

CERTIFICATE OF ANALYSIS

Page 1 of 4

Product Description: V-PLEX® and V-PLEX Proinflammatory Panel 1 (human) Kit

Kit Catalog Numbers: K15049-Series, K15052-Series, K15053-Series

Kit Lot Number: K0081048

Expiration Date: 30 Nov 2019

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	hu IFN- γ	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	hu IL-1 β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
3	hu IL-2	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	hu IL-4	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
5	hu IL-6	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
6	hu IL-8	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
7	hu IL-10	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
8	hu IL-12p70	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
9	hu IL-13	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
10	hu TNF- α	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu IFN- γ	1.76	80% - 120%	< 20%	938	80% - 120%	< 20%
2	hu IL-1 β	0.646	80% - 120%	< 20%	375	80% - 120%	< 20%
3	hu IL-2	0.890	80% - 120%	< 20%	938	80% - 120%	< 20%
4	hu IL-4	0.218	80% - 120%	< 20%	158	80% - 120%	< 20%
5	hu IL-6	0.633	80% - 120%	< 20%	488	80% - 120%	< 20%
6	hu IL-8	0.591	80% - 120%	< 20%	375	80% - 120%	< 20%
7	hu IL-10	0.298	80% - 120%	< 20%	233	80% - 120%	< 20%
8	hu IL-12p70	1.22	80% - 120%	< 20%	315	80% - 120%	< 20%
9	hu IL-13	4.21	80% - 120%	< 20%	353	80% - 120%	< 20%
10	hu TNF- α	0.690	80% - 120%	< 20%	248	80% - 120%	< 20%

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	hu IFN- γ	< 1.0%
2	hu IL-1 β	< 1.0%
3	hu IL-2	< 1.0%
4	hu IL-4	< 1.0%
5	hu IL-6	< 1.0%
6	hu IL-8	< 1.0%
7	hu IL-10	< 1.0%
8	hu IL-12p70	< 1.0%
9	hu IL-13	< 1.0%
10	hu TNF- α	< 1.0%

CERTIFICATE OF ANALYSIS

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IFN- γ	Calibrator	1430	pg/mL
		Proinflammatory Panel 1(human) Control 1	582	pg/mL
		Proinflammatory Panel 1(human) Control 2	164	pg/mL
		Proinflammatory Panel 1(human) Control 3	40.9	pg/mL
2	hu IL-1 β	Calibrator	561	pg/mL
		Proinflammatory Panel 1(human) Control 1	208	pg/mL
		Proinflammatory Panel 1(human) Control 2	52.7	pg/mL
		Proinflammatory Panel 1(human) Control 3	10.6	pg/mL
3	hu IL-2	Calibrator	1430	pg/mL
		Proinflammatory Panel 1(human) Control 1	407	pg/mL
		Proinflammatory Panel 1(human) Control 2	42.1	pg/mL
		Proinflammatory Panel 1(human) Control 3	3.54	pg/mL
4	hu IL-4	Calibrator	227	pg/mL
		Proinflammatory Panel 1(human) Control 1	101	pg/mL
		Proinflammatory Panel 1(human) Control 2	20.6	pg/mL
		Proinflammatory Panel 1(human) Control 3	2.95	pg/mL
5	hu IL-6	Calibrator	750	pg/mL
		Proinflammatory Panel 1(human) Control 1	271	pg/mL
		Proinflammatory Panel 1(human) Control 2	43.5	pg/mL
		Proinflammatory Panel 1(human) Control 3	8.03	pg/mL
6	hu IL-8	Calibrator	616	pg/mL
		Proinflammatory Panel 1(human) Control 1	233	pg/mL
		Proinflammatory Panel 1(human) Control 2	37.4	pg/mL
		Proinflammatory Panel 1(human) Control 3	6.13	pg/mL
7	hu IL-10	Calibrator	337	pg/mL
		Proinflammatory Panel 1(human) Control 1	125	pg/mL
		Proinflammatory Panel 1(human) Control 2	24.2	pg/mL
		Proinflammatory Panel 1(human) Control 3	3.49	pg/mL
8	hu IL-12p70	Calibrator	550	pg/mL
		Proinflammatory Panel 1(human) Control 1	194	pg/mL
		Proinflammatory Panel 1(human) Control 2	36.4	pg/mL
		Proinflammatory Panel 1(human) Control 3	7.29	pg/mL
9	hu IL-13	Calibrator	513	pg/mL
		Proinflammatory Panel 1(human) Control 1	168	pg/mL
		Proinflammatory Panel 1(human) Control 2	48.8	pg/mL
		Proinflammatory Panel 1(human) Control 3	13.4	pg/mL
10	hu TNF- α	Calibrator	346	pg/mL
		Proinflammatory Panel 1(human) Control 1	89.5	pg/mL
		Proinflammatory Panel 1(human) Control 2	17.4	pg/mL
		Proinflammatory Panel 1(human) Control 3	2.21	pg/mL

Controls tested may not be the same lot as shipped.

