

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-12/IL-23p40 Kit  
**Kit Lot Number:** K0081808  
**Expiration Date:** 30 Apr 2023  
**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	3270	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12/IL-23p40 Plate	Z0047577	2-8°C	31 Jan 2024
Anti-hu IL-12/IL-23p40 Antibody	D0081418	2-8°C	30 Apr 2023
Cytokine Panel 1 (human) Calibrator Blend	A0080283	2-8°C	31 Mar 2026
Diluent 43	M0430045	≤ -10°C	30 Nov 2023
Diluent 3	M0030108	≤ -10°C	30 Nov 2023

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	Karen Hamilla	Quality		24 JAN 2022