



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human Eotaxin-3 Kit
Lot Number: K00S0285
Expiration Date: 31 DEC 2026
Catalog Number: K151Q3S-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Eotaxin-3 Antibody	B00T0265	2-8°C	31 DEC 2028
TURBO-BOOST Human Eotaxin-3 Antibody	D00T0265	2-8°C	30 NOV 2028
Human Eotaxin-3 Calibrator	A00S0132	2-8°C	31 MAY 2030
Blocker S1	R0140015	≤ -10°C	31 OCT 2027
Diluent 57	M0570016	≤ -10°C	31 DEC 2026
S-PLEX Enhance E1	Y0550025	≤ -10°C	30 NOV 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	Eotaxin-3	80% - 120%	< 20%

LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
Eotaxin-3	35.3 - 564	70.5	fg/mL	80% - 120%	< 20%	17.8	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Eotaxin-3	1X Calibrator	655000	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		01 OCT 2025