



CERTIFICATE OF ANALYSIS

Page 1 of 4

COA-04154

Product Description: Chemokine Panel 1 (human) Kit V-PLEX™ and V-PLEX Plus
Kit Catalog Numbers: K15047D-1, K15047D-2, K15047D-4, K15047G-1, K15047G-2, K15047G-4
Kit Lot Number: K0080125
Expiration Date: 31Oct2015

Acceptance Data:

			Assay				
Metric		Specifications	hu Eotaxin	hu MIP-1β	hu Eotaxin-3	hu TARC	hu IP-10
% Recovery Concentration	Calibrators	80% - 120%	Pass	Pass	Pass	Pass	Pass
	Controls	80% - 120%	Pass	Pass	Pass	1	Pass
	LLOQ	80% - 120%	Pass	Pass	Pass	Pass	Pass
	ULOQ	80% - 120%	Pass	Pass	Pass	Pass	Pass
% Recovery CV	Calibrators	<20%	Pass	Pass	Pass	Pass	Pass
	Controls	<20%	Pass	Pass	Pass	Pass	Pass
	LLOQ	<20%	Pass	Pass	Pass	Pass	Pass
	ULOQ	<20%	Pass	Pass	Pass	Pass	Pass
Non-specific binding		<1%	Pass	Pass	Pass	Pass	Pass

			Assay			
Metric		Specifications	hu MIP-1α	hu MCP-1	hu MDC	hu MCP-4
% Recovery Concentration	Calibrators	80% - 120%	Pass	Pass	Pass	Pass
	Controls	80% - 120%	Pass	2	3	Pass
	LLOQ	80% - 120%	Pass	Pass	Pass	Pass
	ULOQ	80% - 120%	Pass	Pass	Pass	Pass
% Recovery CV	Calibrators	<20%	Pass	Pass	Pass	Pass
	Controls	<20%	Pass	Pass	Pass	Pass
	LLOQ	<20%	Pass	Pass	Pass	Pass
	ULOQ	<20%	Pass	Pass	Pass	Pass
Non-specific binding		<1%	Pass	Pass	Pass	Pass

Metric		Specifications	hu IL-8
% Recovery Concentration	Calibrators	80% - 120%	Pass
	Controls	80% - 120%	Pass
	LLOQ	75% - 125%	Pass
	ULOQ	75% - 125%	Pass
% Recovery CV	Calibrators	<20%	Pass
	Controls	<20%	Pass
	LLOQ	<20%	Pass
	ULOQ	<20%	Pass
Non-specific binding		<1%	Pass

1. Pass (Control 1, Control 2); Control 3 = 78.5%
2. Pass (Control 1, Control 2); Control 3 = 77.4%
3. Pass (Control 1, Control 2); Control 3 = 78.1%



CERTIFICATE OF ANALYSIS

Page 2 of 4

COA-04154

Concentration of Materials Tested:

Units = pg/mL	Assay									
	hu Eotaxin	hu MIP-1 β	hu Eotaxin-3	hu TARC	hu IP-10	hu MIP-1 α	hu IL-8	hu MCP-1	hu MDC	hu MCP-4
Reconstituted Calibrator	1490	1080	4930	1470	2610	1030	82 000	486	9780	625
Reconstituted Control 1 (A00C0190)	1150	678	3710	1050	1570	773	14 8000	412	6720	516
Reconstituted Control 2 (A00C0191)	297	175	914	260	385	189	52 400	98.9	1400	135
Reconstituted Control 3 (A00C0192)	68.5	36.3	208	61.1	94.8	48.2	N/A	25.7	349	31.8
LLOQ	12.2	2.23	10.2	3.04	1.37	2.13	673	1.01	20.2	5.13
ULOQ	1150	833	3800	1130	500	795	63300	375	7550	482



