



## CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-05088

**Product Description:** NHP IL-12/IL-23 p40 Kit V-PLEX™ and V-PLEX Plus

**Kit Catalog Numbers:** K156RJD-1, K156RJD-2, K156RJD-4, K156RJG-1, K156RJG-2, K156RJG-4

**Kit Lot Number:** K0080040

**Expiration Date:** 31 AUG 2015

### Acceptance Data:

Metric		Specifications	Assay
			NHP IL-12/23 p40
% 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 IL-12/23 p40*
Reconstituted Calibrator	2910
Reconstituted Control 1 (A00C0193)	907
Reconstituted Control 2 (A00C0194)	246
Reconstituted Control 3 (A00C0195)	59.3
ULOQ	2620
LLOQ	1.42

\*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-05088

#### Standard Curve Data from Kit QC:

NHP IL-12/23 p40		
Conc., (pg/mL)	Average Signal	%CV
2910	219 881	4.0
728	51 586	4.5
182	12 563	3.5
45.5	3033	6.1
11.4	808	9.2
2.84	245	7.3
0.710	124	5.4
0	74	9.7

NHP IL-12/IL-23 p40 Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12/IL-23 p40 Plate	Z0044651	2-8°C	30 SEP 2015
Anti-hu IL-12/23 p40 Antibody	D0080089	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), 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	Caitlin Hansen	Quality Control		19 JAN 2015