



**CERTIFICATE OF ANALYSIS**

**Product Description:** V-PLEX® and V-PLEX Plus Angiogenesis Panel 1 (human) Kit  
**Lot Number:** K0082492  
**Expiration Date:** 30 SEP 2025  
**Catalog Number:** K151RZ-Series, K15190-Series, K151KW-Series, K151LT-Series, K151LU-Series, K151MD-Series, K151ME-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Angiogenesis Panel 1 (human) Plate	Z0057396	2-8°C	31 JAN 2026
Anti-hu FGF(basic) Antibody	D0082125	2-8°C	31 JAN 2027
Anti-hu PIGF Antibody	D0082124	2-8°C	31 JAN 2027
Anti-hu Tie-2 Antibody	D0082122	2-8°C	31 JAN 2027
Anti-hu VEGF Antibody	D0082126	2-8°C	31 JAN 2027
Anti-hu Flt-1 Antibody	D0082123	2-8°C	31 JAN 2027
Human VEGF-C Antibody	D0082127	2-8°C	28 FEB 2027
Human VEGF-D Antibody	D0082121	2-8°C	31 JAN 2027
Diluent 11	M0110066	≤ -10°C	30 SEP 2026
Diluent 7	M0070075	≤ -10°C	30 APR 2026
Angiogenesis Panel 1 (human) Calibrator Blend	A0080361	≤ -70°C	30 SEP 2025

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

**Precision, Accuracy, and Sensitivity Specifications:**

Spot	Name	Calibrator		Angiogenesis Control 1		Angiogenesis Control 2		Angiogenesis Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	VEGF-A	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
2	VEGF-C	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
3	VEGF-D	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
4	Tie-2	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
5	Flt-1	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
6	PIGF	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
7	FGF(basic)	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
2	VEGF-C	146	80% - 120%	< 20%	17500	80% - 120%	< 20%
3	VEGF-D	67.1	80% - 120%	< 20%	18800	80% - 120%	< 20%
4	Tie-2	396	80% - 120%	< 20%	63400	80% - 120%	< 20%
5	Flt-1	10	80% - 120%	< 25%	6410	80% - 120%	< 20%
6	PIGF	1.5	80% - 120%	< 25%	800	80% - 120%	< 25%
7	FGF(basic)	2.6	80% - 120%	< 25%	1780	80% - 120%	< 20%



## CERTIFICATE OF ANALYSIS

## Non-Specific Binding Specifications:


Spot	Name	VEGF-A	VEGF-C	VEGF-D	Tie-2	Flt-1	PlGF	FGF (basic)
1	VEGF-A	NA	< 0.5%	< 0.5%	< 0.5%	< 10.0%	< 0.5%	< 0.5%
2	VEGF-C	< 0.5%	NA	< 5.0%	< 0.5%	< 5.0%	< 0.5%	< 0.5%
3	VEGF-D	< 0.5%	< 0.5%	NA	< 0.5%	< 0.5%	< 0.5%	< 0.5%
4	Tie-2	< 0.5%	< 0.5%	< 0.5%	NA	< 0.5%	< 0.5%	< 0.5%
5	Flt-1	< 10.0%	< 0.5%	< 0.5%	< 0.5%	NA	< 0.5%	< 0.5%
6	PlGF	< 0.5%	< 0.5%	< 0.5%	< 0.5%	< 0.5%	NA	< 0.5%
7	FGF (basic)	< 0.5%	< 0.5%	< 0.5%	< 0.5%	< 0.5%	< 0.5%	NA

## Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	VEGF-A	40400	pg/mL
2	VEGF-C	784000	pg/mL
3	VEGF-D	568000	pg/mL
4	Tie-2	1938000	pg/mL
5	Flt-1	208000	pg/mL
6	PlGF	24200	pg/mL
7	FGF(basic)	52200	pg/mL

Dilution to top of Curve: 20X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		09 JUL 2024