

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX Proinflammatory Panel 1 (human) Kit  
**Lot Number:** K00S0267  
**Expiration Date:** 31 MAY 2027  
**Catalog Number:** K15396S-Series; K151AKV-Series

**Lot-Specific Components:**

| Description                                             | Lot Number | Storage Temperature        | Expiration Date |
|---------------------------------------------------------|------------|----------------------------|-----------------|
| S-PLEX Proinflammatory Panel 1 (human) Calibrator Blend | A00S0127   | 2-8°C                      | 30 JUN 2030     |
| TURBO-BOOST Human IFN- $\gamma$ Antibody                | D0TB0027   | 2-8°C                      | 31 OCT 2028     |
| TURBO-BOOST Human IL-10 Antibody                        | D0TB0021   | 2-8°C                      | 31 OCT 2028     |
| TURBO-BOOST Human IL-12p70 Antibody                     | D0TB0022   | 2-8°C                      | 31 OCT 2028     |
| TURBO-BOOST Human IL-17A Antibody                       | D0TB0028   | 2-8°C                      | 31 OCT 2028     |
| TURBO-BOOST Human IL-1 $\beta$ Antibody                 | D0TB0026   | 2-8°C                      | 30 NOV 2028     |
| TURBO-BOOST Human IL-2 Antibody                         | D0TB0024   | 2-8°C                      | 31 OCT 2028     |
| TURBO-BOOST Human IL-4 Antibody                         | D0TB0023   | 2-8°C                      | 31 OCT 2028     |
| TURBO-BOOST Human IL-6 Antibody                         | D0TB0020   | 2-8°C                      | 31 OCT 2028     |
| TURBO-BOOST Human TNF- $\alpha$ Antibody                | D0TB0025   | 2-8°C                      | 31 OCT 2028     |
| Blocker S1                                              | R0140015   | $\leq -10^{\circ}\text{C}$ | 31 OCT 2027     |
| Blocker S2                                              | 521327     | $\leq -10^{\circ}\text{C}$ | 31 MAR 2030     |
| Diluent 58                                              | M0580014   | $\leq -10^{\circ}\text{C}$ | 31 MAY 2027     |
| S-PLEX Enhance E1                                       | Y0550023   | $\leq -10^{\circ}\text{C}$ | 31 AUG 2027     |
| S-PLEX Enhance E2                                       | Y0560023   | $\leq -10^{\circ}\text{C}$ | 31 OCT 2027     |
| S-PLEX Detect D1 with TURBO-TAG <sup>TM</sup> Reagent   | Y0580032   | $\leq -70^{\circ}\text{C}$ | 31 AUG 2027     |
| S-PLEX Detect D2                                        | Y0590025   | $\leq -70^{\circ}\text{C}$ | 30 NOV 2027     |
| S-PLEX Enhance E3                                       | Y0570025   | $\leq -70^{\circ}\text{C}$ | 31 AUG 2027     |
| S-PLEX Proinflammatory Panel 1 (human) Coating Solution | 527691     | $\leq -70^{\circ}\text{C}$ | 31 AUG 2028     |

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

**Precision, Accuracy, and Sensitivity Specifications:**

| Spot Number | Name          | Calibrator |       |
|-------------|---------------|------------|-------|
|             |               | Recovery   | CV    |
| 1           | IL-6          | 80% - 120% | < 20% |
| 2           | IL-10         | 80% - 120% | < 20% |
| 3           | IL-12p70      | 80% - 120% | < 20% |
| 4           | IL-4          | 80% - 120% | < 20% |
| 5           | TNF- $\alpha$ | 80% - 120% | < 20% |
| 6           | IL-2          | 80% - 120% | < 20% |
| 8           | IL-1 $\beta$  | 75% - 125% | < 30% |
| 9           | IFN- $\gamma$ | 80% - 120% | < 20% |
| 10          | IL-17A        | 80% - 120% | < 20% |



## CERTIFICATE OF ANALYSIS

| Name          | LLOQ         |          |       |            |       | LLOD     |       |
|---------------|--------------|----------|-------|------------|-------|----------|-------|
|               | Range Tested | Observed | Units | Recovery   | CV    | Observed | Units |
| IL-6          | 3.42 - 814   | 18.9     | fg/mL | 80% - 120% | < 25% | 6.34     | fg/mL |
| IL-10         | 12.5 - 2980  | 140      | fg/mL | 80% - 120% | < 25% | 15.6     | fg/mL |
| IL-12p70      | 39.2 - 9320  | 128      | fg/mL | 80% - 120% | < 25% | 53.2     | fg/mL |
| IL-4          | 5.11 - 1220  | 16.0     | fg/mL | 80% - 120% | < 25% | 6.26     | fg/mL |
| TNF-a         | 5.39 - 1280  | 28.8     | fg/mL | 80% - 120% | < 25% | 10.5     | fg/mL |
| IL-2          | 8.75 - 2080  | 37.9     | fg/mL | 80% - 120% | < 25% | 9.10     | fg/mL |
| IL-1 $\beta$  | 16.0 - 3810  | 274      | fg/mL | 75% - 125% | < 30% | 88.4     | fg/mL |
| IFN- $\gamma$ | 3.75 - 890   | 21.2     | fg/mL | 80% - 120% | < 25% | 9.14     | fg/mL |
| IL-17A        | 19.2 - 4570  | 166      | fg/mL | 80% - 120% | < 25% | 46.4     | fg/mL |

## Stock Concentration of Calibrator:

| Name          | Item          | Concentration | Units |
|---------------|---------------|---------------|-------|
| IL-6          | 1X Calibrator | 37800         | fg/mL |
| IL-10         | 1X Calibrator | 140000        | fg/mL |
| IL-12p70      | 1X Calibrator | 512000        | fg/mL |
| IL-4          | 1X Calibrator | 64100         | fg/mL |
| TNF-a         | 1X Calibrator | 57600         | fg/mL |
| IL-2          | 1X Calibrator | 75800         | fg/mL |
| IL-1 $\beta$  | 1X Calibrator | 137000        | fg/mL |
| IFN- $\gamma$ | 1X Calibrator | 42300         | fg/mL |
| IL-17A        | 1X Calibrator | 332000        | fg/mL |

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

|                 | Name          | Function | Signature                                                                            | Date        |
|-----------------|---------------|----------|--------------------------------------------------------------------------------------|-------------|
| Review/Approval | Karen Hamilla | Quality  |  | 13 AUG 2025 |