



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX® Human IL-3 Kit

Catalog Number: K151ADQS-Series

Lot Number: K00S0037

Expiration Date: 31 Jan 2023

Precision, Accuracy, and Sensitivity Specifications:

Calibrator		
Name	Recovery	CV
IL-3	80% - 120%	< 20%

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IL-3	1.33 – 21.2	21.2	fg/mL	80% - 120%	< 20%

LLOD		
Name	Observed	Units
IL-3	4.56	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-3	1X Calibrator	38000	fg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-3 Antibody	D00T0043A	2-8°C	31 May 2024
Biotin Human IL-3 Antibody	B00T0043A	2-8°C	30 Apr 2024
Human IL-3 Calibrator	A00S0033	2-8°C	31 Aug 2025
Diluent 58	M0580001	≤ -10°C	28 Feb 2023
S-PLEX Coating Reagent C1	329998	≤ -70°C	31 Aug 2023
Blocker S1	R0140004	≤ -10°C	30 Apr 2023
S-PLEX Enhance E1	Y0550004	≤ -10°C	28 Feb 2023
S-PLEX Enhance E2	Y0560004	≤ -10°C	31 Jan 2023
S-PLEX Enhance E3	Y0570004	≤ -70°C	31 Jan 2023
S-PLEX Detect D1	Y0580006	≤ -70°C	31 Mar 2023
S-PLEX Detect D2	Y0590004	≤ -70°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 FEB 2021