

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX Human IL-5 Kit  
**Catalog Number:** K151J3S-Series; K156J3S-Series  
**Lot Number:** K00S0166  
**Expiration Date:** 31 AUG 2025

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-5 Antibody	B00T0175	2-8°C	28 FEB 2027
Human IL-5 Calibrator	A00S0092	2-8°C	31 OCT 2028
TURBO-BOOST™ Human IL-5 Antibody	D00T0175	2-8°C	31 MAR 2027
Blocker S1	R0140007	≤ -10°C	30 JUN 2026
Diluent 3	M0030123	≤ -10°C	31 JAN 2026
Diluent 43	M0430053	≤ -10°C	31 JAN 2026
S-PLEX Enhance E1	Y0550016	≤ -10°C	31 JUL 2026
S-PLEX Enhance E2	Y0560017	≤ -10°C	31 JUL 2026
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580023	≤ -70°C	31 JUL 2026
S-PLEX Detect D2	Y0590017	≤ -70°C	30 JUN 2026
S-PLEX Enhance E3	Y0570018	≤ -70°C	30 JUN 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%	2.18	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
<b>Review/Approval</b>	Karen Hamilla	Quality		03 APR 2024