

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-12/IL-23p40 Kit

**Kit Catalog Numbers:** K151RJ-Series

**Kit Lot Number:** K0080896

**Expiration Date:** 30 Jun 2019

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu IL-12/IL-23p40	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu IL-12/IL-23p40	5.68	80% - 120%	< 20%	2250	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	hu IL-12/IL-23p40	Calibrator	3400	pg/mL
		Cytokine Panel 1 (human) Control 1	1580	pg/mL
		Cytokine Panel 1 (human) Control 2	236	pg/mL
		Cytokine Panel 1 (human) Control 3	39.4	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

hu IL-12/IL-23p40 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
3400	333,000	1.7
850	89,300	1.6
213	22,600	1.6
53.1	5,770	1.3
13.3	1,500	2.4
3.32	434	1.9
0.830	167	2.8
0.00	82	3.4

**CERTIFICATE OF ANALYSIS**

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12/IL-23p40 Plate	Z0046655	2-8°C	31 Mar 2020
Anti-hu IL-12/23p40 Antibody	D0080889	2-8°C	31 Jul 2019
Cytokine Panel 1 (human) Calibrator Blend	A0080121	2-8°C	28 Feb 2022
Diluent 43	M0430020	≤ -10°C	30 Jun 2019
Diluent 3	M0030074	≤ -10°C	30 Jun 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), 100mL	Not Kit Specific
Cytokine Panel 1 (human) Control 1	Not Kit Specific
Cytokine Panel 1 (human) Control 2	Not Kit Specific
Cytokine 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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	30 OCT 2017