



CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-04434

Product Description: NHP Eotaxin-3 Kit V-PLEX™ and V-PLEX Plus

Kit Catalog Numbers: K156NUD-1, K156NUD-2, K156NUD-4, K156NUG-1, K156NUG-2, K156NUG-4

Kit Lot Number: K0080149

Expiration Date: 30 SEP 2015

Acceptance Data:

Metric		Specifications	Assay
			NHP Eotaxin-3
% Recovery Concentration	Calibrators	80% - 120%	Pass
	Controls	70% - 130%	Pass
	LLOQ	80% - 120%	Pass
	ULOQ	80% - 120%	Pass
% Recovery CV	Calibrators	< 20%	Pass
	Controls	< 20%	Pass
	LLOQ	< 20%	Pass
	ULOQ	< 20%	Pass

Concentration of Materials Tested:

Units = pg/mL	Assay
	NHP Eotaxin-3*
Reconstituted Calibrator	4620
Reconstituted Control 1 (A00C0190)	3710
Reconstituted Control 2 (A00C0191)	914
Reconstituted Control 3 (A00C0192)	208
ULOQ	3750
LLOQ	10.2

*The calibrator and controls for this assay contain a human recombinant chemokine calibrator sequence that is highly homologous to the non-human primate chemokine. 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 Eotaxin-3		
Conc., (pg/mL)	Average Signal	%CV
4620	79 692	3.5
1155	22 463	2.8
289	5998	2.4
72.2	1637	2.7
18.0	488	3.8
4.51	167	3.8
1.13	82	3.6
0	55	4.6




CERTIFICATE OF ANALYSIS

Page 2 of 2

COA-04434

NHP Eotaxin-3 Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Human Eotaxin-3 Plate	Z0045176	2-8°C	31 JUL 2016
Anti-hu Eotaxin-3 Antibody	D0080323	2-8°C	30 SEP 2016
Chemokine Panel 1 (human) Calibrator Blend	A0080017	2-8°C	31 OCT 2015
Diluent 43	M0430007	≤ -10°C	30 APR 2016
Diluent 3	M0030055	≤ -10°C	30 SEP 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
Chemokine Panel 1 (human) Control 1	Not Kit Specific; See Separate Control Pack COA		
Chemokine Panel 1 (human) Control 2			
Chemokine 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		14 MAY 2014