

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human VEGF Kit
Kit Lot Number: K0082059
Expiration Date: 31 May 2024
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	Z0057197	2-8°C	30 Nov 2024
Anti-hu VEGF Antibody	D0081648	2-8°C	30 Sep 2024
Cytokine Panel 1 (human) Calibrator Blend	A0080307	2-8°C	30 Sep 2026
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		05 AUG 2022