

CERTIFICATE OF ANALYSIS

Page 1 of 4

Product Description: V-PLEX® and V-PLEX Plus TH17 Panel 1 (human) Kit
Kit Catalog Numbers: K15085-Series
Kit Lot Number: K0081422
Expiration Date: 28 Feb 2022

Acceptance Data/ Percent Recovery 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%

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	6.05*	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%

*See additional comments

Acceptance Data/ 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 Material Tested :

Spot	Name	Item	Conc	Units
1	IL-17A Gen B.	Calibrator	3140	pg/mL
		TH17 Panel 1 (human) Control 1	1980	pg/mL
		TH17 Panel 1 (human) Control 2	293	pg/mL
		TH17 Panel 1 (human) Control 3	42.5	pg/mL
4	IL-21	Calibrator	1180	pg/mL
		TH17 Panel 1 (human) Control 1	872	pg/mL
		TH17 Panel 1 (human) Control 2	161	pg/mL
		TH17 Panel 1 (human) Control 3	26.8	pg/mL
5	IL-31	Calibrator	964	pg/mL
		TH17 Panel 1 (human) Control 1	1080	pg/mL
		TH17 Panel 1 (human) Control 2	176	pg/mL
		TH17 Panel 1 (human) Control 3	23.9	pg/mL
6	IL-27	Calibrator	18200	pg/mL
		TH17 Panel 1 (human) Control 1	31000	pg/mL
		TH17 Panel 1 (human) Control 2	6560	pg/mL
		TH17 Panel 1 (human) Control 3	1380	pg/mL
7	IL-23	Calibrator	5540	pg/mL
		TH17 Panel 1 (human) Control 1	5360	pg/mL
		TH17 Panel 1 (human) Control 2	602	pg/mL
		TH17 Panel 1 (human) Control 3	69.8	pg/mL
8	IL-22	Calibrator	542	pg/mL
		TH17 Panel 1 (human) Control 1	526	pg/mL
		TH17 Panel 1 (human) Control 2	83.3	pg/mL
		TH17 Panel 1 (human) Control 3	15.4	pg/mL
10	MIP-3 α	Calibrator	508	pg/mL
		TH17 Panel 1 (human) Control 1	428	pg/mL
		TH17 Panel 1 (human) Control 2	175	pg/mL
		TH17 Panel 1 (human) Control 3	64.6	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-17A Gen B. - Spot 1			IL-21 - Spot 4			IL-31 - Spot 5		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
3140	539,000	3.0	1180	165,000	6.9	964	548,000	2.2
785	133,000	3.0	295	36,200	4.6	241	169,000	2.0
196	35,400	4.1	73.8	9,800	6.1	60.3	48,300	2.3
49.1	9,580	3.4	18.4	3,340	4.0	15.1	12,900	2.2
12.3	3,140	3.3	4.61	1,640	4.0	3.77	3,480	1.9
3.07	1,520	3.2	1.15	1,240	4.7	0.941	1,100	2.9
0.767	1,090	4.0	0.288	1,160	6.6	0.235	461	6.9
0.00	940	5.6	0.00	1,120	6.4	0.00	253	7.1

CERTIFICATE OF ANALYSIS

IL-27 - Spot 6		
Conc. pg/mL	ECL Signal	%CV
18200	176,000	3.3
4550	54,800	3.7
1140	16,500	2.8
284	5,590	3.1
71.1	2,930	2.9
17.8	2,150	3.1
4.44	1,970	2.9
0.00	1,890	2.4

IL-23 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
5540	431,000	6.9
1390	120,000	5.0
346	32,600	4.5
86.6	8,620	4.8
21.6	2,620	3.8
5.41	978	3.8
1.35	591	5.1
0.00	445	8.0

IL-22 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
542	394,000	2.8
136	114,000	2.6
33.9	32,000	4.7
8.47	8,710	5.9
2.12	2,900	4.0
0.529	1,410	7.6
0.132	1,070	20.0
0.00	757	9.7

MIP-3 α - Spot 10		
Conc. pg/mL	ECL Signal	%CV
508	333,000	5.5
127	77,000	4.3
31.8	19,500	6.2
7.94	5,230	4.7
1.98	1,830	3.6
0.496	914	6.9
0.124	684	9.4
0.00	580	5.5

Components:

Description	Lot Number	Storage Temperature	Expiration Date
TH17 Panel 1 (human) Plate	Z0047268	2-8°C	31 Jul 2022
Anti-hu IL-17A Antibody	D0081135	2-8°C	30 Apr 2022
Anti-hu IL-21 Antibody	D0081427	2-8°C	31 Jul 2023
Anti-hu IL-31 Antibody	D0081140	2-8°C	28 Feb 2022
Anti-hu IL-27 Antibody	D0081429	2-8°C	31 Jul 2023
Anti-hu IL-23 Antibody	D0081138	2-8°C	31 Mar 2022
Anti-hu IL-22 Antibody	D0081137	2-8°C	28 Feb 2022
Anti-hu MIP-3 α Antibody	D0081428	2-8°C	31 Jul 2023
TH17 Panel 1 (human) Calibrator Blend	A0080241	2-8°C	28 Feb 2025
Diluent 43	M0430036	≤ -10°C	31 May 2022
Diluent 3	M0030096	≤ -10°C	30 Jun 2022
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20X), 100mL	Not Kit Specific
TH17 Panel 1 (human) Control 1	Not Kit Specific
TH17 Panel 1 (human) Control 2	Not Kit Specific
TH17 Panel 1 (human) Control 3	Not Kit Specific

CERTIFICATE OF ANALYSIS


Page 4 of 4

Additional Comments:

	Assay	Spot	Specification	Comment
1	IL-21	4	LLOQ Conc.	Reported LLOQ 6.05 pg/mL

Statement:

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		27 MAR 2020