

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: K0080885
Expiration Date: 31 Jul 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	2450	pg/ml
		Cytokine Panel 1 (Human) Control 1	1160	pg/ml
		Cytokine Panel 1 (Human) Control 2	328	pg/ml
		Cytokine Panel 1 (Human) Control 3	126	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
2450	405,000	2.6
613	86,600	3.8
153	17,300	4.4
38.3	3,890	4.7
9.57	1,090	4.4
2.39	322	2.2
0.598	135	3.5
0.00	73	5.1

CERTIFICATE OF ANALYSIS

Components:


Description	Lot Number	Storage Temperature	Expiration Date
Human IL-16 Plate	Z0046635	2-8°C	28 Feb 2020
Anti-hu IL-16 Antibody	D0080891	2-8°C	31 Jul 2019
Cytokine Panel 1 (human) Calibrator Blend	A0080063	2-8°C	31 Jul 2020
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		09 OCT 2017