

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX® Human IL-22 Kit
Lot Number: K00S0222
Expiration Date: 30 APR 2026
Catalog Number: K151H3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-22 Antibody	B00T0229	2-8°C	31 JAN 2028
Human IL-22 Calibrator	A00S0119	2-8°C	31 OCT 2029
TURBO-BOOST Human IL-22 Antibody	D00T0229	2-8°C	31 DEC 2027
S-PLEX Coating Reagent C1	476750	≤ -70°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580028	≤ -70°C	31 JAN 2027
S-PLEX Detect D2	Y0590021	≤ -70°C	31 MAR 2027
S-PLEX Enhance E3	Y0570021	≤ -70°C	30 NOV 2026
Blocker S1	R0140008	≤ -10°C	31 JUL 2026
Diluent 3	M0030128	≤ -10°C	31 OCT 2026
Diluent 43	M0430055	≤ -10°C	30 APR 2026
S-PLEX Enhance E1	Y0550019	≤ -10°C	31 MAR 2027
S-PLEX Enhance E2	Y0560019	≤ -10°C	30 NOV 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-22	80% - 120%	< 20%

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
IL-22	3.75 - 60.0	15.0	fg/mL	80% - 120%	< 20%	4.04	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-22	15X Calibrator	605000	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 DEC 2024