



CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-04878

Product Description: NHP IL-15 Kit V-PLEX™ and V-PLEX Plus

Kit Catalog Numbers: K156RDD-1, K156RDD-2, K156RDD-4, K156RDG-1, K156RDG-2, K156RDG-4

Kit Lot Number: K0080232

Expiration Date: 30 SEP 2016

Acceptance Data:

Metric		Specifications	Assay
			NHP IL-15
% 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-15*
Reconstituted Calibrator	683
Reconstituted Control 1 (A00C0251)	191
Reconstituted Control 2 (A00C0252)	48.3
Reconstituted Control 3 (A00C0263)	16.2
ULOQ	525
LLOQ	1.40

*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-04878

Standard Curve Data from Kit QC:

NHP IL-15		
Conc., (pg/mL)	Average Signal	%CV
683	157 434	4.3
171	28 411	3.4
42.7	6689	6.0
10.7	1690	4.2
2.67	469	5.7
0.667	152	5.3
0.167	73	5.4
0	47	7.2

NHP IL-15 Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Human IL-15 Plate	Z0045466	2-8°C	31 MAR 2017
Anti-hu IL-15 Antibody	D0080318	2-8°C	30 SEP 2016
Cytokine Panel 1 (human) Calibrator Blend	A0080041	2-8°C	30 JUN 2019
Diluent 43	M0430011	≤ -10°C	28 FEB 2017
Diluent 3	M0030062	≤ -10°C	31 OCT 2016
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	Caitlin Hansen	Quality Control		21 OCT 2014