



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®]Human IL-4 Kit

Catalog Number: K151A3S - Series

Lot Number: K00S0140

Expiration Date: 31 Jul 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-4 Antibody	B00T0149	2-8°C	30 Sep 2026
TURBO-BOOST [™] Human IL-4 Antibody	D00T0149	2-8°C	30 Sep 2026
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 Aug 2025
Blocker S1	R0010831	≤ -10°C	31 Jul 2025
S-PLEX Enhance E1	Y0550011	≤ -10°C	31 Jul 2025
S-PLEX Enhance E2	Y0560013	≤ -10°C	31 Aug 2025
S-PLEX Enhance E3	Y0570013	≤ -70°C	31 Jul 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580017	≤ -70°C	31 Aug 2025
S-PLEX Detect D2	Y0590013	≤ -70°C	31 Jul 2025
Human IL-4 Calibrator	A00S0083	2-8°C	31 Mar 2028
Diluent 2	M0020126	≤ -10°C	31 Aug 2025
Diluent 3	M0030120	≤ -10°C	31 Jul 2025

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

Name	Calibrator	
	Recovery	CV
IL-4	80% - 120%	< 20%

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
IL-4	0.306 - 4.90	1.23	fg/mL	80% - 120%	< 20%	0.387	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-4	20X Calibrator	117000	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		28 JUL 2023