

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human Eotaxin-3 Kit
Lot Number: K00S0306
Expiration Date: 31 DEC 2027
Catalog Number: K151Q3S-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Eotaxin-3 Antibody	B00T0277	2-8°C	31 JUL 2029
Human Eotaxin-3 Calibrator	A00S0132	2-8°C	31 MAY 2030
TURBO-BOOST Human Eotaxin-3 Antibody	D00T0277	2-8°C	31 AUG 2029
Blocker S1	R0140016	≤ -10°C	31 DEC 2027
Diluent 57	M0570021	≤ -10°C	29 FEB 2028
S-PLEX Enhance E1	Y0550027	≤ -10°C	29 FEB 2028
S-PLEX Enhance E2	Y0560025	≤ -10°C	29 FEB 2028
S-PLEX Coating Reagent C1	521330	≤ -70°C	30 JUN 2028
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580036	≤ -70°C	30 APR 2028
S-PLEX Detect D2	Y0590028	≤ -70°C	29 FEB 2028
S-PLEX Enhance E3	Y0570029	≤ -70°C	29 FEB 2028

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	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
Eotaxin-3	35.3 - 564	99.5	fg/mL	80% - 120%	< 20%	26.8	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Eotaxin-3	1X Calibrator	655000	fg/mL

This lot meets MSD's non-specific binding specifications.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		18 MAR 2026