

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: K0081491
Expiration Date: 31 Oct 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	1560	pg/mL
		Proinflammatory Panel 1 (human) Control 1	612	pg/mL
		Proinflammatory Panel 1 (human) Control 2	164	pg/mL
		Proinflammatory Panel 1 (human) