

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Mouse IL-17A/F Kit  
**Kit Catalog Numbers:** K152WN-Series  
**Kit Lot Number:** K0081038  
**Expiration Date:** 31 Aug 2019

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		Cytokine Panel 1 (mouse) Control 1		Cytokine Panel 1 (mouse) Control 2		Cytokine Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	ms IL-17A/F	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-17A/F	1.39	75% - 125%	< 20%	1620	75% - 125%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	ms IL-17A/F	Calibrator	2460	pg/mL
		Cytokine Panel 1 (mouse) Control 1	2220	pg/mL
		Cytokine Panel 1 (mouse) Control 2	212	pg/mL
		Cytokine Panel 1 (mouse) Control 3	18.6	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

ms IL-17A/F - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2460	281,000	7.4
615	62,800	6.1
154	15,400	6.8
38.4	3,680	9.6
9.61	1,010	10.0
2.40	307	9.8
0.601	155	6.7
0.00	92	3.2

## CERTIFICATE OF ANALYSIS


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-17A/F Plate	Z0046390	2-8°C	31 Aug 2019
Anti-ms IL-17A/F Antibody	D0080975	2-8°C	30 Nov 2019
Cytokine Panel 1 (mouse) Calibrator Blend	A0080125	2-8°C	31 Mar 2022
Diluent 41	M0410016	≤ -10°C	31 Aug 2019
Diluent 45	M0450016	≤ -10°C	30 Sep 2019
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Cytokine Panel 1 (mouse) Control 1	Not Kit Specific
Cytokine Panel 1 (mouse) Control 2	Not Kit Specific
Cytokine 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		23 AUG 2018