

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX<sup>®</sup> Human IL-13 Kit  
**Catalog Number:** K151Y9S-Series  
**Lot Number:** K00S0175  
**Expiration Date:** 28 FEB 2025

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-13 Antibody	B00T0111	2-8°C	31 MAR 2026
Human IL-13 Calibrator	A00S0076	2-8°C	31 OCT 2027
TURBO-BOOST™ Human IL-13 Antibody	D00T0111	2-8°C	31 MAY 2026
Blocker S1	R0010811	≤ -10°C	28 FEB 2025
Diluent 2	M0020124	≤ -10°C	28 FEB 2025
S-PLEX Enhance E1	Y0550010	≤ -10°C	31 MAY 2025
S-PLEX Enhance E2	Y0560011	≤ -10°C	31 MAR 2025
S-PLEX Coating Reagent C1	446866	≤ -70°C	31 MAR 2026
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580015	≤ -70°C	30 APR 2025
S-PLEX Detect D2	Y0590011	≤ -70°C	31 MAR 2025
S-PLEX Enhance E3	Y0570011	≤ -70°C	31 MAR 2025

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-13	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-13	5.86 - 93.8	11.7	fg/mL	80% - 120%	< 20%	10.3	fg/mL

**Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
IL-13	1X Calibrator	126000	fg/mL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Michelle Likiam	Quality		11 DEC 2023