

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Proinflammatory Panel 1 (human) Kit
Catalog Number: K15396S-Series
Lot Number: K00S0138
Expiration Date: 31 OCT 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	D00T0127	2-8°C	31 AUG 2026
TURBO-BOOST™ Human IL-10 Antibody	D00T0128	2-8°C	31 JUL 2026
TURBO-BOOST™ Human IL-12p70 Antibody	D00T0129	2-8°C	31 AUG 2026
TURBO-BOOST™ Human IL-17A Antibody	D00T0130	2-8°C	30 SEP 2026
TURBO-BOOST™ Human IL-1β Antibody	D00T0131	2-8°C	31 JUL 2026
TURBO-BOOST™ Human IL-2 Antibody	D00T0132	2-8°C	31 OCT 2026
TURBO-BOOST™ Human IL-4 Antibody	D00T0133	2-8°C	31 JUL 2026
TURBO-BOOST™ Human IL-6 Antibody	D00T0134	2-8°C	31 JUL 2026
TURBO-BOOST™ Human TNF-α Antibody	D00T0135	2-8°C	30 SEP 2026
Blocker S1	R0010832	≤ -10°C	31 OCT 2025
Blocker S2	438076	≤ -10°C	31 MAY 2028
Diluent 58	M0580009	≤ -10°C	30 NOV 2025
S-PLEX Enhance E1	Y0550013	≤ -10°C	30 NOV 2025
S-PLEX Enhance E2	Y0560014	≤ -10°C	31 OCT 2025
S-PLEX Proinflammatory Panel 1 (human) Coating Solution	457686	≤ -70°C	31 DEC 2026
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580018	≤ -70°C	31 OCT 2025
S-PLEX Detect D2	Y0590014	≤ -70°C	31 OCT 2025
S-PLEX Enhance E3	Y0570014	≤ -70°C	31 OCT 2025

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	39.0	fg/mL	80% - 120%	< 25%	17.0	fg/mL
IL-10	12.5 - 2980	83.5	fg/mL	80% - 120%	< 25%	23.9	fg/mL
IL-12p70	39.2 - 9320	505	fg/mL	80% - 120%	< 25%	74.9	fg/mL
IL-4	5.11 - 1220	27.5	fg/mL	80% - 120%	< 25%	15.2	fg/mL
TNF-a	5.39 - 1280	24.6	fg/mL	80% - 120%	< 25%	8.78	fg/mL
IL-2	8.75 - 2080	39.3	fg/mL	80% - 120%	< 25%	14.2	fg/mL
IL-1 β	16.0 - 3810	160	fg/mL	75% - 125%	< 30%	69.4	fg/mL
IFN- γ	3.75 - 890	20.3	fg/mL	80% - 120%	< 25%	6.46	fg/mL
IL-17A	19.2 - 4570	108	fg/mL	80% - 120%	< 25%	31.9	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-1 β	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		23 FEB 2024