



CERTIFICATE OF ANALYSIS

Page 1 of 4

COA-04431

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: K0080135
Expiration Date: 30 SEP 2015

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	Pass	Pass
	ULOQ	80% - 120%	Pass	Pass	Pass	Pass	Pass
	LLOQ	80% - 120%	Pass	Pass	Pass	Pass	Pass
% Recovery CV	Calibrators	< 20%	Pass	Pass	Pass	Pass	Pass
	Controls	< 20%	Pass	Pass	Pass	Pass	Pass
	ULOQ	< 20%	Pass	Pass	Pass	Pass	Pass
	LLOQ	< 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	Pass	Pass	Pass
	ULOQ	80% - 120%	Pass	Pass	Pass	Pass
	LLOQ	80% - 120%	Pass	Pass	Pass	Pass
% Recovery CV	Calibrators	< 20%	Pass	Pass	Pass	Pass
	Controls	< 20%	Pass	Pass	Pass	Pass
	ULOQ	< 20%	Pass	Pass	Pass	Pass
	LLOQ	< 20%	Pass	Pass	Pass	Pass
Non-specific binding		< 1%	Pass	Pass	Pass	Pass

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



CERTIFICATE OF ANALYSIS

Page 2 of 4

COA-04431

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	1380	1020	4370	1430	2380	979	73 900	471	9500	588
Reconstituted Control 1 (A00C0190)	1150	678	3710	1050	1570	773	148 000	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
ULOQ	1180	875	3750	1230	500	840	63 400	404	8150	505
LLOQ	12.0	2.27	9.73	3.18	1.37	8.54	645	1.05	21.1	5.13



CERTIFICATE OF ANALYSIS

COA-04431

Standard Curve Data from Kit QC:

hu Eotaxin		
Conc., (pg/mL)	Average Signal	%CV
1380	569 783	3.5
345	132 957	3.3
86.3	18 773	5.2
21.6	2378	4.4
5.39	584	6.9
1.35	391	9.7
0.337	360	13.2
0	356	10.8

hu MIP-1β		
Conc., (pg/mL)	Average Signal	%CV
1020	1 132 142	4.8
255	286 242	2.8
63.8	46 315	3.2
15.9	6613	3.7
3.98	1199	5.3
0.996	327	7.7
0.249	156	12.1
0	98	15.2

hu Eotaxin-3		
Conc., (pg/mL)	Average Signal	%CV
4370	203 563	5.5
1093	47 430	5.1
273	12 497	3.6
68.3	3350	6.0
17.1	1048	5.8
4.27	390	10.1
1.07	201	12.7
0	127	13.9

hu TARC		
Conc., (pg/mL)	Average Signal	%CV
1430	540 600	2.4
358	172 493	2.8
89.4	41 616	3.2
22.3	11 333	3.7
5.59	3349	3.6
1.40	1323	4.5
0.349	754	7.2
0	558	5.9

hu IP-10		
Conc., (pg/mL)	Average Signal	%CV
2380	724 517	9.7
595	414 340	3.1
149	134 847	3.7
37.2	38 733	4.5
9.30	11 080	4.0
2.32	3035	4.4
0.581	917	5.0
0	100	12.7

hu MIP-1α		
Conc., (pg/mL)	Average Signal	%CV
979	681 662	2.5
245	125 976	2.9
61.2	15 684	2.7
15.3	2070	2.9
3.82	597	3.3
0.956	353	5.3
0.239	307	4.7
0	282	4.7

hu IL-8		
Conc., (pg/mL)	Average Signal	%CV
73 900	786 396	3.9
18 475	148 781	8.5
4619	15 999	6.6
1155	2660	5.6
289	754	7.1
72.2	284	10.9
18.0	171	11.8
0	143	24.1

hu MCP-1		
Conc., (pg/mL)	Average Signal	%CV
471	522 536	8.2
118	121 664	8.0
29.4	27 068	13.1
7.36	6953	9.3
1.84	1920	9.8
0.460	634	12.6
0.115	285	11.5
0	149	18.9

hu MDC		
Conc., (pg/mL)	Average Signal	%CV
9500	1 224 357	3.8
2375	317 668	4.7
594	46 800	4.5
148	7278	3.8
37.1	1651	4.5
9.28	510	4.6
2.32	251	10.2
0	158	12.1

hu MCP-4		
Conc., (pg/mL)	Average Signal	%CV
588	973 764	2.5
147	163 860	4.6
36.8	19 322	4.5
9.19	1974	4.7
2.30	303	6.8
0.574	137	27.2
0.144	117	21.2
0	126	24.5



CERTIFICATE OF ANALYSIS

COA-04431

Chemokine Panel 1 (human) Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0045237	2-8°C	30 SEP 2016
Anti-hu Eotaxin Antibody	D0080302	2-8°C	31 AUG 2016
Anti-hu MIP-1β Antibody	D0080303	2-8°C	31 AUG 2016
Anti-hu Eotaxin-3 Antibody	D0080304	2-8°C	31 AUG 2016
Anti-hu TARC Antibody	D0080305	2-8°C	31 AUG 2016
Anti-hu IP-10 Antibody	D0080306	2-8°C	31 AUG 2016
Anti-hu MIP-1α Antibody	D0080307	2-8°C	31 AUG 2016
Anti-hu IL-8 (HA) Antibody	D0080308	2-8°C	31 AUG 2016
Anti-hu MCP-1 Antibody	D0080309	2-8°C	31 AUG 2016
Anti-hu MDC Antibody	D0080310	2-8°C	31 AUG 2016
Anti-hu MCP-4 Antibody	D0080311	2-8°C	31 AUG 2016
Chemokine Panel 1 (human) Calibrator Blend	A0080037	2-8°C	30 SEP 2016
Diluent 43	M0430007	≤ -10°C	30 APR 2016
Diluent 3	M0030055	≤ -10°C	30 SEP 2015
MSD Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A
Additional Components Included in V-PLEX Plus			
Wash Buffer (20X), 100 mL	Not Kit Specific	Room Temperature	N/A
Plate Seals	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		19 JUN 2014