

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-12/IL-23p40 Kit

Kit Lot Number: K0081747

Expiration Date: 28 Feb 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	LLO			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	3650	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12/IL-23 p40 Plate	Z0047534	2-8°C	31 Oct 2023
Anti-hu IL12/23p40 Antibody	D0081362	2-8°C	28 Feb 2023
Cytokine Panel 1 (human) Calibrator Blend	A0080253	2-8°C	30 Apr 2025
Diluent 43	M0430044	≤ -10°C	30 Sep 2023
Diluent 3	M0030107	≤ -10°C	30 Sep 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		15 JUL 2021