

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: K0081218
Expiration Date: 30 Jun 2021

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

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-23	Calibrator	8650	pg/mL
		TH17 Panel 1 (mouse) Control 1	6450	pg/mL
		TH17 Panel 1 (mouse) Control 2	1330	pg/mL
		TH17 Panel 1 (mouse) Control 3	199	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-23 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
8650	516,000	1.4
2160	154,000	2.6
541	39,500	1.6
135	9,960	1.9
33.8	2,520	1.7
8.45	676	3.3
2.11	223	2.7
0.00	75	8.0

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-23 SECTOR Plate	Z0047056	2-8°C	31 Aug 2021
Anti-ms IL-23 Antibody	D0081276	2-8°C	30 Sep 2022
TH17 Panel 1 (mouse) Calibrator Blend	A0080157	2-8°C	31 Oct 2022
Diluent 41	M0410026	≤ -10°C	30 Jun 2021
Diluent 45	M0450026	≤ -10°C	30 Jun 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	Karen Hamilla	Quality		04 JUN 2019