



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX and V-PLEX Plus Human IL-12p70 Kit**Lot Number:** K0082763**Expiration Date:** 31 JAN 2027**Catalog Number:** K151QV-Series**Lot-Specific Components:**

| Description                                      | Lot Number | Storage Temperature | Expiration Date |
|--------------------------------------------------|------------|---------------------|-----------------|
| Anti-hu IL-12p70 Antibody                        | D0082119   | 2-8°C               | 31 JAN 2027     |
| Human IL-12 p70, 96-Well Small Spot SECTOR Plate | Z0048154   | 2-8°C               | 31 DEC 2027     |
| Proinflammatory Panel 1 (human) Calibrator Blend | A0080414   | 2-8°C               | 31 JAN 2030     |
| Diluent 2                                        | M0020134   | ≤ -10°C             | 30 SEP 2027     |
| Diluent 3                                        | M0030140   | ≤ -10°C             | 30 SEP 2027     |

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-12p70 | 80% - 120% | < 20% | 80% - 120%                                | < 20% | 80% - 120%                                | < 20% | 80% - 120%                                | < 20% |

|      |          | LLOQ  |            |       | ULOQ  |            |       |
|------|----------|-------|------------|-------|-------|------------|-------|
| Spot | Name     | Value | Recovery   | CV    | Value | Recovery   | CV    |
| 1    | IL-12p70 | 1.22  | 80% - 120% | < 20% | 315   | 80% - 120% | < 20% |

**Stock Concentration of Calibrator:**

| Spot | Name     | Conc. | Units |
|------|----------|-------|-------|
| 1    | IL-12p70 | 497   | 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  |  | 14 AUG 2025 |