

CERTIFICATE OF ANALYSIS

Page 1 of 4

Product Description: V-PLEX® and V-PLEX Plus Angiogenesis Panel 1 (human) Kit
Kit Catalog Numbers: K15190-Series
Kit Lot Number: K0081029
Expiration Date: 31 Mar 2019

Acceptance Data/ Percent Recovery Specifications:

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

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu VEGF	5.00	80% - 120%	< 20%	1510	80% - 120%	< 20%
2	hu VEGF-C	146	80% - 120%	< 20%	17500	80% - 120%	< 20%
3	hu VEGF-D	67.1	80% - 120%	< 20%	18800	80% - 120%	< 20%
4	hu Tie-2	396	80% - 120%	< 20%	63400	80% - 120%	< 20%
5	hu Flt-1	10.0	80% - 120%	< 25%	6410	80% - 120%	< 20%
6	hu PIGF	1.50	80% - 120%	< 25%	800	80% - 120%	< 25%
7	hu bFGF	2.60	80% - 120%	< 25%	1780	80% - 120%	< 20%

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	hu VEGF	hu VEGF-C	hu VEGF-D	hu Tie-2	hu Flt-1
1	hu VEGF	NA	< 0.5%	< 0.5%	< 0.5%	< 10.0%
2	hu VEGF-C	< 0.5%	NA	< 0.5%	< 0.5%	< 0.5%
3	hu VEGF-D	< 0.5%	< 5.0%	NA	< 0.5%	< 0.5%
4	hu Tie-2	< 0.5%	< 0.5%	< 0.5%	NA	< 0.5%
5	hu Flt-1	< 10.0%	< 5.0%	< 0.5%	< 0.5%	NA
6	hu PIGF	< 0.5%	< 0.5%	< 0.5%	< 0.5%	< 0.5%
7	hu bFGF	< 0.5%	< 0.5%	< 0.5%	< 0.5%	< 0.5%

Spot	Name	hu PIGF	hu bFGF
1	hu VEGF	< 0.5%	< 0.5%
2	hu VEGF-C	< 0.5%	< 0.5%
3	hu VEGF-D	< 0.5%	< 0.5%
4	hu Tie-2	< 0.5%	< 0.5%
5	hu Flt-1	< 0.5%	< 0.5%
6	hu PIGF	NA	< 0.5%
7	hu bFGF	< 0.5%	NA

CERTIFICATE OF ANALYSIS

Page 2 of 4

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu VEGF	20X Calibrator	43000	pg/mL
		Angiogenesis Control 1	437	pg/mL
		Angiogenesis Control 2	88.9	pg/mL
		Angiogenesis Control 3	16	pg/mL
2	hu VEGF-C	20X Calibrator	414000	pg/mL
		Angiogenesis Control 1	5710	pg/mL
		Angiogenesis Control 2	1580	pg/mL
		Angiogenesis Control 3	366	pg/mL
3	hu VEGF-D	20X Calibrator	476000	pg/mL
		Angiogenesis Control 1	7070	pg/mL
		Angiogenesis Control 2	990	pg/mL
		Angiogenesis Control 3	94.4	pg/mL
4	hu Tie-2	20X Calibrator	1414000	pg/mL
		Angiogenesis Control 1	18700	pg/mL
		Angiogenesis Control 2	6320	pg/mL
		Angiogenesis Control 3	427	pg/mL
5	hu Flt-1	20X Calibrator	162800	pg/mL
		Angiogenesis Control 1	2320	pg/mL
		Angiogenesis Control 2	298	pg/mL
		Angiogenesis Control 3	26.3	pg/mL
6	hu PlGF	20X Calibrator	24000	pg/mL
		Angiogenesis Control 1	253	pg/mL
		Angiogenesis Control 2	61	pg/mL
		Angiogenesis Control 3	12.2	pg/mL
7	hu bFGF	20X Calibrator	37000	pg/mL
		Angiogenesis Control 1	534	pg/mL
		Angiogenesis Control 2	60.9	pg/mL
		Angiogenesis Control 3	5.5	pg/mL

Controls tested may not be the same lot as shipped.

Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentrations, please refer to the Standard Curve Data

CERTIFICATE OF ANALYSIS

Standard Curve Data:

hu VEGF - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2150	658,000	3.1
538	122,000	3.1
134	23,700	1.7
33.6	5,400	3.7
8.40	1,520	3.0
2.10	523	2.7
0.525	284	3.3
0.00	205	2.6

hu VEGF-C - Spot 2		
Conc. pg/mL	ECL Signal	%CV
20700	580,000	3.8
5180	101,000	5.7
1290	18,200	5.4
323	4,030	6.4
80.9	1,280	6.2
20.2	555	4.4
5.05	368	3.9
0.00	311	5.6

hu VEGF-D - Spot 3		
Conc. pg/mL	ECL Signal	%CV
23800	558,000	3.7
5950	175,000	3.4
1490	35,700	5.0
372	7,070	7.3
93.0	1,830	7.9
23.2	632	6.8
5.81	363	20.0
0.00	230	8.2

hu Tie-2 - Spot 4		
Conc. pg/mL	ECL Signal	%CV
70700	139,000	4.6
17700	34,300	4.8
4420	8,540	3.5
1100	2,340	2.9
276	833	3.3
69.0	448	4.4
17.3	363	7.6
0.00	328	6.7

hu Flt-1 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
8140	475,000	2.4
2040	121,000	3.0
509	29,500	2.0
127	7,350	3.0
31.8	2,040	2.4
7.95	645	3.4
1.99	291	4.7
0.00	179	6.1

hu PIGF - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1200	729,000	2.4
300	184,000	3.8
75.0	42,800	1.7
18.8	10,500	1.7
4.69	3,110	1.4
1.17	1,110	3.4
0.293	591	5.3
0.00	443	8.2

hu bFGF - Spot 7		
Conc. pg/mL	ECL Signal	%CV
1850	739,000	4.2
463	199,000	4.0
116	48,500	4.0
28.9	12,100	3.8
7.23	3,340	4.8
1.81	924	4.8
0.452	333	5.8
0.00	147	9.8

CERTIFICATE OF ANALYSIS

Page 4 of 4

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Angiogenesis Panel 1 (human) Plate	Z0056282	2-8°C	30 Nov 2020
Anti-hu VEGF Antibody	D0080912	2-8°C	31 Jul 2019
Anti-hu VEGF-C Antibody	D0081073	2-8°C	30 Sep 2021
Anti-hu VEGF-D Antibody	D0081069	2-8°C	30 Sep 2021
Anti-hu Tie-2 Antibody	D0081070	2-8°C	30 Sep 2021
Anti-hu Flt-1 Antibody	D0081071	2-8°C	31 Jul 2020
Anti-hu PIGF Antibody	D0081101	2-8°C	31 Jan 2022
Anti-hu bFGF Antibody	D0081072	2-8°C	31 Jul 2020
Angiogenesis Panel 1 (human) Calibrator Blend	A0080104	≤ -70°C	31 Mar 2019
Blocker A	Not Kit Specific	Not Kit Specific	N/A
Diluent 7	M0070053	≤ -10°C	31 Jul 2020
Diluent 11	M0110035	≤ -10°C	31 Dec 2020
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20x), 100 mL	Not Kit Specific
Angiogenesis Control 1	Not Kit Specific
Angiogenesis Control 2	Not Kit Specific
Angiogenesis Control 3	Not Kit Specific

Statement:

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	13 SEP 2018