



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX® Human IL-1 β Kit
Catalog Number: K151ADS-Series; K156ADS-Series
Lot Number: K00S0040
Expiration Date: 31 JAN 2023

Precision, Accuracy, and Sensitivity Specifications:

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

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IL-1 β	48.8 - 780	97.5	fg/mL	80% - 120%	< 20%

LLOD		
Name	Observed	Units
IL-1 β	19.3	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-1 β	1X Calibrator	145000	fg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-1 β Detection Antibody	D00T0041A	2-8°C	30 Apr 2024
Biotin Human IL-1 β Capture Antibody	B00T0041A	2-8°C	30 Apr 2024
Human IL-1 β Calibrator	A00S0038	2-8°C	30 Sep 2025
Diluent 58	M0580001	$\leq -10^{\circ}\text{C}$	28 Feb 2023
S-PLEX Coating Reagent C1	329998	$\leq -70^{\circ}\text{C}$	31 Aug 2023
Blocker S1	R0140004	$\leq -10^{\circ}\text{C}$	30 Apr 2023
S-PLEX Enhance E1	Y0550004	$\leq -10^{\circ}\text{C}$	28 Feb 2023
S-PLEX Enhance E2	Y0560004	$\leq -10^{\circ}\text{C}$	31 Jan 2023
S-PLEX Enhance E3	Y0570004	$\leq -70^{\circ}\text{C}$	31 Jan 2023
S-PLEX Detect D1	Y0580006	$\leq -70^{\circ}\text{C}$	31 Mar 2023
S-PLEX Detect D2	Y0590004	$\leq -70^{\circ}\text{C}$	31 Jan 2023

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		22 DEC 2020