



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: K0081771

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081771 has been extended to 31 AUG 2023.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 43	M0430037	31 AUG 2022	31 AUG 2023
Human Eotaxin-3 Plate	Z0047464	31 JUL 2023	31 AUG 2023
Anti-hu Eotaxin-3 Antibody	D0081386	31 MAR 2023	31 AUG 2023
Diluent 3	M0030103	30 APR 2023	31 AUG 2023

Function	Name	Signature	Date
Quality	Karen Hamilla		22 JUL 2022

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: K0081771
Expiration Date: 31 Aug 2022
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	5550	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	D0081386	2-8°C	31 Mar 2023
Chemokine Panel 1 (human) Calibrator Blend	A0080280	2-8°C	28 Feb 2026
Diluent 43	M0430037	≤ -10°C	31 Aug 2022
Diluent 3	M0030103	≤ -10°C	30 Apr 2023

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		01 SEP 2021