



CERTIFICATE OF ANALYSIS

COA-5858

Page 1 of 3

Product Description: V-PLEX® and V-PLEX Plus NHP IL-8 Kit
Kit Catalog Numbers: K156RAD-1, K156RAD-2, K156RAD-4, K156RAG-1, K156RAG-2, K156RAG-4
Kit Lot Number: K0080391
Expiration Date: 30 Nov 2016

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	NHP IL-8	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	NHP IL-8	1.13	80% - 120%	< 20%	375	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP IL-8	Calibrator	529	pg/mL
		Proinflammatory Panel 1 (human) Control 1	129	pg/mL
		Proinflammatory Panel 1 (human) Control 2	37.6	pg/mL
		Proinflammatory Panel 1 (human) Control 3	5.19	pg/mL

Controls tested may not be the same lot as shipped.

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-5858

Page 2 of 3

Standard Curve Data:

NHP IL-8 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
529	238,501	4.7
132	61,893	3.2
33.1	14,784	5.2
8.27	3,703	7.0
2.07	1,019	2.7
0.517	291	3.8
0.129	116	9.4
0	54	6.4

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-8 Plate	Z0044973	2-8°C	29 Feb 2016
Anti-hu IL-8 Antibody	D0080238	2-8°C	30 Apr 2016
Proinflammatory Panel 1 (human) Calibrator Blend	A0080031	2-8°C	31 Mar 2016
Diluent 2	M0020079	≤ -10°C	30 Nov 2015
Diluent 3	M0030059	≤ -10°C	31 Mar 2016
MSD Read Buffer T (4x)	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20x), 100mL	Not Kit Specific
Proinflammatory Panel 1 (human) Control 1	Not Kit Specific
Proinflammatory Panel 1 (human) Control 2	Not Kit Specific
Proinflammatory Panel 1 (human) Control 3	Not Kit Specific

Statement:

The above kit 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.



CERTIFICATE OF ANALYSIS

COA-5858

Page 3 of 3

	Name	Function	Signature	Date
Review/Approval	Monika Jain	Quality Assurance	<i>Monika Jain</i>	13 Aug 2015




Meso Scale Diagnostics, L.L.C.

Description: V-PLEX® and V-PLEX Plus Human IL-8 Kit

Catalog Number(s): K151RAD/G – 1/2/4

Lot #: K0080391

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of Lot # K0080391 has been extended to November 30, 2016.

Function	Name	Signature	Date
Quality Assurance	Michael Grossman		13 AUG 2015

This memo supersedes all other statements of date of expiration for this kit, Lot # K0080391.