

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: K0080724
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	2230	pg/mL
		Cytokine Panel 1 (mouse) Control 2	211	pg/mL
		Cytokine Panel 1 (mouse) Control 3	19.1	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	332,000	5.0
615	80,400	7.2
154	20,900	3.1
38.4	4,750	7.0
9.61	1,330	6.7
2.40	367	4.3
0.601	150	5.2
0.00	74	4.6

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	D0080936	2-8°C	31 Aug 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

Plate Seals	Not Kit Specific
Wash Buffer (20x), 100mL	Not Kit Specific
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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	28 JUL 2017