

CERTIFICATE OF ANALYSIS

Page 1 of 4

Product Description: V-PLEX® and V-PLEX Plus Proinflammatory Panel 1 (human) Kit
Kit Catalog Numbers: K151QR-Series; K151OD-Series; K15049-Series; K15052-Series; K15053-Series; K15056-Series; K15348-Series, K15346-Series, K15349-Series, K15345-Series
Kit Lot Number: K0081459
Expiration Date: 30 Apr 2022

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

Acceptance Data/ Non Specific Binding Specifications:

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

CERTIFICATE OF ANALYSIS

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IFN- γ	Calibrator	1500	pg/mL
		Proinflammatory Panel 1 (human) Control 1	546	pg/mL
		Proinflammatory Panel 1 (human) Control 2	167	pg/mL
		Proinflammatory Panel 1 (human) Control 3	32.2	pg/mL
2	IL-1 β	Calibrator	610	pg/mL
		Proinflammatory Panel 1 (human) Control 1	209	pg/mL
		Proinflammatory Panel 1 (human) Control 2	49.7	pg/mL
		Proinflammatory Panel 1 (human) Control 3	10	pg/mL
3	IL-2	Calibrator	1360	pg/mL
		Proinflammatory Panel 1 (human) Control 1	384	pg/mL
		Proinflammatory Panel 1 (human) Control 2	37.7	pg/mL
		Proinflammatory Panel 1 (human) Control 3	2.96	pg/mL
4	IL-4	Calibrator	211	pg/mL
		Proinflammatory Panel 1 (human) Control 1	109	pg/mL
		Proinflammatory Panel 1 (human) Control 2	20.2	pg/mL
		Proinflammatory Panel 1 (human) Control 3	3.11	pg/mL
5	IL-6	Calibrator	690	pg/mL
		Proinflammatory Panel 1 (human) Control 1	289	pg/mL
		Proinflammatory Panel 1 (human) Control 2	49.1	pg/mL
		Proinflammatory Panel 1 (human) Control 3	7.89	pg/mL
6	IL-8	Calibrator	612	pg/mL
		Proinflammatory Panel 1 (human) Control 1	220	pg/mL
		Proinflammatory Panel 1 (human) Control 2	33.2	pg/mL
		Proinflammatory Panel 1 (human) Control 3	5.42	pg/mL
7	IL-10	Calibrator	356	pg/mL
		Proinflammatory Panel 1 (human) Control 1	135	pg/mL
		Proinflammatory Panel 1 (human) Control 2	24.8	pg/mL
		Proinflammatory Panel 1 (human) Control 3	3.62	pg/mL
8	IL-12p70	Calibrator	464	pg/mL
		Proinflammatory Panel 1 (human) Control 1	158	pg/mL
		Proinflammatory Panel 1 (human) Control 2	31.4	pg/mL
		Proinflammatory Panel 1 (human) Control 3	5.92	pg/mL
9	IL-13	Calibrator	509	pg/mL
		Proinflammatory Panel 1 (human) Control 1	121	pg/mL
		Proinflammatory Panel 1 (human) Control 2	40.3	pg/mL
		Proinflammatory Panel 1 (human) Control 3	13.2	pg/mL
10	TNF- α	Calibrator	367	pg/mL
		Proinflammatory Panel 1 (human) Control 1	110	pg/mL
		Proinflammatory Panel 1 (human) Control 2	16.4	pg/mL
		Proinflammatory Panel 1 (human) Control 3	2.37	pg/mL

Controls tested may not be the same lot as shipped.

CERTIFICATE OF ANALYSIS

Standard Curve Data:

IFN- γ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1500	412,000	2.1
375	112,000	2.9
93.8	27,200	5.3
23.4	6,870	5.5
5.86	2,150	2.5
1.46	870	4.9
0.366	537	12.0
0.00	419	14.0

IL-1 β - Spot 2		
Conc. pg/mL	ECL Signal	%CV
610	1,630,000	4.9
153	469,000	3.5
38.1	119,000	4.0
9.53	30,800	5.0
2.38	8,840	3.0
0.596	3,100	2.8
0.149	1,580	3.3
0.00	1,140	5.0

IL-2 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
1360	1,300,000	3.6
340	369,000	2.2
85.0	100,000	2.7
21.3	24,900	2.5
5.31	6,930	3.2
1.33	1,980	2.7
0.332	799	4.0
0.00	387	6.6

IL-4 - Spot 4		
Conc. pg/mL	ECL Signal	%CV
211	949,000	3.1
52.8	298,000	2.3
13.2	79,100	3.0
3.30	20,300	2.6
0.824	5,420	3.0
0.206	1,540	2.9
0.0515	549	5.9
0.00	219	12.0

IL-6 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
690	1,020,000	2.7
173	231,000	3.2
43.1	47,300	4.5
10.8	10,700	3.4
2.70	3,000	2.8
0.674	930	4.1
0.168	403	6.8
0.00	230	8.9

IL-8 - Spot 6		
Conc. pg/mL	ECL Signal	%CV
612	1,010,000	2.1
153	258,000	2.0
38.3	63,900	3.4
9.56	15,900	2.8
2.39	4,330	3.5
0.598	1,250	2.6
0.149	472	4.2
0.00	224	8.6

IL-10 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
356	857,000	1.7
89.0	252,000	1.1
22.3	65,700	2.1
5.56	16,500	1.9
1.39	4,500	2.1
0.348	1,350	2.5
0.0869	548	4.6
0.00	289	8.0

IL-12p70 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
464	341,000	5.7
116	107,000	3.4
29.0	27,400	6.3
7.25	7,080	3.4
1.81	2,090	5.1
0.453	739	5.2
0.113	371	8.4
0.00	281	15.0

IL-13 - Spot 9		
Conc. pg/mL	ECL Signal	%CV
509	956,000	3.0
127	163,000	2.6
31.8	19,200	4.1
7.95	2,810	4.5
1.99	831	4.4
0.497	356	7.8
0.124	238	10.0
0.00	212	10.0

CERTIFICATE OF ANALYSIS

TNF- α - Spot 10		
Conc. pg/mL	ECL Signal	%CV
367	414,000	3.4
91.8	102,000	2.0
22.9	24,900	2.8
5.73	6,030	4.9
1.43	1,900	3.3
0.358	728	3.9
0.0896	431	4.8
0.00	330	6.5


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Plate	Z0047299	2-8°C	31 Aug 2022
Anti-hu IFN- γ Antibody	D0081426	2-8°C	30 Jun 2023
Anti-hu IL-1 β Antibody	D0081182	2-8°C	30 Apr 2022
Anti-hu IL-2 Antibody	D0081320	2-8°C	31 Dec 2022
Anti-hu IL-4 Antibody	D0081342	2-8°C	31 May 2023
Anti-hu IL-6 Antibody	D0081398	2-8°C	30 Apr 2023
Anti-hu IL-8 Antibody	D0081453	2-8°C	31 Jul 2023
Anti-hu IL-10 Antibody	D0081437	2-8°C	30 Jun 2023
Anti-hu IL-12p70 Antibody	D0081307	2-8°C	31 Oct 2022
Anti-hu IL-13 Antibody	D0081381	2-8°C	31 Mar 2023
Anti-hu TNF- α Antibody	D0081486	2-8°C	30 Sep 2023
Proinflammatory Panel 1 (human) Calibrator Blend	A0080242	2-8°C	31 Mar 2025
Diluent 2	M0020112	\leq -10°C	31 Jul 2022
Diluent 3	M0030097	\leq -10°C	31 Aug 2022
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	Karen Hamilla	Quality		25 AUG 2020