



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IL-17A Kit
Lot Number: K00S0264
Expiration Date: 28 FEB 2027
Catalog Number: K151C3S-Series; K156C3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-17A Antibody	B00T0249	2-8°C	31 AUG 2028
Human IL-17A Calibrator	A00S0110	2-8°C	31 JUL 2029
TURBO-BOOST Human IL-17A Antibody	D00T0249	2-8°C	31 AUG 2028
Blocker S1	R0140013	≤ -10°C	31 JUL 2027
Diluent 3	M0030138	≤ -10°C	31 AUG 2027
Diluent 43	M0430058	≤ -10°C	28 FEB 2027
S-PLEX Enhance E1	Y0550022	≤ -10°C	31 JUL 2027
S-PLEX Enhance E2	Y0560021	≤ -10°C	30 APR 2027
S-PLEX Coating Reagent C1	521329	≤ -70°C	31 DEC 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580032	≤ -70°C	31 AUG 2027
S-PLEX Detect D2	Y0590024	≤ -70°C	31 AUG 2027
S-PLEX Enhance E3	Y0570025	≤ -70°C	31 AUG 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-17A	75% - 125%	< 25%

LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-17A	7.50 - 120	30.0	fg/mL	75% - 125%	< 25%	18.4	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		12 JUN 2025