

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX<sup>®</sup> Proinflammatory Panel 1 (human) Kit  
**Lot Number:** K00S0225  
**Expiration Date:** 30 SEP 2026  
**Catalog Number:** K15396S-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
S-PLEX Proinflammatory Panel 1 (human) Calibrator Blend	A00S0107	2-8°C	31 JUL 2029
TURBO-BOOST Human IFN- $\gamma$ Antibody	D0TB0006	2-8°C	30 NOV 2027
TURBO-BOOST Human IL-10 Antibody	D0TB0001	2-8°C	30 NOV 2027
TURBO-BOOST Human IL-12p70 Antibody	D0TB0002	2-8°C	30 NOV 2027
TURBO-BOOST Human IL-17A Antibody	D0TB0007	2-8°C	31 DEC 2027
TURBO-BOOST Human IL-1 $\beta$ Antibody	D0TB0005	2-8°C	30 NOV 2027
TURBO-BOOST Human IL-2 Antibody	D00T0213	2-8°C	31 OCT 2027
TURBO-BOOST Human IL-4 Antibody	D0TB0003	2-8°C	30 NOV 2027
TURBO-BOOST Human IL-6 Antibody	D00T0209	2-8°C	31 OCT 2027
TURBO-BOOST Human TNF- $\alpha$ Antibody	D0TB0004	2-8°C	30 NOV 2027
Blocker S1	R0140010	$\leq -10^{\circ}\text{C}$	31 DEC 2026
Blocker S2	438078	$\leq -10^{\circ}\text{C}$	30 NOV 2028
Diluent 58	M0580012	$\leq -10^{\circ}\text{C}$	31 OCT 2026
S-PLEX Enhance E1	Y0550018	$\leq -10^{\circ}\text{C}$	30 NOV 2027
S-PLEX Enhance E2	Y0560019	$\leq -10^{\circ}\text{C}$	30 NOV 2026
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> Reagent	Y0580027	$\leq -70^{\circ}\text{C}$	31 OCT 2026
S-PLEX Detect D2	Y0590019	$\leq -70^{\circ}\text{C}$	30 SEP 2026
S-PLEX Enhance E3	Y0570020	$\leq -70^{\circ}\text{C}$	31 OCT 2026
S-PLEX Proinflammatory Panel 1 (human) Coating Solution	496112	$\leq -70^{\circ}\text{C}$	30 NOV 2027

Additional components may be provided to execute the assay. Additional components are not lot-specific.

**Precision, Accuracy, and Sensitivity Specifications:**

Spot Number	Name	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-1 $\beta$	75% - 125%	< 30%
9	IFN- $\gamma$	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	18.0	fg/mL	80% - 120%	< 25%	6.00	fg/mL
IL-10	12.5 - 2980	93.0	fg/mL	80% - 120%	< 25%	22.5	fg/mL
IL-12p70	39.2 - 9320	125	fg/mL	80% - 120%	< 25%	33.2	fg/mL
IL-4	5.11 - 1220	31.0	fg/mL	80% - 120%	< 25%	5.50	fg/mL
TNF-a	5.39 - 1280	23.1	fg/mL	80% - 120%	< 25%	7.54	fg/mL
IL-2	8.75 - 2080	43.1	fg/mL	80% - 120%	< 25%	14.8	fg/mL
IL-1β	16.0 - 3810	88.0	fg/mL	75% - 125%	< 30%	32.1	fg/mL
IFN-γ	3.75 - 890	11.5	fg/mL	80% - 120%	< 25%	4.51	fg/mL
IL-17A	19.2 - 4570	29.9	fg/mL	80% - 120%	< 25%	14.7	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-6	1X Calibrator	36000	fg/mL
IL-10	1X Calibrator	186000	fg/mL
IL-12p70	1X Calibrator	501000	fg/mL
IL-4	1X Calibrator	62000	fg/mL
TNF-a	1X Calibrator	46200	fg/mL
IL-2	1X Calibrator	86200	fg/mL
IL-1β	1X Calibrator	176000	fg/mL
IFN-γ	1X Calibrator	45900	fg/mL
IL-17A	1X Calibrator	239000	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		27 SEP 2024