

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: K0081202
Expiration Date: 30 Jun 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%	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
Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Eotaxin	Calibrator	1960	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	1240	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	5960	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	1800	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	2480	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	72100	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	537	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	10200	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	751	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
1960	583,000	7.4
490	108,000	7.5
123	13,000	11.0
30.6	1,350	10.0
7.66	293	8.9
1.91	186	13.0
0.479	162	14.0
0.00	169	15.0

MIP-1 β - Spot 2		
Conc. pg/mL	ECL Signal	%CV
1240	1,450,000	4.8
310	346,000	4.5
77.5	56,700	2.6
19.4	7,760	4.9
4.84	1,400	3.5
1.21	399	5.7
0.303	205	11.0
0.00	158	17.0

Eotaxin-3 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
5960	203,000	5.6
1490	64,700	4.5
373	17,700	3.9
93.1	4,410	5.7
23.3	1,250	5.0
5.82	407	7.2
1.46	205	11.0
0.00	130	15.0

TARC - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1800	691,000	4.4
450	218,000	3.6
113	51,100	5.7
28.1	12,700	4.1
7.03	3,430	3.5
1.76	1,260	4.4
0.439	641	7.0
0.00	443	8.6

IP-10 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
2480	569,000	5.1
620	329,000	4.2
155	101,000	4.2
38.8	26,800	5.3
9.69	7,060	3.5
2.42	1,870	4.8
0.605	571	7.5
0.00	126	21.0

MIP-1 α - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1090	499,000	5.3
273	76,000	6.3
68.1	7,980	7.1
17.0	810	7.2
4.26	279	9.8
1.06	222	18.0
0.266	216	23.0
0.00	207	17.0

IL-8 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
72100	654,000	8.3
18000	81,600	11.0
4510	6,550	9.9
1130	797	12.0
282	326	8.0
70.4	246	25.0
17.6	224	32.0
0.00	198	19.0

MCP-1 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
537	1,030,000	5.1
134	254,000	6.0
33.6	57,700	6.1
8.39	13,000	7.6
2.10	3,610	6.0
0.524	1,160	6.4
0.131	547	12.0
0.00	338	16.0

MDC - Spot 9		
Conc. pg/mL	ECL Signal	%CV
10200	1,680,000	5.4
2550	232,000	4.8
638	28,900	5.4
159	4,950	4.7
39.8	1,290	5.3
9.96	504	12.0
2.49	320	26.0
0.00	243	20.0

CERTIFICATE OF ANALYSIS

MCP-4 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
751	974,000	4.1
188	124,000	5.6
46.9	12,500	7.4
11.7	1,220	12.0
2.93	286	10.0
0.733	240	29.0
0.183	234	39.0
0.00	207	22.0


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0047039	2-8°C	31 Aug 2021
Anti-hu Eotaxin Antibody	D0081259	2-8°C	31 Aug 2022
Anti-hu MIP-1β Antibody	D0081162	2-8°C	28 Feb 2022
Anti-hu Eotaxin-3 Antibody	D0081122	2-8°C	31 Jan 2022
Anti-hu TARC Antibody	D0081261	2-8°C	31 Aug 2022
Anti-hu IP-10 Antibody	D0081227	2-8°C	31 May 2022
Anti-hu MIP-1α Antibody	D0081123	2-8°C	31 Jan 2022
Anti-hu IL-8 (HA) Antibody	D0081165	2-8°C	28 Feb 2022
Anti-hu MCP-1 Antibody	D0081124	2-8°C	31 Jan 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	A0080183	2-8°C	31 Oct 2023
Diluent 43	M0430030	≤ -10°C	30 Jun 2021
Diluent 3	M0030087	≤ -10°C	30 Jun 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	Carissa Hubbard	Quality		18 JUN 2019