

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Proinflammatory Panel 1 (human) Kit
Catalog Number: K15396S-Series
Lot Number: K00S0200
Expiration Date: 28 FEB 2026

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
S-PLEX Proinflammatory Panel 1 (human) Calibrator Blend	A00S0093	2-8°C	31 OCT 2028
TURBO-BOOST™ Human IFN-γ Antibody	D00T0157	2-8°C	30 APR 2027
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	D00T0161	2-8°C	31 JAN 2027
TURBO-BOOST™ Human IL-2 Antibody	D00T0144	2-8°C	31 OCT 2026
TURBO-BOOST™ Human IL-4 Antibody	D00T0163	2-8°C	31 MAR 2027
TURBO-BOOST™ Human IL-6 Antibody	D00T0146	2-8°C	31 OCT 2026
TURBO-BOOST™ Human TNF-α Antibody	D00T0165	2-8°C	31 JAN 2027
Blocker S1	R0010871	≤ -10°C	31 MAR 2026
Blocker S2	438077	≤ -10°C	30 SEP 2028
Diluent 58	M0580011	≤ -10°C	31 JUL 2026
S-PLEX Enhance E1	Y0550014	≤ -10°C	28 FEB 2026
S-PLEX Enhance E2	Y0560015	≤ -10°C	28 FEB 2026
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580021	≤ -70°C	30 APR 2026
S-PLEX Detect D2	Y0590016	≤ -70°C	30 APR 2026
S-PLEX Enhance E3	Y0570017	≤ -70°C	31 MAR 2026
S-PLEX Proinflammatory Panel 1(human) Coating Solution	468937	≤ -70°C	31 MAR 2027

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	16.4	fg/mL	80% - 120%	< 25%	5.60	fg/mL
IL-10	12.5 - 2980	340	fg/mL	80% - 120%	< 25%	28.6	fg/mL
IL-12p70	39.2 - 9320	4610	fg/mL	80% - 120%	< 25%	109	fg/mL
IL-4	5.11 - 1220	14.0	fg/mL	80% - 120%	< 25%	6.76	fg/mL
TNF-a	5.39 - 1280	13.7	fg/mL	80% - 120%	< 25%	9.66	fg/mL
IL-2	8.75 - 2080	41.5	fg/mL	80% - 120%	< 25%	22.4	fg/mL
IL-1 β	16.0 - 3810	40.9	fg/mL	75% - 125%	< 30%	31.9	fg/mL
IFN- γ	3.75 - 890	8.36	fg/mL	80% - 120%	< 25%	4.53	fg/mL
IL-17A	19.2 - 4570	37.1	fg/mL	80% - 120%	< 25%	20.3	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-6	1X Calibrator	32800	fg/mL
IL-10	1X Calibrator	170000	fg/mL
IL-12p70	1X Calibrator	576000	fg/mL
IL-4	1X Calibrator	56000	fg/mL
TNF-a	1X Calibrator	54850	fg/mL
IL-2	1X Calibrator	83050	fg/mL
IL-1 β	1X Calibrator	163500	fg/mL
IFN- γ	1X Calibrator	66900	fg/mL
IL-17A	1X Calibrator	297000	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		19 APR 2024