

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, K156NO-Series, K156NP-Series, K151NP-Series, K151NS-Series
Kit Lot Number: K0081355
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%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 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%	7500	80% - 120%	< 20%
10	MCP-4	5.13	80% - 120%	< 20%	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	< 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	1790	pg/mL
		Chemokine Panel 1 (human) Control 1	1650	pg/mL
		Chemokine Panel 1 (human) Control 2	596	pg/mL
		Chemokine Panel 1 (human) Control 3	184	pg/mL
2	MIP-1 β	Calibrator	1280	pg/mL
		Chemokine Panel 1 (human) Control 1	1030	pg/mL
		Chemokine Panel 1 (human) Control 2	176	pg/mL
		Chemokine Panel 1 (human) Control 3	27.8	pg/mL
3	Eotaxin-3	Calibrator	5710	pg/mL
		Chemokine Panel 1 (human) Control 1	4250	pg/mL
		Chemokine Panel 1 (human) Control 2	593	pg/mL
		Chemokine Panel 1 (human) Control 3	94.6	pg/mL
4	TARC	Calibrator	1870	pg/mL
		Chemokine Panel 1 (human) Control 1	1300	pg/mL
		Chemokine Panel 1 (human) Control 2	216	pg/mL
		Chemokine Panel 1 (human) Control 3	37	pg/mL
5	IP-10	Calibrator	2370	pg/mL
		Chemokine Panel 1 (human) Control 1	589	pg/mL
		Chemokine Panel 1 (human) Control 2	86.8	pg/mL
		Chemokine Panel 1 (human) Control 3	14.3	pg/mL
6	MIP-1 α	Calibrator	1090	pg/mL
		Chemokine Panel 1 (human) Control 1	920	pg/mL
		Chemokine Panel 1 (human) Control 2	403	pg/mL
		Chemokine Panel 1 (human) Control 3	169	pg/mL
7	IL-8	Calibrator	66700	pg/mL
		Chemokine Panel 1 (human) Control 1	65000	pg/mL
		Chemokine Panel 1 (human) Control 2	22200	pg/mL
		Chemokine Panel 1 (human) Control 3	8800	pg/mL
8	MCP-1	Calibrator	544	pg/mL
		Chemokine Panel 1 (human) Control 1	468	pg/mL
		Chemokine Panel 1 (human) Control 2	71.4	pg/mL
		Chemokine Panel 1 (human) Control 3	12.1	pg/mL
9	MDC	Calibrator	9350	pg/mL
		Chemokine Panel 1 (human) Control 1	9830	pg/mL
		Chemokine Panel 1 (human) Control 2	2780	pg/mL
		Chemokine Panel 1 (human) Control 3	664	pg/mL
10	MCP-4	Calibrator	683	pg/mL
		Chemokine Panel 1 (human) Control 1	704	pg/mL
		Chemokine Panel 1 (human) Control 2	246	pg/mL
		Chemokine Panel 1 (human) Control 3	71.8	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
1790	762,000	4.8
448	106,000	6.2
112	10,800	7.9
28.0	907	13.0
6.99	209	11.0
1.75	140	16.0
0.437	125	15.0
0.00	126	22.0

MIP-1 β - Spot 2		
Conc. pg/mL	ECL Signal	%CV
1280	1,730,000	3.1
320	381,000	2.3
80.0	55,000	2.7
20.0	6,610	3.3
5.00	1,080	4.4
1.25	299	7.1
0.313	153	11.0
0.00	115	20.0

Eotaxin-3 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
5710	260,000	3.9
1430	71,000	2.6
357	18,000	3.3
89.2	4,420	3.0
22.3	1,240	3.6
5.58	373	5.6
1.39	164	13.0
0.00	90	29.0

TARC - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1870	1,010,000	4.5
468	287,000	4.7
117	70,000	4.7
29.2	16,700	7.4
7.30	4,720	6.7
1.83	1,490	12.0
0.457	718	7.7
0.00	389	30.0

IP-10 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
2370	800,000	3.2
593	306,000	2.4
148	82,600	3.6
37.0	21,700	3.5
9.26	5,640	3.2
2.31	1,500	4.0
0.579	440	4.8
0.00	103	17.0

MIP-1 α - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1090	691,000	3.8
273	109,000	3.4
68.1	10,700	6.8
17.0	852	6.0
4.26	203	7.8
1.06	153	17.0
0.266	140	12.0
0.00	130	18.0

IL-8 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
66700	805,000	12.0
16700	149,000	15.0
4170	13,200	16.0
1040	1,380	18.0
261	411	12.0
65.1	190	12.0
16.3	146	15.0
0.00	129	24.0

MCP-1 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
544	1,020,000	4.9
136	262,000	3.4
34.0	60,400	2.8
8.50	14,500	4.9
2.13	3,810	3.6
0.531	1,100	5.1
0.133	386	7.7
0.00	142	23.0

MDC - Spot 9		
Conc. pg/mL	ECL Signal	%CV
9350	2,070,000	4.3
2340	275,000	3.7
584	34,400	5.1
146	5,980	5.5
36.5	1,550	4.5
9.13	495	7.5
2.28	235	7.4
0.00	147	13.0

CERTIFICATE OF ANALYSIS

Page 4 of 4

MCP-4 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
683	611,000	3.5
171	75,000	5.0
42.7	6,640	7.1
10.7	530	16.0
2.67	211	9.2
0.667	168	14.0
0.167	164	13.0
0.00	157	19.0


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0047206	2-8°C	31 Mar 2022
Anti-hu Eotaxin Antibody	D0081287	2-8°C	31 Dec 2022
Anti-hu MIP-1 β Antibody	D0081181	2-8°C	31 Jul 2022
Anti-hu Eotaxin-3 Antibody	D0081260	2-8°C	31 Oct 2022
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	D0081263	2-8°C	31 Aug 2022
Anti-hu IL-8 (HA) Antibody	D0081165	2-8°C	28 Feb 2022
Anti-hu MCP-1 Antibody	D0081265	2-8°C	31 Aug 2022
Anti-hu MDC Antibody	D0081166	2-8°C	28 Feb 2022
Anti-hu MCP-4 Antibody	D0081125	2-8°C	31 Mar 2022
Chemokine Panel 1 (human) Calibrator Blend	A0080204	2-8°C	30 Apr 2024
Diluent 43	M0430035	\leq -10°C	31 Mar 2022
Diluent 3	M0030092	\leq -10°C	31 Mar 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
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		18 JUN 2020