



## RFE MEMO

Page 1 of 1

**Product Description(s):** S-PLEX® Human IL-1 $\beta$  Kit


**Catalog Number(s) / Size(s):** K151ADSS-Series; K156ADSS-Series

**Lot Number:** K00S0072

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K00S0072 has been extended to 29 FEB 2024.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 58	M0580001	28 Feb 2023	29 Feb 2024
Blocker S1	R0010757	31 Jan 2024	29 Feb 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		17 NOV 2022

This memo supersedes all other statements regarding date of expiration for this product.



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human IL-1 $\beta$  Kit  
**Catalog Number:** K151ADSS-Series; K156ADSS-Series  
**Lot Number:** K00S0072  
**Expiration Date:** 28 FEB 2023

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST <sup>™</sup> Human IL-1 $\beta$ Antibody	D00T0080	2-8°C	31 MAY 2025
Biotin Human IL-1 $\beta$ Antibody	B00T0080	2-8°C	31 MAY 2025
Human IL-1 $\beta$ Calibrator	A00S0059	2-8°C	31 DEC 2026
Diluent 58	M0580001	$\leq -10^{\circ}\text{C}$	28 FEB 2023
S-PLEX Coating Reagent C1	351987	$\leq -70^{\circ}\text{C}$	30 APR 2024
Blocker S1	R0010757	$\leq -10^{\circ}\text{C}$	31 JAN 2024
S-PLEX Enhance E1	Y0550007	$\leq -10^{\circ}\text{C}$	31 AUG 2024
S-PLEX Enhance E2	Y0560010	$\leq -10^{\circ}\text{C}$	31 JAN 2025
S-PLEX Enhance E3	Y0570009	$\leq -70^{\circ}\text{C}$	30 SEP 2024
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> reagent	Y0580012	$\leq -70^{\circ}\text{C}$	31 OCT 2024
S-PLEX Detect D2	Y0590008	$\leq -70^{\circ}\text{C}$	31 JUL 2024

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 $\beta$	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-1 $\beta$	48.7 - 780	48.8	fg/mL	80% - 120%	< 20%	18.2	fg/mL

### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-1 $\beta$	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		11 OCT 2022