

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: K0080886
Expiration Date: 31 Jan 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		TH17 Panel 1 (human) Control 1		TH17 Panel 1 (human) Control 2		TH17 Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu 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	hu IL-23	4.60	80% - 120%	< 20%	3250	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-23	Calibrator	4350	pg/mL
		TH17 Panel 1 (human) Control 1	4173	pg/mL
		TH17 Panel 1 (human) Control 2	464	pg/mL
		TH17 Panel 1 (human) Control 3	53	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-23 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
4350	209,000	3.5
1090	61,400	3.8
272	16,000	3.0
68.0	4,090	2.7
17.0	1,130	2.9
4.25	356	2.9
1.06	166	2.8
0.00	101	3.6

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-23 Plate	Z0046637	2-8°C	28 Feb 2020
Anti-hu IL-23 Antibody	D0080759	2-8°C	31 Jan 2019
TH17 Panel 1 (human) Calibrator Blend	A0080111	2-8°C	30 Nov 2021
Diluent 43	M0430023	≤ -10°C	31 Dec 2019
Diluent 3	M0030078	≤ -10°C	31 Jan 2020
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), 100mL	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>	13 OCT 2017