



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Chemokine Panel 1 (human) Kit
Catalog Number(s) / Size(s): K15047-Series; K15055-Series; K156NO-Series; K151NO-Series
Lot Number: K0081091

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081091 has been extended to 31 OCT 2020.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Anti-hu Eotaxin Antibody	D0080990	30 APR 2020	31 OCT 2020
Anti-hu IP-10 Antibody	D0080994	31 Dec 2019	31 OCT 2020
Anti-hu MIP-1 α Antibody	D0080995	31 Oct 2019	31 OCT 2020
Anti-hu IL-8 (HA) Antibody	D0080996	31 Oct 2019	31 OCT 2020
Anti-hu MCP-1 Antibody	D0080997	31 Oct 2019	31 OCT 2020
Anti-hu MDC Antibody	D0080998	31 Oct 2019	31 OCT 2020
Anti-hu MCP-4 Antibody	D0081068	31 Jul 2020	31 OCT 2020

Function	Name	Signature	Date
Quality	Karen Hamilla		10 JUN 2019

This memo supersedes all other statements regarding date of expiration for this product.

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; K15055-Series; K156NO-Series; K151NO-Series
Kit Lot Number: K0081091
Expiration Date: 31 Oct 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	Eotaxin	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	MIP-1 β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
3	Eotaxin-3	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	TARC	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
5	IP-10	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
6	MIP-1 α	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
7	IL-8	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
8	MCP-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
9	MDC	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
10	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	Eotaxin	12.3	80% - 120%	< 20%	1120	80% - 120%	< 20%
2	MIP-1 β	1.02	80% - 120%	< 20%	750	80% - 120%	< 20%
3	Eotaxin-3	10.2	80% - 120%	< 20%	3750	80% - 120%	< 20%
4	TARC	3.32	80% - 120%	< 20%	1120	80% - 120%	< 20%
5	IP-10	1.37	80% - 120%	< 20%	500	80% - 120%	< 20%
6	MIP-1 α	13.8	80% - 120%	< 20%	743	80% - 120%	< 20%
7	IL-8	713	75% - 125%	< 20%	43400	75% - 125%	< 20%
8	MCP-1	1.09	80% - 120%	< 20%	375	80% - 120%	< 20%
9	MDC	88.3	80% - 120%	< 20%	7500	80% - 120%*	< 20%
10	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	Eotaxin	< 1.0%
2	MIP-1 β	< 1.0%
3	Eotaxin-3	< 1.0%
4	TARC	< 1.0%
5	IP-10	< 1.0%
6	MIP-1 α	< 1.0%
7	IL-8	< 1.0%
8	MCP-1	< 1.0%
9	MDC	< 1.0%
10	MCP-4	< 1.0%

CERTIFICATE OF ANALYSIS

Page 2 of 5

Stock Concentration of Material Tested:

Spot	Name	Item	Conc	Units
1	Eotaxin	Calibrator	1790	pg/mL
		Chemokine Panel 1 (human) Control 1	1650	pg/mL
		Chemokine Panel 1 (human) Control 2	596	pg/mL
		Chemokine Panel 1 (human) Control 3	184	pg/mL
2	MIP-1 β	Calibrator	1180	pg/mL
		Chemokine Panel 1 (human) Control 1	1030	pg/mL
		Chemokine Panel 1 (human) Control 2	176	pg/mL
		Chemokine Panel 1 (human) Control 3	27.8	pg/mL
3	Eotaxin-3	Calibrator	5180	pg/mL
		Chemokine Panel 1 (human) Control 1	4250	pg/mL
		Chemokine Panel 1 (human) Control 2	593	pg/mL
		Chemokine Panel 1 (human) Control 3	94.6	pg/mL
4	TARC	Calibrator	1700	pg/mL
		Chemokine Panel 1 (human) Control 1	1300	pg/mL
		Chemokine Panel 1 (human) Control 2	216	pg/mL
		Chemokine Panel 1 (human) Control 3	37	pg/mL
5	IP-10	Calibrator	2090	pg/mL
		Chemokine Panel 1 (human) Control 1	589	pg/mL
		Chemokine Panel 1 (human) Control 2	86.8	pg/mL
		Chemokine Panel 1 (human) Control 3	14.3	pg/mL
6	MIP-1 α	Calibrator	1060	pg/mL
		Chemokine Panel 1 (human) Control 1	920	pg/mL
		Chemokine Panel 1 (human) Control 2	403	pg/mL
		Chemokine Panel 1 (human) Control 3	169	pg/mL
7	IL-8	Calibrator	62200	pg/mL
		Chemokine Panel 1 (human) Control 1	65000	pg/mL
		Chemokine Panel 1 (human) Control 2	22200	pg/mL
		Chemokine Panel 1 (human) Control 3	8800	pg/mL
8	MCP-1	Calibrator	540	pg/mL
		Chemokine Panel 1 (human) Control 1	468	pg/mL
		Chemokine Panel 1 (human) Control 2	71.4	pg/mL
		Chemokine Panel 1 (human) Control 3	12.1	pg/mL
9	MDC	Calibrator	9980	pg/mL
		Chemokine Panel 1 (human) Control 1	9830	pg/mL
		Chemokine Panel 1 (human) Control 2	2780	pg/mL
		Chemokine Panel 1 (human) Control 3	664	pg/mL
10	MCP-4	Calibrator	714	pg/mL
		Chemokine Panel 1 (human) Control 1	704	pg/mL
		Chemokine Panel 1 (human) Control 2	246	pg/mL
		Chemokine Panel 1 (human) Control 3	71.8	pg/mL

