



CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-05206

Product Description: NHP IL-6 Kit V-PLEX™ and V-PLEX Plus

Kit Catalog Numbers: K156QXD-1, K156QXD-2, K156QXD-4, K156QXG-1, K156QXG-2, K156QXG-4

Kit Lot Number: K0080271

Expiration Date: 30 APR 2017

Acceptance Data:

Metric		Specifications	Assay
			NHP IL-6
% 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-6*
Reconstituted Calibrator	749
Reconstituted Control 1 (A00C0273)	128
Reconstituted Control 2 (A00C0274)	58.3
Reconstituted Control 3 (A00C0293)	8.27
ULOQ	488
LLOQ	1.58

*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-05206

Standard Curve Data from Kit QC:

NHP IL-6		
Conc., (pg/mL)	Average Signal	%CV
749	471 116	4.6
187	103 571	4.7
46.8	23 394	5.4
11.7	5776	5.6
2.93	1528	5.2
0.731	429	4.3
0.183	152	3.3
0	61	4.9

NHP IL-6 Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Human IL-6 Plate	Z0045549	2-8°C	30 JUN 2017
Anti-hu IL-6 Antibody	D0080492	2-8°C	31 MAY 2017
Proinflammatory Panel 1 (human) Calibrator Blend	A0080045	2-8°C	30 NOV 2019
Diluent 2	M0020085	≤ -10°C	31 JAN 2017
Diluent 3	M0030063	≤ -10°C	30 APR 2017
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
Proinflammatory Panel 1 (human) Control 1	Not Kit Specific; See Separate Control Pack COA		
Proinflammatory Panel 1 (human) Control 2			
Proinflammatory 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		11 FEB 2015