

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Human IL-21 Kit
Lot Number: K00S0192
Expiration Date: 31 MAR 2026
Catalog Number: K151ADNS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-21 Antibody	B00T0201	2-8°C	31 OCT 2027
Human IL-21 Calibrator	A00S0072	2-8°C	31 AUG 2027
TURBO-BOOST Human IL-21 Antibody	D00T0201	2-8°C	31 OCT 2027
Blocker S1	R0140009	≤ -10°C	31 OCT 2026
Diluent 2	M0020128	≤ -10°C	31 MAR 2026
S-PLEX Enhance E1	Y0550017	≤ -10°C	31 AUG 2026
S-PLEX Enhance E2	Y0560018	≤ -10°C	31 AUG 2026
S-PLEX Coating Reagent C1	476750	≤ -70°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG [™] Reagent	Y0580025	≤ -70°C	30 SEP 2026
S-PLEX Detect D2	Y0590019	≤ -70°C	30 SEP 2026
S-PLEX Enhance E3	Y0570020	≤ -70°C	31 OCT 2026

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-21	80% - 120%	< 20%

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
IL-21	18.8 - 300	37.5	fg/mL	80% - 120%	< 20%	9.45	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-21	1X Calibrator	86700	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		25 JUL 2024