

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® AND V-PLEX Plus Human IL-12/IL-23p40 Kit
Kit Catalog Numbers: K156RJ-Series, K151RJ-Series
Kit Lot Number: K0081461
Expiration Date: 30 Sep 2021

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	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	IL-12/IL-23p40	1.32	80% - 120%	< 20%	2250	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-12/IL-23p40	Calibrator	3900	pg/mL
		Cytokine Panel 1 (human) Control 1	1490	pg/mL
		Cytokine Panel 1 (human) Control 2	225	pg/mL
		Cytokine Panel 1 (human) Control 3	39.8	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-12/IL-23p40 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
3900	324,000	1.9
975	84,500	3.7
244	21,300	1.9
60.9	5,330	2.3
15.2	1,370	2.2
3.81	398	3.0
0.952	163	3.3
0.00	84	4.3

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12/IL-23 p40 Plate	Z0047301	2-8°C	30 Sep 2022
Anti-hu IL-12/23p40 Antibody	D0081080	2-8°C	30 Sep 2021
Cytokine Panel 1 (human) Calibrator Blend	A0080227	2-8°C	31 Oct 2024
Dilent 43	M0430035	≤ -10°C	31 Mar 2022
Dileunt 3	M0030092	≤ -10°C	31 Mar 2022
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
MSD Wash Buffer (20X), 100mL	Not Kit Specific
Plate Seals	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
Review/Approval	Karen Hamilla	Quality		21 APR 2020