



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX® Human IL-5 Kit
Catalog Number: K151J3S-Series; K156J3S-Series
Lot Number: K00S0141
Expiration Date: 31 Mar 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-5 Antibody	B00T0150	2-8°C	30 Sep 2026
TURBO-BOOST™ Human IL-5 Antibody	D00T0150	2-8°C	31 Oct 2026
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 Aug 2025
Human IL-5 Calibrator	A00S0084	2-8°C	31 Mar 2028
Diluent 43	M0430050	≤ -10°C	31 Mar 2025
Diluent 3	M0030122	≤ -10°C	30 Nov 2025
Blocker S1	R0010831	≤ -10°C	31 Jul 2025
S-PLEX Enhance E1	Y0550012	≤ -10°C	31 Aug 2025
S-PLEX Enhance E2	Y0560013	≤ -10°C	31 Aug 2025
S-PLEX Enhance E3	Y0570013	≤ -70°C	31 Jul 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580017	≤ -70°C	31 Aug 2025
S-PLEX Detect D2	Y0590013	≤ -70°C	31 Jul 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-5	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-5	3.00 – 48.0	12.0	fg/mL	80% - 120%	< 20%	1.90	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-5	30X Calibrator	1080000	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		14 JUL 2023