



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX® Human IL-1 $\beta$  Kit  
**Catalog Number:** K151ADSS-Series; K156ADSS - Series  
**Lot Number:** K00S0134  
**Expiration Date:** 31 Jul 2025

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-1 $\beta$ Antibody	B00T0148	2-8°C	31 Aug 2026
TURBO-BOOST™ Human IL-1 $\beta$ Antibody	D00T0148	2-8°C	31 Oct 2026
S-PLEX Coating Reagent C1	417723	$\leq -70^{\circ}\text{C}$	31 Aug 2025
Blocker S1	R0010831	$\leq -10^{\circ}\text{C}$	31 Jul 2025
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 Enhance E3	Y0570013	$\leq -70^{\circ}\text{C}$	31 Jul 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580017	$\leq -70^{\circ}\text{C}$	31 Aug 2025
S-PLEX Detect D2	Y0590014	$\leq -70^{\circ}\text{C}$	31 Oct 2025
Human IL-1 $\beta$ Calibrator	A00S0086	2-8°C	30 Apr 2028
Diluent 58	M0580009	$\leq -10^{\circ}\text{C}$	30 Nov 2025

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

### Precision, Accuracy, and Sensitivity Specifications:

Name	Calibrator		LLOQ				LLOD		
	Recovery	CV	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-1 $\beta$	80% - 120%	< 20%	48.7 - 780	97.5	fg/mL	80% - 120%	< 20%	18.9	fg/mL

### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-1 $\beta$	1X Calibrator	148000	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		25 JUL 2023