

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human IL-1 β Kit
Lot Number: K00S0251
Expiration Date: 31 OCT 2026
Catalog Number: K151ADSS-Series; K156ADSS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-1 β Antibody	B00T0221	2-8°C	31 OCT 2027
Human IL-1 β Calibrator	A00S0086	2-8°C	30 APR 2028
TURBO-BOOST Human IL-1 β Antibody	D00T0221	2-8°C	31 OCT 2027
Blocker S1	R0140009	$\leq -10^{\circ}\text{C}$	31 OCT 2026
Diluent 58	M0580012	$\leq -10^{\circ}\text{C}$	31 OCT 2026
S-PLEX Enhance E1	Y0550020	$\leq -10^{\circ}\text{C}$	30 APR 2027
S-PLEX Enhance E2	Y0560020	$\leq -10^{\circ}\text{C}$	31 MAR 2027
S-PLEX Coating Reagent C1	476750	$\leq -70^{\circ}\text{C}$	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG TM Reagent	Y0580029	$\leq -70^{\circ}\text{C}$	30 APR 2027
S-PLEX Detect D2	Y0590022	$\leq -70^{\circ}\text{C}$	30 APR 2027
S-PLEX Enhance E3	Y0570023	$\leq -70^{\circ}\text{C}$	30 APR 2027

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	780	fg/mL	80% - 120%	< 20%	11.8	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		23 JAN 2025