



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IL-1 β Kit
Lot Number: K00S0293
Expiration Date: 31 JUL 2027
Catalog Number: K151ADSS-Series, K156ADSS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-1 β Antibody	B00T0271	2-8°C	31 MAR 2029
Human IL-1 β Calibrator	A00S0129	2-8°C	31 MAR 2030
TURBO-BOOST Human IL-1 β Antibody	D00T0271	2-8°C	30 APR 2029
Blocker S1	R0140014	$\leq -10^{\circ}\text{C}$	30 SEP 2027
Diluent 58	M0580015	$\leq -10^{\circ}\text{C}$	31 JUL 2027
S-PLEX Enhance E1	Y0550025	$\leq -10^{\circ}\text{C}$	30 NOV 2027
S-PLEX Enhance E2	Y0560025	$\leq -10^{\circ}\text{C}$	29 FEB 2028
S-PLEX Coating Reagent C1	521329	$\leq -70^{\circ}\text{C}$	31 DEC 2027
S-PLEX Detect D1 with TURBO-TAG TM Reagent	Y0580035	$\leq -70^{\circ}\text{C}$	31 JAN 2028
S-PLEX Detect D2	Y0590027	$\leq -70^{\circ}\text{C}$	31 JAN 2028
S-PLEX Enhance E3	Y0570028	$\leq -70^{\circ}\text{C}$	31 JAN 2028

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	390	fg/mL	80% - 120%	< 20%	37.7	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-1 β	1X Calibrator	115000	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		17 DEC 2025