

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-12/IL-23p40 Kit  
**Kit Lot Number:** K0081567  
**Expiration Date:** 30 Apr 2022  
**Kit Catalog Numbers:** K151RJ-Series; K156RJ-Series

**Precision and Accuracy 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 Calibrator:**


Spot	Name	Conc.	Units
1	IL-12/IL-23p40	3380	pg/ml

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12/IL-23p40 plate	Z0047407	2-8°C	31 Mar 2023
Anti-hu IL-12/23p40 Antibody	D0081186	2-8°C	30 Apr 2022
Cytokine Panel 1(human) Calibrator Blend	A0080240	2-8°C	30 Nov 2024
Diluent 43	M0430039	≤ -10°C	30 Sep 2022
Diluent 3	M0030097	≤ -10°C	31 Aug 2022

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Carissa B. Hubbard	Quality		16 NOV 2020