

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Human IL-22 Kit
Catalog Number: K151H3S-Series
Lot Number: K00S0193
Expiration Date: 31 JUL 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-22 Antibody	B00T0202	2-8°C	31 AUG 2027
Human IL-22 Calibrator	A00S0025	2-8°C	31 JUL 2025
TURBO-BOOST™ Human IL-22 Antibody	D00T0202	2-8°C	31 AUG 2027
Blocker S1	R0140007	≤ -10°C	30 JUN 2026
Diluent 3	M0030126	≤ -10°C	31 JUL 2026
Diluent 43	M0430055	≤ -10°C	30 APR 2026
S-PLEX Enhance E1	Y0550016	≤ -10°C	31 JUL 2026
S-PLEX Enhance E2	Y0560017	≤ -10°C	31 JUL 2026
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580023	≤ -70°C	31 JUL 2026
S-PLEX Detect D2	Y0590017	≤ -70°C	30 JUN 2026
S-PLEX Enhance E3	Y0570018	≤ -70°C	30 JUN 2026

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-22	80% - 120%	< 20%

		LLOQ				LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-22	3.75 - 60.0	7.50	fg/mL	80% - 120%	< 20%	3.11	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-22	15X Calibrator	537000	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		10 APR 2024