

**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:** K0081007  
**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 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 bFGF	2.60	80% - 120%	< 20%	1780	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	hu bFGF	20X Calibrator	37000	pg/mL
		Angiogenesis Control 1	544	pg/mL
		Angiogenesis Control 2	62.6	pg/mL
		Angiogenesis Control 3	7.37	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:**

hu bFGF - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1850	285,000	3.1
463	77,100	1.5
116	19,100	1.7
28.9	5,020	2.4
7.23	1,320	2.0
1.81	391	2.7
0.452	155	3.9
0.00	76	3.9

**Components:**

**CERTIFICATE OF ANALYSIS**

Page 2 of 2

Description	Lot Number	Storage Temperature	Expiration Date
Human bFGF Plate	Z0056225	2-8°C	31 Oct 2020
Anti-hu bFGF Antibody	D0080911	2-8°C	30 Sep 2019
Angiogenesis Panel 1 (human) Calibrator Blend	A0080104	≤ -70°C	31 Mar 2019
Blocker A	Not Kit Specific	Not Kit Specific	N/A
Diluent 7	M0070051	≤ -10°C	31 Mar 2019
Diluent 11	M0110031	≤ -10°C	31 Aug 2019
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), 1000mL	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>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	13 JUN 2018