



EXTENSION MEMO


Page 1 of 1

Product Description(s): S-PLEX® Human IL-5 Kit
Catalog Number(s) / Size(s): K151J3S-Series; K156J3S-Series
Lot Number: K00S0108

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K00S0108 has been extended to 30 APR 2025.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 43	M0430050	31 MAR 2025	30 APR 2025
S-PLEX Coating Reagent C1	351987	30 APR 2024	30 APR 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		07 JUN 2024

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX® Human IL-5 Kit
Catalog Number: K151J3S-Series; K156J3S-Series
Lot Number: K00S0108
Expiration Date: 30 Apr 2024

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-5 Antibody	D00T0113	2-8°C	30 Jun 2026
Biotin Human IL-5 Antibody	B00T0113	2-8°C	31 Mar 2026
Human IL-5 Calibrator	A00S0057	2-8°C	30 Nov 2026
Diluent 43	M0430050	≤ -10°C	31 Mar 2025
Diluent 3	M0030119	≤ -10°C	30 Apr 2025
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 Apr 2024
Blocker S1	R0010824	≤ -10°C	30 Jun 2025
S-PLEX Enhance E1	Y0550011	≤ -10°C	31 Jul 2025
S-PLEX Enhance E2	Y0560012	≤ -10°C	31 May 2025
S-PLEX Enhance E3	Y0570012	≤ -70°C	31 May 2025
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580016	≤ -70°C	31 May 2025
S-PLEX Detect D2	Y0590012	≤ -70°C	30 Apr 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-5	80% - 120%	< 20%

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
IL-5	3.00 – 48.0	3.00	fg/mL	80% - 120%	< 20%	1.68	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		08 MAR 2023