

CERTIFICATE OF ANALYSIS

Page 1 of 5

Product Description: V-PLEX® and V-PLEX Plus Chemokine Panel 1 (human) Kit

Kit Catalog Numbers: K15047-Series

Kit Lot Number: K0080919

Expiration Date: 31 Jul 2019

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	hu Eotaxin	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	hu MIP-1β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
3	hu Eotaxin-3	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	hu TARC	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
5	hu IP-10	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
6	hu MIP-1α	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
7	hu IL-8	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
8	hu MCP-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
9	hu MDC	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
10	hu 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	hu Eotaxin	12.3	80% - 120%	< 20%	1120	80% - 120%	< 20%
2	hu MIP-1β	2.27	80% - 120%	< 20%	750	80% - 120%	< 20%
3	hu Eotaxin-3	10.2	80% - 120%	< 20%	3750	80% - 120%	< 20%
4	hu TARC	3.32	80% - 120%	< 20%	1120	80% - 120%	< 20%
5	hu IP-10	1.37	80% - 120%	< 20%	500	80% - 120%	< 20%
6	hu MIP-1α	13.8	80% - 120%	< 20%	743	80% - 120%	< 20%
7	hu IL-8	713	75% - 125%	< 20%	43400	75% - 125%	< 20%
8	hu MCP-1	1.09	80% - 120%	< 20%	375	80% - 120%	< 20%
9	hu MDC	88.3	80% - 120%	< 20%	7500	80% - 120% *	< 20%
10	hu MCP-4	5.13	80% - 120%	< 20%	469	80% - 120%	< 20%

* See Additional Comments.

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	hu Eotaxin	< 1.0%
2	hu MIP-1β	< 1.0%
3	hu Eotaxin-3	< 1.0%
4	hu TARC	< 1.0%
5	hu IP-10	< 1.0%
6	hu MIP-1α	< 1.0%
7	hu IL-8	< 1.0%
8	hu MCP-1	< 1.0%
9	hu MDC	< 1.0%
10	hu MCP-4	< 1.0%

CERTIFICATE OF ANALYSIS

Page 2 of 5

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu Eotaxin	Calibrator	1860	pg/mL
		Chemokine Panel 1 (human) Control 1	1660	pg/mL
		Chemokine Panel 1 (human) Control 2	330	pg/mL
		Chemokine Panel 1 (human) Control 3	180	pg/mL
2	hu MIP-1 β	Calibrator	1050	pg/mL
		Chemokine Panel 1 (human) Control 1	980	pg/mL
		Chemokine Panel 1 (human) Control 2	170	pg/mL
		Chemokine Panel 1 (human) Control 3	30	pg/mL
3	hu Eotaxin-3	Calibrator	5410	pg/mL
		Chemokine Panel 1 (human) Control 1	4050	pg/mL
		Chemokine Panel 1 (human) Control 2	580	pg/mL
		Chemokine Panel 1 (human) Control 3	90	pg/mL
4	hu TARC	Calibrator	1630	pg/mL
		Chemokine Panel 1 (human) Control 1	1460	pg/mL
		Chemokine Panel 1 (human) Control 2	220	pg/mL
		Chemokine Panel 1 (human) Control 3	40	pg/mL
5	hu IP-10	Calibrator	2420	pg/mL
		Chemokine Panel 1 (human) Control 1	590	pg/mL
		Chemokine Panel 1 (human) Control 2	90	pg/mL
		Chemokine Panel 1 (human) Control 3	16	pg/mL
6	hu MIP-1 α	Calibrator	1030	pg/mL
		Chemokine Panel 1 (human) Control 1	910	pg/mL
		Chemokine Panel 1 (human) Control 2	410	pg/mL
		Chemokine Panel 1 (human) Control 3	160	pg/mL
7	hu IL-8	Calibrator	86200	pg/mL
		Chemokine Panel 1 (human) Control 1	66500	pg/mL
		Chemokine Panel 1 (human) Control 2	22500	pg/mL
		Chemokine Panel 1 (human) Control 3	8800	pg/mL
8	hu MCP-1	Calibrator	686	pg/mL
		Chemokine Panel 1 (human) Control 1	610	pg/mL
		Chemokine Panel 1 (human) Control 2	160	pg/mL
		Chemokine Panel 1 (human) Control 3	16	pg/mL
9	hu MDC	Calibrator	12200	pg/mL
		Chemokine Panel 1 (human) Control 1	11600	pg/mL
		Chemokine Panel 1 (human) Control 2	3300	pg/mL
		Chemokine Panel 1 (human) Control 3	1010	pg/mL
10	hu MCP-4	Calibrator	787	pg/mL
		Chemokine Panel 1 (human) Control 1	710	pg/mL
		Chemokine Panel 1 (human) Control 2	270	pg/mL
		Chemokine Panel 1 (human) Control 3	82	pg/mL

Controls tested may not be the same lot as shipped.

