

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus TH17 Panel 1 (human) Kit  
**Kit Lot Number:** K0081636  
**Expiration Date:** 30 Jun 2023  
**Kit Catalog Numbers:** K15085-Series; K151XD-Series; K151XA-Series; K151WU-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 Gen. B	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%

\* See Additional Comments.

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-17A Gen. B	5.86	80% - 120%	< 20%	1950	80% - 120%	< 20%
4	IL-21	21.0	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 Gen. B	< 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 Gen. B	3110	pg/mL
4	IL-21	932	pg/mL
5	IL-31	970	pg/mL
6	IL-27	19900	pg/mL
7	IL-23	5110	pg/mL
8	IL-22	521	pg/mL
10	MIP-3 $\alpha$	506	pg/mL

**Lot-Specific Components:**


Description	Lot Number	Storage Temperature	Expiration Date
TH17 Panel 1 (human) Plate	Z0047457	2-8°C	30 Jun 2023
Anti-hu IL-17A Antibody	D0081613	2-8°C	31 Jul 2024
Anti-hu IL-21 Antibody	D0081614	2-8°C	31 Jul 2024
Anti-hu IL-31 Antibody	D0081631	2-8°C	31 Jul 2024
Anti-hu IL-27 Antibody	D0081616	2-8°C	31 Jul 2024
Anti-hu IL-23 Antibody	D0081617	2-8°C	30 Jun 2024
Anti-hu IL-22 Antibody	D0081618	2-8°C	31 Jul 2024
Anti-hu MIP-3 $\alpha$ Antibody	D0081537	2-8°C	31 Mar 2024
TH17 Panel 1 (human) Calibrator Blend	A0080263	2-8°C	30 Sep 2025
Diluent 43	M0430043	$\leq$ -10°C	31 Jul 2023
Diluent 3	M0030104	$\leq$ -10°C	30 Jun 2023

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

**Additional Comments:**

Assay	Spot	Specification	Comment
IL-21	4	Control Recovery = 75% - 125%	Control 1 = 64.7% Control 2 = 64.8% Control 3 = 53.3%

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

Name	Function	Signature	Date
Review/Approval Karen Hamilla	Quality		12 MAY 2021