

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: K0081216
Expiration Date: 31 Jan 2021

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	49800	pg/mL
		Angiogenesis Control 1	559	pg/mL
		Angiogenesis Control 2	62.7	pg/mL
		Angiogenesis Control 3	5.88	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
2490	397,000	2.2
623	108,000	1.6
156	27,700	5.8
38.9	6,960	2.9
9.73	1,840	4.9
2.43	500	3.4
0.608	178	3.2
0.00	73	4.5

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human bFGF Plate	Z0056490	2-8°C	31 Aug 2021
Anti-hu bFGF Antibody	D0081194	2-8°C	31 May 2022
Angiogenesis Panel 1 (human) Calibrator Blend	A0080185	≤ -70°C	31 Mar 2021
Blocker A Solution	Not Kit Specific	Not Kit Specific	N/A
Diluent 7	M0070056	≤ -70°C	31 Jan 2021
Diluent 11	M0110040	≤ -70°C	30 Sep 2021
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		20 JAN 2020