

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Mouse IL-17A Kit  
**Kit Catalog Numbers:** K152RF-Series  
**Kit Lot Number:** K0081237  
**Expiration Date:** 30 Jun 2021

### Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		TH17 Panel 1 (mouse) Control 1		TH17 Panel 1 (mouse) Control 2		TH17 Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-17A	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-17A	0.255	75% - 125%	< 20%	360	75% - 125%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-17A	Calibrator	638	pg/mL
		TH17 Panel 1 (mouse) Control 1	494	pg/mL
		TH17 Panel 1 (mouse) Control 2	85.2	pg/mL
		TH17 Panel 1 (mouse) Control 3	14.8	pg/mL

Controls tested may not be the same lot as shipped.

### Standard Curve Data:

IL-17A - Spot 1		
Conc. pg/mL	ECL Signal	%CV
638	259,000	0.8
160	67,000	2.5
39.9	16,700	1.3
9.97	4,180	1.1
2.49	1,080	1.0
0.623	316	2.1
0.156	126	3.8
0.00	64	6.1

**CERTIFICATE OF ANALYSIS**

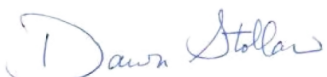
**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-17A Plate	Z0047078	2-8°C	30 Sep 2021
Anti-ms IL-17A Antibody	D0081190	2-8°C	31 Jul 2022
TH17 Panel 1 (mouse) Calibrator Blend	A0080157	2-8°C	31 Oct 2022
Diluent 41	M0410027	≤ -10°C	31 Aug 2021
Diluent 45	M0450026	≤ -10°C	30 Jun 2021
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
TH17 Panel 1 (mouse) Control 1	Not Kit Specific
TH17 Panel 1 (mouse) Control 2	Not Kit Specific
TH17 Panel 1 (mouse) 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		11 JUN 2019