

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: K0081263
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%	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

Standard Curve Data:

Eotaxin - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1790	660,000	4.5
448	97,600	5.8
112	11,000	6.9
28.0	1,130	6.3
6.99	252	10.0
1.75	157	17.0
0.437	140	18.0
0.00	133	15.0

MIP-1β - Spot 2		
Conc. pg/mL	ECL Signal	%CV
1280	1,570,000	5.6
320	366,000	4.5
80.0	57,800	2.8
20.0	8,540	4.1
5.00	1,530	4.7
1.25	430	4.6
0.313	188	9.2
0.00	130	17.0

Eotaxin-3 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
5710	220,000	4.3
1430	59,700	4.3
357	15,200	3.4
89.2	3,860	4.7
22.3	1,060	4.6
5.58	356	7.0
1.39	164	16.0
0.00	105	25.0

TARC - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1870	979,000	3.9
468	282,000	4.7
117	71,100	5.0
29.2	18,100	3.9
7.30	5,110	5.8
1.83	1,690	5.8
0.457	853	6.2
0.00	564	13.0

IP-10 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
2370	637,000	2.9
593	251,000	3.3
148	67,800	4.3
37.0	17,500	3.9
9.26	4,550	5.6
2.31	1,210	5.3
0.579	367	6.7
0.00	94	19.0

MIP-1α - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1090	621,000	3.0
273	105,000	3.9
68.1	11,000	3.7
17.0	1,020	4.8
4.26	246	8.4
1.06	169	10.0
0.266	159	9.8
0.00	161	14.0

IL-8 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
66700	829,000	6.7
16700	167,000	12.0
4170	16,700	14.0
1040	1,650	9.0
261	463	11.0
65.1	233	13.0
16.3	190	12.0
0.00	179	16.0

MCP-1 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
544	947,000	5.0
136	231,000	2.4
34.0	52,300	3.3
8.50	12,900	4.2
2.13	3,180	5.5
0.531	943	4.8
0.133	319	7.7
0.00	132	15.0

MDC - Spot 9		
Conc. pg/mL	ECL Signal	%CV
9350	1,600,000	2.9
2340	216,000	3.6
584	27,600	5.2
146	5,150	5.0
36.5	1,330	5.1
9.13	473	4.5
2.28	253	6.5
0.00	189	11.0

MCP-4 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
683	565,000	3.2
171	74,500	4.5
42.7	7,320	6.6
10.7	727	7.9
2.67	218	11.0
0.667	151	17.0
0.167	139	18.0
0.00	148	16.0

CERTIFICATE OF ANALYSIS

Page 4 of 4


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0047111	2-8°C	30 Nov 2021
Anti-hu Eotaxin Antibody	D0081259	2-8°C	31 Aug 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	D0081261	2-8°C	31 Aug 2022
Anti-hu IP-10 Antibody	D0081227	2-8°C	31 May 2022
Anti-hu MIP-1 α 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	$\leq -10^{\circ}\text{C}$	30 Sep 2021
Diluent 3	M0030090	$\leq -10^{\circ}\text{C}$	31 Dec 2021
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
MSD Wash Buffer (20X), 100 mL	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		29 AUG 2019