



CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-04841

Product Description: NHP TARC Kit V-PLEX™ and V-PLEX Plus

Kit Catalog Numbers: K156NTD-1, K156NTD-2, K156NTD-4, K156NTG-1, K156NTG-2, K156NTG-4

Kit Lot Number: K0080222

Expiration Date: 31 OCT 2015

Acceptance Data:

Metric		Specifications	Assay
			NHP TARC
% 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 TARC*
Reconstituted Calibrator	1450
Reconstituted Control 1 (A00C0267)	971
Reconstituted Control 2 (A00C0268)	235
Reconstituted Control 3 (A00C0269)	63.4
ULOQ	1120
LLOQ	3.32

*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.



CERTIFICATE OF ANALYSIS

COA-04841

Standard Curve Data from Kit QC:

NHP TARC		
Conc., (pg/mL)	Average Signal	%CV
1450	317 563	4.3
363	95 137	3.4
90.6	23 655	6.1
22.7	6201	3.3
5.66	1697	2.9
1.42	475	3.9
0.354	158	3.0
0	49	7.6

NHP TARC Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Human TARC Plate	Z0045439	2-8°C	28 FEB 2017
Anti-hu TARC Antibody	D0080190	2-8°C	30 APR 2016
Chemokine Panel 1 (human) Calibrator Blend	A0080017	2-8°C	31 OCT 2015
Diluent 43	M0430009	≤ -10°C	30 NOV 2016
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
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	Caitlin Hansen	Quality Control		14 OCT 2014