Controls tested may not be the same lot as shipped.

CERTIFICATE OF ANALYSIS

Standard Curve Data:

Eotaxin - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1790	903,000	3.5
448	146,000	6.3
112	17,100	8.4
28.0	1,670	9.8
6.99	267	7.7
1.75	122	12.0
0.437	103	18.0
0.00	106	17.0

MIP-1 β - Spot 2		
Conc. pg/mL	ECL Signal	%CV
1180	1,590,000	3.2
295	358,000	2.4
73.8	54,500	3.6
18.4	7,950	4.7
4.61	1,550	3.9
1.15	427	3.8
0.288	184	12.0
0.00	112	20.0

Eotaxin-3 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
5180	262,000	4.0
1300	74,200	6.3
324	19,200	4.4
80.9	4,930	4.4
20.2	1,370	4.9
5.06	403	7.6
1.26	163	12.0
0.00	80	21.0

TARC - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1700	946,000	2.9
425	263,000	5.2
106	63,000	4.1
26.6	16,400	4.7
6.64	4,170	4.4
1.66	1,350	4.0
0.415	593	4.2
0.00	322	9.9

IP-10 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
2090	780,000	3.4
523	320,000	3.0
131	92,500	3.1
32.7	24,800	3.1
8.16	6,420	4.7
2.04	1,670	3.4
0.510	494	5.1
0.00	90	22.0

MIP-1 α - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1060	627,000	1.9
265	96,500	1.6
66.3	9,730	3.7
16.6	997	5.9
4.14	290	9.1
1.04	195	6.7
0.259	188	7.2
0.00	181	14.0

IL-8 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
62200	801,000	4.3
15600	149,000	7.9
3890	14,600	13.0
972	2,020	5.3
243	541	7.9
60.7	256	8.5
15.2	198	12.0
0.00	179	18.0

MCP-1 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
540	837,000	5.2
135	219,000	3.1
33.8	53,200	4.8
8.44	13,600	5.2
2.11	3,450	4.7
0.527	1,010	4.1
0.132	347	7.6
0.00	162	18.0

MDC - Spot 9		
Conc. pg/mL	ECL Signal	%CV
9980	2,280,000	3.4
2500	323,000	7.0
624	39,000	6.9
156	7,040	7.0
39.0	1,710	4.8
9.75	522	5.9
2.44	240	8.5
0.00	183	24.0

CERTIFICATE OF ANALYSIS

MCP-4 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
714	595,000	5.5
179	73,700	7.0
44.6	6,930	11.0
11.2	678	19.0
2.79	180	12.0
0.697	134	15.0
0.174	121	15.0
0.00	157	27.0

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0046916	2-8°C	28 Feb 2021
Anti-hu Eotaxin Antibody	D0080990	2-8°C	30 Apr 2020
Anti-hu MIP-1β Antibody	D0081121	2-8°C	31 Jan 2022
Anti-hu Eotaxin-3 Antibody	D0081122	2-8°C	31 Jan 2022
Anti-hu TARC Antibody	D0081090	2-8°C	31 Oct 2021
Anti-hu IP-10 Antibody	D0080994	2-8°C	31 Dec 2019
Anti-hu MIP-1α Antibody	D0080995	2-8°C	31 Oct 2019
Anti-hu IL-8 (HA) Antibody	D0080996	2-8°C	31 Oct 2019
Anti-hu MCP-1 Antibody	D0080997	2-8°C	31 Oct 2019
Anti-hu MDC Antibody	D0080998	2-8°C	31 Oct 2019
Anti-hu MCP-4 Antibody	D0081068	2-8°C	31 Jul 2020
Chemokine Panel 1 (human) Calibrator Blend	A0080171	2-8°C	30 Jun 2023
Diluent 43	M0430027	≤ -10°C	31 Dec 2020
Diluent 3	M0030084	≤ -10°C	31 Jan 2021
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), 100 mL	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	MDC	9	ULOQ Conc.	ULOQ Recovery 121.2% (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	Karen Hamilla	Quality		10 JUN 2019