



## CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-04594

**Product Description:** NHP MDC Kit V-PLEX™ and V-PLEX Plus

**Kit Catalog Numbers:** K156NPD-1, K156NPD-2, K156NPD-4, K156NPG-1, K156NPG-2, K156NPG-4

**Kit Lot Number:** K0080184

**Expiration Date:** 31 DEC 2015

### Acceptance Data:

Metric		Specifications	Assay
			NHP MDC
% Recovery Concentration	Calibrators	80% - 120%	Pass
	Controls	75% - 125%	Pass
	ULOQ	80% - 120%	Pass
	LLOQ	80% - 120%	Pass
% Recovery CV	Calibrators	< 20%	Pass
	Controls	< 20%	Pass
	ULOQ	< 20%	Pass
	LLOQ	< 20%	Pass

### Concentration of Materials Tested:

Units = pg/mL	Assay
	NHP MDC*
Reconstituted Calibrator	9220
Reconstituted Control 1 (A00C0190)	6720
Reconstituted Control 2 (A00C0191)	1400
Reconstituted Control 3 (A00C0192)	349
ULOQ	7500
LLOQ	88.3

\*The calibrator and controls for this assay contain a human recombinant chemokine calibrator sequence that is highly homologous to the non-human primate chemokine. This assay has been verified to cross-react with homologs of rhesus and cynomolgus monkeys. See Product Insert for experimental data.

### Standard Curve Data from Kit QC:

NHP MDC		
Conc., (pg/mL)	Average Signal	%CV
9220	442 831	9.9
2305	117 941	3.9
576	18 209	6.2
144	3140	4.5
36.0	763	3.9
9.00	236	2.2
2.25	110	2.4
0	70	5.8



## CERTIFICATE OF ANALYSIS

Page 2 of 2

COA-04594

NHP MDC Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Human MDC Plate	Z0045319	2-8°C	30 NOV 2016
Anti-hu MDC Antibody	D0080119	2-8°C	31 DEC 2015
Chemokine Panel 1 (human) Calibrator Blend	A0080038	2-8°C	30 SEP 2016
Diluent 43	M0430009	≤ -10°C	30 NOV 2016
Diluent 3	M0030061	≤ -10°C	31 AUG 2016
MSD Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A
Additional Components Included in V-PLEX Plus			
Wash Buffer (20X), 100 mL	Not Kit Specific	Room Temperature	N/A
Plate Seals	Not Kit Specific	Room Temperature	N/A
Chemokine Panel 1 (human) Control 1	Not Kit Specific; See Separate Control Pack COA		
Chemokine Panel 1 (human) Control 2			
Chemokine Panel 1 (human) Control 3			

**Statement:** The above kit lot was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Caitlin Hansen	Quality Control		06 AUG 2014