



## CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-04480

**Product Description:** NHP TNF- $\beta$  Kit V-PLEX™ and V-PLEX Plus

**Kit Catalog Numbers:** K156RGD-1, K156RGD-2, K156RGD-4, K156RGG-1, K156RGG-2, K156RGG-4

**Kit Lot Number:** K0080041

**Expiration Date:** 31 AUG 2015

### Acceptance Data:

Metric		Specifications	Assay
			NHP TNF- $\beta$
% Recovery Concentration	Calibrators	80% - 120%	Pass
	Controls	80% - 120%	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 TNF- $\beta$ *
Reconstituted Calibrator	589
Reconstituted Control 1 (A00C0193)	165
Reconstituted Control 2 (A00C0194)	43.6
Reconstituted Control 3 (A00C0195)	10.0
ULOQ	531
LLOQ	0.290

\*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-04480

#### Standard Curve Data from Kit QC:

NHP TNF-β		
Conc., (pg/mL)	Average Signal	%CV
589	414 030	2.2
147	98 089	1.8
37	24 341	2.4
9.20	6113	1.9
2.30	1739	1.6
0.575	515	2.7
0.144	218	4.1
0	123	5.3

NHP TNF-β V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Human TNF-β Plate	Z0044652	2-8°C	30 SEP 2015
Anti-hu TNF-β Antibody	D0080090	2-8°C	30 NOV 2015
Cytokine Panel 1 (human) Calibrator Blend	A0080006	2-8°C	31 AUG 2015
Diluent 43	M0430003	≤ -10°C	30 SEP 2015
Diluent 3	M0030055	≤ -10°C	30 SEP 2015
Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A
Additional Components Included in V-PLEX Plus			
Wash Buffer (20X), 100mL	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		30 MAY 2014