

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX<sup>®</sup> Proinflammatory Panel 1 (human) Kit  
**Catalog Number:** K15396S-Series  
**Lot Number:** K00S0133  
**Expiration Date:** 30 APR 2025

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
S-PLEX Proinflammatory Panel 1 (human) Calibrator Blend	A00S0078	2-8°C	31 DEC 2027
TURBO-BOOST™ Human IFN-γ Antibody	D00T0139	2-8°C	30 SEP 2026
TURBO-BOOST™ Human IL-10 Antibody	D00T0140	2-8°C	31 OCT 2026
TURBO-BOOST™ Human IL-12p70 Antibody	D00T0141	2-8°C	31 AUG 2026
TURBO-BOOST™ Human IL-17A Antibody	D00T0142	2-8°C	30 SEP 2026
TURBO-BOOST™ Human IL-1β Antibody	D00T0143	2-8°C	31 AUG 2026
TURBO-BOOST™ Human IL-2 Antibody	D00T0144	2-8°C	31 OCT 2026
TURBO-BOOST™ Human IL-4 Antibody	D00T0145	2-8°C	30 SEP 2026
TURBO-BOOST™ Human IL-6 Antibody	D00T0146	2-8°C	31 OCT 2026
TURBO-BOOST™ Human TNF-α Antibody	D00T0147	2-8°C	31 OCT 2026
Blocker S1	R0010824	≤ -10°C	30 JUN 2025
Diluent 58 - 10 mL	M0580007	≤ -10°C	30 APR 2025
S-PLEX Enhance E1	Y0550010	≤ -10°C	31 MAY 2025
S-PLEX Enhance E2	Y0560012	≤ -10°C	31 MAY 2025
Blocker S2	437541	≤ -70°C	31 DEC 2027
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580016	≤ -70°C	31 MAY 2025
S-PLEX Detect D2	Y0590012	≤ -70°C	30 APR 2025
S-PLEX Enhance E3	Y0570012	≤ -70°C	31 MAY 2025
S-PLEX Proinflammatory Panel 1 (human) Coating Solution	430391	≤ -70°C	31 MAY 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	78.0	fg/mL	80% - 120%	< 25%	25.4	fg/mL
IL-10	12.5 - 2980	1340	fg/mL	80% - 120%	< 25%	69.4	fg/mL
IL-12p70	39.2 - 9320	1010	fg/mL	80% - 120%	< 25%	169	fg/mL
IL-4	5.11 - 1220	55.0	fg/mL	80% - 120%	< 25%	22.3	fg/mL
TNF-a	5.39 - 1280	49.2	fg/mL	80% - 120%	< 25%	22.9	fg/mL
IL-2	8.75 - 2080	629	fg/mL	80% - 120%	< 25%	34.8	fg/mL
IL-1β	16.0 - 3810	3810	fg/mL	75% - 125%	< 30%	125	fg/mL
IFN-γ	3.75 - 890	324	fg/mL	80% - 120%	< 25%	15.4	fg/mL
IL-17A	19.2 - 4570	216	fg/mL	80% - 120%	< 25%	109	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-6	1X Calibrator	39000	fg/mL
IL-10	1X Calibrator	167000	fg/mL
IL-12p70	1X Calibrator	505000	fg/mL
IL-4	1X Calibrator	55000	fg/mL
TNF-a	1X Calibrator	49200	fg/mL
IL-2	1X Calibrator	78600	fg/mL
IL-1β	1X Calibrator	160000	fg/mL
IFN-γ	1X Calibrator	40500	fg/mL
IL-17A	1X Calibrator	216000	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		22 FEB 2024