



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IL-4 Kit
Lot Number: K00S0300
Expiration Date: 30 JUN 2027
Catalog Number: K151A3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-4 Antibody	B00T0276	2-8°C	31 MAY 2029
Human IL-4 Calibrator	A00S0111	2-8°C	31 JAN 2030
TURBO-BOOST Human IL-4 Antibody	D00T0276	2-8°C	30 JUN 2029
Blocker S1	R0140016	≤ -10°C	31 DEC 2027
Diluent 2	M0020133	≤ -10°C	30 JUN 2027
Diluent 3	M0030145	≤ -10°C	29 FEB 2028
S-PLEX Enhance E1	Y0550026	≤ -10°C	31 JAN 2028
S-PLEX Enhance E2	Y0560025	≤ -10°C	29 FEB 2028
S-PLEX Coating Reagent C1	521330	≤ -70°C	30 JUN 2028
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580035	≤ -70°C	31 JAN 2028
S-PLEX Detect D2	Y0590027	≤ -70°C	31 JAN 2028
S-PLEX Enhance E3	Y0570028	≤ -70°C	31 JAN 2028

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%	0.948	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-4	20X Calibrator	125000	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		22 JAN 2026