

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: K0081444

Expiration Date: 30 Jun 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%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
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	1.88	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	6.27	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	1870	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	1330	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	5970	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	1830	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	2360	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	1060	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	64500	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	552	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	724	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
1870	718,000	3.9
468	103,000	4.6
117	10,600	7.7
29.2	1,040	7.7
7.30	256	9.0
1.83	173	13.0
0.457	150	20.0
0.00	157	19.0

MIP-1 β - Spot 2		
Conc. pg/mL	ECL Signal	%CV
1330	1,890,000	3.5
333	418,000	4.0
83.1	64,600	3.8
20.8	8,970	3.2
5.20	1,790	3.9
1.30	488	4.7
0.325	240	17.0
0.00	143	19.0

Eotaxin-3 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
5970	292,000	4.1
1490	75,900	3.6
373	18,900	4.2
93.3	4,610	3.3
23.3	1,270	3.4
5.83	406	7.3
1.46	213	21.0
0.00	128	25.0

TARC - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1830	1,060,000	4.2
458	311,000	2.8
114	75,800	4.1
28.6	19,500	3.0
7.15	5,270	5.4
1.79	1,760	2.9
0.447	805	8.0
0.00	479	7.3

IP-10 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
2360	803,000	2.3
590	317,000	2.9
148	87,200	4.4
36.9	22,100	3.4
9.22	5,900	3.4
2.30	1,530	4.5
0.576	503	12.0
0.00	138	29.0

MIP-1 α - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1060	743,000	1.8
265	121,000	3.1
66.3	12,600	3.5
16.6	1,170	2.3
4.14	274	6.0
1.04	186	10.0
0.259	214	33.0
0.00	227	32.0

IL-8 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
64500	849,000	4.6
16100	142,000	10.0
4030	12,900	9.4
1010	1,630	6.6
252	485	7.7
63.0	233	10.0
15.7	221	31.0
0.00	269	46.0

MCP-1 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
552	1,000,000	4.4
138	268,000	4.4
34.5	61,700	3.6
8.63	15,100	4.2
2.16	4,050	4.4
0.539	1,290	2.1
0.135	629	11.0
0.00	444	20.0

MDC - Spot 9		
Conc. pg/mL	ECL Signal	%CV
10200	1,880,000	3.1
2550	242,000	3.4
638	33,300	4.0
159	6,370	5.1
39.8	1,730	6.8
9.96	588	4.8
2.49	370	31.0
0.00	302	43.0

CERTIFICATE OF ANALYSIS

MCP-4 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
724	560,000	3.3
181	66,100	3.3
45.3	5,900	7.6
11.3	637	8.8
2.83	284	16.0
0.707	228	9.5
0.177	363	54.0
0.00	373	45.0


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0047276	2-8°C	31 Jul 2022
Anti-hu Eotaxin Antibody	D0081259	2-8°C	31 Aug 2022
Anti-hu MIP-1β Antibody	D0081312	2-8°C	31 Jan 2023
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-1α Antibody	D0081387	2-8°C	31 Mar 2023
Anti-hu IL-8 (HA) Antibody	D0081264	2-8°C	31 Aug 2022
Anti-hu MCP-1 Antibody	D0081388	2-8°C	31 Mar 2023
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)	A0080228	2-8°C	31 Aug 2024
Diluent 43	M0430037	≤ -10°C	31 Aug 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
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		22 MAY 2020