



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0396-2
Lot Number: A00S0160
Expiration Date: 31 JUL 2031
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-6	37400	fg/mL
IL-10	146000	fg/mL
IL-12p70	447000	fg/mL
IL-4	62800	fg/mL
TNF- α	52600	fg/mL
IL-2	71400	fg/mL
IL-1 β	147000	fg/mL
IFN- γ	44300	fg/mL
IL-17A	226000	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		24 AUG 2026