



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX Human Eotaxin-3 Gen. B Kit
Lot Number: K0082672
Expiration Date: 31 MAY 2027
Catalog Number: K151ANJ-Series, K156ANJ-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Gen. B Calibrator Blend	A0080416	2-8°C	31 DEC 2029
Human Eotaxin-3 Antibody	D0082257	2-8°C	30 NOV 2027
Human Eotaxin-3 Gen. B 96-Well Small Spot SECTOR Plate	Z0048097	2-8°C	31 AUG 2027
Diluent 3	M0030137	≤ -10°C	31 JUL 2027
Diluent 57	M0570018	≤ -10°C	31 MAY 2027

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

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator		Chemokine Panel 1 (human) Gen. B Control 1		Chemokine Panel 1 (human) Gen. B Control 2		Chemokine Panel 1 (human) Gen. B Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Eotaxin-3	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	Eotaxin-3	11.1	80% - 120%	< 20%	11100	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Eotaxin-3	21500	pg/mL

Dilution to top of Curve: 1X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		29 APR 2025