

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

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

**Kit Catalog Numbers:** K151RH-Series; K156RH-Series

**Kit Lot Number:** K0081359

**Expiration Date:** 31 Dec 2021

**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	VEGF	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

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

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	VEGF	Calibrator	1080	pg/mL
		Cytokine Panel 1 (human) Control 1	472	pg/mL
		Cytokine Panel 1 (human) Control 2	143	pg/mL
		Cytokine Panel 1 (human) Control 3	49.6	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

VEGF - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1080	292,000	1.5
270	66,700	1.8
67.5	14,500	2.2
16.9	3,580	2.2
4.22	957	2.6
1.05	289	2.8
0.264	115	4.0
0.00	57	6.8

## CERTIFICATE OF ANALYSIS

Page 2 of 2


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human VEGF Plate	Z0056556	2-8°C	31 Dec 2021
Anti-hu VEGF Antibody	D0081393	2-8°C	30 Apr 2023
Cytokine Panel 1 (human) Calibrator Blend	A0080188	2-8°C	30 Nov 2023
Diluent 43	M0430033	≤ -10°C	31 Jan 2022
Diluent 3	M0030090	≤ -10°C	31 Dec 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
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
<b>Review/Approval</b>	Karen Hamilla	Quality		11 MAR 2020