



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IL-13 Kit

Catalog Number: K151Y9S-Series

Lot Number: K00S0092

Expiration Date: 31 OCT 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human IL-13 Antibody	D00T0092	2-8°C	31 DEC 2025
Biotin Human IL-13 Antibody	B00T0092	2-8°C	28 FEB 2026
Human IL-13 Calibrator	A00S0027	2-8°C	31 AUG 2025
Diluent 2	M0020115	≤ -10°C	31 OCT 2023
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 APR 2024
Blocker S1	R010768	≤ -10°C	31 MAY 2024
S-PLEX Enhance E1	Y0550008	≤ -10°C	30 NOV 2024
S-PLEX Enhance E2	Y0560010	≤ -10°C	31 JAN 2025
S-PLEX Enhance E3	Y0570010	≤ -70°C	31 JAN 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580013	≤ -70°C	31 DEC 2024
S-PLEX Detect D2	Y0590010	≤ -70°C	29 FEB 2025

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-13	80% - 120%	< 20%

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-13	5.86 - 93.8	23.5	fg/mL	80% - 120%	< 20%	8.43	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-13	1X Calibrator	128000	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		10 OCT 2022