

CERTIFICATE OF ANALYSIS

Page 1 of 4

Product Description: V-PLEX® and V-PLEX Plus Chemokine Panel 1 (NHP) Kit

Kit Catalog Numbers: K15055-Series

Kit Lot Number: K0080872

Expiration Date: 30 Jun 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
2	NHP MIP-1 β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
3	NHP Eotaxin-3	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	NHP TARC	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
5	NHP IP-10	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
6	NHP MIP-1 α	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
7	NHP IL-8	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
8	NHP MCP-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
9	NHP MDC	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
10	NHP MCP-4	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
2	NHP MIP-1 β	2.27	80% - 120%	< 20%	750	80% - 120%	< 20%
3	NHP Eotaxin-3	10.2	80% - 120%	< 20%	3750	80% - 120%	< 20%
4	NHP TARC	3.32	80% - 120%	< 20%	1120	80% - 120%	< 20%
5	NHP IP-10	1.37	80% - 120%	< 20%	500	80% - 120%	< 20%
6	NHP MIP-1 α	13.8	80% - 120%	< 20%	743	80% - 120%	< 20%
7	NHP IL-8	713	75% - 125%	< 20%	43400	75% - 125%	< 20%
8	NHP MCP-1	1.09	80% - 120%	< 20%	375	80% - 120%	< 20%
9	NHP MDC	88.3	80% - 120%	< 20%	7500	80% - 120%*	< 20%
10	NHP MCP-4	5.13	80% - 120%	< 20%	469	80% - 120%	< 20%

* See Additional Comments.

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
2	NHP MIP-1 β	< 1.0%
3	NHP Eotaxin-3	< 1.0%
4	NHP TARC	< 1.0%
5	NHP IP-10	< 1.0%
6	NHP MIP-1 α	< 1.0%
7	NHP IL-8	< 1.0%
8	NHP MCP-1	< 1.0%
9	NHP MDC	< 1.0%
10	NHP MCP-4	< 1.0%

CERTIFICATE OF ANALYSIS
Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
2	NHP MIP-1 β	Calibrator	1050	pg/mL
		Chemokine Panel 1 (human) Control 1	797	pg/mL
		Chemokine Panel 1 (human) Control 2	130	pg/mL
		Chemokine Panel 1 (human) Control 3	22.4	pg/mL
3	NHP Eotaxin-3	Calibrator	5410	pg/mL
		Chemokine Panel 1 (human) Control 1	4760	pg/mL
		Chemokine Panel 1 (human) Control 2	755	pg/mL
		Chemokine Panel 1 (human) Control 3	118	pg/mL
4	NHP TARC	Calibrator	1630	pg/mL
		Chemokine Panel 1 (human) Control 1	1130	pg/mL
		Chemokine Panel 1 (human) Control 2	197	pg/mL
		Chemokine Panel 1 (human) Control 3	34.2	pg/mL
5	NHP IP-10	Calibrator	2420	pg/mL
		Chemokine Panel 1 (human) Control 1	348	pg/mL
		Chemokine Panel 1 (human) Control 2	57.3	pg/mL
		Chemokine Panel 1 (human) Control 3	10.3	pg/mL
6	NHP MIP-1 α	Calibrator	1030	pg/mL
		Chemokine Panel 1 (human) Control 1	824	pg/mL
		Chemokine Panel 1 (human) Control 2	371	pg/mL
		Chemokine Panel 1 (human) Control 3	152	pg/mL
7	NHP IL-8	Calibrator	86200	pg/mL
		Chemokine Panel 1 (human) Control 1	171000	pg/mL
		Chemokine Panel 1 (human) Control 2	58100	pg/mL
		Chemokine Panel 1 (human) Control 3	22500	pg/mL
8	NHP MCP-1	Calibrator	686	pg/mL
		Chemokine Panel 1 (human) Control 1	420	pg/mL
		Chemokine Panel 1 (human) Control 2	66.6	pg/mL
		Chemokine Panel 1 (human) Control 3	11.1	pg/mL
9	NHP MDC	Calibrator	12200	pg/mL
		Chemokine Panel 1 (human) Control 1	11100	pg/mL
		Chemokine Panel 1 (human) Control 2	3460	pg/mL
		Chemokine Panel 1 (human) Control 3	907	pg/mL
10	NHP MCP-4	Calibrator	787	pg/mL
		Chemokine Panel 1 (human) Control 1	635	pg/mL
		Chemokine Panel 1 (human) Control 2	232	pg/mL
		Chemokine Panel 1 (human) Control 3	82.2	pg/mL

Controls tested may not be the same lot as shipped.

