

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human Eotaxin-3 Kit
Kit Lot Number: K0082281
Expiration Date: 30 Sep 2024
Kit Catalog Numbers: K151NU-Series; K156NU-Series

Precision and Accuracy 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	6760	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human Eotaxin-3 Plate	Z0047857	2-8°C	31 Oct 2025
Anti-hu Eotaxin-3 Antibody	D0081677	2-8°C	30 Sep 2024
Chemokine Panel 1 (human) Calibrator Blend	A0080338	2-8°C	31 Aug 2027
Diluent 3	M0030117	≤ -10°C	31 Jan 2025
Diluent 43	M0430050	≤ -10°C	31 Mar 2025

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	Xiomara Talavera	Quality		22 MAY 2023