



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX[®] Human IL-5 Kit
Lot Number: K00S0223
Expiration Date: 31 AUG 2026
Catalog Number: K151J3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-5 Antibody	B00T0230	2-8°C	31 DEC 2027
Human IL-5 Calibrator	A00S0092	2-8°C	31 OCT 2028
TURBO-BOOST Human IL-5 Antibody	D00T0230	2-8°C	31 JAN 2028
Blocker S1	R0140009	≤ -10°C	31 OCT 2026
Diluent 3	M0030129	≤ -10°C	30 NOV 2026
Diluent 43	M0430056	≤ -10°C	30 SEP 2026
S-PLEX Enhance E1	Y0550017	≤ -10°C	31 AUG 2026
S-PLEX Enhance E2	Y0560018	≤ -10°C	31 AUG 2026
S-PLEX Coating Reagent C1	476750	≤ -70°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG [™] Reagent	Y0580025	≤ -70°C	30 SEP 2026
S-PLEX Detect D2	Y0590019	≤ -70°C	30 SEP 2026
S-PLEX Enhance E3	Y0570020	≤ -70°C	31 OCT 2026

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

Precision, Accuracy, and Sensitivity Specifications:

Spot Number	Name	Calibrator	
		Recovery	CV
1	IL-5	80% - 120%	< 20%

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
IL-5	3.00 - 48.0	6.00	fg/mL	80% - 120%	< 20%	1.95	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-5	30X Calibrator	965000	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		19 AUG 2024