

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-16 Kit

Kit Catalog Numbers: K152RE-Series

Kit Lot Number: K0080784

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-16	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-16	15.0	75% - 125%	< 25%	1800	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IL-16	Calibrator	2420	pg/mL
		TH17 Panel 1 (mouse) Control 1	1670	pg/mL
		TH17 Panel 1 (mouse) Control 2	924	pg/mL
		TH17 Panel 1 (mouse) Control 3	519	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms IL-16 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2420	217,000	2.8
605	34,600	1.6
151	6,470	3.2
37.8	1,840	5.1
9.45	928	5.4
2.36	669	11.0
0.591	614	8.3
0.00	618	11.0

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-16 Plate	Z0046475	2-8°C	30 Sep 2019
Anti-ms IL-16 Antibody	D0080956	2-8°C	31 Oct 2019
TH17 Panel 1 (mouse) Calibrator Blend	A0080146	2-8°C	31 Aug 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