


**Extension Memo**

Product Description: V-PLEX and V-PLEX Plus Human VEGF Kit
Lot Number: K0082483
New Expiration Date: 31 DEC 2026
Catalog Number(s): K151RH-Series, K156RH-Series

Based on additional testing, the original date of expiration for this kit has been extended by 12 months.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Human VEGF-A 96-Well Small Spot Plate	Z0057472	30 NOV 2026	31 DEC 2026
Anti-hu VEGF Antibody	D0081882	31 DEC 2025	31 DEC 2026
Diluent 3	M0030127	30 SEP 2026	31 DEC 2026
Diluent 43	M0430056	30 SEP 2026	31 DEC 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		19 Dec 2025



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human VEGF Kit
Lot Number: K0082483
Expiration Date: 31 DEC 2025
Catalog Number: K151RH-Series; K156RH-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu VEGF Antibody	D0081882	2-8°C	31 DEC 2025
Cytokine Panel 1 (human) Calibrator Blend	A0080374	2-8°C	31 JUL 2028
Human VEGF Plate	Z0057472	2-8°C	30 NOV 2026
Diluent 3	M0030127	≤ -10°C	30 SEP 2026
Diluent 43	M0430056	≤ -10°C	30 SEP 2026

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	VEGF	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%


		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	VEGF	7.7	80% - 120%	< 20%	562	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	VEGF	1660	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		24 JUN 2024