



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-1 $\beta$  Kit  
**Lot Number:** K0082956  
**Expiration Date:** 29 FEB 2028  
**Catalog Number:** K151QP-Series, K156QP-Series

**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

| Description                                        | Lot Number | Storage Temperature        | Expiration Date |
|----------------------------------------------------|------------|----------------------------|-----------------|
| Anti-hu IL-1 $\beta$ Antibody                      | D0082366   | 2-8°C                      | 31 MAR 2028     |
| Human IL-1 $\beta$ 96-Well Small Spot SECTOR Plate | Z0048264   | 2-8°C                      | 30 NOV 2028     |
| Proinflammatory Panel 1 (human) Calibrator Blend   | A0080461   | 2-8°C                      | 31 MAR 2031     |
| Diluent 2                                          | M0020137   | $\leq -10^{\circ}\text{C}$ | 29 FEB 2028     |
| Diluent 3                                          | M0030147   | $\leq -10^{\circ}\text{C}$ | 31 MAY 2028     |

Additional components may be provided to execute the assay. Additional components are not lot-specific.

**Precision, Accuracy, and Sensitivity 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-1 $\beta$ | 80% - 120% | < 20% | 80% - 120%                                | < 20% | 80% - 120%                                | < 20% | 80% - 120%                                | < 20% |

|      |              | LLOQ  |            |       | ULOQ  |            |       |
|------|--------------|-------|------------|-------|-------|------------|-------|
| Spot | Name         | Value | Recovery   | CV    | Value | Recovery   | CV    |
| 1    | IL-1 $\beta$ | 2.14  | 80% - 120% | < 20% | 375   | 80% - 120% | < 20% |

**Stock Concentration of Calibrator:**

| Spot | Name         | Conc. | Units |
|------|--------------|-------|-------|
| 1    | IL-1 $\beta$ | 600   | pg/mL |

Dilution to top of Curve: 1X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

|                 | Name          | Function | Signature                                                                            | Date        |
|-----------------|---------------|----------|--------------------------------------------------------------------------------------|-------------|
| Review/Approval | Karen Hamilla | Quality  |  | 16 JUL 2026 |