

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: K0080783
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-17A	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	0.255	75% - 125%	< 20%	360	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IL-17A	Calibrator	585	pg/mL
		TH17 Panel 1 (mouse) Control 1	508	pg/mL
		TH17 Panel 1 (mouse) Control 2	101	pg/mL
		TH17 Panel 1 (mouse) Control 3	14.6	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms IL-17A - Spot 1		
Conc. pg/mL	ECL Signal	%CV
585	286,000	0.8
146	73,700	1.3
36.6	18,300	2.4
9.14	4,540	3.2
2.29	1,190	1.8
0.571	334	3.5
0.143	128	2.9
0.00	60	5.0

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-17A Plate	Z0046474	2-8°C	30 Sep 2019
Anti-ms IL-17A Antibody	D0080957	2-8°C	30 Nov 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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	02 JAN 2018