



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K151A3S - Series

Lot Number: K00S0165

Expiration Date: 31 JAN 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-4 Antibody	B00T0174	2-8°C	31 JAN 2027
TURBO-BOOST™ Human IL-4 Antibody	D00T0174	2-8°C	28 FEB 2027
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
Blocker S1	R0010871	≤ -10°C	31 MAR 2026
S-PLEX Enhance E1	Y0550014	≤ -10°C	28 FEB 2026
S-PLEX Enhance E2	Y0560010	≤ -10°C	31 JAN 2025
S-PLEX Enhance E3	Y0570015	≤ -70°C	30 NOV 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580018	≤ -70°C	31 OCT 2025
S-PLEX Detect D2	Y0590015	≤ -70°C	31 JAN 2026
Human IL-4 Calibrator	A00S0083	2-8°C	31 MAR 2028
Diluent 2	M0020127	≤ -10°C	30 NOV 2025
Diluent 3	M0030123	≤ -10°C	31 JAN 2026

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	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
IL-4	0.306 - 4.90	4.90	fg/mL	80% - 120%	< 20%	0.709	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		15 NOV 2023