

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX<sup>®</sup> Proinflammatory Panel 1 (human) Kit  
**Catalog Number:** K15396S-Series  
**Lot Number:** K00S0197  
**Expiration Date:** 31 JUL 2026

**Lot-Specific Components:**

| Description   | Lot Number | Storage Temperature | Expiration Date |
|---|------------|---------------------|-----------------|
| S-PLEX Proinflammatory Panel 1 (human) Calibrator Blend | A00S0093   | 2-8°C               | 31 OCT 2028     |
| TURBO-BOOST™ Human IFN-γ Antibody                       | D00T0188   | 2-8°C               | 30 JUN 2027     |
| TURBO-BOOST™ Human IL-10 Antibody                       | D00T0189   | 2-8°C               | 31 JUL 2027     |
| TURBO-BOOST™ Human IL-12p70 Antibody                    | D00T0190   | 2-8°C               | 31 JUL 2027     |
| TURBO-BOOST™ Human IL-17A Antibody                      | D00T0191   | 2-8°C               | 30 SEP 2027     |
| TURBO-BOOST™ Human IL-1β Antibody                       | D00T0192   | 2-8°C               | 30 SEP 2027     |
| TURBO-BOOST™ Human IL-2 Antibody                        | D00T0193   | 2-8°C               | 31 OCT 2027     |
| TURBO-BOOST™ Human IL-4 Antibody                        | D00T0194   | 2-8°C               | 31 JUL 2027     |
| TURBO-BOOST™ Human IL-6 Antibody                        | D00T0195   | 2-8°C               | 31 JUL 2027     |
| TURBO-BOOST™ Human TNF-α Antibody                       | D00T0196   | 2-8°C               | 31 JUL 2027     |
| Blocker S1  | R0140008   | ≤ -10°C             | 31 JUL 2026     |
| Blocker S2  | 438078     | ≤ -10°C             | 30 NOV 2028     |
| Diluent 58  | M0580011   | ≤ -10°C             | 31 JUL 2026     |
| S-PLEX Enhance E1                                       | Y0550017   | ≤ -10°C             | 31 AUG 2026     |
| S-PLEX Enhance E2                                       | Y0560017   | ≤ -10°C             | 31 JUL 2026     |
| S-PLEX Proinflammatory Panel 1 (human) Coating Solution | 477701     | ≤ -70°C             | 30 JUN 2027     |
| S-PLEX Detect D1 with TURBO-TAG™ reagent                | Y0580025   | ≤ -70°C             | 30 SEP 2026     |
| S-PLEX Detect D2  | Y0590018   | ≤ -70°C             | 31 JUL 2026     |
| S-PLEX Enhance E3                                       | Y0570019   | ≤ -70°C             | 31 JUL 2026     |

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-a    | 80% - 120% | < 20% |
| 6           | IL-2     | 80% - 120% | < 20% |
| 8           | IL-1β    | 75% - 125% | < 30% |
| 9           | IFN-γ    | 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   | 16.4     | fg/mL | 80% - 120% | < 25% | 7.72     | fg/mL |
| IL-10         | 12.5 - 2980  | 85.0     | fg/mL | 80% - 120% | < 25% | 23.4     | fg/mL |
| IL-12p70      | 39.2 - 9320  | 288      | fg/mL | 80% - 120% | < 25% | 84.4     | fg/mL |
| IL-4          | 5.11 - 1220  | 28.0     | fg/mL | 80% - 120% | < 25% | 8.50     | fg/mL |
| TNF-a         | 5.39 - 1280  | 27.4     | fg/mL | 80% - 120% | < 25% | 12.5     | fg/mL |
| IL-2          | 8.75 - 2080  | 41.5     | fg/mL | 80% - 120% | < 25% | 19.4     | fg/mL |
| IL-1 $\beta$  | 16.0 - 3810  | 81.8     | fg/mL | 75% - 125% | < 30% | 33.9     | fg/mL |
| IFN- $\gamma$ | 3.75 - 890   | 134      | fg/mL | 80% - 120% | < 25% | 9.48     | fg/mL |
| IL-17A        | 19.2 - 4570  | 149      | fg/mL | 80% - 120% | < 25% | 23.5     | fg/mL |

## Stock Concentration of Calibrator:

| Name          | Item          | Concentration | Units |
|---------------|---------------|---------------|-------|
| IL-6          | 1X Calibrator | 32800         | fg/mL |
| IL-10         | 1X Calibrator | 170000        | fg/mL |
| IL-12p70      | 1X Calibrator | 576000        | fg/mL |
| IL-4          | 1X Calibrator | 56000         | fg/mL |
| TNF-a         | 1X Calibrator | 54850         | fg/mL |
| IL-2          | 1X Calibrator | 83050         | fg/mL |
| IL-1 $\beta$  | 1X Calibrator | 163500        | fg/mL |
| IFN- $\gamma$ | 1X Calibrator | 66900         | fg/mL |
| IL-17A        | 1X Calibrator | 297000        | 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  |  | 17 MAY 2024 |