



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX[®] Human IL-4 Kit
Lot Number: K00S0094
Expiration Date: 31 AUG 2023
Catalog Number: K151A3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-4 Capture Antibody	B00T0094	2-8°C	28 FEB 2026
Human IL-4 Calibrator	A00S0041	2-8°C	30 NOV 2025
TURBO-BOOST [™] Human IL-4 Detection Antibody	D00T0094	2-8°C	31 JAN 2026
Blocker S1	R0010757	≤ -10°C	31 JAN 2024
Diluent 2	M0020122	≤ -10°C	31 OCT 2024
Diluent 3	M0030114	≤ -10°C	30 SEP 2024
S-PLEX Enhance E1	Y0550009	≤ -10°C	28 FEB 2025
S-PLEX Enhance E2	Y0560011	≤ -10°C	31 MAR 2025
S-PLEX Coating Reagent C1	329998	≤ -70°C	31 AUG 2023
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580014	≤ -70°C	28 FEB 2025
S-PLEX Detect D2	Y0590010	≤ -70°C	28 FEB 2025
S-PLEX Enhance E3	Y0570008	≤ -70°C	30 JUN 2024

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

Precision, Accuracy, and Sensitivity Specifications:

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

LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-4	0.306 - 4.90	4.90	fg/mL	80% - 120%	< 20%	0.670	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-4	20X Calibrator	114000	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		02 SEP 2025