



## CERTIFICATE OF ANALYSIS

**Product Description:** S-PLEX Human IL-5 Kit  
**Lot Number:** K00S0229  
**Expiration Date:** 30 SEP 2026  
**Catalog Number:** K151J3S-Series; K156J3S-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-5 Antibody	B00T0232	2-8°C	31 MAR 2028
Human IL-5 Calibrator	A00S0057	2-8°C	30 NOV 2026
TURBO-BOOST Human IL-5 Antibody	D00T0232	2-8°C	31 MAR 2028
Blocker S1	R0140011	≤ -10°C	28 FEB 2027
Diluent 3	M0030130	≤ -10°C	31 JAN 2027
Diluent 43	M0430056	≤ -10°C	30 SEP 2026
S-PLEX Enhance E1	Y0550018	≤ -10°C	30 NOV 2026
S-PLEX Enhance E2	Y0560019	≤ -10°C	30 NOV 2026
S-PLEX Coating Reagent C1	516221	≤ -70°C	31 OCT 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580027	≤ -70°C	31 OCT 2026
S-PLEX Detect D2	Y0590020	≤ -70°C	30 NOV 2026
S-PLEX Enhance E3	Y0570021	≤ -70°C	30 NOV 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	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
IL-5	3.00 - 48.0	6.00	fg/mL	80% - 120%	< 20%	2.19	fg/mL

**Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
IL-5	30X Calibrator	909000	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		10 DEC 2024