

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: K0081210
Expiration Date: 30 Jun 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	1120	pg/mL
		Cytokine Panel 1 (human) Control 1	384	pg/mL
		Cytokine Panel 1 (human) Control 2	114	pg/mL
		Cytokine Panel 1 (human) Control 3	37.5	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

VEGF - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1120	227,000	2.2
280	39,300	4.1
70.0	7,850	2.4
17.5	1,870	2.3
4.38	535	2.8
1.09	178	4.1
0.273	92	3.8
0.00	61	7.4

CERTIFICATE OF ANALYSIS

Components:


Description	Lot Number	Storage Temperature	Expiration Date
Human VEGF Plate	Z0056486	2-8°C	31 Aug 2021
Anti-hu VEGF Antibody	D0081148	2-8°C	28 Feb 2022
Cytokine Panel 1 (human) Calibrator Blend	A0080170	2-8°C	30 Jun 2023
Diluent 43	M0430030	≤ -10°C	30 Jun 2021
Diluent 3	M0030087	≤ -10°C	30 Jun 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) 100 mL	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	Karen Hamilla	Quality		05 JUN 2019