The calibrator and controls for this assay contain a human recombinant cytokine calibrator sequence that is highly homologous to the non-human primate cytokine. This assay has been verified to cross-react with homologs of rhesus and cynomolgus monkeys. See Product Insert for experimental data.

CERTIFICATE OF ANALYSIS

Standard Curve Data:

NHP MIP-1 β - Spot 2		
Conc. pg/mL	ECL Signal	%CV
1050	1,800,000	3.0
263	473,000	1.8
65.6	89,600	2.2
16.4	16,400	3.2
4.10	3,620	4.3
1.03	926	4.3
0.256	310	7.0
0.00	116	20.0

NHP Eotaxin-3 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
5410	144,000	3.1
1350	44,400	3.1
338	12,200	3.6
84.5	3,020	3.7
21.1	893	3.9
5.28	299	6.8
1.32	151	14.0
0.00	101	30.0

NHP TARC - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1630	853,000	2.3
408	262,000	3.1
102	65,600	3.8
25.5	17,400	4.3
6.37	4,870	3.2
1.59	1,740	4.1
0.398	871	5.4
0.00	599	7.5

NHP IP-10 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
2420	557,000	4.2
605	336,000	4.2
151	117,000	4.1
37.8	31,400	5.0
9.45	8,650	2.9
2.36	2,250	2.5
0.591	624	4.6
0.00	90	26.0

NHP MIP-1 α - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1030	594,000	2.0
258	93,700	2.5
64.4	9,410	2.2
16.1	818	4.2
4.02	244	8.9
1.01	165	9.8
0.251	158	13.0
0.00	157	14.0

NHP IL-8 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
86200	880,000	3.8
21600	157,000	5.7
5390	16,700	6.5
1350	2,200	7.3
337	639	6.3
84.2	278	12.0
21.0	219	15.0
0.00	181	20.0

NHP MCP-1 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
686	956,000	5.1
172	251,000	4.5
42.9	58,800	4.9
10.7	13,700	6.8
2.68	3,480	5.8
0.670	984	5.8
0.167	342	10.0
0.00	143	17.0

NHP MDC - Spot 9		
Conc. pg/mL	ECL Signal	%CV
12200	2,190,000	3.5
3050	449,000	3.2
763	51,300	3.6
191	7,870	2.7
47.7	1,880	2.9
11.9	580	5.2
2.98	295	10.0
0.00	197	17.0

NHP MCP-4 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
787	1,210,000	2.9
197	217,000	3.3
49.2	27,800	5.0
12.3	2,700	6.4
3.07	374	7.6
0.769	163	15.0
0.192	148	22.0
0.00	151	24.0

CERTIFICATE OF ANALYSIS

Page 4 of 4

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0046612	2-8°C	31 Jan 2020
Anti-hu MIP-1β Antibody	D0080861	2-8°C	30 Jun 2019
Anti-hu Eotaxin-3 Antibody	D0080899	2-8°C	31 Jul 2019
Anti-hu TARC Antibody	D0080900	2-8°C	31 Jul 2019
Anti-hu IP-10 Antibody	D0080863	2-8°C	30 Jun 2019
Anti-hu MIP-1α Antibody	D0080902	2-8°C	31 Jul 2019
Anti-hu IL-8 (HA) Antibody	D0080864	2-8°C	30 Jun 2019
Anti-hu MCP-1 Antibody	D0080904	2-8°C	31 Jul 2019
Anti-hu MDC Antibody	D0080866	2-8°C	30 Jun 2019
Anti-hu MCP-4 Antibody	D0080867	2-8°C	30 Jun 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080131	2-8°C	31 Jul 2022
Diluent 43	M0430023	≤ -10°C	31 Dec 2019
Diluent 3	M0030078	≤ -10°C	31 Jan 2020
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)	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

Additional Comments:

	Assay	Spot	Specification	Comment
1	NHP MDC	9	ULOQ Conc.	ULOQ = 121.3%

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 Assurance		05 OCT 2017