CERTIFICATE OF ANALYSIS

COA-04154

Standard Curve Data from Kit QC:

hu Eotaxin		
Conc., (pg/ml)	Average Signal	%CV
1490	663 497	3.0
373	136 035	3.3
93.1	19 768	5.3
23.3	2351	4.6
5.82	364	5.1
1.46	153	11.1
0.364	131	24.9
0	118	16.0

hu MIP-1 β		
Conc., (pg/ml)	Average Signal	%CV
1080	1 304 876	3.0
270	276 302	3.8
67.5	38 484	4.2
16.9	5018	3.6
4.22	895	4.2
1.05	260	9.7
0.264	159	25.2
0	105	20.7

hu Eotaxin-3		
Conc., (pg/ml)	Average Signal	%CV
4930	319 412	4.3
1233	55 027	4.8
308	14 400	6.1
77.0	4026	6.0
19.3	1178	4.1
4.81	367	8.6
1.20	184	22.3
0	99	26.9

hu TARC		
Conc., (pg/ml)	Average Signal	%CV
1470	621 339	2.7
368	180 857	4.6
91.9	42 664	4.3
23.0	11 209	5.3
5.74	3124	5.9
1.44	1031	6.0
0.359	436	9.2
0	257	12.8

hu IP-10		
Conc., (pg/ml)	Average Signal	%CV
2610	559 394	7.4
653	331 056	5.9
163	106 145	6.3
40.8	31 197	5.4
10.2	8852	4.2
2.55	2435	5.1
0.637	699	4.3
0	87	16.4

hu MIP-1 α		
Conc., (pg/ml)	Average Signal	%CV
1030	747 895	2.4
258	120 821	4.1
64.4	14 208	4.5
16.1	1768	4.3
4.02	455	6.1
1.01	245	6.3
0.251	210	12.0
0	197	8.7

hu IL-8		
Conc., (pg/ml)	Average Signal	%CV
82 000	721 364	2.9
20 500	134 073	10.9
5125	13 285	6.7
1281	2204	6.4
320	631	4.9
80.1	270	9.5
20.0	188	15.9
0	160	17.2

hu MCP-1		
Conc., (pg/ml)	Average Signal	%CV
486	747 383	4.1
122	172 876	5.7
30.4	38 421	7.1
7.59	9972	4.9
1.90	2646	6.7
0.475	803	8.3
0.119	302	6.9
0	141	16.2

hu MDC		
Conc., (pg/ml)	Average Signal	%CV
9780	1 107 106	4.4
2445	277 253	4.8
611	42 801	6.7
153	6645	4.6
38.2	1571	3.0
9.55	482	5.2
2.39	231	8.7
0	148	14.6

hu MCP-4		
Conc., (pg/ml)	Average Signal	%CV
625	727 208	4.0
156	115 183	7.8
39.1	14 458	7.8
9.77	1 475	5.7
2.44	242	9.3
0.610	114	19.9
0.153	97	17.7
0	100	25.4



CERTIFICATE OF ANALYSIS

COA-04154

Chemokine Panel 1 (human) Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0045125	2-8°C	30Jun2016
Anti-hu Eotaxin Antibody	D0080187	2-8°C	30Apr2016
Anti-hu MIP-1β Antibody	D0080188	2-8°C	30Apr2016
Anti-hu Eotaxin-3 Antibody	D0080189	2-8°C	30Apr2016
Anti-hu TARC Antibody	D0080190	2-8°C	30Apr2016
Anti-hu IP-10 Antibody	D0080191	2-8°C	30Apr2016
Anti-hu MIP-1α Antibody	D0080192	2-8°C	30Apr2016
Anti-hu IL-8 (HA) Antibody	D0080193	2-8°C	30Apr2016
Anti-hu MCP-1 Antibody	D0080194	2-8°C	30Apr2016
Anti-hu MDC Antibody	D0080195	2-8°C	30Apr2016
Anti-hu MCP-4 Antibody	D0080196	2-8°C	30Apr2016
Chemokine Panel 1 (human) Calibrator Blend	A0080016	2-8°C	31Oct2015
Diluent 43	M0430005	≤-10°C	31Mar2016
Diluent 3	M0030057	≤-10°C	31Oct2015
MSD Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A
Additional Components Included in V-PLEX Plus			
Wash Buffer (20X), 100mL	Not Kit Specific	Room Temperature	N/A
Plate Seals, pack of 3	Not Kit Specific	Room Temperature	N/A
Chemokine Panel 1 (human) Control 1	Not Kit Specific; See Separate Control Pack CoA		
Chemokine Panel 1 (human) Control 2			
Chemokine Panel 1 (human) Control 3			

Statement: The above kit lot was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Caitlin Hansen	Quality Control		11Feb2014