

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human VEGF Kit  
**Kit Lot Number:** K0082183  
**Expiration Date:** 31 Mar 2025  
**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%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 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	1140	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human VEGF Plate	Z0057296	2-8°C	30 Apr 2025
Anti-hu VEGF Antibody	D0081772	2-8°C	30 Sep 2025
Cytokine Panel 1(human) Calibrator Blend	A0080307	2-8°C	30 Sep 2026
Diluent 3	M0030118	≤ -10°C	31 Mar 2025
Diluent 43	M0430050	≤ -10°C	31 Mar 2025

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	Xiomara Talavera	Quality		19 DEC 2022