

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Proinflammatory Panel 1 (human) Kit
Lot Number: K00S0299
Expiration Date: 31 MAY 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	A00S0127	2-8°C	30 JUN 2030
TURBO-BOOST Human IFN- γ Antibody	D0TB0039	2-8°C	30 NOV 2028
TURBO-BOOST Human IL-10 Antibody	D0TB0033	2-8°C	28 FEB 2029
TURBO-BOOST Human IL-12p70 Antibody	D0TB0034	2-8°C	30 NOV 2028
TURBO-BOOST Human IL-17A Antibody	D0TB0040	2-8°C	31 MAR 2029
TURBO-BOOST Human IL-1 β Antibody	D0TB0038	2-8°C	31 JAN 2029
TURBO-BOOST Human IL-2 Antibody	D0TB0036	2-8°C	30 NOV 2028
TURBO-BOOST Human IL-4 Antibody	D0TB0035	2-8°C	31 JAN 2029
TURBO-BOOST Human IL-6 Antibody	D0TB0032	2-8°C	30 NOV 2028
TURBO-BOOST Human TNF- α Antibody	D0TB0037	2-8°C	30 NOV 2028
Blocker S1	R0140015	$\leq -10^{\circ}\text{C}$	31 OCT 2027
Blocker S2	521326	$\leq -10^{\circ}\text{C}$	31 DEC 2029
Diluent 58	M0580014	$\leq -10^{\circ}\text{C}$	31 MAY 2027
S-PLEX Enhance E1	Y0550023	$\leq -10^{\circ}\text{C}$	31 AUG 2027
S-PLEX Enhance E2	Y0560025	$\leq -10^{\circ}\text{C}$	29 FEB 2028
S-PLEX Detect D1 with TURBO-TAG TM Reagent	Y0580033	$\leq -70^{\circ}\text{C}$	30 NOV 2027
S-PLEX Detect D2	Y0590027	$\leq -70^{\circ}\text{C}$	31 JAN 2028
S-PLEX Enhance E3	Y0570026	$\leq -70^{\circ}\text{C}$	31 OCT 2027
S-PLEX Proinflammatory Panel 1 (human) Coating Solution	528464	$\leq -70^{\circ}\text{C}$	31 OCT 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	18.9	fg/mL	80% - 120%	< 25%	5.19	fg/mL
IL-10	12.5 - 2980	70.0	fg/mL	80% - 120%	< 25%	13.3	fg/mL
IL-12p70	39.2 - 9320	128	fg/mL	80% - 120%	< 25%	32.6	fg/mL
IL-4	5.11 - 1220	16.0	fg/mL	80% - 120%	< 25%	5.10	fg/mL
TNF-a	5.39 - 1280	14.4	fg/mL	80% - 120%	< 25%	5.68	fg/mL
IL-2	8.75 - 2080	37.9	fg/mL	80% - 120%	< 25%	8.10	fg/mL
IL-1 β	16.0 - 3810	137	fg/mL	75% - 125%	< 30%	18.5	fg/mL
IFN- γ	3.75 - 890	10.6	fg/mL	80% - 120%	< 25%	2.79	fg/mL
IL-17A	19.2 - 4570	83.0	fg/mL	80% - 120%	< 25%	17.8	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-6	1X Calibrator	37800	fg/mL
IL-10	1X Calibrator	140000	fg/mL
IL-12p70	1X Calibrator	512000	fg/mL
IL-4	1X Calibrator	64100	fg/mL
TNF-a	1X Calibrator	57600	fg/mL
IL-2	1X Calibrator	75800	fg/mL
IL-1 β	1X Calibrator	137000	fg/mL
IFN- γ	1X Calibrator	42300	fg/mL
IL-17A	1X Calibrator	332000	fg/mL

This lot meets MSD's non-specific binding specifications.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		21 NOV 2025