



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human IL-5 Kit  
**Catalog Number:** K151J3S-Series; K156J3S-Series  
**Lot Number:** K00S0073  
**Expiration Date:** 31 JAN 2024

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST <sup>™</sup> Human IL-5 Antibody	D00T0081	2-8°C	31 OCT 2025
Biotin Human IL-5 Antibody	B00T0081	2-8°C	31 OCT 2025
Human IL-5 Calibrator	A00S0057	2-8°C	30 NOV 2026
Diluent 43	M0430047	≤ -10°C	31 MAY 2024
Diluent 3	M0030116	≤ -10°C	30 NOV 2024
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 APR 2024
Blocker S1	R0010757	≤ -10°C	31 JAN 2024
S-PLEX Enhance E1	Y0550007	≤ -10°C	31 AUG 2024
S-PLEX Enhance E2	Y0560010	≤ -10°C	31 JAN 2025
S-PLEX Enhance E3	Y0570009	≤ -70°C	30 SEP 2024
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> Reagent	Y0580013	≤ -70°C	31 DEC 2024
S-PLEX Detect D2	Y0590009	≤ -70°C	30 NOV 2024

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-5	80% - 120%	< 20%

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-5	3.00 - 48.0	6.00	fg/mL	80% - 120%	< 20%	1.21	fg/mL

### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-5	30X Calibrator	909000	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		30 SEP 2022