

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: K0080993

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-1β	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-1α	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-1β	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-1α	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-1β	< 1.0%
3	NHP Eotaxin-3	< 1.0%
4	NHP TARC	< 1.0%
5	NHP IP-10	< 1.0%
6	NHP MIP-1α	< 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	1040	pg/mL
		Chemokine Panel 1 (human) Control 1	940	pg/mL
		Chemokine Panel 1 (human) Control 2	158	pg/mL
		Chemokine Panel 1 (human) Control 3	26.9	pg/mL
3	NHP Eotaxin-3	Calibrator	4550	pg/mL
		Chemokine Panel 1 (human) Control 1	3920	pg/mL
		Chemokine Panel 1 (human) Control 2	553	pg/mL
		Chemokine Panel 1 (human) Control 3	86.3	pg/mL
4	NHP TARC	Calibrator	1600	pg/mL
		Chemokine Panel 1 (human) Control 1	1390	pg/mL
		Chemokine Panel 1 (human) Control 2	213	pg/mL
		Chemokine Panel 1 (human) Control 3	37.8	pg/mL
5	NHP IP-10	Calibrator	2380	pg/mL
		Chemokine Panel 1 (human) Control 1	574	pg/mL
		Chemokine Panel 1 (human) Control 2	87.5	pg/mL
		Chemokine Panel 1 (human) Control 3	15.4	pg/mL
6	NHP MIP-1 α	Calibrator	1020	pg/mL
		Chemokine Panel 1 (human) Control 1	887	pg/mL
		Chemokine Panel 1 (human) Control 2	393	pg/mL
		Chemokine Panel 1 (human) Control 3	152	pg/mL
7	NHP IL-8	Calibrator	70900	pg/mL
		Chemokine Panel 1 (human) Control 1	70100	pg/mL
		Chemokine Panel 1 (human) Control 2	23500	pg/mL
		Chemokine Panel 1 (human) Control 3	9330	pg/mL
8	NHP MCP-1	Calibrator	671	pg/mL
		Chemokine Panel 1 (human) Control 1	595	pg/mL
		Chemokine Panel 1 (human) Control 2	155	pg/mL
		Chemokine Panel 1 (human) Control 3	15.2	pg/mL
9	NHP MDC	Calibrator	11600	pg/mL
		Chemokine Panel 1 (human) Control 1	10800	pg/mL
		Chemokine Panel 1 (human) Control 2	3120	pg/mL
		Chemokine Panel 1 (human) Control 3	982	pg/mL
10	NHP MCP-4	Calibrator	769	pg/mL
		Chemokine Panel 1 (human) Control 1	688	pg/mL
		Chemokine Panel 1 (human) Control 2	258	pg/mL
		Chemokine Panel 1 (human) Control 3	78.2	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
1040	1,680,000	3.7
260	469,000	4.5
65.0	88,900	3.8
16.3	15,600	3.4
4.06	2,910	5.7
1.02	722	4.9
0.254	245	8.8
0.00	110	14.0

NHP Eotaxin-3 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
4550	120,000	4.7
1140	36,300	6.3
284	9,010	4.0
71.1	2,280	5.6
17.8	668	6.3
4.44	246	10.0
1.11	139	12.0
0.00	93	25.0

NHP TARC - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1600	983,000	3.4
400	291,000	6.6
100	68,800	4.2
25.0	18,200	5.7
6.25	5,060	5.6
1.56	1,850	7.4
0.391	1,000	8.0
0.00	737	8.9

NHP IP-10 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
2380	714,000	3.2
595	341,000	3.5
149	103,000	2.9
37.2	28,200	2.6
9.30	7,280	2.8
2.32	1,910	2.5
0.581	558	4.2
0.00	97	17.0

NHP MIP-1 α - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1020	703,000	2.0
255	130,000	4.6
63.8	13,700	2.4
15.9	1,170	4.1
3.98	237	6.2
0.996	153	11.0
0.249	138	16.0
0.00	132	8.9

NHP IL-8 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
70900	931,000	4.7
17700	217,000	8.8
4430	18,500	8.5
1110	1,810	7.9
277	463	5.9
69.2	209	11.0
17.3	175	15.0
0.00	153	16.0

NHP MCP-1 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
671	1,150,000	3.9
168	315,000	3.7
41.9	71,700	2.4
10.5	17,000	3.8
2.62	4,290	3.2
0.655	1,200	3.4
0.164	391	6.1
0.00	130	18.0

NHP MDC - Spot 9		
Conc. pg/mL	ECL Signal	%CV
11600	2,000,000	3.3
2900	291,000	6.4
725	32,800	5.0
181	5,700	4.2
45.3	1,420	4.8
11.3	454	5.3
2.83	233	6.0
0.00	149	14.0

NHP MCP-4 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
769	877,000	2.3
192	140,000	4.7
48.1	15,300	4.9
12.0	1,400	7.9
3.00	243	8.6
0.751	139	11.0
0.188	126	20.0
0.00	120	20.0

CERTIFICATE OF ANALYSIS

Page 4 of 4


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0046801	2-8°C	30 Sep 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	D0080996	2-8°C	31 Oct 2019
Anti-hu MCP-1 Antibody	D0080997	2-8°C	31 Oct 2019
Anti-hu MDC Antibody	D0080905	2-8°C	31 Jul 2019
Anti-hu MCP-4 Antibody	D0081068	2-8°C	31 Jul 2020
Chemokine Panel 1 (human) Calibrator Blend	A0080159	2-8°C	31 Dec 2022
Diluent 43	M0430024	≤ -10°C	31 Jan 2020
Diluent 3	M0030080	≤ -10°C	31 May 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), 1000mL	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		02 MAY 2018