



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Human Eotaxin-3 Kit


Catalog Number(s) / Size(s): K151NU-Series; K156NU-Series

Lot Number: K0082251

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0082251 has been extended to 31 JUL 2024.

The expiration date for the following kit component has been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Human Eotaxin-3 Plate	Z0047464	31 JUL 2023	31 JUL 2024
Diluent 43	M0430047	31 MAY 2024	31 JUL 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		17 MAR 2023

This memo supersedes all other statements regarding date of expiration for this product.

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human Eotaxin-3 Kit
Kit Lot Number: K0082251
Expiration Date: 31 Jul 2023
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	6120	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human Eotaxin-3 Plate	Z0047464	2-8°C	31 Jul 2023
Anti-hu Eotaxin-3 Antibody	D0081652	2-8°C	30 Sep 2024
Chemokine Panel 1 (human) Calibrator Blend	A0080347	2-8°C	31 Oct 2027
Diluent 3	M0030114	≤ -10°C	30 Sep 2024
Diluent 43	M0430047	≤ -10°C	31 May 2024

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	Karen Hamilla	Quality		17 MAR 20232