

## 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:** K0080947  
**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
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	671	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

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
671	239,000	1.9
168	62,600	2.6
41.9	16,400	3.9
10.5	4,370	3.6
2.62	1,160	3.3
0.655	339	4.3
0.164	129	4.0
0.00	60	4.7

### CERTIFICATE OF ANALYSIS

**Components:**


Description	Lot Number	Storage Temperature	Expiration Date
Human MCP-1 Plate	Z0046742	2-8°C	30 Jun 2020
Anti-hu MCP-1 Antibody	D0080904	2-8°C	31 Jul 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080159	2-8°C	31 Dec 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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		14 FEB 2018