

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus TH17 Panel 1 (human) Kit
Kit Lot Number: K0082041
Expiration Date: 31 May 2024
Kit Catalog Numbers: K15085-Series; K151XD-Series; K151WU-Series; K151XA-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-17A	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	IL-21	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
5	IL-31	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
6	IL-27	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
7	IL-23	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
8	IL-22	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
10	MIP-3α	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-17A	5.86	80% - 120%	< 20%	1950	80% - 120%	< 20%
4	IL-21	6.12	75% - 125%	< 25%	650	75% - 125%	< 25%
5	IL-31	4.22	80% - 120%	< 20%	650	80% - 120%	< 20%
6	IL-27	38.7	80% - 120%	< 20%	13000	80% - 120%	< 20%
7	IL-23	4.60	80% - 120%	< 20%	3250	80% - 120%	< 20%
8	IL-22	2.78	80% - 120%	< 20%	325	80% - 120%	< 20%
10	MIP-3α	0.750	80% - 120%	< 20%	325	80% - 120%	< 20%

Non-Specific Binding Specifications:

Spot	Name	NSB Spec
1	IL-17A	< 1.0%
4	IL-21	< 1.0%
5	IL-31	< 1.0%
6	IL-27	< 1.0%
7	IL-23	< 1.0%
8	IL-22	< 1.0%
10	MIP-3α	< 1.0%

CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	IL-17A	2580	pg/mL
4	IL-21	970	pg/mL
5	IL-31	948	pg/mL
6	IL-27	17700	pg/mL
7	IL-23	4910	pg/mL
8	IL-22	503	pg/mL
10	MIP-3 α	473	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TH17 Panel 1 (human) Plate	Z0047706	2-8°C	30 Nov 2024
Anti-hu IL-17A Antibody	D0081862	2-8°C	30 Nov 2025
Anti-hu IL-21 Antibody	D0081863	2-8°C	31 Oct 2025
Anti-hu IL-31 Antibody	D0081864	2-8°C	31 Oct 2025
Anti-hu IL-27 Antibody	D0081865	2-8°C	31 Oct 2025
Anti-hu IL-23 Antibody	D0081866	2-8°C	30 Nov 2025
Anti-hu IL-22 Antibody	D0081867	2-8°C	31 Dec 2025
Anti-hu MIP-3 α Antibody	D0081868	2-8°C	31 Oct 2025
TH17 Panel 1 (human) Calibrator Blend	A0080324	2-8°C	31 Mar 2027
Diluent 3	M0030114	$\leq -10^{\circ}\text{C}$	30 Sep 2024
Diluent 43	M0430047	$\leq -10^{\circ}\text{C}$	31 May 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		08 SEP 2022