

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Mouse IL-17F Kit

**Kit Catalog Numbers:** K152WS-Series

**Kit Lot Number:** K0080782

**Expiration Date:** 30 Sep 2019

**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	ms IL-17F	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms IL-17F	320	75% - 125%	< 20%	52000	75% - 125%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	ms IL-17F	Calibrator	87800	pg/mL
		TH17 Panel 1 (mouse) Control 1	51900	pg/mL
		TH17 Panel 1 (mouse) Control 2	10800	pg/mL
		TH17 Panel 1 (mouse) Control 3	2370	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

ms IL-17F - Spot 1		
Conc. pg/mL	ECL Signal	%CV
87800	233,000	3.4
22000	67,100	3.8
5490	17,300	2.9
1370	4,180	4.8
343	1,090	3.8
85.7	312	4.7
21.4	122	3.9
0.00	59	6.5

## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-17F Plate	Z0046472	2-8°C	30 Sep 2019
Anti-ms IL-17F Antibody	D0080955	2-8°C	31 Oct 2019
TH17 Panel 1 (mouse) Calibrator Blend	A0080154	2-8°C	31 Oct 2022
Diluent 41	M0410018	≤ -10°C	30 Nov 2019
Diluent 45	M0450019	≤ -10°C	31 Dec 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
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 Assurance	<i>Dawn Stollar</i>	02 JAN 2018