

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX Plus Human IL-12/IL-23p40 Kit
Lot Number: K0082600
Expiration Date: 31 JUL 2026
Catalog Number: K151RJ-Series, K156RJ-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-12/23p40 Antibody	D0082041	2-8°C	30 SEP 2026
Cytokine Panel 1 (human) Calibrator Blend	A0080300	2-8°C	31 JUL 2026
Human IL-12/IL-23 p40 SECTOR Plate	Z0048047	2-8°C	31 MAY 2027
Diluent 3	M0030129	≤ -10°C	30 NOV 2026
Diluent 43	M0430056	≤ -10°C	30 SEP 2026

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

Precision, Accuracy, and Sensitivity 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	3350	pg/mL

Dilution to top of Curve: 1X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		17 JAN 2025