

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: K0080918

Expiration Date: 31 Jul 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%

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

Page 2 of 4

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	980	pg/mL
		Chemokine Panel 1 (human) Control 2	170	pg/mL
		Chemokine Panel 1 (human) Control 3	30	pg/mL
3	NHP Eotaxin-3	Calibrator	5410	pg/mL
		Chemokine Panel 1 (human) Control 1	4050	pg/mL
		Chemokine Panel 1 (human) Control 2	580	pg/mL
		Chemokine Panel 1 (human) Control 3	90	pg/mL
4	NHP TARC	Calibrator	1630	pg/mL
		Chemokine Panel 1 (human) Control 1	1460	pg/mL
		Chemokine Panel 1 (human) Control 2	220	pg/mL
		Chemokine Panel 1 (human) Control 3	40	pg/mL
5	NHP IP-10	Calibrator	2420	pg/mL
		Chemokine Panel 1 (human) Control 1	590	pg/mL
		Chemokine Panel 1 (human) Control 2	90	pg/mL
		Chemokine Panel 1 (human) Control 3	16	pg/mL
6	NHP MIP-1 α	Calibrator	1030	pg/mL
		Chemokine Panel 1 (human) Control 1	910	pg/mL
		Chemokine Panel 1 (human) Control 2	410	pg/mL
		Chemokine Panel 1 (human) Control 3	160	pg/mL
7	NHP IL-8	Calibrator	86200	pg/mL
		Chemokine Panel 1 (human) Control 1	66500	pg/mL
		Chemokine Panel 1 (human) Control 2	22500	pg/mL
		Chemokine Panel 1 (human) Control 3	8800	pg/mL
8	NHP MCP-1	Calibrator	686	pg/mL
		Chemokine Panel 1 (human) Control 1	610	pg/mL
		Chemokine Panel 1 (human) Control 2	160	pg/mL
		Chemokine Panel 1 (human) Control 3	16	pg/mL
9	NHP MDC	Calibrator	12200	pg/mL
		Chemokine Panel 1 (human) Control 1	11600	pg/mL
		Chemokine Panel 1 (human) Control 2	3300	pg/mL
		Chemokine Panel 1 (human) Control 3	1010	pg/mL
10	NHP MCP-4	Calibrator	787	pg/mL
		Chemokine Panel 1 (human) Control 1	710	pg/mL
		Chemokine Panel 1 (human) Control 2	270	pg/mL
		Chemokine Panel 1 (human) Control 3	82	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	1,570,000	3.3
263	426,000	2.6
65.6	83,800	3.2
16.4	15,200	2.9
4.10	3,090	4.4
1.03	791	3.9
0.256	273	6.7
0.00	110	20.0

NHP Eotaxin-3 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
5410	121,000	2.9
1350	33,300	2.7
338	8,620	4.4
84.5	2,160	4.2
21.1	641	6.5
5.28	240	10.0
1.32	133	19.0
0.00	98	25.0

NHP TARC - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1630	841,000	3.6
408	263,000	4.1
102	65,500	3.1
25.5	17,500	4.1
6.37	4,780	4.6
1.59	1,650	5.7
0.398	804	5.1
0.00	525	10.0

NHP IP-10 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
2420	500,000	4.8
605	291,000	4.2
151	85,000	3.8
37.8	22,800	2.6
9.45	5,800	3.0
2.36	1,560	4.4
0.591	461	5.1
0.00	93	25.0

NHP MIP-1 α - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1030	642,000	3.0
258	106,000	3.0
64.4	11,100	3.6
16.1	1,000	7.3
4.02	237	8.9
1.01	154	13.0
0.251	155	14.0
0.00	140	13.0

NHP IL-8 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
86200	694,000	3.6
21600	96,500	7.5
5390	8,260	7.3
1350	1,080	9.7
337	379	9.2
84.2	226	13.0
21.0	193	15.0
0.00	174	17.0

NHP MCP-1 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
686	841,000	5.9
172	223,000	4.1
42.9	50,800	4.2
10.7	11,800	4.0
2.68	2,990	5.0
0.670	888	6.5
0.167	329	10.0
0.00	164	24.0

NHP MDC - Spot 9		
Conc. pg/mL	ECL Signal	%CV
12200	2,130,000	3.7
3050	279,000	4.6
763	33,700	5.1
191	5,860	4.9
47.7	1,540	4.9
11.9	504	6.5
2.98	240	11.0
0.00	152	16.0

NHP MCP-4 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
787	1,100,000	2.9
197	140,000	3.1
49.2	14,100	7.0
12.3	1,260	7.4
3.07	245	10.0
0.769	147	18.0
0.192	123	20.0
0.00	129	21.0

CERTIFICATE OF ANALYSIS

Page 4 of 4


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0046693	2-8°C	30 Apr 2020
Anti-hu MIP-1β Antibody	D0080898	2-8°C	30 Sep 2019
Anti-hu Eotaxin-3 Antibody	D0080899	2-8°C	31 Jul 2019
Anti-hu TARC Antibody	D0080993	2-8°C	31 Oct 2019
Anti-hu IP-10 Antibody	D0080901	2-8°C	31 Jul 2019
Anti-hu MIP-1α Antibody	D0080902	2-8°C	31 Jul 2019
Anti-hu IL-8 (HA) Antibody	D0080903	2-8°C	31 Jul 2019
Anti-hu MCP-1 Antibody	D0080904	2-8°C	31 Jul 2019
Anti-hu MDC Antibody	D0080905	2-8°C	31 Jul 2019
Anti-hu MCP-4 Antibody	D0080906	2-8°C	31 Aug 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080131	2-8°C	31 Jul 2022
Diluent 43	M0430024	≤ -10°C	31 Jan 2020
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), 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 Assurance		12 FEB 2018