



RFE MEMO

Page 1 of 1

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

Catalog Number(s) / Size(s): K15396S-Series

Lot Number: K00S0100

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K00S0100 has been extended to 29 FEB 2024.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 58	M0580001	28 Feb 2023	29 Feb 2024
Blocker S1	R0010757	31 Jan 2024	29 Feb 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		08 NOV 2022

This memo supersedes all other statements regarding date of expiration for this product.

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Proinflammatory Panel 1 (human) Kit
Catalog Number: K15396S-Series
Lot Number: K00S0100
Expiration Date: 28 FEB 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
S-PLEX Proinflammatory Panel 1 (human) Calibrator Blend	A00S0065	2-8°C	30 JUN 2027
TURBO-BOOST™ Human IFN- γ Antibody	D00T0107	2-8°C	31 MAR 2026
TURBO-BOOST™ Human IL-10 Antibody	D00T0101	2-8°C	31 MAR 2026
TURBO-BOOST™ Human IL-12p70 Antibody	D00T0102	2-8°C	31 MAR 2026
TURBO-BOOST™ Human IL-17A Antibody	D00T0108	2-8°C	31 MAR 2026
TURBO-BOOST™ Human IL-1 β Antibody	D00T0106	2-8°C	31 MAR 2026
TURBO-BOOST™ Human IL-2 Antibody	D00T0104	2-8°C	31 MAR 2026
TURBO-BOOST™ Human IL-4 Antibody	D00T0103	2-8°C	31 MAR 2026
TURBO-BOOST™ Human IL-6 Antibody	D00T0100	2-8°C	31 MAR 2026
TURBO-BOOST™ Human TNF- α Antibody	D00T0105	2-8°C	31 MAR 2026
Blocker S1	R0010757	$\leq -10^{\circ}\text{C}$	31 JAN 2024
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	Y0550009	$\leq -10^{\circ}\text{C}$	28 FEB 2025
S-PLEX Enhance E2	Y0560011	$\leq -10^{\circ}\text{C}$	31 MAR 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580014	$\leq -70^{\circ}\text{C}$	28 FEB 2025
S-PLEX Detect D2	Y0590008	$\leq -70^{\circ}\text{C}$	31 JUL 2024
S-PLEX Enhance E3	Y0570009	$\leq -70^{\circ}\text{C}$	30 SEP 2024
S-PLEX Proinflammatory Panel 1(human) Coating Solution	426479	$\leq -70^{\circ}\text{C}$	31 MAR 2026

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	158	fg/mL	80% - 120%	< 25%	26.7	fg/mL
IL-10	12.5 - 2980	148	fg/mL	80% - 120%	< 25%	50.4	fg/mL
IL-12p70	39.2 - 9320	484	fg/mL	80% - 120%	< 25%	117	fg/mL
IL-4	5.11 - 1220	55.3	fg/mL	80% - 120%	< 25%	13.9	fg/mL
TNF- α	5.39 - 1280	107	fg/mL	80% - 120%	< 25%	15.1	fg/mL
IL-2	8.75 - 2080	85.8	fg/mL	80% - 120%	< 25%	12.9	fg/mL
IL-1 β	16.0 - 3810	318	fg/mL	75% - 125%	< 30%	110	fg/mL
IFN- γ	3.75 - 890	57.6	fg/mL	80% - 120%	< 25%	12.2	fg/mL
IL-17A	19.2 - 4570	210	fg/mL	80% - 120%	< 25%	98.5	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-6	1X Calibrator	39500	fg/mL
IL-10	1X Calibrator	148000	fg/mL
IL-12p70	1X Calibrator	484000	fg/mL
IL-4	1X Calibrator	55300	fg/mL
TNF- α	1X Calibrator	53400	fg/mL
IL-2	1X Calibrator	85800	fg/mL
IL-1 β	1X Calibrator	159000	fg/mL
IFN- γ	1X Calibrator	28800	fg/mL
IL-17A	1X Calibrator	210000	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