

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX PLUS TH17 Panel 1 (human) Kit
Kit Lot Number: K0082272
Expiration Date: 31 Oct 2024
Kit Catalog Numbers: K151WU-Series; K151XA-Series; K151XD-Series; K15085-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-3 α	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-3 α	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-3 α	< 1.0%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-17A	2850	pg/mL
4	IL-21	1030	pg/mL
5	IL-31	1070	pg/mL
6	IL-27	17800	pg/mL
7	IL-23	5160	pg/mL
8	IL-22	494	pg/mL
10	MIP-3 α	495	pg/mL


CERTIFICATE OF ANALYSIS

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TH17 Panel 1 (human) SECTOR Plate	Z0047849	2-8°C	30 Sep 2025
Anti-hu IL-17A Antibody	D0082048	2-8°C	30 Sep 2026
Anti-hu IL-21 Antibody	D0082049	2-8°C	30 Sep 2026
Anti-hu IL-31 Antibody	D0082050	2-8°C	31 Oct 2026
Anti-hu IL-27 Antibody	D0082051	2-8°C	30 Sep 2026
Anti-hu IL-23 Antibody	D0082052	2-8°C	30 Sep 2026
Anti-hu IL-22 Antibody	D0082053	2-8°C	31 Oct 2026
Anti-hu MIP-3α Antibody	D0082054	2-8°C	30 Sep 2026
TH17 Panel 1 (human) Calibrator Blend	A0080354	2-8°C	31 Dec 2027
Diluent 3	M0030120	≤ -10°C	31 Jul 2025
Diluent 43	M0430049	≤ -10°C	31 Oct 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		01 JUN 2023