

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX Proinflammatory Panel 1 (human) Kit  
**Lot Number:** K00S0303  
**Expiration Date:** 31 JUL 2027  
**Catalog Number:** K15396S-Series, K151AKV-Series

**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
S-PLEX Proinflammatory Panel 1 (human) Calibrator Blend	A00S0141	2-8°C	28 FEB 2031
TURBO-BOOST Human IFN- $\gamma$ Antibody	D0TB0027	2-8°C	31 OCT 2028
TURBO-BOOST Human IL-10 Antibody	D0TB0021	2-8°C	31 OCT 2028
TURBO-BOOST Human IL-12p70 Antibody	D0TB0013	2-8°C	31 MAY 2028
TURBO-BOOST Human IL-17A Antibody	D0TB0028	2-8°C	31 OCT 2028
TURBO-BOOST Human IL-1 $\beta$ Antibody	D0TB0026	2-8°C	30 NOV 2028
TURBO-BOOST Human IL-2 Antibody	D0TB0024	2-8°C	31 OCT 2028
TURBO-BOOST Human IL-4 Antibody	D0TB0014	2-8°C	31 MAY 2028
TURBO-BOOST Human IL-6 Antibody	D0TB0020	2-8°C	31 OCT 2028
TURBO-BOOST Human TNF- $\alpha$ Antibody	D0TB0025	2-8°C	31 OCT 2028
Blocker S1	R0140018	$\leq -10^{\circ}\text{C}$	30 APR 2028
Diluent 58	M0580015	$\leq -10^{\circ}\text{C}$	31 JUL 2027
S-PLEX Enhance E1	Y0550026	$\leq -10^{\circ}\text{C}$	31 JAN 2028
S-PLEX Enhance E2	Y0560026	$\leq -10^{\circ}\text{C}$	29 FEB 2028
Blocker S2	562164	$\leq -10^{\circ}\text{C}$	31 OCT 2030
S-PLEX Detect D1 with TURBO-TAG <sup>TM</sup> Reagent	Y0580035	$\leq -70^{\circ}\text{C}$	31 JAN 2028
S-PLEX Detect D2	Y0590028	$\leq -70^{\circ}\text{C}$	29 FEB 2028
S-PLEX Enhance E3	Y0570028	$\leq -70^{\circ}\text{C}$	31 JAN 2028
S-PLEX Proinflammatory Panel 1 (human) Coating Solution	568801	$\leq -70^{\circ}\text{C}$	30 JUN 2029

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

**CERTIFICATE OF ANALYSIS****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-a	80% - 120%	< 20%
6	IL-2	80% - 120%	< 20%
8	IL-1 $\beta$	75% - 125%	< 30%
9	IFN- $\gamma$	80% - 120%	< 20%
10	IL-17A	80% - 120%	< 20%

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
IL-6	3.42 - 814	25.2	fg/mL	80% - 120%	< 25%	12.5	fg/mL
IL-10	12.5 - 2980	65.7	fg/mL	80% - 120%	< 25%	19.2	fg/mL
IL-12p70	39.2 - 9320	385	fg/mL	80% - 120%	< 25%	130	fg/mL
IL-4	5.11 - 1220	30.8	fg/mL	80% - 120%	< 25%	8.92	fg/mL
TNF-a	5.39 - 1280	68.2	fg/mL	80% - 120%	< 25%	14.0	fg/mL
IL-2	8.75 - 2080	57.8	fg/mL	80% - 120%	< 25%	18.9	fg/mL
IL-1 $\beta$	16.0 - 3810	94.9	fg/mL	75% - 125%	< 30%	37.8	fg/mL
IFN- $\gamma$	3.75 - 890	21.1	fg/mL	80% - 120%	< 25%	5.91	fg/mL
IL-17A	19.2 - 4570	237	fg/mL	80% - 120%	< 25%	37.9	fg/mL

**Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
IL-6	1X Calibrator	39600	fg/mL
IL-10	1X Calibrator	150000	fg/mL
IL-12p70	1X Calibrator	440000	fg/mL
IL-4	1X Calibrator	64000	fg/mL
TNF-a	1X Calibrator	53700	fg/mL
IL-2	1X Calibrator	75500	fg/mL
IL-1 $\beta$	1X Calibrator	137000	fg/mL
IFN- $\gamma$	1X Calibrator	43600	fg/mL
IL-17A	1X Calibrator	252000	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
<b>Review/Approval</b>	Karen Hamilla	Quality		11 JUN 2026