



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IL-17A Kit
Catalog Number: K151C3S-Series; K156C3S-Series
Lot Number: K00S0162
Expiration Date: 31 AUG 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human IL-17A Antibody	D00T0171	2-8°C	31 MAR 2027
Biotin Human IL-17A Antibody	B00T0171	2-8°C	31 JAN 2027
Human IL-17A Calibrator	A00S0045	2-8°C	31 DEC 2025
Diluent 43	M0430053	≤ -10°C	31 JAN 2026
Diluent 3	M0030123	≤ -10°C	31 JAN 2026
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
Blocker S1	R0010832	≤ -10°C	31 OCT 2025
S-PLEX Enhance E1	Y0550013	≤ -10°C	30 NOV 2025
S-PLEX Enhance E2	Y0560015	≤ -10°C	28 FEB 2026
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

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-17A	75% - 125%	< 25%

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-17A	7.50 - 120	30.0	fg/mL	75% - 125%	< 25%	14.8	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-17A	10X Calibrator	2400000	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		31 OCT 2023