

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Angiogenesis Panel 1 (human) Kit**Lot Number:** K0082746**Expiration Date:** 31 MAR 2027**Catalog Number:** K15190-Series, K151KW-Series, K151LT-Series, K151LU-Series, K151MD-Series, K151ME-Series, K151RZ-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Angiogenesis Panel 1 (human) 96-Well 7-Spot SECTOR Plate	Z0057532	2-8°C	31 MAY 2027
Anti-hu bFGF Antibody	D0082382	2-8°C	31 MAY 2028
Anti-hu Flt-1 Antibody	D0082380	2-8°C	31 MAY 2028
Anti-hu PlGF Antibody, 1 Plate	D0082381	2-8°C	31 MAY 2028
Anti-hu Tie-2 Antibody	D0082379	2-8°C	31 MAY 2028
Anti-hu VEGF Antibody	D0082383	2-8°C	31 MAY 2028
Human VEGF-C Antibody, 1 Plate	D0082384	2-8°C	31 MAY 2028
Human VEGF-D Antibody, 1 Plate	D0082378	2-8°C	31 MAY 2028
Diluent 11	M0110072	≤ -10°C	30 NOV 2027
Diluent 7	M0070077	≤ -10°C	31 MAR 2027
Angiogenesis Panel 1 (human) Calibrator Blend	A0080415	≤ -70°C	30 SEP 2027

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

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator		Angiogenesis Control 1		Angiogenesis Control 2		Angiogenesis Control 3	
Spot	Name	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	PlGF	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%



CERTIFICATE OF ANALYSIS

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	VEGF-A	5.00	80% - 120%	< 20%	1510	80% - 120%	< 20%
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	10.0	80% - 120%	< 25%	6410	80% - 120%	< 20%
6	PlGF	1.50	80% - 120%	< 25%	800	80% - 120%	< 25%
7	FGF (basic)	2.60	80% - 120%	< 25%	1780	80% - 120%	< 20%

Non-Specific Binding Specifications:

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

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	VEGF-A	39400	pg/mL
2	VEGF-C	686000	pg/mL
3	VEGF-D	582000	pg/mL
4	Tie-2	1864000	pg/mL
5	Flt-1	204000	pg/mL
6	PlGF	24400	pg/mL
7	FGF (basic)	51200	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		01 AUG 2025