CERTIFICATE OF ANALYSIS

Standard Curve Data:

hu IFN- γ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1430	227,000	2.1
358	60,800	2.7
89.4	16,400	10.0
22.3	4,360	10.0
5.59	1,350	3.5
1.40	712	4.5
0.349	584	14.0
0.00	522	13.0

hu IL-1 β - Spot 2		
Conc. pg/mL	ECL Signal	%CV
561	1,440,000	2.0
140	401,000	1.7
35.1	109,000	7.5
8.77	27,900	7.2
2.19	7,210	4.7
0.548	2,450	3.4
0.137	1,310	11.0
0.00	877	11.0

hu IL-2 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
1430	1,110,000	2.2
358	331,000	1.7
89.4	93,500	8.6
22.3	24,600	8.2
5.59	6,100	4.5
1.40	1,980	4.0
0.349	939	12.0
0.00	538	11.0

hu IL-4 - Spot 4		
Conc. pg/mL	ECL Signal	%CV
227	763,000	2.0
56.8	241,000	3.3
14.2	68,700	8.6
3.55	17,900	9.6
0.887	4,260	2.8
0.222	1,250	4.1
0.0554	502	15.0
0.00	211	12.0

hu IL-6 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
750	1,360,000	2.1
188	338,000	1.9
46.9	77,200	6.8
11.7	17,800	9.9
2.93	4,300	2.3
0.732	1,230	2.5
0.183	515	11.0
0.00	232	14.0

hu IL-8 - Spot 6		
Conc. pg/mL	ECL Signal	%CV
616	874,000	2.7
154	233,000	3.2
38.5	62,500	6.9
9.63	15,800	7.5
2.41	4,000	3.4
0.602	1,300	5.4
0.150	659	11.0
0.00	392	13.0

hu IL-10 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
337	726,000	1.1
84.3	215,000	1.5
21.1	60,200	8.3
5.27	15,400	8.5
1.32	4,000	2.9
0.329	1,350	3.5
0.0823	756	11.0
0.00	476	9.3

hu IL-12p70 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
550	396,000	4.1
138	124,000	2.0
34.4	34,600	6.6
8.59	8,870	9.6
2.15	2,290	3.2
0.537	829	7.8
0.134	489	14.0
0.00	343	18.0

hu IL-13 - Spot 9		
Conc. pg/mL	ECL Signal	%CV
513	1,170,000	2.6
128	213,000	2.5
32.1	29,100	7.4
8.02	4,550	9.1
2.00	1,130	4.4
0.501	428	5.3
0.125	271	16.0
0.00	192	14.0

CERTIFICATE OF ANALYSIS

Page 4 of 4

hu TNF- α - Spot 10		
Conc. pg/mL	ECL Signal	%CV
346	338,000	3.6
86.5	84,600	2.5
21.6	22,700	7.3
5.41	5,680	9.0
1.35	1,750	3.0
0.338	780	5.2
0.0845	593	11.0
0.00	494	15.0


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Plate	Z0046868	2-8°C	31 Dec 2020
Anti-hu IFN- γ Antibody	D0081094	2-8°C	30 Nov 2021
Anti-hu IL-1 β Antibody	D0081095	2-8°C	31 Dec 2021
Anti-hu IL-2 Antibody	D0081036	2-8°C	29 Feb 2020
Anti-hu IL-4 Antibody	D0081017	2-8°C	31 Jan 2020
Anti-hu IL-6 Antibody	D0081078	2-8°C	31 Oct 2021
Anti-hu IL-8 Antibody	D0081097	2-8°C	30 Nov 2021
Anti-hu IL-10 Antibody	D0081098	2-8°C	31 Jan 2022
Anti-hu IL-12p70 Antibody	D0081008	2-8°C	30 Nov 2019
Anti-hu IL-13 Antibody	D0081092	2-8°C	31 Oct 2021
Anti-hu TNF- α Antibody	D0081099	2-8°C	30 Nov 2021
Proinflammatory Panel 1(human) Calibrator Blend	A0080172	2-8°C	30 Jun 2023
Diluent 2	M0020099	\leq -10°C	31 Mar 2020
Diluent 3	M0030082	\leq -10°C	31 Oct 2020
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
MSD Wash Buffer 20X, 100mL	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
Review/Approval	Dawn Stollar	Quality Assurance		19 SEP 2018