

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human bFGF Kit  
**Kit Catalog Numbers:** K151MD-Series  
**Kit Lot Number:** K0081293  
**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	bFGF	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	bFGF	2.60	80% - 120%	< 20%	1780	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	bFGF	20X Calibrator	50200	pg/mL
		Angiogenesis Control 1	545	pg/mL
		Angiogenesis Control 2	62	pg/mL
		Angiogenesis Control 3	7.46	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:

bFGF - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2510	329,000	1.9
628	87,100	3.0
157	21,000	5.2
39.2	5,510	1.7
9.80	1,450	3.8
2.45	410	2.7
0.613	147	2.8
0.00	63	7.9

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human bFGF Plate	Z0056561	2-8°C	31 Jan 2022
Anti-hu bFGF Antibody	D0081235	2-8°C	30 Jun 2022
Angiogenesis Panel 1 (human) Calibrator Blend	A0080213	≤ -70°C	31 May 2022
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
<b>Review/Approval</b>	Karen Hamilla	Quality		10 FEB 2020