



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX® Human IL-4 Kit
Lot Number: K00S0254
Expiration Date: 30 NOV 2025
Catalog Number: K151A3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-4 Antibody	B00T0187	2-8°C	31 JUL 2027
Human IL-4 Calibrator	A00S0083	2-8°C	31 MAR 2028
TURBO-BOOST Human IL-4 Antibody	D00T0187	2-8°C	30 JUN 2027
S-PLEX Coating Reagent C1	476750	≤ -70°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580030	≤ -70°C	30 APR 2027
S-PLEX Detect D2	Y0590022	≤ -70°C	30 APR 2027
S-PLEX Enhance E3	Y0570023	≤ -70°C	30 APR 2027
Blocker S1	R0140009	≤ -10°C	31 OCT 2026
Diluent 2	M0020127	≤ -10°C	30 NOV 2025
Diluent 3	M0030128	≤ -10°C	31 OCT 2026
S-PLEX Enhance E1	Y0550021	≤ -10°C	30 JUN 2027
S-PLEX Enhance E2	Y0560021	≤ -10°C	30 APR 2027

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	2.45	fg/mL	80% - 120%	< 20%	0.720	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-4	20X Calibrator	117000	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 FEB 2025