

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: K0081161

Expiration Date: 31 Jan 2021

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	1.02	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	1180	pg/mL
		Chemokine Panel 1 (human) Control 1	1000	pg/mL
		Chemokine Panel 1 (human) Control 2	176	pg/mL
		Chemokine Panel 1 (human) Control 3	25.8	pg/mL
3	NHP Eotaxin-3	Calibrator	6240	pg/mL
		Chemokine Panel 1 (human) Control 1	4450	pg/mL
		Chemokine Panel 1 (human) Control 2	661	pg/mL
		Chemokine Panel 1 (human) Control 3	100	pg/mL
4	NHP TARC	Calibrator	1780	pg/mL
		Chemokine Panel 1 (human) Control 1	1350	pg/mL
		Chemokine Panel 1 (human) Control 2	213	pg/mL
		Chemokine Panel 1 (human) Control 3	36	pg/mL
5	NHP IP-10	Calibrator	2360	pg/mL
		Chemokine Panel 1 (human) Control 1	601	pg/mL
		Chemokine Panel 1 (human) Control 2	92.1	pg/mL
		Chemokine Panel 1 (human) Control 3	15.5	pg/mL
6	NHP MIP-1 α	Calibrator	1140	pg/mL
		Chemokine Panel 1 (human) Control 1	929	pg/mL
		Chemokine Panel 1 (human) Control 2	411	pg/mL
		Chemokine Panel 1 (human) Control 3	167	pg/mL
7	NHP IL-8	Calibrator	79700	pg/mL
		Chemokine Panel 1 (human) Control 1	66700	pg/mL
		Chemokine Panel 1 (human) Control 2	23000	pg/mL
		Chemokine Panel 1 (human) Control 3	9290	pg/mL
8	NHP MCP-1	Calibrator	562	pg/mL
		Chemokine Panel 1 (human) Control 1	478	pg/mL
		Chemokine Panel 1 (human) Control 2	73.3	pg/mL
		Chemokine Panel 1 (human) Control 3	12.8	pg/mL
9	NHP MDC	Calibrator	11800	pg/mL
		Chemokine Panel 1 (human) Control 1	9580	pg/mL
		Chemokine Panel 1 (human) Control 2	2840	pg/mL
		Chemokine Panel 1 (human) Control 3	948	pg/mL
10	NHP MCP-4	Calibrator	724	pg/mL
		Chemokine Panel 1 (human) Control 1	673	pg/mL
		Chemokine Panel 1 (human) Control 2	251	pg/mL
		Chemokine Panel 1 (human) Control 3	83.1	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
1180	1,800,000	4.5
295	457,000	3.8
73.8	71,800	2.3
18.4	9,150	4.1
4.61	1,240	3.8
1.15	341	5.6
0.288	193	11.0
0.00	166	8.8

NHP Eotaxin-3 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
6240	235,000	4.4
1560	67,400	3.8
390	16,900	4.3
97.5	4,250	6.7
24.4	1,120	3.8
6.09	385	5.9
1.52	191	11.0
0.00	134	16.0

NHP TARC - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1780	857,000	8.9
445	255,000	8.5
111	55,000	11.0
27.8	14,500	8.2
6.95	4,110	4.1
1.74	1,780	7.0
0.435	983	7.2
0.00	922	8.0

NHP IP-10 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
2360	805,000	9.7
590	387,000	5.1
148	111,000	8.0
36.9	29,700	4.9
9.22	7,660	3.1
2.30	2,000	3.6
0.576	588	5.6
0.00	136	14.0

NHP MIP-1 α - Spot 6		
Conc. pg/mL	ECL Signal	%CV
1140	724,000	10.0
285	130,000	5.7
71.3	14,200	4.1
17.8	1,440	7.5
4.45	372	5.2
1.11	255	6.4
0.278	250	7.3
0.00	253	6.8

NHP IL-8 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
79700	779,000	16.0
19900	137,000	25.0
4980	11,200	20.0
1250	1,330	13.0
311	452	7.1
77.8	304	7.5
19.5	301	15.0
0.00	266	11.0

NHP MCP-1 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
562	1,130,000	4.0
141	302,000	4.6
35.1	66,400	5.2
8.78	14,900	8.5
2.20	3,840	6.4
0.549	1,200	7.1
0.137	552	6.3
0.00	335	5.5

NHP MDC - Spot 9		
Conc. pg/mL	ECL Signal	%CV
11800	2,690,000	3.8
2950	415,000	3.7
738	45,800	7.4
184	7,270	6.0
46.1	1,830	2.6
11.5	667	4.0
2.88	393	6.3
0.00	305	10.0

NHP MCP-4 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
724	806,000	6.4
181	122,000	5.9
45.3	13,900	11.0
11.3	1,380	8.6
2.83	324	9.1
0.707	217	11.0
0.177	216	18.0
0.00	208	11.0

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Plate	Z0046943	2-8°C	31 Mar 2021
Anti-hu MIP-1β Antibody	D0081121	2-8°C	31 Jan 2022
Anti-hu Eotaxin-3 Antibody	D0081122	2-8°C	31 Jan 2022
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	A0080175	2-8°C	31 Jul 2023
Diluent 43	M0430028	≤ -10°C	31 Jan 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), 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		06 FEB 2019