

## 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:** K0081339

**Expiration Date:** 31 Dec 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	566	pg/mL
		TH17 Panel 1 (mouse) Control 1	508	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
566	168,000	20.0
142	47,500	2.9
35.4	11,500	2.9
8.84	2,850	1.9
2.21	739	2.8
0.553	222	3.2
0.138	95	3.1
0.00	58	4.5

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-17A SECTOR Plate	Z0047189	2-8°C	31 Mar 2022
Anti-ms IL-17A Antibody	D0081190	2-8°C	31 Jul 2022
TH17 Panel 1 (mouse) Calibrator Blend	A0080192	2-8°C	31 May 2024
Diluent 41	M0410029	≤ -10°C	31 Dec 2021
Diluent 45	M0450030	≤ -10°C	28 Feb 2022
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 (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>	Karen Hamilla	Quality		18 NOV 2019