



CERTIFICATE OF ANALYSIS

COA-5852

Page 1 of 3

Product Description: V-PLEX® and V-PLEX Plus NHP IFN- γ Kit
Kit Catalog Numbers: K156QOD-1, K156QOD-2, K156QOD-4, K156QOG-1, K156QOG-2, K156QOG-4
Kit Lot Number: K0080375
Expiration Date: 31 Jul 2017

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 IFN- γ	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	NHP IFN- γ	7.47	80% - 120%	< 20%	938	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP IFN- γ	Calibrator	1270	pg/mL
		Proinflammatory Panel 1 (human) Control 1	345	pg/mL
		Proinflammatory Panel 1 (human) Control 2	123	pg/mL
		Proinflammatory Panel 1 (human) Control 3	37.9	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-5852

Standard Curve Data:

NHP IFN- γ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1270	106,210	1.8
318	27,273	1.5
79.4	6,866	1.7
19.8	1,824	1.4
4.96	523	3.2
1.24	211	5.1
0.310	128	7.1
0	96	7.4

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IFN- γ Plate	Z0045802	2-8°C	31 Dec 2017
Anti-hu IFN- γ Antibody	D0080510	2-8°C	31 Jul 2017
Proinflammatory Panel 1 (human) Calibrator Blend	A0080062	2-8°C	31 Jul 2020
Diluent 2	M0020087	\leq -10°C	31 Oct 2017
Diluent 3	M0030067	\leq -10°C	30 Nov 2017
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-5852

Page 3 of 3

	Name	Function	Signature	Date
Review/Approval	Monika Jain	Quality Assurance	<i>Monika Jain</i>	06Aug2015