



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX Plus Human MIP-1 $\beta$  Kit**Lot Number:** K0082759**Expiration Date:** 28 FEB 2027**Catalog Number:** K151NR-Series, K156NR-Series**Lot-Specific Components:**

| Description                                         | Lot Number | Storage Temperature        | Expiration Date |
|-----------------------------------------------------|------------|----------------------------|-----------------|
| Anti-hu MIP-1 $\beta$ Antibody                      | D0082473   | 2-8°C                      | 30 SEP 2028     |
| Chemokine Panel 1 (human) Calibrator Blend          | A0080418   | 2-8°C                      | 28 FEB 2030     |
| Human MIP-1 $\beta$ 96-Well Small Spot SECTOR Plate | Z0048153   | 2-8°C                      | 31 JAN 2028     |
| Diluent 3                                           | M0030140   | $\leq -10^{\circ}\text{C}$ | 30 SEP 2027     |
| Diluent 43                                          | M0430058   | $\leq -10^{\circ}\text{C}$ | 28 FEB 2027     |

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

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

| Spot | Name          | LLOQ  |             |       | ULOQ  |            |       |
|------|---------------|-------|-------------|-------|-------|------------|-------|
|      |               | Value | Recovery    | CV    | Value | Recovery   | CV    |
| 1    | MIP-1 $\beta$ | 1.88  | 80% - 120%* | < 20% | 520   | 80% - 120% | < 20% |

\*See Additional Comments

**Stock Concentration of Calibrator:**

| Spot | Name          | Conc. | Units |
|------|---------------|-------|-------|
| 1    | MIP-1 $\beta$ | 954   | pg/mL |

Dilution to top of Curve: 1X

**Additional Comments:**

| Spot | Name          | Specification | Comment                                                                                              |
|------|---------------|---------------|------------------------------------------------------------------------------------------------------|
| 1    | MIP-1 $\beta$ | LLOQ Conc.    | % LLOQ measured 139% (specification 80-120%) at target 1.88pg/mL, was fully conforming at 7.52pg/mL. |

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

|                 | Name          | Function | Signature                                                                            | Date        |
|-----------------|---------------|----------|--------------------------------------------------------------------------------------|-------------|
| Review/Approval | Karen Hamilla | Quality  |  | 04 NOV 2025 |