



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Human Eotaxin-3 Kit
Lot Number: K0082786
Expiration Date: 30 NOV 2026
Catalog Number: K151NU-Series; K156NU-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu Eotaxin-3 Antibody	D0082077	2-8°C	30 NOV 2026
Chemokine Panel 1 (human) Calibrator Blend	A0080369	2-8°C	30 JUN 2028
Human Eotaxin-3, 96-Well Small Spot SECTOR Plate	Z0048172	2-8°C	29 FEB 2028
Diluent 3	M0030140	≤ -10°C	30 SEP 2027
Diluent 43	M0430058	≤ -10°C	28 FEB 2027

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

Precision, Accuracy, and Sensitivity Specifications:

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

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Eotaxin-3	10.2	80% - 120%	< 20%	3750	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Eotaxin-3	6230	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		05 NOV 2025