



**RFE MEMO**


Page 1 of 1

**Product Description(s):** Anti-hu Eotaxin-3 Antibody

**Catalog Number(s) / Size(s):** D21NU-2

**Lot Number:** D0080992

Based on additional testing, the original date of expiration for this detection antibody has been extended by 12 months. The expiration date of D0080992 has been extended to 31 OCT 2020.

<b>Function</b>	<b>Name</b>	<b>Signature</b>	<b>Date</b>
<b>Quality</b>	Dawn Stollar		06 JUN 2019

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

## 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, K15055-Series; K156NO-Series; K151NO-Series  
**Kit Lot Number:** K0081148  
**Expiration Date:** 30 Apr 2020

### 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-1β	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-1α	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-1β	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-1α	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%	5830	80% - 120%	< 20%
10	MCP-4	5.13	80% - 120%	< 20%	469	80% - 120%	< 20%

### Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	Eotaxin	< 1.0%
2	MIP-1β	< 1.0%
3	Eotaxin-3	< 1.0%
4	TARC	< 1.0%
5	IP-10	< 1.0%
6	MIP-1α	< 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 4

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Eotaxin	Calibrator	1960	pg/mL
		Chemokine Panel 1 (human) Control 1	1550	pg/mL
		Chemokine Panel 1 (human) Control 2	594	pg/mL
		Chemokine Panel 1 (human) Control 3	182	pg/mL
2	MIP-1 $\beta$	Calibrator	1240	pg/mL
		Chemokine Panel 1 (human) Control 1	991	pg/mL
		Chemokine Panel 1 (human) Control 2	172	pg/mL
		Chemokine Panel 1 (human) Control 3	25.5	pg/mL
3	Eotaxin-3	Calibrator	5960	pg/mL
		Chemokine Panel 1 (human) Control 1	4400	pg/mL
		Chemokine Panel 1 (human) Control 2	654	pg/mL
		Chemokine Panel 1 (human) Control 3	99.6	pg/mL
4	TARC	Calibrator	1800	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	35.4	pg/mL
5	IP-10	Calibrator	2480	pg/mL
		Chemokine Panel 1 (human) Control 1	544	pg/mL
		Chemokine Panel 1 (human) Control 2	86.4	pg/mL
		Chemokine Panel 1 (human) Control 3	14.2	pg/mL
6	MIP-1 $\alpha$	Calibrator	1090	pg/mL
		Chemokine Panel 1 (human) Control 1	926	pg/mL
		Chemokine Panel 1 (human) Control 2	404	pg/mL
		Chemokine Panel 1 (human) Control 3	162	pg/mL
7	IL-8	Calibrator	72100	pg/mL
		Chemokine Panel 1 (human) Control 1	63500	pg/mL
		Chemokine Panel 1 (human) Control 2	21900	pg/mL
		Chemokine Panel 1 (human) Control 3	8970	pg/mL
8	MCP-1	Calibrator	537	pg/mL
		Chemokine Panel 1 (human) Control 1	465	pg/mL
		Chemokine Panel 1 (human) Control 2	71.7	pg/mL
		Chemokine Panel 1 (human) Control 3	12.1	pg/mL
9	MDC	Calibrator	10200	pg/mL
		Chemokine Panel 1 (human) Control 1	8970	pg/mL
		Chemokine Panel 1 (human) Control 2	2710	pg/mL
		Chemokine Panel 1 (human) Control 3	907	pg/mL
10	MCP-4	Calibrator	751	pg/mL
		Chemokine Panel 1 (human) Control 1	650	pg/mL
		Chemokine Panel 1 (human) Control 2	246	pg/mL
		Chemokine Panel 1 (human) Control 3	79.2	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
1960	929,000	4.2
490	179,000	4.0
123	23,000	7.3
30.6	2,130	6.8
7.66	292	8.8
1.91	126	15.0
0.479	102	18.0
0.00	101	19.0

MIP-1 $\beta$ - Spot 2		
Conc. pg/mL	ECL Signal	%CV
1240	1,310,000	7.3
310	353,000	5.7
77.5	59,200	5.6
19.4	9,010	4.9
4.84	1,660	6.4
1.21	440	6.3
0.303	175	8.9
0.00	101	19.0

Eotaxin-3 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
5960	149,000	11.0
1490	41,500	11.0
373	11,200	8.0
93.1	2,920	6.0
23.3	817	8.0
5.82	266	10.0
1.46	133	16.0
0.00	92	17.0

TARC - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1800	746,000	5.8
450	217,000	6.9
113	55,800	4.7
28.1	13,400	8.1
7.03	4,050	6.4
1.76	1,320	13.0
0.439	674	10.0
0.00	463	13.0

IP-10 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
2480	535,000	3.6
620	321,000	4.1
155	100,000	2.9
38.8	27,000	3.0
9.69	7,060	3.6
2.42	1,870	5.3
0.605	543	6.1
0.00	104	16.0

MIP-1 $\alpha$ - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1090	883,000	2.1
273	160,000	2.2
68.1	17,100	3.3
17.0	1,580	4.8
4.26	278	5.8
1.06	148	13.0
0.266	128	13.0
0.00	128	17.0

IL-8 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
72100	837,000	6.4
18000	113,000	8.6
4510	9,440	8.7
1130	1,250	7.6
282	407	5.5
70.4	193	8.2
17.6	141	16.0
0.00	129	27.0

MCP-1 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
537	930,000	5.1
134	235,000	5.2
33.6	52,900	4.5
8.39	12,600	7.0
2.10	3,100	4.3
0.524	910	6.7
0.131	319	7.6
0.00	143	21.0

MDC - Spot 9		
Conc. pg/mL	ECL Signal	%CV
10200	1,790,000	5.0
2550	263,000	3.9
638	34,200	6.2
159	6,250	4.0
39.8	1,580	6.0
9.96	508	5.8
2.49	233	8.8
0.00	169	23.0

# CERTIFICATE OF ANALYSIS

MCP-4 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
751	900,000	3.8
188	110,000	5.0
46.9	10,700	8.6
11.7	922	11.0
2.93	214	8.9
0.733	131	19.0
0.183	117	18.0
0.00	123	26.0


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0046956	2-8°C	31 May 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	D0080992	2-8°C	31 Oct 2020
Anti-hu TARC Antibody	D0081163	2-8°C	28 Feb 2022
Anti-hu IP-10 Antibody	D0081227	2-8°C	31 May 2022
Anti-hu MIP-1α Antibody	D0081123	2-8°C	31 Jan 2022
Anti-hu IL-8 (HA) Antibody	D0081165	2-8°C	28 Feb 2022
Anti-hu MCP-1 Antibody	D0081124	2-8°C	31 Jan 2022
Anti-hu MDC Antibody	D0081166	2-8°C	28 Feb 2022
Anti-hu MCP-4 Antibody	D0081125	2-8°C	31 Mar 2022
Chemokine Panel 1 (human) Calibrator Blend	A0080183	2-8°C	31 Oct 2023
Diluent 43	M0430029	≤ -10°C	30 Apr 2021
Diluent 3	M0030085	≤ -10°C	28 Feb 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), 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

### 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		06 JUN 2019