

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX and V-PLEX Plus Angiogenesis Panel 1 (human) Kit, SECTOR
Lot Number: K0082843
Expiration Date: 31 MAR 2027
Catalog Number: K15190-Series, K151KW-Series, K151LT-Series, K151LU-Series, K151MD-Series, K151ME-Series, K151RZ-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Angiogenesis Panel 1 (human) 96-Well 7-Spot SECTOR Plate	Z0057608	2-8°C	29 FEB 2028
Anti-hu bFGF Antibody	D0082521	2-8°C	28 FEB 2029
Anti-hu Flt-1 Antibody	D0082519	2-8°C	28 FEB 2029
Anti-hu PlGF Antibody	D0082520	2-8°C	28 FEB 2029
Anti-hu Tie-2 Antibody	D0082518	2-8°C	31 MAR 2029
Anti-hu VEGF-A Antibody	D0082522	2-8°C	28 FEB 2029
Human VEGF-C Antibody	D0082523	2-8°C	28 FEB 2029
Human VEGF-D Antibody	D0082517	2-8°C	28 FEB 2029
Diluent 11	M0110073	≤ -10°C	29 FEB 2028
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	VEGFR-1/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%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	VEGF-A	5	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	VEGFR-1/Flt-1	10	80% - 120%	< 25%	6410	80% - 120%	< 20%
6	PlGF	1.5	80% - 120%	< 25%	800	80% - 120%	< 25%
7	FGF (basic)	2.6	80% - 120%	< 25%	1780	80% - 120%	< 20%

*See Additional Comments



CERTIFICATE OF ANALYSIS

Non-Specific Binding Specifications:

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

Dilution to top of Curve: 20X

Additional Comments:

Spot	Name	Specification	Comment
4	Tie-2	LLOQ Conc.	Specification for LLOQ is 80-120%, actual value 79.5% at target of 396 pg/mL, was passing at 550 pg/mL.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		13 FEB 2026