



## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** S-PLEX® Proinflammatory Panel 1 (human) Kit

**Catalog Number:** K15396S-Series

**Lot Number:** K00S0062

**Expiration Date:** 28 FEB 2023

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-6 Antibody	D00T0051A	2-8°C	31 Aug 2024
TURBO-BOOST™ Human IL-10 Antibody	D00T0052A	2-8°C	31 Oct 2024
TURBO-BOOST™ Human IL-12p70 Antibody	D00T0053A	2-8°C	31 Aug 2024
TURBO-BOOST™ Human IL-4 Antibody	D00T0054A	2-8°C	31 Aug 2024
TURBO-BOOST™ Human TNF-α Antibody	D00T0069A	2-8°C	31 Oct 2024
TURBO-BOOST™ Human IL-2 Antibody	D00T0055A	2-8°C	31 Aug 2024
TURBO-BOOST™ Human IL-1β Antibody	D00T0056A	2-8°C	31 Aug 2024
TURBO-BOOST™ Human IFN-γ Antibody	D00T0057A	2-8°C	31 Oct 2024
TURBO-BOOST™ Human IL-17A Antibody	D00T0058A	2-8°C	31 Aug 2024
S-PLEX Proinflammatory Panel 1 (human) Coating Solution	370063	≤ -70°C	30 Nov 2024
S-PLEX Proinflammatory Panel 1 (human) Calibrator Blend	A00S0042	2-8°C	31 May 2026
Diluent 58	M0580001	≤ -10°C	28 Feb 2023
Blocker S1	R0140004	≤ -10°C	30 Apr 2023
Blocker S2	338741	≤ -10°C	31 Dec 2025
S-PLEX Enhance E1	Y0550006	≤ -10°C	31 Oct 2023
S-PLEX Enhance E2	Y0560007	≤ -10°C	31 Oct 2023
S-PLEX Enhance E3	Y0570006	≤ -70°C	31 Oct 2023
S-PLEX Detect D1	Y0580008	≤ -70°C	30 Nov 2023
S-PLEX Detect D2	Y0590006	≤ -70°C	31 Oct 2023

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-α	80% - 120%	< 20%
6	IL-2	80% - 120%	< 20%
8	IL-1β	80% - 120%	< 20%
9	IFN-γ	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	53.9	fg/mL	80% - 120%	< 25%	16.1	fg/mL
IL-10	12.5 - 2980	174	fg/mL	80% - 120%	< 25%	43.8	fg/mL
IL-12p70	39.2 - 9320	486	fg/mL	80% - 120%	< 25%	122	fg/mL
IL-4	5.11 - 1220	27.1	fg/mL	80% - 120%	< 25%	11.3	fg/mL
TNF- $\alpha$	5.39 - 1280	52.6	fg/mL	80% - 120%	< 25%	27.4	fg/mL
IL-2	8.75 - 2080	80.2	fg/mL	80% - 120%	< 25%	26.2	fg/mL
IL-1 $\beta$	16.0 - 3810	168	fg/mL	80% - 120%	< 25%	111	fg/mL
IFN- $\gamma$	3.75 - 890	127	fg/mL	80% - 120%	< 25%	12.3	fg/mL
IL-17A	19.2 - 4570	130	fg/mL	80% - 120%	< 25%	46.7	fg/mL

#### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-6	1X Calibrator	53900	fg/mL
IL-10	1X Calibrator	174000	fg/mL
IL-12p70	1X Calibrator	486000	fg/mL
IL-4	1X Calibrator	54200	fg/mL
TNF- $\alpha$	1X Calibrator	52600	fg/mL
IL-2	1X Calibrator	80200	fg/mL
IL-1 $\beta$	1X Calibrator	168000	fg/mL
IFN- $\gamma$	1X Calibrator	31700	fg/mL
IL-17A	1X Calibrator	259000	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		19 NOV 2021