



EXTENSION MEMO

Page 1 of 1

Product Description(s): S-PLEX® Human Eotaxin-3 Kit


Catalog Number(s) / Size(s): K151Q3S-Series

Lot Number: K00S0185

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K00S0185 has been extended to 31 JAN 2025.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Biotin Human Eotaxin-3 Antibody	B00T0036A	29 FEB 2024	31 JAN 2025
TURBO-BOOST™ Human Eotaxin-3 Antibody	D00T0036A	31 JAN 2024	31 JAN 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		29 FEB 2024

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

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Human Eotaxin-3 Kit
Catalog Number: K151Q3S-Series
Lot Number: K00S0185
Expiration Date: 31 JAN 2024

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Eotaxin-3 Antibody	B00T0036A	2-8°C	29 FEB 2024
Human Eotaxin-3 Calibrator	A00S0028	2-8°C	31 AUG 2025
TURBO-BOOST™ Human Eotaxin-3 Antibody	D00T0036A	2-8°C	31 JAN 2024
Blocker S1	R0010871	≤ -10°C	31 MAR 2026
Diluent 57	M0570011	≤ -10°C	31 JUL 2025
S-PLEX Enhance E1	Y0550014	≤ -10°C	28 FEB 2026
S-PLEX Enhance E2	Y0560016	≤ -10°C	31 MAY 2026
S-PLEX Coating Reagent C1	446866	≤ -70°C	31 MAR 2026
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580022	≤ -70°C	31 MAY 2026
S-PLEX Detect D2	Y0590016	≤ -70°C	30 APR 2026
S-PLEX Enhance E3	Y0570017	≤ -70°C	31 MAR 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	Eotaxin-3	80% - 120%	< 20%

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
Eotaxin-3	35.3 - 564	282	fg/mL	80% - 120%	< 20%	51.7	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Eotaxin-3	1X Calibrator	609000	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		29 FEB 2024