



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human Eotaxin-3 Kit
Catalog Number: K151Q3S-Series
Lot Number: K00S0034
Expiration Date: 31 MAR 2022

Precision, Accuracy, and Sensitivity Specifications:

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

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

LLOD		
Name	Observed	Units
Eotaxin-3	58.9	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Eotaxin-3	1X Calibrator	609000	fg/mL

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	Y0560003	≤ -10°C	31 MAR 2022
S-PLEX Enhance E3	Y0570003	≤ -70°C	31 DEC 2022
S-PLEX Detect D1	Y0580005	≤ -70°C	31MAR 2023
S-PLEX Detect D2	Y0590003	≤ -70°C	31AUG 2022

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		04 FEB 2021