

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-Plex® and V-Plex Plus Human IL-17C Kit  
**Kit Catalog Numbers:** K151WP-Series  
**Kit Lot Number:** K0080683  
**Expiration Date:** 30 Apr 2019

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		Cytokine Panel 2 (human) Control 1		Cytokine Panel 2 (human) Control 2		Cytokine Panel 2 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu IL-17C	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu IL-17C	3.84	75% - 125%	< 20%	1620	75% - 125%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	hu IL-17C	Calibrator	2500	pg/ml
		Cytokine Panel 2 (human) Control 1	1041	pg/ml
		Cytokine Panel 2 (human) Control 2	165	pg/ml
		Cytokine Panel 2 (human) Control 3	19.2	pg/ml

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

hu IL-17C - Spot 1		
Conc. pg/ml	ECL Signal	%CV
2500	129,000	3.2
625	33,100	3.5
156	7,670	3.3
39.1	1,870	2.9
9.77	482	2.4
2.44	187	3.9
0.610	114	6.8
0.00	88	9.3

**CERTIFICATE OF ANALYSIS**

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-17C Plate	Z0046316	2-8°C	31 May 2019
Anti-hu IL-17C Antibody	D0080768	2-8°C	30 Apr 2019
Cytokine Panel 2 (human) Calibrator Blend	A0080116	2-8°C	31 Jan 2022
Diluent 43	M0430020	≤ -10°C	30 Jun 2019
Diluent 3	M0030074	≤ -10°C	30 Jun 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), 100 mL	Not Kit Specific
Cytokine Panel 2 (human) Control 1	Not Kit Specific
Cytokine Panel 2 (human) Control 2	Not Kit Specific
Cytokine Panel 2 (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>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	27 JUL 2017