



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IL-5 Kit
Lot Number: K00S0274
Expiration Date: 28 FEB 2027
Catalog Number: K151J3S-Series, K156J3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-5 Antibody	B00T0255	2-8°C	31 OCT 2028
Human IL-5 Calibrator	A00S0131	2-8°C	31 MAR 2030
TURBO-BOOST Human IL-5 Antibody	D00T0255	2-8°C	30 NOV 2028
Blocker S1	R0140016	≤ -10°C	31 DEC 2027
Diluent 3	M0030143	≤ -10°C	31 JAN 2028
Diluent 43	M0430058	≤ -10°C	28 FEB 2027
S-PLEX Enhance E1	Y0550024	≤ -10°C	31 OCT 2027
S-PLEX Enhance E2	Y0560024	≤ -10°C	31 DEC 2027
S-PLEX Coating Reagent C1	521330	≤ -70°C	30 JUN 2028
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580034	≤ -70°C	31 DEC 2027
S-PLEX Detect D2	Y0590026	≤ -70°C	31 DEC 2027
S-PLEX Enhance E3	Y0570027	≤ -70°C	31 DEC 2027

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-5	80% - 120%	< 20%

LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-5	3.00 - 48.0	6.00	fg/mL	80% - 120%	< 20%	1.73	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-5	30X Calibrator	964000	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		31 OCT 2025