



## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus NHP MCP-1 Kit  
**Kit Catalog Numbers:** K156NN - Series  
**Kit Lot Number:** K0080660  
**Expiration Date:** 31 Jan 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
1	NHP MCP-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	NHP MCP-1	1.09	80% - 120%	< 20%	375	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP MCP-1	Calibrator	530	pg/mL
		Chemokine Panel 1 (human) Control 1	437	pg/mL
		Chemokine Panel 1 (human) Control 2	71.5	pg/mL
		Chemokine Panel 1 (human) Control 3	11.6	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.

### Standard Curve Data:

NHP MCP-1 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
530	218,000	1.5
133	51,900	2.3
33.1	13,000	2.5
8.28	3,390	2.1
2.07	881	2.9
0.518	273	2.8
0.129	110	3.9
0.00	59	4.5



## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MCP-1 Plate	Z0046298	2-8°C	31 May 2019
Anti-hu MCP-1 Antibody	D0080865	2-8°C	30 Jun 2019
Chemokine Panel 1 (human) Calibrator	A0080106	2-8°C	30 Nov 2021
Diluent 43	M0430018	≤ -10°C	31 Jan 2019
Diluent 3	M0030073	≤ -10°C	31 Mar 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	<i>Dawn Stollar</i>	27 JAN 2017