

CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human IL-17A Gen. B Kit
Kit Catalog Numbers: K151WM-Series
Kit Lot Number: K0080589
Expiration Date: 31 Jan 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		TH17 Panel 1 (human) Control 1		TH17 Panel 1 (human) Control 2		TH17 Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-17A	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-17A	5.86	80% - 120%	< 20%	1950	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-17A	Calibrator	2750	pg/ml
		TH17 Panel 1 (human) Control 1	1360	pg/ml
		TH17 Panel 1 (human) Control 2	214	pg/ml
		TH17 Panel 1 (human) Control 3	39.2	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
2750	177,000	2.1
688	41,000	3.4
172	9,920	2.4
43.0	2,600	4.7
10.7	672	4.0
2.69	226	3.3
0.671	105	4.5
0.00	72	5.5


CERTIFICATE OF ANALYSIS
Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-17A Gen. B Plate	Z0046169	2-8°C	31 Jan 2019
Anti-hu IL-17A Antibody	D0080754	2-8°C	31 Jan 2019
TH17 Panel 1 (human) Calibrator Blend	A0080101	2-8°C	30 Sep 2021
Diluent 43	M0430018	≤ -10°C	31 Jan 2019
Diluent 3	M0030073	≤ -10°C	31 Mar 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
TH17 Panel 1 (human) Control 1	Not Kit Specific
TH17 Panel 1 (human) Control 2	Not Kit Specific
TH17 Panel 1 (human) Control 3	Not Kit Specific

Statement:

The above kit was tested using MSD procedures and it conforms to MSD's specifications. 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		15 FEB 2017