

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Proinflammatory Panel 1 (human) Kit
Catalog Number: K15396S-Series
Lot Number: K00S0075
Expiration Date: 28 FEB 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
S-PLEX Proinflammatory Panel 1 (human) Calibrator Blend	A00S0042	2-8°C	31 MAY 2026
TURBO-BOOST [™] Human IFN- γ Antibody	D00T0057A	2-8°C	31 OCT 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-17A Antibody	D00T0058A	2-8°C	31 AUG 2024
TURBO-BOOST [™] Human IL-1 β Antibody	D00T0056A	2-8°C	31 AUG 2024
TURBO-BOOST [™] Human IL-2 Antibody	D00T0055A	2-8°C	31 AUG 2024
TURBO-BOOST [™] Human IL-4 Antibody	D00T0054A	2-8°C	31 AUG 2024
TURBO-BOOST [™] Human IL-6 Antibody	D00T0051A	2-8°C	31 AUG 2024
TURBO-BOOST [™] Human TNF- α Antibody	D00T0069A	2-8°C	31 OCT 2024
Blocker S1	R0140004	$\leq -10^{\circ}\text{C}$	30 APR 2023
Blocker S2	338741	$\leq -10^{\circ}\text{C}$	31 DEC 2025
Diluent 58	M0580001	$\leq -10^{\circ}\text{C}$	28 FEB 2023
S-PLEX Enhance E1	Y0550006	$\leq -10^{\circ}\text{C}$	31 OCT 2023
S-PLEX Enhance E2	Y0560008	$\leq -10^{\circ}\text{C}$	30 APR 2024
S-PLEX Coating Reagent C1	371318	$\leq -70^{\circ}\text{C}$	30 JUN 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580008	$\leq -70^{\circ}\text{C}$	30 NOV 2023
S-PLEX Detect D2	Y0590006	$\leq -70^{\circ}\text{C}$	31 OCT 2023
S-PLEX Enhance E3	Y0570006	$\leq -70^{\circ}\text{C}$	31 OCT 2023

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

Precision, Accuracy, and Sensitivity Specifications:

Spot 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	27.0	fg/mL	80% - 120%	< 25%	12.8	fg/mL
IL-10	12.5 - 2980	43.5	fg/mL	80% - 120%	< 25%	25.1	fg/mL
IL-12p70	39.2 - 9320	122	fg/mL	80% - 120%	< 25%	67.3	fg/mL
IL-4	5.11 - 1220	13.6	fg/mL	80% - 120%	< 25%	6.84	fg/mL
TNF- α	5.39 - 1280	52.6	fg/mL	80% - 120%	< 25%	12.7	fg/mL
IL-2	8.75 - 2080	80.2	fg/mL	80% - 120%	< 25%	13.6	fg/mL
IL-1 β	16.0 - 3810	168	fg/mL	75% - 125%	< 30%	79.1	fg/mL
IFN- γ	3.75 - 890	31.7	fg/mL	80% - 120%	< 25%	8.39	fg/mL
IL-17A	19.2 - 4570	130	fg/mL	80% - 120%	< 25%	39.9	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- α	1X Calibrator	52600	fg/mL
IL-2	1X Calibrator	80200	fg/mL
IL-1 β	1X Calibrator	168000	fg/mL
IFN- γ	1X Calibrator	31700	fg/mL
IL-17A	1X Calibrator	259000	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		21 FEB 2024