

CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Proinflammatory Panel 1 (NHP) Kit
Kit Catalog Numbers: K15056-series
Kit Lot Number: K0081035
Expiration Date: 31 Jan 2020

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%
2	NHP IL-1 β	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- γ	1.76	80% - 120%	< 20%	938	80% - 120%	< 20%
2	NHP IL-1 β	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- γ	< 1.0%
2	NHP IL-1 β	< 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

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP IFN- γ	Calibrator	1430	pg/mL
		Proinflammatory Panel 1 (human) Control 1	574	pg/mL
		Proinflammatory Panel 1 (human) Control 2	155	pg/mL
		Proinflammatory Panel 1 (human) Control 3	41.8	pg/mL
2	NHP IL-1 β	Calibrator	561	pg/mL
		Proinflammatory Panel 1 (human) Control 1	202	pg/mL
		Proinflammatory Panel 1 (human) Control 2	50.4	pg/mL
		Proinflammatory Panel 1 (human) Control 3	10.7	pg/mL
3	NHP IL-2	Calibrator	1430	pg/mL
		Proinflammatory Panel 1 (human) Control 1	400	pg/mL
		Proinflammatory Panel 1 (human) Control 2	41.2	pg/mL
		Proinflammatory Panel 1 (human) Control 3	3.48	pg/mL
5	NHP IL-6	Calibrator	750	pg/mL
		Proinflammatory Panel 1 (human) Control 1	264	pg/mL
		Proinflammatory Panel 1 (human) Control 2	41.2	pg/mL
		Proinflammatory Panel 1 (human) Control 3	8.24	pg/mL
6	NHP IL-8	Calibrator	616	pg/mL
		Proinflammatory Panel 1 (human) Control 1	233	pg/mL
		Proinflammatory Panel 1 (human) Control 2	36.9	pg/mL
		Proinflammatory Panel 1 (human) Control 3	5.93	pg/mL
7	NHP IL-10	Calibrator	337	pg/mL
		Proinflammatory Panel 1 (human) Control 1	125	pg/mL
		Proinflammatory Panel 1 (human) Control 2	23.3	pg/mL
		Proinflammatory Panel 1 (human) Control 3	3.47	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			NHP IL-1 β - 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
1430	214,000	1.5	561	1,290,000	2.9	1430	1,220,000	1.8
358	56,800	1.6	140	352,000	1.9	358	343,000	1.2
89.4	14,100	2.7	35.1	88,700	1.7	89.4	93,300	1.8
22.3	3,590	2.8	8.77	22,200	2.0	22.3	23,600	2.7
5.59	1,080	2.8	2.19	6,020	2.7	5.59	6,460	5.7
1.40	469	5.1	0.548	1,880	3.5	1.40	1,790	1.7
0.349	302	9.2	0.137	861	3.5	0.349	727	4.9
0.00	247	9.0	0.00	507	9.1	0.00	338	6.5

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
750	1,350,000	2.8	616	931,000	4.4	337	577,000	1.5
188	311,000	1.8	154	229,000	3.9	84.3	169,000	1.8
46.9	69,200	2.3	38.5	56,500	2.7	21.1	44,400	2.0
11.7	15,500	2.8	9.63	14,200	3.1	5.27	11,100	2.2
2.93	4,010	2.4	2.41	3,770	2.5	1.32	3,090	1.7
0.732	1,070	3.2	0.602	1,110	3.0	0.329	984	3.0
0.183	388	4.3	0.150	447	5.4	0.0823	453	5.5
0.00	143	9.9	0.00	234	6.8	0.00	290	6.1

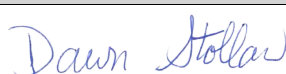
Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Plate	Z0046854	2-8°C	30 Nov 2020
Anti-hu IFN-γ Antibody	D0081074	2-8°C	31 Aug 2021
Anti-hu IL-1β 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	D0081078	2-8°C	31 Oct 2021
Anti-hu IL-8 Antibody	D0081037	2-8°C	29 Feb 2020
Anti-hu IL-10 Antibody	D0081024	2-8°C	31 Jan 2020
Proinflammatory Panel 1 (human) Calibrator Blend	A0080172	2-8°C	30 Jun 2023
Diluent 2	M0020099	≤ -10°C	31 Mar 2020
Diluent 3	M0030082	≤ -10°C	31 Oct 2020
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 (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
Review/Approval	Dawn Stollar	Quality Assurance		06 AUG 2018