



CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-04380

Product Description: NHP VEGF Kit V-PLEX™ and V-PLEX Plus

Kit Catalog Numbers: K156RHD-1, K156RHD-2, K156RHD-4, K156RHG-1, K156RHG-2, K156RHG-4

Kit Lot Number: K0080127

Expiration Date: 31 OCT 2015

Acceptance Data:

Metric		Specifications	Assay
			NHP VEGF
% Recovery Concentration	Calibrators	80% - 120%	Pass
	Controls	70% - 130%	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 VEGF*
Reconstituted Calibrator	894
Reconstituted Control 1 (A00C0209)	279
Reconstituted Control 2 (A00C0210)	88.1
Reconstituted Control 3 (A00C0211)	23.5
ULOQ	562
LLOQ	7.70

*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 from Kit QC:

NHP VEGF		
Conc., (pg/mL)	Average Signal	%CV
894	203 517	5.7
224	35 527	4.8
55.9	8090	3.9
14.0	1796	5.3
3.49	548	3.4
0.873	173	2.4
0.218	86	6.1
0	52	7.1



CERTIFICATE OF ANALYSIS

Page 2 of 2

COA-04380

NHP VEGF Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Human VEGF Plate	Z0054868	2-8°C	30 JUN 2016
Anti-hu VEGF Antibody	D0080291	2-8°C	31 AUG 2016
Cytokine Panel 1 (human) Calibrator Blend	A0080033	2-8°C	31 MAR 2016
Diluent 43	M0430007	≤ -10°C	30 APR 2016
Diluent 3	M0030057	≤ -10°C	31 OCT 2015
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
Cytokine Panel 1 (human) Control 1	Not Kit Specific; See Separate Control Pack COA		
Cytokine Panel 1 (human) Control 2			
Cytokine 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	Dana Short	Quality Control		09 MAY 2014