

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human VEGF Kit

Kit Catalog Numbers: K151RH-Series

Kit Lot Number: K0081032

Expiration Date: 31 May 2020

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu VEGF	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu VEGF	7.70	80% - 120%	< 20%	562	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu VEGF	Calibrator	1490	pg/mL
		Cytokine Panel 1 (human) Control 1	410	pg/mL
		Cytokine Panel 1 (human) Control 2	140	pg/mL
		Cytokine Panel 1 (human) Control 3	50	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu VEGF - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1490	331,000	1.5
373	62,800	1.5
93.1	13,000	3.1
23.3	3,100	3.7
5.82	819	2.0
1.46	244	3.3
0.364	106	3.2
0.00	59	6.8

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human VEGF Plate	Z0056286	2-8°C	30 Nov 2020
Anti-hu VEGF Antibody	D0081058	2-8°C	31 May 2020
Cytokine Panel 1 (human) Calibrator Blend	A0080144	2-8°C	31 Aug 2022
Diluent 43	M0430026	≤ -10°C	31 Aug 2020
Diluent 3	M0030082	≤ -10°C	31 Oct 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), 100mL	Not Kit Specific
Cytokine Panel 1 (human) Control 1	Not Kit Specific
Cytokine Panel 1 (human) Control 2	Not Kit Specific
Cytokine Panel 1 (human) 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		12 JUL 2018