

**CERTIFICATE OF ANALYSIS**

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-17A Gen. B Kit  
**Kit Catalog Numbers:** K151WM-Series  
**Kit Lot Number:** K0080720  
**Expiration Date:** 31 Jan 2019

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		TH17 Panel 1 (human) Control 1		TH17 Panel 1 (human) Control 2		TH17 Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu IL-17A	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-17A	5.86	80% - 120%	< 20%	1950	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	hu IL-17A	Calibrator	2730	pg/mL
		TH17 Panel 1 (human) Control 1	2725	pg/mL
		TH17 Panel 1 (human) Control 2	413	pg/mL
		TH17 Panel 1 (human) Control 3	57	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

hu IL-17A - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2730	180,000	1.3
683	42,200	2.0
171	9,810	2.8
42.7	2,560	1.6
10.7	698	2.3
2.67	223	2.2
0.667	101	3.3
0.00	62	5.7

## CERTIFICATE OF ANALYSIS

Page 2 of 2

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-17A Gen. B Plate	Z0046395	2-8°C	31 Jul 2019
Anti-hu IL-17A Antibody	D0080754	2-8°C	31 Jan 2019
TH17 Panel 1 (human) Calibrator Blend	A0080101	2-8°C	30 Sep 2021
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), 100mL	Not Kit Specific
TH17 Panel 1 (human) Control 1	Not Kit Specific
TH17 Panel 1 (human) Control 2	Not Kit Specific
TH17 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>	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	03MAY2017