

CERTIFICATE OF ANALYSIS

Page 1 of 4

Product Description: V-PLEX® and V-PLEX Plus Chemokine Panel 1 (human) Kit
Kit Catalog Numbers: K15047-Series; K15055-Series; K151NO-Series; K151NP-Series; K151NS-Series; K156NO-Series; K156NP-Series
Kit Lot Number: K0081512
Expiration Date: 28 Feb 2022

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Chemokine Panel 1 (human) Control 1		Chemokine Panel 1 (human) Control 2		Chemokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Eotaxin	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	MIP-1β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
3	Eotaxin-3	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	TARC	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
5	IP-10	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
6	MIP-1α	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
7	IL-8	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
8	MCP-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
9	MDC	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
10	MCP-4	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Eotaxin	12.3	80% - 120%	< 20%	1120	80% - 120%	< 20%
2	MIP-1β	1.02	80% - 120%	< 20%	750	80% - 120%	< 20%
3	Eotaxin-3	10.2	80% - 120%	< 20%	3750	80% - 120%	< 20%
4	TARC	3.32	80% - 120%	< 20%	1120	80% - 120%	< 20%
5	IP-10	1.37	80% - 120%	< 20%	500	80% - 120%	< 20%
6	MIP-1α	13.8	80% - 120%	< 20%	743	80% - 120%	< 20%
7	IL-8	713	75% - 125%	< 20%	43400	75% - 125%	< 20%
8	MCP-1	1.09	80% - 120%	< 20%	375	80% - 120%	< 20%
9	MDC	88.3	80% - 120%	< 20%	6590	80% - 120%	< 20%
10	MCP-4	8.85	75% - 125%	< 25%	469	80% - 120%	< 20%

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	Eotaxin	< 1.0%
2	MIP-1β	< 1.0%
3	Eotaxin-3	< 1.0%
4	TARC	< 1.0%
5	IP-10	< 1.0%
6	MIP-1α	< 1.0%
7	IL-8 (HA)	< 1.0%
8	MCP-1	< 1.0%
9	MDC	< 1.0%
10	MCP-4	< 1.0%

CERTIFICATE OF ANALYSIS

Page 2 of 4

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Eotaxin	Calibrator	1880	pg/mL
		Chemokine Panel 1 (human) Control 1	1610	pg/mL
		Chemokine Panel 1 (human) Control 2	590	pg/mL
		Chemokine Panel 1 (human) Control 3	182	pg/mL
2	MIP-1 β	Calibrator	1270	pg/mL
		Chemokine Panel 1 (human) Control 1	1040	pg/mL
		Chemokine Panel 1 (human) Control 2	178	pg/mL
		Chemokine Panel 1 (human) Control 3	27.1	pg/mL
3	Eotaxin-3	Calibrator	4990	pg/mL
		Chemokine Panel 1 (human) Control 1	4300	pg/mL
		Chemokine Panel 1 (human) Control 2	615	pg/mL
		Chemokine Panel 1 (human) Control 3	95.7	pg/mL
4	TARC	Calibrator	1800	pg/mL
		Chemokine Panel 1 (human) Control 1	1340	pg/mL
		Chemokine Panel 1 (human) Control 2	223	pg/mL
		Chemokine Panel 1 (human) Control 3	36.9	pg/mL
5	IP-10	Calibrator	2340	pg/mL
		Chemokine Panel 1 (human) Control 1	561	pg/mL
		Chemokine Panel 1 (human) Control 2	85.6	pg/mL
		Chemokine Panel 1 (human) Control 3	14	pg/mL
6	MIP-1 α	Calibrator	1100	pg/mL
		Chemokine Panel 1 (human) Control 1	911	pg/mL
		Chemokine Panel 1 (human) Control 2	401	pg/mL
		Chemokine Panel 1 (human) Control 3	166	pg/mL
7	IL-8	Calibrator	78600	pg/mL
		Chemokine Panel 1 (human) Control 1	61100	pg/mL
		Chemokine Panel 1 (human) Control 2	21300	pg/mL
		Chemokine Panel 1 (human) Control 3	8370	pg/mL
8	MCP-1	Calibrator	553	pg/mL
		Chemokine Panel 1 (human) Control 1	465	pg/mL
		Chemokine Panel 1 (human) Control 2	71.3	pg/mL
		Chemokine Panel 1 (human) Control 3	11.5	pg/mL
9	MDC	Calibrator	9730	pg/mL
		Chemokine Panel 1 (human) Control 1	8960	pg/mL
		Chemokine Panel 1 (human) Control 2	2630	pg/mL
		Chemokine Panel 1 (human) Control 3	645	pg/mL
10	MCP-4	Calibrator	705	pg/mL
		Chemokine Panel 1 (human) Control 1	660	pg/mL
		Chemokine Panel 1 (human) Control 2	239	pg/mL
		Chemokine Panel 1 (human) Control 3	71.2	pg/mL

Controls tested may not be the same lot as shipped.

