

**CERTIFICATE OF ANALYSIS**

**Product Description:** V-PLEX® and V-PLEX Plus Human Eotaxin-3 Kit  
**Lot Number:** K0082839  
**Expiration Date:** 30 JUN 2027  
**Catalog Number:** K151NU-Series, K156NU-Series

**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu Eotaxin-3 Antibody	D0082173	2-8°C	30 JUN 2027
Chemokine Panel 1 (human) Calibrator Blend	A0080418	2-8°C	28 FEB 2030
Human Eotaxin-3, 96-Well Small Spot SECTOR Plate	Z0048200	2-8°C	30 APR 2028
Diluent 3	M0030141	≤ -10°C	30 NOV 2027
Diluent 43	M0430059	≤ -10°C	30 SEP 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) Control 1		Chemokine Panel 1 (human) Control 2		Chemokine Panel 1 (human) 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	10.2	80% - 120%	< 20%	3750	80% - 120%	< 20%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	Eotaxin-3	5770	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		15 JAN 2026