

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-23 Kit
Kit Lot Number: K0081948
Expiration Date: 30 Nov 2023
Kit Catalog Numbers: K151WW-Series

Precision and Accuracy Specifications:

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

Spot	Name	LLOQ			ULOQ		
		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	5220	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-23 Plate	Z0047651	2-8°C	30 Jun 2024
Anti-hu IL-23Antibody	D0081521	2-8°C	30 Nov 2023
TH17 Panel 1 (human) Calibrator Blend	A0080303	2-8°C	31 Aug 2026
Diluent 43	M0430046	≤ -10°C	29 Feb 2024
Diluent 3	M0030110	≤ -10°C	31 Mar 2024

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		17 MAY 2022