

**CERTIFICATE OF ANALYSIS**

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-17A Kit  
**Kit Catalog Numbers:** K151RF-series  
**Kit Lot Number:** K0080751  
**Expiration Date:** 30 Jun 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-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	9.32	80% - 120%	< 20%	3650	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	hu IL-17A	Calibrator	6560	pg/mL
		Cytokine Panel 1 (human) Control 1	2590	pg/mL
		Cytokine Panel 1 (human) Control 2	327	pg/mL
		Cytokine Panel 1 (human) Control 3	57.5	pg/mL

Controls tested may not be the same lot as shipped.

**CERTIFICATE OF ANALYSIS**

**Standard Curve Data:**

hu IL-17A - Spot 1		
Conc. pg/mL	ECL Signal	%CV
6560	823,000	1.5
1640	220,000	1.9
410	55,900	2.2
103	13,700	3.4
25.6	3,490	2.5
6.41	932	3.3
1.60	302	2.3
0.00	99	6.0


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-17A	Z0046421	2-8°C	31 Aug 2019
Anti-hu IL-17A Antibody	D0080892	2-8°C	31 Jul 2019
Cytokine Panel 1 (human) Calibrator Blend	A0080073	2-8°C	31 Jan 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
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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		27 APR 2017