



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IL-5 Kit
Catalog Number: K151J3S-Series; K156J3S-Series
Lot Number: K00S0053
Expiration Date: 30 APR 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-5 Antibody	D00T0062	2-8°C	31 JUL 2025
Biotin Human IL-5 Antibody	B00T0062	2-8°C	31 JUL 2025
Human IL-5 Calibrator	A00S0054	2-8°C	31 JUL 2026
Diluent 43	M0430046	≤ -10°C	29 FEB 2024
Diluent 3	M0030111	≤ -10°C	31 MAY 2024
S-PLEX Coating Reagent C1	329998	≤ -70°C	31 AUG 2023
Blocker S1	R0140004	≤ -10°C	30 APR 2023
S-PLEX Enhance E1	R0140006	≤ -10°C	31 JAN 2024
S-PLEX Enhance E2	Y0560008	≤ -10°C	30 APR 2024
S-PLEX Enhance E3	Y0570007	≤ -70°C	29 FEB 2024
S-PLEX Detect D1	Y0580009	≤ -70°C	29 FEB 2024
S-PLEX Detect D2	Y0590007	≤ -70°C	29 FEB 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%	1.57	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-5	30X Calibrator	1040000	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		16 MAR 2022