



Extension Memo

Page 1 of 1

Product Description: S-PLEX Human IL-1 β Kit
Catalog Number(s): K151ADSS-Series; K156ADSS-Series
Lot Number: K00S0163
New Expiration Date: 31 Aug 2026

Based on additional testing, the original date of expiration for this kit has been extended by 12 months.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 58	M0580011	31 Jul 2026	31 Aug 2026
S-PLEX Coating Reagent C1	417723	31 Aug 2025	31 Aug 2026
Blocker S1	R0140007	30 Jun 2026	31 Aug 2026
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580023	31 Jul 2026	31 Aug 2026
S-PLEX Detect D2	Y0590017	30 Jun 2026	31 Aug 2026
S-PLEX Enhance E1	Y0550016	31 Jul 2026	31 Aug 2026
S-PLEX Enhance E2	Y0560017	31 Jul 2026	31 Aug 2026
S-PLEX Enhance E3	Y0570018	30 Jun 2026	31 Aug 2026

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 Jul 2025



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IL-1 β Kit
Catalog Number: K151ADSS-Series; K156ADSS-Series
Lot Number: K00S0163
Expiration Date: 31 AUG 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-1 β Antibody	B00T0172	2-8°C	31 MAR 2027
Human IL-1 β Calibrator	A00S0090	2-8°C	31 OCT 2028
TURBO-BOOST™ Human IL-1 β Antibody	D00T0172	2-8°C	28 FEB 2027
Blocker S1	R0140007	≤ -10°C	30 JUN 2026
Diluent 58	M0580011	≤ -10°C	31 JUL 2026
S-PLEX Enhance E1	Y0550016	≤ -10°C	31 JUL 2026
S-PLEX Enhance E2	Y0560017	≤ -10°C	31 JUL 2026
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580023	≤ -70°C	31 JUL 2026
S-PLEX Detect D2	Y0590017	≤ -70°C	30 JUN 2026
S-PLEX Enhance E3	Y0570018	≤ -70°C	30 JUN 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-1 β	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-1 β	48.7 - 780	97.5	fg/mL	80% - 120%	< 20%	9.76	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-1 β	1X Calibrator	135000	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		03 APR 2024