

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human Eotaxin-3 Gen. B Kit
Kit Lot Number: K0082355
Expiration Date: 31 Jul 2025
Kit Catalog Numbers: K151ANJ-Series, K156ANJ-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human Eotaxin-3 Gen. B SECTOR Plate	Z0047910	2-8°C	28 Feb 2026
Human Eotaxin-3 Antibody	D0082135	2-8°C	31 Jan 2027
Chemokine Panel 1 (hu) Gen. B Calibrator	A0080377	2-8°C	31 Aug 2028
Diluent 3	M0030123	≤ -10°C	31 Jan 2026
Diluent 57	M0570011	≤ -10°C	31 Jul 2025

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

Precision and Accuracy Specifications:

Spot	Name	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	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Eotaxin-3	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%


Spot	Name	LLOQ			ULOQ		
		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	17900	pg/mL

Dilution to Top of Curve = 1X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		19 Dec 2023