

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human IL-9 Kit
Catalog Number: K151R3S-Series
Lot Number: K00S0177
Expiration Date: 31 AUG 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-9 Antibody	B00T0182	2-8°C	31 JUL 2027
Human IL-9 Calibrator	A00S0100	2-8°C	31 DEC 2028
TURBO-BOOST™ Human IL-9 Antibody	D00T0182	2-8°C	31 JUL 2027
Blocker S1	R0140007	≤ -10°C	30 JUN 2026
Diluent 3	M0030125	≤ -10°C	31 MAY 2026
Diluent 43	M0430055	≤ -10°C	30 APR 2026
S-PLEX Enhance E1	Y0550015	≤ -10°C	31 MAY 2026
S-PLEX Enhance E2	Y0560016	≤ -10°C	31 MAY 2026
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580022	≤ -70°C	31 MAY 2026
S-PLEX Detect D2	Y0590017	≤ -70°C	30 JUN 2026
S-PLEX Enhance E3	Y0570017	≤ -70°C	31 MAR 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	39.0	fg/mL	75% - 125%	< 25%	28.4	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
Review/Approval	Karen Hamilla	Quality		05 MAR 2024