



Extension Memo

Page 1 of 1

Product Description: S-PLEX Human IL-3 Kit
Catalog Number(s): K151ADQS-Series
Lot Number: K00S0235
New Expiration Date: 30 Apr 2025

Based on additional testing, the original date of expiration for this kit has been extended by 12 months.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Biotin Human IL-3 Antibody	B00T0043	30 Apr 2024	30 Apr 2025
TURBO-TAG Human IL-3 Antibody	D00T0043	31 May 2024	30 Apr 2025

Function	Name	Signature	Date
Quality	Michelle Likiam		27 Nov 2024

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human IL-3 Kit
Lot Number: K00S0235
Expiration Date: 30 APR 2024
Catalog Number: K151ADQS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-3 Calibrator	A00S0033	2-8°C	31 AUG 2025
Biotin Human IL-3 Antibody	B00T0043	2-8°C	30 APR 2024
TURBO-TAG Human IL-3 Antibody	D00T0043	2-8°C	31 MAY 2024
Blocker S1	R0140010	≤ -10°C	31 DEC 2026
Diluent 58	M0580012	≤ -10°C	31 OCT 2026
S-PLEX Enhance E1	Y0550019	≤ -10°C	31 MAR 2027
S-PLEX Enhance E2	Y0560019	≤ -10°C	30 NOV 2026
S-PLEX Coating Reagent C1	476750	≤ -70°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG reagent	Y0580028	≤ -70°C	31 JAN 2027
S-PLEX Detect D2	Y0590020	≤ -70°C	30 NOV 2026
S-PLEX Enhance E3	Y0570021	≤ -70°C	30 NOV 2026

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

Spot Number	Name	Calibrator	
		Recovery	CV
1	IL-3	80% - 120%	< 20%

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
IL-3	10.6 - 170	42.5	fg/mL	80% - 120%	< 20%	6.54	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-3	1X Calibrator	38000	fg/mL

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		27 NOV 2024