

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus NHP IFN- γ Kit
Kit Catalog Numbers: K156QO-Series
Kit Lot Number: K0080670
Expiration Date: 28 Feb 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (mouse) control 1		Proinflammatory Panel 1 (mouse) control 2		Proinflammatory Panel 1 (mouse) control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	NHP IFN- γ	80% - 120%	< 20%	80% - 130%	< 25%	80% - 130%	< 25%	80% - 130%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	NHP IFN- γ	0.390	75% - 125%	< 25%	570	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP IFN- γ	Calibrator	933	pg/mL
		Proinflammatory Panel 1 (mouse) control 1	479	pg/mL
		Proinflammatory Panel 1 (mouse) control 2	25.7	pg/mL
		Proinflammatory Panel 1 (mouse) control 3	1.53	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.

Standard Curve Data:

NHP IFN- γ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
933	404,000	1.9
233	102,000	2.2
58.3	25,700	1.5
14.6	6,820	1.2
3.64	1,800	1.3
0.911	539	1.8
0.228	204	2.3
0.00	81	7.1

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IFN- γ Plate	Z0046313	2-8°C	31 May 2019
Anti-ms IFN- γ Antibody	D0080785	2-8°C	31 Mar 2019
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080110	2-8°C	31 Dec 2021
Diluent 41	M0410015	$\leq -10^{\circ}\text{C}$	30 Jun 2019
Diluent 45	M0450012	$\leq -10^{\circ}\text{C}$	28 Feb 2019
MSD Read Buffer T	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 (mouse) Control 1	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 2	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 3	Not Kit Specific

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance		11 JAN 2017