

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX Human IL-9 Kit  
**Lot Number:** K00S0210  
**Expiration Date:** 30 SEP 2026  
**Catalog Number:** K151R3S-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-9 Antibody	B00T0223	2-8°C	31 JAN 2028
Human IL-9 Calibrator	A00S0100	2-8°C	31 DEC 2028
TURBO-BOOST Human IL-9 Antibody	D00T0223	2-8°C	31 OCT 2027
Blocker S1	R0140010	≤ -10°C	31 DEC 2026
Diluent 3	M0030129	≤ -10°C	30 NOV 2026
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	476750	≤ -70°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580027	≤ -70°C	31 OCT 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-9	75% - 125%	< 25%

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
IL-9	9.75 - 156	19.5	fg/mL	75% - 125%	< 25%	12.1	fg/mL

**Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
IL-9	15X Calibrator	1550000	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		30 OCT 2024