CERTIFICATE OF ANALYSIS

Standard Curve Data:

hu Eotaxin - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1860	845,000	3.0
465	122,000	4.6
116	13,900	4.4
29.1	1,350	7.0
7.27	262	7.7
1.82	144	12.0
0.454	112	19.0
0.00	117	19.0

hu MIP-1β - Spot 2		
Conc. pg/mL	ECL Signal	%CV
1050	1,530,000	3.4
263	404,000	3.5
65.6	77,300	4.0
16.4	12,800	3.7
4.10	2,410	5.1
1.03	610	4.3
0.256	228	8.9
0.00	111	16.0

hu Eotaxin-3 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
5410	118,000	3.7
1350	34,600	3.9
338	9,300	5.2
84.5	2,370	5.4
21.1	704	6.0
5.28	255	6.6
1.32	135	17.0
0.00	100	24.0

hu TARC - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1630	876,000	3.8
408	269,000	2.7
102	67,400	2.3
25.5	17,600	3.7
6.37	4,860	3.6
1.59	1,740	5.2
0.398	857	5.8
0.00	585	8.9

hu IP-10 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
2420	516,000	4.2
605	299,000	3.6
151	92,500	3.8
37.8	24,400	3.2
9.45	6,320	3.3
2.36	1,700	4.3
0.591	489	4.5
0.00	93	24.0

hu MIP-1α - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1030	658,000	2.5
258	116,000	2.5
64.4	13,100	2.3
16.1	1,160	3.5
4.02	238	6.1
1.01	151	10.0
0.251	138	12.0
0.00	132	13.0

hu IL-8 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
86200	759,000	5.0
21600	119,000	9.3
5390	11,400	8.4
1350	1,500	9.0
337	469	7.4
84.2	249	11.0
21.0	206	11.0
0.00	176	9.5

hu MCP-1 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
686	888,000	5.7
172	232,000	4.9
42.9	53,200	4.1
10.7	12,300	6.3
2.68	3,050	6.5
0.670	991	9.6
0.167	365	18.0
0.00	184	30.0

hu MDC - Spot 9		
Conc. pg/mL	ECL Signal	%CV
12200	2,040,000	2.7
3050	273,000	4.4
763	33,600	3.7
191	5,780	4.1
47.7	1,470	2.6
11.9	484	6.4
2.98	232	7.7
0.00	149	13.0

CERTIFICATE OF ANALYSIS

Page 4 of 5

hu MCP-4 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
787	965,000	3.0
197	133,000	4.2
49.2	15,000	4.6
12.3	1,360	7.1
3.07	249	7.8
0.769	148	13.0
0.192	132	15.0
0.00	126	20.0

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0046694	2-8°C	30 Apr 2020
Anti-hu Eotaxin Antibody	D0080959	2-8°C	30 Sep 2019
Anti-hu MIP-1 β Antibody	D0080991	2-8°C	31 Dec 2019
Anti-hu Eotaxin-3 Antibody	D0080899	2-8°C	31 Jul 2019
Anti-hu TARC Antibody	D0080993	2-8°C	31 Oct 2019
Anti-hu IP-10 Antibody	D0080901	2-8°C	31 Jul 2019
Anti-hu MIP-1 α Antibody	D0080995	2-8°C	31 Oct 2019
Anti-hu IL-8 (HA) Antibody	D0080903	2-8°C	31 Jul 2019
Anti-hu MCP-1 Antibody	D0080904	2-8°C	31 Jul 2019
Anti-hu MDC Antibody	D0080905	2-8°C	31 Jul 2019
Anti-hu MCP-4 Antibody	D0080999	2-8°C	31 May 2020
Chemokine Panel 1 (human) Calibrator Blend	A0080131	2-8°C	31 Jul 2022
Diluent 43	M0430024	\leq -10°C	31 Jan 2020
Diluent 3	M0030078	\leq -10°C	31 Jan 2020
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

CERTIFICATE OF ANALYSIS

Page 5 of 5

Additional Comments:

	Assay	Spot	Specification	Comment
1	hu MDC	9	ULOQ Conc.	ULOQ recovery at 120.9% (Specifications 80-120%)

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	10 MAY 2018