

CERTIFICATE OF ANALYSIS

Page 1 of 4

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

Kit Catalog Numbers: K15055D-Series

Kit Lot Number: K0080860

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%

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	1,440,000	4.4
263	422,000	3.5
65.6	87,600	3.5
16.4	15,600	4.0
4.10	3,370	5.3
1.03	870	4.9
0.256	318	8.7
0.00	124	13.0

NHP Eotaxin-3 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
5410	136,000	3.5
1350	41,600	4.2
338	11,700	4.9
84.5	3,000	5.4
21.1	925	5.0
5.28	328	8.4
1.32	181	7.7
0.00	122	18.0

NHP TARC - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1630	715,000	3.4
408	235,000	4.8
102	56,700	4.4
25.5	15,400	4.1
6.37	4,380	4.7
1.59	1,680	5.4
0.398	933	6.6
0.00	681	9.5

NHP IP-10 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
2420	495,000	4.3
605	335,000	3.2
151	116,000	3.9
37.8	32,000	6.1
9.45	8,930	3.7
2.36	2,410	2.7
0.591	690	3.6
0.00	107	16.0

NHP MIP-1 α - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1030	612,000	4.5
258	102,000	3.5
64.4	11,400	3.0
16.1	1,000	6.1
4.02	280	5.8
1.01	196	11.0
0.251	187	10.0
0.00	182	12.0

NHP IL-8 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
86200	770,000	4.5
21600	116,000	7.1
5390	11,800	8.6
1350	1,640	9.2
337	581	6.9
84.2	305	8.7
21.0	270	7.2
0.00	242	9.3

NHP MCP-1 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
686	969,000	6.3
172	259,000	4.6
42.9	60,700	5.2
10.7	14,100	7.1
2.68	3,630	5.2
0.670	1,050	3.6
0.167	420	7.3
0.00	197	10.0

NHP MDC - Spot 9		
Conc. pg/mL	ECL Signal	%CV
12200	1,670,000	4.0
3050	259,000	6.1
763	31,400	4.7
191	4,970	6.2
47.7	1,320	5.3
11.9	461	6.1
2.98	278	6.0
0.00	205	11.0

NHP MCP-4 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
787	1,400,000	2.4
197	257,000	4.1
49.2	35,800	4.1
12.3	3,670	6.6
3.07	496	5.6
0.769	209	15.0
0.192	177	15.0
0.00	163	16.0

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0046507	2-8°C	31 Oct 2019
Anti-hu MIP-1 β 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-1 α Antibody	D0080902	2-8°C	31 Jul 2019

CERTIFICATE OF ANALYSIS


Page 4 of 4

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	D0080867	2-8°C	30 Jun 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080131	2-8°C	31 Jul 2022
Diluent 43	M0430021	≤ -10°C	30 Sep 2019
Diluent 3	M0030077	≤ -10°C	30 Nov 2019
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		09 AUG 2017