CERTIFICATE OF ANALYSIS

Page 3 of 4

Standard Curve Data:

Eotaxin - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1880	571,000	3.0
470	74,800	4.4
118	7,780	4.5
29.4	744	5.1
7.34	239	6.6
1.84	162	15.0
0.459	159	17.0
0.00	162	19.0

MIP-1 β - Spot 2		
Conc. pg/mL	ECL Signal	%CV
1270	1,380,000	3.5
318	309,000	3.3
79.4	51,500	2.0
19.8	7,500	3.6
4.96	1,440	3.8
1.24	393	6.1
0.310	190	8.4
0.00	142	22.0

Eotaxin-3 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
4990	200,000	3.0
1250	51,600	1.4
312	13,500	1.5
78.0	3,390	1.8
19.5	978	3.2
4.87	327	7.2
1.22	177	12.0
0.00	120	23.0

TARC - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1800	1,020,000	2.6
450	304,000	2.7
113	76,400	3.2
28.1	20,700	3.7
7.03	5,500	3.3
1.76	1,840	3.7
0.439	838	3.7
0.00	517	5.9

IP-10 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
2340	606,000	2.5
585	186,000	2.5
146	47,700	3.0
36.6	12,000	3.5
9.14	3,230	4.1
2.29	897	3.7
0.571	309	8.7
0.00	114	17.0

MIP-1 α - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1100	579,000	2.6
275	91,700	3.8
68.8	9,550	4.0
17.2	908	3.8
4.30	273	6.2
1.07	201	6.0
0.269	200	7.5
0.00	189	9.0

IL-8 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
78600	897,000	3.3
19700	218,000	6.1
4910	21,300	5.9
1230	2,050	4.0
307	468	6.3
76.8	224	13.0
19.2	165	14.0
0.00	155	20.0

MCP-1 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
553	791,000	4.1
138	203,000	4.1
34.6	49,900	3.2
8.64	12,700	2.9
2.16	3,400	5.9
0.540	1,010	4.1
0.135	376	9.0
0.00	165	18.0

MDC - Spot 9		
Conc. pg/mL	ECL Signal	%CV
9730	1,500,000	3.5
2430	172,000	3.0
608	26,300	5.5
152	5,410	4.3
38.0	1,510	4.8
9.50	500	4.6
2.38	242	15.0
0.00	163	14.0

CERTIFICATE OF ANALYSIS

Page 4 of 4

MCP-4 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
705	315,000	4.3
176	33,100	4.5
44.1	2,830	6.6
11.0	239	21.0
2.75	157	12.0
0.688	129	21.0
0.172	115	27.0
0.00	117	20.0


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0047359	2-8°C	31 Dec 2022
Anti-hu Eotaxin Antibody	D0081287	2-8°C	31 Dec 2022
Anti-hu MIP-1 β Antibody	D0081312	2-8°C	31 Jan 2023
Anti-hu Eotaxin-3 Antibody	D0081386	2-8°C	31 Mar 2023
Anti-hu TARC Antibody	D0081313	2-8°C	30 Nov 2022
Anti-hu IP-10 Antibody	D0081262	2-8°C	31 Oct 2022
Anti-hu MIP-1 α Antibody	D0081508	2-8°C	30 Sep 2023
Anti-hu IL-8 (HA) Antibody	D0081165	2-8°C	28 Feb 2022
Anti-hu MCP-1 Antibody	D0081388	2-8°C	31 Mar 2023
Anti-hu MDC Antibody	D0081507	2-8°C	30 Sep 2023
Anti-hu MCP-4 Antibody	D0081506	2-8°C	30 Sep 2023
Chemokine Panel 1 (human) Calibrator Blend	A0080259	2-8°C	31 May 2025
Diluent 43	M0430040	$\leq -10^{\circ}\text{C}$	30 Nov 2022
Diluent 3	M0030100	$\leq -10^{\circ}\text{C}$	31 Oct 2022
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

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

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		11 SEP 2020