



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IL-1 β Kit
Catalog Number: K151ADS-Series; K156ADS-Series
Lot Number: K00S0107
Expiration Date: 30 APR 2024

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human IL-1 β Antibody	D00T0112	2-8°C	31 MAR 2026
Biotin Human IL-1 β Antibody	B00T0112	2-8°C	31 MAR 2026
Human IL-1 β Calibrator	A00S0059	2-8°C	31 DEC 2026
Diluent 58	M0580007	$\leq -10^{\circ}\text{C}$	30 APR 2025
S-PLEX Coating Reagent C1	351987	$\leq -70^{\circ}\text{C}$	30 APR 2024
Blocker S1	R0010811	$\leq -10^{\circ}\text{C}$	28 FEB 2025
S-PLEX Enhance E1	Y0550009	$\leq -10^{\circ}\text{C}$	28 FEB 2025
S-PLEX Enhance E2	Y0560011	$\leq -10^{\circ}\text{C}$	31 MAR 2025
S-PLEX Enhance E3	Y0570011	$\leq -70^{\circ}\text{C}$	31 MAR 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580015	$\leq -70^{\circ}\text{C}$	30 APR 2025
S-PLEX Detect D2	Y0590011	$\leq -70^{\circ}\text{C}$	31 MAR 2025

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

Precision, Accuracy, and Sensitivity Specifications:

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

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-1 β	48.7 - 780	48.8	fg/mL	80% - 120%	< 20%	20.8	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-1 β	1X Calibrator	136000	fg/mL

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		09 JAN 2023