

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
Kit Lot Number: K0081223
Expiration Date: 30 Sep 2021

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%	5850	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

Standard Curve Data:

Eotaxin - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1790	640,000	4.3
448	103,000	5.3
112	11,900	4.7
28.0	1,300	5.3
6.99	293	8.9
1.75	193	15.0
0.437	160	16.0
0.00	160	13.0

MIP-1β - Spot 2		
Conc. pg/mL	ECL Signal	%CV
1280	1,470,000	5.0
320	358,000	3.1
80.0	60,500	2.9
20.0	9,360	3.4
5.00	1,830	3.7
1.25	520	4.7
0.313	226	5.7
0.00	141	17.0

Eotaxin-3 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
5710	232,000	9.6
1430	71,000	5.6
357	17,800	6.3
89.2	4,660	4.6
22.3	1,360	4.3
5.58	454	7.2
1.39	208	10.0
0.00	141	22.0

TARC - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1870	820,000	4.3
468	252,000	4.5
117	61,700	4.2
29.2	16,700	5.4
7.30	4,840	4.9
1.83	1,840	3.8
0.457	1,040	7.4
0.00	806	7.3

IP-10 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
2370	614,000	3.6
593	272,000	3.3
148	81,600	3.5
37.0	22,000	3.0
9.26	5,870	2.5
2.31	1,540	3.1
0.579	484	5.0
0.00	117	14.0

MIP-1α - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1090	610,000	3.3
273	101,000	3.3
68.1	10,800	3.8
17.0	1,040	4.6
4.26	263	6.8
1.06	195	10.0
0.266	176	9.8
0.00	180	7.9

IL-8 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
66700	789,000	4.3
16700	164,000	6.4
4170	14,300	8.4
1040	1,580	6.4
261	449	7.5
65.1	263	9.0
16.3	212	9.9
0.00	209	13.0

MCP-1 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
544	838,000	3.8
136	213,000	4.8
34.0	51,100	3.7
8.50	12,600	4.3
2.13	3,400	3.0
0.531	1,010	4.6
0.133	411	5.6
0.00	193	12.0

MDC - Spot 9		
Conc. pg/mL	ECL Signal	%CV
9350	1,730,000	2.7
2340	240,000	4.7
584	31,000	2.5
146	5,700	4.4
36.5	1,540	5.1
9.13	552	5.6
2.28	313	7.1
0.00	233	8.9

MCP-4 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
683	609,000	6.2
171	83,700	6.9
42.7	8,380	7.5
10.7	854	7.0
2.67	262	8.0
0.667	187	8.9
0.167	183	12.0
0.00	173	16.0

CERTIFICATE OF ANALYSIS

Page 4 of 4


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0047063	2-8°C	30 Sep 2021
Anti-hu Eotaxin Antibody	D0081259	2-8°C	31 Aug 2022
Anti-hu MIP-1β Antibody	D0081181	2-8°C	31 Jul 2022
Anti-hu Eotaxin-3 Antibody	D0081260	2-8°C	31 Oct 2022
Anti-hu TARC Antibody	D0081261	2-8°C	31 Aug 2022
Anti-hu IP-10 Antibody	D0081262	2-8°C	31 Oct 2022
Anti-hu MIP-1α Antibody	D0081263	2-8°C	31 Aug 2022
Anti-hu IL-8 (HA) Antibody	D0081264	2-8°C	31 Aug 2022
Anti-hu MCP-1 Antibody	D0081265	2-8°C	31 Aug 2022
Anti-hu MDC Antibody	D0081266	2-8°C	31 Aug 2022
Anti-hu MCP-4 Antibody	D0081267	2-8°C	31 Dec 2022
Chemokine Panel 1 (human) Calibrator Blend	A0080204	2-8°C	30 Apr 2024
Diluent 43	M0430032	≤ -10°C	30 Sep 2021
Diluent 3	M0030089	≤ -10°C	30 Sep 2021
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		28 AUG 2019