

**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