

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human IL-17A Kit
Lot Number: K00S0248
Expiration Date: 30 APR 2026
Catalog Number: K151C3S-Series; K156C3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-17A Antibody	B00T0220	2-8°C	30 NOV 2027
Human IL-17A Calibrator	A00S0110	2-8°C	31 JUL 2029
TURBO-BOOST Human IL-17A Antibody	D00T0220	2-8°C	30 NOV 2027
S-PLEX Coating Reagent C1	476750	≤ -70°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580029	≤ -70°C	30 APR 2027
S-PLEX Detect D2	Y0590021	≤ -70°C	31 MAR 2027
S-PLEX Enhance E3	Y0570022	≤ -70°C	30 APR 2027
Blocker S1	R0140009	≤ -10°C	31 OCT 2026
Diluent 3	M0030128	≤ -10°C	31 OCT 2026
Diluent 43	M0430055	≤ -10°C	30 APR 2026
S-PLEX Enhance E1	Y0550020	≤ -10°C	30 APR 2027
S-PLEX Enhance E2	Y0560020	≤ -10°C	31 MAR 2027

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

Precision, Accuracy, and Sensitivity Specifications:

Spot Number	Name	Calibrator	
		Recovery	CV
1	IL-17A	75% - 125%	< 25%

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
IL-17A	7.50 - 120	60.0	fg/mL	75% - 125%	< 25%	26.3	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-17A	10X Calibrator	3120000	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		26 DEC 2024