



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Angiogenesis Panel 1 (human) Kit, SECTOR
Lot Number: K0082872
Expiration Date: 30 SEP 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	Z0057651	2-8°C	30 JUN 2028
Anti-hu bFGF Antibody	D0082576	2-8°C	30 JUN 2029
Anti-hu Flt-1 Antibody	D0082574	2-8°C	30 JUN 2029
Anti-hu PIGF Antibody	D0082575	2-8°C	31 JUL 2029
Anti-hu Tie-2 Antibody	D0082573	2-8°C	30 JUN 2029
Anti-hu VEGF-A Antibody	D0082577	2-8°C	30 JUN 2029
Human VEGF-C Antibody	D0082578	2-8°C	31 JUL 2029
Human VEGF-D Antibody	D0082572	2-8°C	30 JUN 2029
Diluent 11	M0110073	≤ -10°C	29 FEB 2028
Diluent 7	M0070078	≤ -10°C	29 FEB 2028
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:

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	VEGFR-1/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
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	VEGFR-1/Flt-1	10.0	80% - 120%	< 25%	6410	80% - 120%	< 20%
6	PIGF	1.50	80% - 120%	< 25%	800	80% - 120%	< 25%
7	FGF(basic)	2.60	80% - 120%	< 25%	1780	80% - 120%	< 20%



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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		23 FEB 2026