



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IL-1 β Kit
Lot Number: K00S0214
Expiration Date: 31 JUL 2025
Catalog Number: K151ADSS-Series; K156ADSS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-1 β Antibody	B00T0148	2-8°C	31 AUG 2026
Human IL-1 β Calibrator	A00S0086	2-8°C	30 APR 2028
TURBO-BOOST Human IL-1 β Antibody	D00T0148	2-8°C	31 OCT 2026
Blocker S1	R0010831	$\leq -10^{\circ}\text{C}$	31 JUL 2025
Diluent 58	M0580012	$\leq -10^{\circ}\text{C}$	31 OCT 2026
S-PLEX Enhance E1	Y0550012	$\leq -10^{\circ}\text{C}$	31 AUG 2025
S-PLEX Enhance E2	Y0560013	$\leq -10^{\circ}\text{C}$	31 AUG 2025
S-PLEX Coating Reagent C1	417723	$\leq -70^{\circ}\text{C}$	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG TM Reagent	Y0580017	$\leq -70^{\circ}\text{C}$	31 AUG 2025
S-PLEX Detect D2	Y0590014	$\leq -70^{\circ}\text{C}$	31 OCT 2025
S-PLEX Enhance E3	Y0570013	$\leq -70^{\circ}\text{C}$	31 JUL 2025

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%	8.15	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-1 β	1X Calibrator	148000	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		01 JUL 2024