

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human PIGF Kit
Kit Catalog Numbers: K151ME-Series
Kit Lot Number: K0081421
Expiration Date: 31 Jan 2022

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	PIGF	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	PIGF	1.50	80% - 120%	< 25%	800	80% - 120%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	PIGF	20X Calibrator	22500	pg/mL
		Angiogenesis Control 1	231	pg/mL
		Angiogenesis Control 2	51.7	pg/mL
		Angiogenesis Control 3	15.3	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

Standard Curve Data:

PIGF - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1130	185,000	5.1
281	37,100	2.5
70.3	7,340	5.6
17.6	1,800	4.4
4.39	546	2.3
1.10	211	2.9
0.275	128	7.2
0.00	104	5.9

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human PIGF Plate	Z0056659	2-8°C	30 Jun 2022
Anti-hu PIGF Antibody	D0081101	2-8°C	31 Jan 2022
Angiogenesis Panel 1 (human) Calibrator Blend	A0080213	≤ -70°C	31 May 2022
MSD Blocker A	Not Kit Specific	Not Kit Specific	N/A
Diluent 7	M0070061	≤ -10°C	30 Apr 2022
Diluent 11	M0110044	≤ -10°C	30 Jun 2022
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), 100mL	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	Karen Hamilla	Quality		02 MAR 2020