



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IL-1 β Kit
Lot Number: K00S0272
Expiration Date: 30 APR 2027
Catalog Number: K151ADSS-Series; K156ADSS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-1 β Antibody	B00T0253	2-8°C	31 AUG 2028
Human IL-1 β Calibrator	A00S0129	2-8°C	31 MAR 2030
TURBO-BOOST Human IL-1 β Antibody	D00T0253	2-8°C	31 AUG 2028
Blocker S1	R0140013	$\leq -10^{\circ}\text{C}$	31 JUL 2027
Diluent 58	M0580014	$\leq -10^{\circ}\text{C}$	31 MAY 2027
S-PLEX Enhance E1	Y0550022	$\leq -10^{\circ}\text{C}$	31 JUL 2027
S-PLEX Enhance E2	Y0560021	$\leq -10^{\circ}\text{C}$	30 APR 2027
S-PLEX Coating Reagent C1	521329	$\leq -70^{\circ}\text{C}$	31 DEC 2027
S-PLEX Detect D1 with TURBO-TAG TM Reagent	Y0580032	$\leq -70^{\circ}\text{C}$	31 AUG 2027
S-PLEX Detect D2	Y0590024	$\leq -70^{\circ}\text{C}$	31 AUG 2027
S-PLEX Enhance E3	Y0570025	$\leq -70^{\circ}\text{C}$	31 AUG 2027

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator	
Spot Number	Name	Recovery	CV
1	IL-1 β	80% - 120%	< 20%

LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-1 β	48.7 - 780	48.8	fg/mL	80% - 120%	< 20%	12.6	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		25 JUN 2025