

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Proinflammatory Panel 1 (human) Kit
Lot Number: K00S0279
Expiration Date: 31 JAN 2027
Catalog Number: K15396S-Series; K151AKV-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- γ Antibody	D0TB0018	2-8°C	30 JUN 2028
TURBO-BOOST Human IL-10 Antibody	D0TB0012	2-8°C	31 MAY 2028
TURBO-BOOST Human IL-12p70 Antibody	D0TB0002	2-8°C	30 NOV 2027
TURBO-BOOST Human IL-17A Antibody	D0TB0019	2-8°C	30 JUN 2028
TURBO-BOOST Human IL-1 β Antibody	D0TB0017	2-8°C	31 MAY 2028
TURBO-BOOST Human IL-2 Antibody	D0TB0015	2-8°C	31 JUL 2028
TURBO-BOOST Human IL-4 Antibody	D0TB0003	2-8°C	30 NOV 2027
TURBO-BOOST Human IL-6 Antibody	D0TB0011	2-8°C	31 JUL 2028
TURBO-BOOST Human TNF- α Antibody	D0TB0016	2-8°C	31 JUL 2028
Blocker S1	R0140012	$\leq -10^{\circ}\text{C}$	30 APR 2027
Blocker S2	476748	$\leq -10^{\circ}\text{C}$	30 JUN 2029
Diluent 58	M0580014	$\leq -10^{\circ}\text{C}$	31 MAY 2027
S-PLEX Enhance E1	Y0550019	$\leq -10^{\circ}\text{C}$	31 MAR 2027
S-PLEX Enhance E2	Y0560020	$\leq -10^{\circ}\text{C}$	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG TM Reagent	Y0580028	$\leq -70^{\circ}\text{C}$	31 JAN 2027
S-PLEX Detect D2	Y0590021	$\leq -70^{\circ}\text{C}$	31 MAR 2027
S-PLEX Enhance E3	Y0570022	$\leq -70^{\circ}\text{C}$	30 APR 2027
S-PLEX Proinflammatory Panel 1 (human) Coating Solution	512758	$\leq -70^{\circ}\text{C}$	31 MAR 2028

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	18.0	fg/mL	80% - 120%	< 25%	6.91	fg/mL
IL-10	12.5 - 2980	372	fg/mL	80% - 120%	< 25%	24.2	fg/mL
IL-12p70	39.2 - 9320	125	fg/mL	80% - 120%	< 25%	59.9	fg/mL
IL-4	5.11 - 1220	15.5	fg/mL	80% - 120%	< 25%	6.89	fg/mL
TNF- α	5.39 - 1280	23.1	fg/mL	80% - 120%	< 25%	10.9	fg/mL
IL-2	8.75 - 2080	43.1	fg/mL	80% - 120%	< 25%	12.7	fg/mL
IL-1 β	16.0 - 3810	352	fg/mL	75% - 125%	< 30%	96.0	fg/mL
IFN- γ	3.75 - 890	45.9	fg/mL	80% - 120%	< 25%	9.29	fg/mL
IL-17A	19.2 - 4570	120	fg/mL	80% - 120%	< 25%	38.7	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- α	1X Calibrator	46200	fg/mL
IL-2	1X Calibrator	86200	fg/mL
IL-1 β	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		23 APR 2025