



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0396-2
Lot Number: A00S0141
Expiration Date: 28 FEB 2031
Storage Temperature: 2-8°C

Concentration Assignment:

| Analyte | Stock Concentration of Reconstituted Calibrator | Unit of Measure |
|---------------|-------------------------------------------------|-----------------|
| IL-6 | 39600 | fg/mL |
| IL-10 | 150000 | fg/mL |
| IL-12p70 | 440000 | fg/mL |
| IL-4 | 64000 | fg/mL |
| TNF- α | 53700 | fg/mL |
| IL-2 | 75500 | fg/mL |
| IL-1 β | 137000 | fg/mL |
| IFN- γ | 43600 | fg/mL |
| IL-17A | 252000 | fg/mL |

Reconstitution Volume: 1000 μ L

This product was manufactured and tested according to MSD documents. This lot meets MSD’s specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|------------------------|---------------|----------|-----------|-------------|
| Review/Approval | Karen Hamilla | Quality | | 29 APR 2026 |