

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX<sup>®</sup> Human IL-22 Kit  
**Catalog Number:** K151H3S-Series  
**Lot Number:** K00S0174  
**Expiration Date:** 31 OCT 2024

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-22 Antibody	B00T0181	2-8°C	31 MAY 2027
Human IL-22 Calibrator	A00S0025	2-8°C	31 JUL 2025
TURBO-BOOST <sup>™</sup> Human IL-22 Antibody	D00T0181	2-8°C	31 MAY 2027
Blocker S1	R0010871	≤ -10°C	31 MAR 2026
Diluent 3	M0030117	≤ -10°C	31 JAN 2025
Diluent 43	M0430049	≤ -10°C	31 OCT 2024
S-PLEX Enhance E1	Y0550015	≤ -10°C	31 MAY 2026
S-PLEX Enhance E2	Y0560016	≤ -10°C	31 MAY 2026
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> reagent	Y0580018	≤ -70°C	31 OCT 2025
S-PLEX Detect D2	Y0590014	≤ -70°C	31 OCT 2025
S-PLEX Enhance E3	Y0570014	≤ -70°C	31 OCT 2025

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	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-22	3.75 - 60.0	15.0	fg/mL	80% - 120%	< 20%	3.14	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
<b>Review/Approval</b>	Karen Hamilla	Quality		22 DEC 2023