



CERTIFICATE OF ANALYSIS

Page 1 of 4

COA-04781

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: K0080216
Expiration Date: 30 APR 2016

Acceptance Data:

Metric		Specifications	Assay				
			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

Metric		Specifications	Assay			
			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

Metric		Specifications	Assay
			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-04781

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	1510	1050	5020	1470	2650	1040	93 700	488	10 100	645
Reconstituted Control 1 (A00C0267)	1040	603	3720	971	1470	709	161 000	382	5790	475
Reconstituted Control 2 (A00C0268)	283	156	897	235	378	188	52 300	88.9	1330	131
Reconstituted Control 3 (A00C0269)	65.6	36.7	226	63.4	93.5	42.8	16 100	22.7	333	32.2
ULOQ	1160	807	3860	1130	500	799	72 000	375	7760	496
LLOQ	11.5	2.13	10.2	2.99	1.37	7.91	713	0.992	20.5	4.91



CERTIFICATE OF ANALYSIS

COA-04781

Standard Curve Data from Kit QC:

hu Eotaxin		
Conc., (pg/ml)	Average Signal	%CV
1510	606 737	3.6
378	126 304	3.1
94.4	17 683	3.9
23.6	2039	3.3
5.90	415	7.3
1.47	252	11.5
0.369	249	10.4
0	244	12.5

hu MIP-1 β		
Conc., (pg/ml)	Average Signal	%CV
1050	1 178 388	3.4
263	255 925	2.4
65.6	36 233	3.9
16.4	4608	3.3
4.10	717	5.1
1.03	220	8.5
0.256	144	12.1
0	118	14.3

hu Eotaxin-3		
Conc., (pg/ml)	Average Signal	%CV
5020	201 746	4.8
1255	48 665	5.3
314	12 817	6.6
78.4	3411	4.3
19.6	973	4.5
4.90	335	8.7
1.23	189	10.4
0	136	18.8

hu TARC		
Conc., (pg/ml)	Average Signal	%CV
1470	649 004	2.7
368	189 199	3.8
91.9	43 239	3.9
23.0	11 007	3.6
5.74	2874	4.1
1.44	940	5.9
0.359	440	8.7
0	264	11.6

hu IP-10		
Conc., (pg/ml)	Average Signal	%CV
2650	527 067	6.6
663	334 406	3.9
166	109 372	2.7
41.4	30 525	2.7
10.4	7955	3.9
2.59	2126	3.5
0.647	619	5.9
0	81	20.0

hu MIP-1 α		
Conc., (pg/ml)	Average Signal	%CV
1040	651 701	2.3
260	108 512	1.8
65.0	11 835	3.8
16.3	1380	2.4
4.06	434	4.9
1.02	311	5.9
0.254	292	6.3
0	286	5.8

hu IL-8		
Conc., (pg/ml)	Average Signal	%CV
93 700	1 432 765	4.9
23 425	205 788	10.1
5856	13 905	9.2
1464	2083	4.1
366	539	7.0
91.5	208	10.0
22.9	140	16.5
0	119	17.9

hu MCP-1		
Conc., (pg/ml)	Average Signal	%CV
488	543 102	6.0
122	123 412	5.1
30.5	27 320	5.8
7.63	6884	5.5
1.91	1750	6.0
0.477	527	9.2
0.119	218	9.7
0	121	20.6

hu MDC		
Conc., (pg/ml)	Average Signal	%CV
10 100	1 460 445	4.1
2525	368 005	6.3
631	51 414	5.1
158	6739	6.0
39.5	1345	3.8
9.86	397	6.2
2.47	200	8.7
0	140	14.5

hu MCP-4		
Conc., (pg/ml)	Average Signal	%CV
645	945 142	2.7
161	137 408	5.1
40.3	15 191	5.4
10.1	1384	4.9
2.52	215	11.0
0.630	118	16.5
0.157	105	21.1
0	101	20.7



CERTIFICATE OF ANALYSIS

COA-04781

Chemokine Panel 1 (human) Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0045260	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	D0080448	2-8°C	31 JAN 2017
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	A0080044	2-8°C	31 JAN 2017
Diluent 43	M0430007	≤ -10°C	30 APR 2016
Diluent 3	M0030062	≤ -10°C	31 OCT 2016
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	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		23 SEP 2014