

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-16 Kit

Kit Catalog Numbers: K151RE - Series

Kit Lot Number: K0081117

Expiration Date: 30 Sep 2019

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 IL-16	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu IL-16	19.1	80% - 120%	< 20%	1870	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-16	Calibrator	2800	pg/mL
		Cytokine Panel 1(human) Control 1	1210	pg/mL
		Cytokine Panel 1(human) Control 2	329	pg/mL
		Cytokine Panel 1(human) Control 3	104	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-16 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2800	404,000	3.3
700	86,400	2.7
175	17,600	2.6
43.8	3,900	3.0
10.9	1,040	2.0
2.73	300	1.4
0.684	131	2.9
0.00	74	3.4

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-16 Plate	Z0046635	2-8°C	28 Feb 2020
Anti-hu IL-16 Antibody	D0081056	2-8°C	30 Apr 2020
Cytokine Panel 1(human) Calibrator Blend	A0080121	2-8°C	28 Feb 2022
Diluent 43	M0430021	≤ -10°C	30 Sep 2019
Diluent 3	M0030077	≤ -10°C	30 Nov 2019
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	<i>Dawn Stollar</i>	07 NOV 2018