

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: K0081105

Expiration Date: 30 Nov 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	638	pg/mL
		TH17 Panel 1 (mouse) Control 1	494	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:

ms IL-17A - Spot 1		
Conc. pg/mL	ECL Signal	%CV
638	328,000	1.9
160	85,800	1.3
39.9	21,600	2.5
9.97	5,420	1.5
2.49	1,420	1.6
0.623	405	2.0
0.156	153	2.3
0.00	74	6.7

CERTIFICATE OF ANALYSIS

Components:


Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-17A Plate	Z0046933	2-8°C	31 Mar 2021
Anti ms IL-17A Antibody	D0080957	2-8°C	30 Nov 2019
TH17 Panel 1 (mouse) Calibrator Blend	A0080157	2-8°C	31 Oct 2022
Diluent 41	M0410024	≤ -10°C	30 Sep 2020
Diluent 45	M0450023	≤ -10°C	31 Jan 2021
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		07 NOV 2018