



## EXTENSION MEMO

Page 1 of 1

**Product Description(s):** V-PLEX® and V-PLEX Plus Human VEGF Kit


**Catalog Number(s) / Size(s):** K151RH-Series; K156RH-Series

**Lot Number:** K0081817

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081817 has been extended to 30 APR 2024.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Human VEGF Plate	Z0056994	31 JAN 2024	30 APR 2024
Anti-hu VEGF Antibody	D0081393	30 APR 2023	30 APR 2024
Diluent 43	M0430045	30 NOV 2023	30 APR 2024
Diluent 3	M0030109	31 JAN 2024	30 APR 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		18 JAN 2024

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 VEGF Kit  
**Kit Lot Number:** K0081817  
**Expiration Date:** 30 Apr 2023  
**Kit Catalog Numbers:** K151RH-Series; K156RH-Series;

**Precision and Accuracy Specifications:**

Spot	Name	Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	VEGF	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	VEGF	7.70	80% - 120%	< 20%	562	80% - 120%	< 20%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	VEGF	1150	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human VEGF Plate	Z0056994	2-8°C	31 Jan 2024
Anti-hu VEGF Antibody	D0081393	2-8°C	30 Apr 2023
Cytokine Panel 1 (human) Calibrator Blend	A0080283	2-8°C	31 Mar 2026
Diluent 43	M0430045	≤ -10°C	30 Nov 2023
Diluent 3	M0030109	≤ -10°C	31 Jan 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		28 OCT 2021