

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-23 Kit

Kit Catalog Numbers: K152WW-Series

Kit Lot Number: K0080777

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-23	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-23	4.19	75% - 125%	< 25%	7600	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IL-23	Calibrator	8690	pg/mL
		TH17 Panel 1 (mouse) Control 1	6450	pg/mL
		TH17 Panel 1 (mouse) Control 2	1270	pg/mL
		TH17 Panel 1 (mouse) Control 3	272	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms IL-23 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
8690	420,000	2.3
2170	129,000	1.0
543	33,600	2.2
136	8,480	1.6
33.9	2,180	1.5
8.49	609	1.7
2.12	204	1.3
0.00	78	6.4

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-23 Plate	Z0046461	2-8°C	30 Sep 2019
Anti-ms IL-23 Antibody	D0080951	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