

**CERTIFICATE OF ANALYSIS**


**Product Description:** S-PLEX Proinflammatory Panel 1 (human) Calibrator Blend  
**Catalog Number:** C0396-2  
**Lot Number:** A00S0127  
**Expiration Date:** 30 JUN 2030  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-6	37800	fg/mL
IL-10	140000	fg/mL
IL-12p70	512000	fg/mL
IL-4	64100	fg/mL
TNF- $\alpha$	57600	fg/mL
IL-2	75800	fg/mL
IL-1 $\beta$	137000	fg/mL
IFN- $\gamma$	42300	fg/mL
IL-17A	332000	fg/mL

**Reconstitution Volume:** 1000  $\mu$ 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
<b>Review/Approval</b>	Karen Hamilla	Quality		25 JUL 2025