

**CERTIFICATE OF ANALYSIS**

**Product Description:** V-PLEX® and V-PLEX Plus Human Eotaxin-3 Gen. B Kit, SECTOR  
**Lot Number:** K0082957  
**Expiration Date:** 30 SEP 2027  
**Catalog Number:** K151ANJ-Series, K156ANJ-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
Chemokine Panel 1 (human) Gen. B Calibrator	A0080416	2-8°C	31 DEC 2029
Human Eotaxin-3 Antibody	D0082318	2-8°C	31 JAN 2028
Human Eotaxin-3 Gen. B 96-Well Small Spot SECTOR Plate	Z0048266	2-8°C	30 NOV 2028
Diluent 3	M0030139	≤ -10°C	30 SEP 2027
Diluent 57	M0570020	≤ -10°C	31 OCT 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) 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%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

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	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		07 JUL 2026