

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Angiogenesis Panel 1 (human) Kit

**Kit Lot Number:** K0081690

**Expiration Date:** 31 May 2022

**Kit Catalog Numbers:** K15190-Series; K151RZ-Series; K151KW-Series; K151MD-Series; K151ME-Series; K151LT-Series; K151LU-Series

**Precision and Accuracy 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	bFGF	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	5.00	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	bFGF	2.60	80% - 120%	< 25%	1780	80% - 120%	< 20%

**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	bFGF	< 0.5%	< 0.5%	< 0.5%	< 0.5%	< 0.5%	< 0.5%	NA

## CERTIFICATE OF ANALYSIS

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	VEGF-A	40200	pg/mL
2	VEGF-C	536000	pg/mL
3	VEGF-D	612000	pg/mL
4	Tie-2	2020000	pg/mL
5	Flt-1	204000	pg/mL
6	PIGF	24000	pg/mL
7	bFGF	54400	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Angiogenesis Panel 1 (human) Plate	Z0056805	2-8°C	31 Mar 2023
Anti-hu VEGF Antibody	D0081494	2-8°C	30 Sep 2023
Anti-hu VEGF-C Antibody	D0081195	2-8°C	31 May 2022
Anti-hu VEGF-D Antibody	D0081576	2-8°C	31 Mar 2024
Anti-hu Tie-2 Antibody	D0081449	2-8°C	30 Sep 2023
Anti-hu Flt-1 Antibody	D0081578	2-8°C	31 Mar 2024
Anti-hu PIGF Antibody	D0081492	2-8°C	31 Oct 2023
Anti-hu bFGF Antibody	D0081493	2-8°C	30 Sep 2023
Angiogenesis Panel 1 (human) Calibrator Blend	A0080260	≤ -70°C	31 Dec 2022
Diluent 7	M0070065	≤ -10°C	31 Jan 2023
Diluent 11	M0110045	≤ -10°C	31 Jan 2023

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		13 MAY 2021