

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

Kit Lot Number: K0080811

Expiration Date: 30 Jun 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

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu Eotaxin	Calibrator	1860	pg/mL
		Chemokine Panel 1 (human) Control 1	1370	pg/mL
		Chemokine Panel 1 (human) Control 2	518	pg/mL
		Chemokine Panel 1 (human) Control 3	160	pg/mL
2	hu MIP-1 β	Calibrator	1050	pg/mL
		Chemokine Panel 1 (human) Control 1	797	pg/mL
		Chemokine Panel 1 (human) Control 2	130	pg/mL
		Chemokine Panel 1 (human) Control 3	22.4	pg/mL
3	hu Eotaxin-3	Calibrator	5410	pg/mL
		Chemokine Panel 1 (human) Control 1	4760	pg/mL
		Chemokine Panel 1 (human) Control 2	755	pg/mL
		Chemokine Panel 1 (human) Control 3	118	pg/mL
4	hu TARC	Calibrator	1630	pg/mL
		Chemokine Panel 1 (human) Control 1	1130	pg/mL
		Chemokine Panel 1 (human) Control 2	197	pg/mL
		Chemokine Panel 1 (human) Control 3	34.2	pg/mL
5	hu IP-10	Calibrator	2420	pg/mL
		Chemokine Panel 1 (human) Control 1	348	pg/mL
		Chemokine Panel 1 (human) Control 2	57.3	pg/mL
		Chemokine Panel 1 (human) Control 3	10.3	pg/mL
6	hu MIP-1 α	Calibrator	1030	pg/mL
		Chemokine Panel 1 (human) Control 1	824	pg/mL
		Chemokine Panel 1 (human) Control 2	371	pg/mL
		Chemokine Panel 1 (human) Control 3	152	pg/mL
7	hu IL-8	Calibrator	86200	pg/mL
		Chemokine Panel 1 (human) Control 1	171000	pg/mL
		Chemokine Panel 1 (human) Control 2	58100	pg/mL
		Chemokine Panel 1 (human) Control 3	22500	pg/mL
8	hu MCP-1	Calibrator	686	pg/mL
		Chemokine Panel 1 (human) Control 1	420	pg/mL
		Chemokine Panel 1 (human) Control 2	66.6	pg/mL
		Chemokine Panel 1 (human) Control 3	11.1	pg/mL
9	hu MDC	Calibrator	12200	pg/mL
		Chemokine Panel 1 (human) Control 1	11100	pg/mL
		Chemokine Panel 1 (human) Control 2	3460	pg/mL
		Chemokine Panel 1 (human) Control 3	907	pg/mL
10	hu MCP-4	Calibrator	787	pg/mL
		Chemokine Panel 1 (human) Control 1	635	pg/mL
		Chemokine Panel 1 (human) Control 2	232	pg/mL
		Chemokine Panel 1 (human) Control 3	82.2	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	947,000	5.2
465	183,000	2.4
116	25,400	4.0
29.1	2,480	4.0
7.27	346	7.1
1.82	148	18.0
0.454	118	21.0
0.00	111	17.0

hu MIP-1β - Spot 2		
Conc. pg/mL	ECL Signal	%CV
1050	1,620,000	3.2
263	433,000	3.1
65.6	85,000	2.1
16.4	14,400	4.0
4.10	3,140	5.7
1.03	797	4.5
0.256	280	7.6
0.00	106	16.0

hu Eotaxin-3 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
5410	142,000	4.8
1350	45,500	3.9
338	12,200	5.1
84.5	3,000	5.5
21.1	901	5.7
5.28	303	7.9
1.32	159	14.0
0.00	96	25.0

hu TARC - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1630	794,000	3.8
408	251,000	2.8
102	59,400	3.1
25.5	15,900	3.9
6.37	4,130	4.9
1.59	1,500	3.6
0.398	721	5.3
0.00	476	8.9

hu IP-10 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
2420	519,000	3.3
605	332,000	3.1
151	111,000	3.1
37.8	29,300	3.8
9.45	7,840	2.7
2.36	2,080	2.5
0.591	586	3.8
0.00	95	21.0

hu MIP-1α - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1030	594,000	3.1
258	91,600	3.2
64.4	9,240	3.2
16.1	712	5.9
4.02	209	7.9
1.01	146	15.0
0.251	137	12.0
0.00	132	13.0

hu IL-8 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
86200	841,000	3.6
21600	123,000	9.0
5390	12,200	12.0
1350	1,500	8.5
337	498	6.0
84.2	238	11.0
21.0	182	13.0
0.00	166	16.0

hu MCP-1 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
686	983,000	4.1
172	257,000	2.6
42.9	59,200	4.9
10.7	14,300	5.0
2.68	3,680	4.8
0.670	1,040	4.7
0.167	357	6.6
0.00	138	20.0

hu MDC - Spot 9		
Conc. pg/mL	ECL Signal	%CV
12200	1,880,000	4.0
3050	352,000	6.2
763	41,900	5.8
191	6,090	4.5
47.7	1,440	3.5
11.9	443	6.0
2.98	209	8.3
0.00	139	13.0

hu MCP-4 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
787	1,260,000	1.9
197	226,000	3.6
49.2	29,900	5.0
12.3	2,660	6.4
3.07	356	6.3
0.769	148	20.0
0.192	125	17.0
0.00	121	18.0

CERTIFICATE OF ANALYSIS

Page 4 of 4

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0046614	2-8°C	31 Oct 2019
Anti-hu Eotaxin Antibody	D0080959	2-8°C	30 Sep 2019
Anti-hu MIP-1β Antibody	D0080898	2-8°C	30 Sep 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-1α Antibody	D0080902	2-8°C	31 Jul 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	D0080867	2-8°C	30 Jun 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080131	2-8°C	31 Jul 2022
Diluent 43	M0430021	≤ -10°C	30 Sep 2019
Diluent 3	M0030077	≤ -10°C	30 Nov 2019
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

Additional Comments:

	Assay	Spot	Specification	Comment
1	hu IP-10	5	ULOQ % Recovery	Observed 123.5% (Specifications 80-120%)
2	hu IP-10	5	ULOQ CV	Observed 20.6% (specifications NMT 20%)

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		23 AUG 2017