

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-23 Kit
Kit Catalog Numbers: K151WW-Series
Kit Lot Number: K0080594
Expiration Date: 31 Jan 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		TH17 Panel 1 (human) Control 1		TH17 Panel 1 (human) Control 3		TH17 Panel 1 (human) Control 2	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-23	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-23	4.60	80% - 120%	< 20%	3250	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-23	Calibrator	4900	pg/mL
		TH17 Panel 1 (human) Control 1	2637	pg/mL
		TH17 Panel 1 (human) Control 2	357	pg/mL
		TH17 Panel 1 (human) Control 3	32.8	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
4900	223,000	4.8
1230	68,800	3.4
306	18,700	3.5
76.6	4,740	2.4
19.1	1,300	3.2
4.79	386	4.4
1.20	154	5.9
0.00	76	4.2

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-23 Plate	Z0046179	2-8°C	28 Feb 2019
Anti-hu IL-23 Antibody	D0080759	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), 100 mL	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 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>	12 JAN 2017