



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®]Human Eotaxin-3 Kit

Catalog Number: K151Q3S-Series

Lot Number: K00S0088

Expiration Date: 28 FEB 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human Eotaxin-3 Antibody	D00T0036A	2-8°C	31 JAN 2024
Biotin Human Eotaxin-3 Antibody	B00T0036A	2-8°C	29 FEB 2024
Human Eotaxin-3 Calibrator	A00S0028	2-8°C	31 AUG 2025
Diluent 57	M0570001	≤ -10°C	28 FEB 2023
S-PLEX Coating Reagent C1	314374	≤ -70°C	28 FEB 2023
Blocker S1	R0140004	≤ -10°C	30 APR 2023
S-PLEX Enhance E1	Y0550004	≤ -10°C	28 FEB 2023
S-PLEX Enhance E2	Y0560007	≤ -10°C	31 OCT 2023
S-PLEX Enhance E3	Y0570008	≤ -70°C	30 JUN 2024
S-PLEX Detect D1	Y0580005	≤ -70°C	31 MAR 2023
S-PLEX Detect D2	Y0590008	≤ -70°C	31 JUL 2024

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

Precision, Accuracy, and Sensitivity Specifications:

Name	Calibrator	
	Recovery	CV
Eotaxin-3	80% - 120%	< 20%

Name	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
Eotaxin-3	35.3 - 564	141	fg/mL	80% - 120%	< 20%	33.7	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Eotaxin-3	1X Calibrator	609000	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		13 MAY 2022