

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-1RA Kit
Kit Lot Number: K0081617
Expiration Date: 30 Nov 2021
Kit Catalog Numbers: K151WT-Series

Precision and Accuracy Specifications:

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

| Spot | Name | LLOQ | | | ULOQ | | |
|------|--------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | IL-1RA | 9.19 | 75% - 125% | < 25% | 650 | 80% - 120% | < 20% |

Stock Concentration of Calibrator:


| Spot | Name | Conc. | Units |
|------|--------|-------|-------|
| 1 | IL-1RA | 905 | pg/mL |

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|---|------------|---------------------|-----------------|
| Human IL-1RA Plate | Z0047106 | 2-8°C | 30 Nov 2021 |
| Anti-hu IL-1RA Antibody | D0081093 | 2-8°C | 31 Dec 2021 |
| Cytokine Panel 2 (human) Calibrator Blend | A0080193 | 2-8°C | 29 Feb 2024 |
| Diluent 43 | M0430042 | ≤ -10°C | 30 Apr 2023 |
| Diluent 3 | M0030101 | ≤ -10°C | 31 Dec 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 | Karen Hamilla | Quality |  | 04 JAN 2021 |