

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human Flt-1 Kit
Kit Catalog Numbers: K151RZ-Series
Kit Lot Number: K0081181
Expiration Date: 31 Jul 2020

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 Flt-1	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu Flt-1	10.0	80% - 120%	< 20%	6410	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu Flt-1	20X Calibrator	196600	pg/mL
		Angiogenesis Control 1	2320	pg/mL
		Angiogenesis Control 2	298	pg/mL
		Angiogenesis Control 3	26.3	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:

hu Flt-1 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
9830	235,000	4.5
2460	60,300	2.0
614	14,600	1.2
154	3,800	2.2
38.4	999	1.4
9.60	304	4.2
2.40	135	3.2
0.00	78	6.6

CERTIFICATE OF ANALYSIS

Page 2 of 2

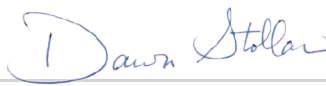
Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human Flt-1 Plate	Z0056457	2-8°C	30 Jun 2021
Anti-hu Flt-1 Antibody	D0081071	2-8°C	31 Jul 2020
Angiogenesis Panel 1 (human) Calibrator Blend	A0080185	≤ -70°C	31 Mar 2021
Blocker A	Not Kit Specific	Not Kit Specific	N/A
Diluent 7	M0070057	≤ -10°C	31 May 2021
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		21 MAR 2019