | 2-8°C | 30 Sep 2024 |
| Diluent 2 | M0020122 | ≤ -10°C | 31 Oct 2024 |
| Diluent 3 | M0030116 | ≤ -10°C | 30 Nov 2024 |

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 |  | 27 FEB 2023 |