



## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** S-PLEX<sup>®</sup> Proinflammatory Panel 1 (human) Kit

**Catalog Number:** K15396S - Series

**Lot Number:** K00S0124

**Expiration Date:** 30 Apr 2025

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST <sup>™</sup> Human IL-6 Antibody	D00T0100	2-8°C	31 Mar 2026
TURBO-BOOST <sup>™</sup> Human IL-10 Antibody	D00T0101	2-8°C	31 Mar 2026
TURBO-BOOST <sup>™</sup> Human IL-12p70 Antibody	D00T0102	2-8°C	31 Mar 2026
TURBO-BOOST <sup>™</sup> Human IL-4 Antibody	D00T0103	2-8°C	31 Mar 2026
TURBO-BOOST <sup>™</sup> Human TNF- $\alpha$ Antibody	D00T0105	2-8°C	31 Mar 2026
TURBO-BOOST <sup>™</sup> Human IL-2 Antibody	D00T0104	2-8°C	31 Mar 2026
TURBO-BOOST <sup>™</sup> Human IL-1 $\beta$ Antibody	D00T0106	2-8°C	31 Mar 2026
TURBO-BOOST <sup>™</sup> Human IFN- $\gamma$ Antibody	D00T0107	2-8°C	31 Mar 2026
TURBO-BOOST <sup>™</sup> Human IL-17A Antibody	D00T0108	2-8°C	31 Mar 2026
S-PLEX Proinflammatory Panel 1 (human) Coating Solution	430391	$\leq -70^{\circ}\text{C}$	31 May 2026
S-PLEX Proinflammatory Panel 1 (human) Calibrator	A00S0065	2-8°C	30 Jun 2027
Diluent 58	M0580007	$\leq -10^{\circ}\text{C}$	30 Apr 2025
Blocker S1	R0010824	$\leq -10^{\circ}\text{C}$	30 Jun 2025
Blocker S2	420803	$\leq -10^{\circ}\text{C}$	30 Sep 2027
S-PLEX Enhance E1	Y0550010	$\leq -10^{\circ}\text{C}$	31 May 2025
S-PLEX Enhance E2	Y0560012	$\leq -10^{\circ}\text{C}$	31 May 2025
S-PLEX Enhance E3	Y0570012	$\leq -70^{\circ}\text{C}$	31 May 2025
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> reagent	Y0580016	$\leq -70^{\circ}\text{C}$	31 May 2025
S-PLEX Detect D2	Y0590012	$\leq -70^{\circ}\text{C}$	30 Apr 2025

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

### Precision, Accuracy, and Sensitivity Specifications:

Spot	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- $\alpha$	80% - 120%	< 20%
6	IL-2	80% - 120%	< 20%
8	IL-1 $\beta$	75% - 125%	< 25%
9	IFN- $\gamma$	80% - 120%	< 20%
10	IL-17A	80% - 120%	< 20%



### CERTIFICATE OF ANALYSIS

Page 2 of 2

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-6	3.42 - 814	79.0	fg/mL	80% - 120%	< 25%	25.1	fg/mL
IL-10	12.5 - 2980	148	fg/mL	80% - 120%	< 25%	30.4	fg/mL
IL-12p70	39.2 - 9320	484	fg/mL	80% - 120%	< 25%	181	fg/mL
IL-4	5.11 - 1220	55.3	fg/mL	80% - 120%	< 25%	9.46	fg/mL
TNF- $\alpha$	5.39 - 1280	53.4	fg/mL	80% - 120%	< 25%	14.0	fg/mL
IL-2	8.75 - 2080	42.9	fg/mL	80% - 120%	< 25%	14.4	fg/mL
IL-1 $\beta$	16.0 - 3810	318	fg/mL	75% - 125%	< 30%	88.4	fg/mL
IFN- $\gamma$	3.75 - 890	28.8	fg/mL	80% - 120%	< 25%	11.1	fg/mL
IL-17A	19.2 - 4570	210	fg/mL	80% - 120%	< 25%	117	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- $\alpha$	1X Calibrator	53400	fg/mL
IL-2	1X Calibrator	85800	fg/mL
IL-1 $\beta$	1X Calibrator	159000	fg/mL
IFN- $\gamma$	1X Calibrator	28800	fg/mL
IL-17A	1X Calibrator	210000	fg/mL

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		25 JAN 2023