

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Proinflammatory Panel 1 (human) Kit
Catalog Number: K15396S-Series
Lot Number: K00S0102
Expiration Date: 28 FEB 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
S-PLEX Proinflammatory Panel 1 (human) Calibrator Blend	A00S0065	2-8°C	30 JUN 2027
TURBO-BOOST™ Human IFN- γ Antibody	D00T0107	2-8°C	31 MAR 2026
TURBO-BOOST™ Human IL-10 Antibody	D00T0101	2-8°C	31 MAR 2026
TURBO-BOOST™ Human IL-12p70 Antibody	D00T0102	2-8°C	31 MAR 2026
TURBO-BOOST™ Human IL-17A Antibody	D00T0108	2-8°C	31 MAR 2026
TURBO-BOOST™ Human IL-1 β Antibody	D00T0106	2-8°C	31 MAR 2026
TURBO-BOOST™ Human IL-2 Antibody	D00T0104	2-8°C	31 MAR 2026
TURBO-BOOST™ Human IL-4 Antibody	D00T0103	2-8°C	31 MAR 2026
TURBO-BOOST™ Human IL-6 Antibody	D00T0100	2-8°C	31 MAR 2026
TURBO-BOOST™ Human TNF- α Antibody	D00T0105	2-8°C	31 MAR 2026
Blocker S1	R0010811	$\leq -10^{\circ}\text{C}$	28 FEB 2025
Blocker S2	420803	$\leq -10^{\circ}\text{C}$	30 SEP 2027
Diluent 58	M0580007	$\leq -10^{\circ}\text{C}$	30 APR 2025
S-PLEX Enhance E1	Y0550009	$\leq -10^{\circ}\text{C}$	28 FEB 2025
S-PLEX Enhance E2	Y0560011	$\leq -10^{\circ}\text{C}$	31 MAR 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580014	$\leq -70^{\circ}\text{C}$	28 FEB 2025
S-PLEX Detect D2	Y0590011	$\leq -70^{\circ}\text{C}$	31 MAR 2025
S-PLEX Enhance E3	Y0570011	$\leq -70^{\circ}\text{C}$	31 MAR 2025
S-PLEX Proinflammatory Panel 1(human) Coating Solution	429069	$\leq -70^{\circ}\text{C}$	30 APR 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- α	80% - 120%	< 20%
6	IL-2	80% - 120%	< 20%
8	IL-1 β	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	79.0	fg/mL	80% - 120%	< 25%	16.1	fg/mL
IL-10	12.5 - 2980	74.0	fg/mL	80% - 120%	< 25%	23.2	fg/mL
IL-12p70	39.2 - 9320	484	fg/mL	80% - 120%	< 25%	96.6	fg/mL
IL-4	5.11 - 1220	27.7	fg/mL	80% - 120%	< 25%	9.94	fg/mL
TNF- α	5.39 - 1280	53.4	fg/mL	80% - 120%	< 25%	14.2	fg/mL
IL-2	8.75 - 2080	21.5	fg/mL	80% - 120%	< 25%	12.3	fg/mL
IL-1 β	16.0 - 3810	159	fg/mL	75% - 125%	< 30%	66.5	fg/mL
IFN- γ	3.75 - 890	14.4	fg/mL	80% - 120%	< 25%	7.81	fg/mL
IL-17A	19.2 - 4570	210	fg/mL	80% - 120%	< 25%	68.2	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-6	1X Calibrator	39500	fg/mL
IL-10	1X Calibrator	148000	fg/mL
IL-12p70	1X Calibrator	484000	fg/mL
IL-4	1X Calibrator	55300	fg/mL
TNF- α	1X Calibrator	53400	fg/mL
IL-2	1X Calibrator	85800	fg/mL
IL-1 β	1X Calibrator	159000	fg/mL
IFN- γ	1X Calibrator	28800	fg/mL
IL-17A	1X Calibrator	210000	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