

## CERTIFICATE OF ANALYSIS

Page 1 of 3

**Product Description:** V-PLEX® and V-PLEX Plus Proinflammatory Panel 1 (NHP) Kit  
**Kit Catalog Numbers:** K15056-Series  
**Kit Lot Number:** K0081072  
**Expiration Date:** 31 Jan 2021

### 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- $\gamma$	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	NHP IL-1 $\beta$	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
3	NHP IL-2	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
5	NHP IL-6	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
6	NHP IL-8	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
7	NHP IL-10	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- $\gamma$	1.76	80% - 120%	< 20%	938	80% - 120%	< 20%
2	NHP IL-1 $\beta$	0.646	80% - 120%	< 20%	375	80% - 120%	< 20%
3	NHP IL-2	0.890	80% - 120%	< 20%	938	80% - 120%	< 20%
5	NHP IL-6	0.633	80% - 120%	< 20%	488	80% - 120%	< 20%
6	NHP IL-8	0.591	80% - 120%	< 20%	375	80% - 120%	< 20%
7	NHP IL-10	0.298	80% - 120%	< 20%	233	80% - 120%	< 20%

### Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	NHP IFN- $\gamma$	< 1.0%
2	NHP IL-1 $\beta$	< 1.0%
3	NHP IL-2	< 1.0%
5	NHP IL-6	< 1.0%
6	NHP IL-8	< 1.0%
7	NHP IL-10	< 1.0%

## CERTIFICATE OF ANALYSIS

Page 2 of 3

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP IFN- $\gamma$	Calibrator	1230	pg/mL
		Proinflammatory Panel 1(human) Control 1	512	pg/mL
		Proinflammatory Panel 1(human) Control 2	132	pg/mL
		Proinflammatory Panel 1(human) Control 3	40.9	pg/mL
2	NHP IL-1 $\beta$	Calibrator	599	pg/mL
		Proinflammatory Panel 1(human) Control 1	203	pg/mL
		Proinflammatory Panel 1(human) Control 2	47.1	pg/mL
		Proinflammatory Panel 1(human) Control 3	10.9	pg/mL
3	NHP IL-2	Calibrator	1450	pg/mL
		Proinflammatory Panel 1(human) Control 1	390	pg/mL
		Proinflammatory Panel 1(human) Control 2	38.3	pg/mL
		Proinflammatory Panel 1(human) Control 3	3.34	pg/mL
5	NHP IL-6	Calibrator	768	pg/mL
		Proinflammatory Panel 1(human) Control 1	264	pg/mL
		Proinflammatory Panel 1(human) Control 2	46.8	pg/mL
		Proinflammatory Panel 1(human) Control 3	8.18	pg/mL
6	NHP IL-8	Calibrator	628	pg/mL
		Proinflammatory Panel 1(human) Control 1	235	pg/mL
		Proinflammatory Panel 1(human) Control 2	36.2	pg/mL
		Proinflammatory Panel 1(human) Control 3	5.93	pg/mL
7	NHP IL-10	Calibrator	356	pg/mL
		Proinflammatory Panel 1(human) Control 1	133	pg/mL
		Proinflammatory Panel 1(human) Control 2	25.3	pg/mL
		Proinflammatory Panel 1(human) Control 3	3.32	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- $\gamma$ - Spot 1			NHP IL-1 $\beta$ - Spot 2			NHP IL-2 - Spot 3		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
1230	193,000	2.2	599	1,570,000	2.4	1450	1,310,000	3.2
308	50,700	2.0	150	417,000	3.0	363	381,000	2.5
76.9	13,200	2.3	37.4	107,000	2.1	90.6	104,000	2.1
19.2	3,540	1.9	9.36	26,700	3.0	22.7	26,100	3.0
4.80	1,300	3.6	2.34	7,640	2.6	5.66	7,320	3.0
1.20	772	4.6	0.585	2,810	2.6	1.42	2,360	3.5
0.300	610	2.9	0.146	1,610	2.9	0.354	1,140	4.3
0.00	562	5.2	0.00	1,190	5.9	0.00	711	6.2

## CERTIFICATE OF ANALYSIS

Page 3 of 3

NHP IL-6 - Spot 5			NHP IL-8 - Spot 6			NHP IL-10 - Spot 7		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
768	1,450,000	2.0	628	793,000	3.0	356	737,000	2.5
192	350,000	1.9	157	205,000	4.3	89.0	217,000	1.9
48.0	77,800	1.5	39.3	51,100	3.4	22.3	57,500	2.4
12.0	18,200	2.6	9.81	12,800	3.5	5.56	14,300	2.4
3.00	4,550	2.1	2.45	3,580	2.8	1.39	3,980	2.5
0.750	1,300	2.7	0.613	1,150	3.2	0.348	1,290	2.1
0.188	502	4.4	0.153	541	4.4	0.0869	625	4.8
0.00	216	8.4	0.00	345	5.0	0.00	388	5.2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1(human) Plate	Z0046894	2-8°C	31 Jan 2021
Anti-hu IFN- $\gamma$ Antibody	D0081094	2-8°C	30 Nov 2021
Anti-hu IL-1 $\beta$ Antibody	D0081095	2-8°C	31 Dec 2021
Anti-hu IL-2 Antibody	D0081096	2-8°C	30 Nov 2021
Anti-hu IL-6 Antibody	D0081130	2-8°C	31 Jan 2022
Anti-hu IL-8 Antibody	D0081097	2-8°C	30 Nov 2021
Anti-hu IL-10 Antibody	D0081098	2-8°C	31 Jan 2022
Proinflammatory Panel 1(human) Calibrator Blend	A0080176	2-8°C	31 Jul 2023
Diluent 2	M0020102	$\leq -10^{\circ}\text{C}$	28 Feb 2021
Diluent 3	M0030084	$\leq -10^{\circ}\text{C}$	31 Jan 2021
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Wash Buffer 20X, 100 mL	Not Kit Specific
Plate Seals	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 product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		08 OCT 2018