



CERTIFICATE OF ANALYSIS

COA-6070

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus NHP MDC Kit
Kit Catalog Numbers: K156NPD-1, K156NPD-2, K156NPD-4, K156NPG-1, K156NPG-2, K156NPG-4
Kit Lot Number: K0080184
Expiration Date: 31 Dec 2016

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 MDC	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	NHP MDC	88.3	80% - 120%	< 20%	7500	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP MDC	Calibrator	9220	pg/mL
		Chemokine panel 1 (human) Control 1	8670	pg/mL
		Chemokine panel 1 (human) Control 2	2660	pg/mL
		Chemokine panel 1 (human) Control 3	703	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

COA-6070

Page 2 of 2

Standard Curve Data:

NHP MDC - Spot 1		
Conc. pg/mL	ECL Signal	%CV
9220	454,000	3.7
2310	120,000	4.6
576	21,600	6.0
144	4,550	3.2
36.0	1,140	3.3
9.00	330	2.7
2.25	126	2.3
0.00	61	7.3

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	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 kit 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	Monika Jain	Quality Assurance		15 OCT 2015




Meso Scale Diagnostics, L.L.C.

Description: V-PLEX® and V-PLEX Plus NHP MDC Kit

Catalog Number(s): K156NP D/G-1/2/4

Lot #: K0080184

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of Lot # K0080184 has been extended to December 31, 2016.

Function	Name	Signature	Date
Quality Assurance	Michael Grossman		14 OCT 2015

This memo supersedes all other statements of date of expiration for this kit, Lot # K0080184.