

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-23 Kit  
**Kit Lot Number:** K0082364  
**Expiration Date:** 30 Jun 2025  
**Kit Catalog Numbers:** K151WW-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-23 SECTOR Plate	Z0047909	2-8°C	28 Feb 2026
Anti-hu IL-23 Antibody	D0081784	2-8°C	30 Jun 2025
TH17 Panel 1 (human) Calibrator Blend	A0080366	2-8°C	31 May 2028
Diluent 3	M0030123	≤ -10°C	31 Jan 2026
Diluent 43	M0430053	≤ -10°C	31 Jan 2026

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

**Precision and Accuracy Specifications:**

		Calibrator		TH17 Panel 1 (human) Control 1		TH17 Panel 1 (human) Control 2		TH17 Panel 1 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-23	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%


		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IL-23	4.60	80% - 120%	< 20%	3250	80% - 120%	< 20%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	IL-23	5270	pg/mL

Dilution to Top of Curve = 1X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		03 NOV 2023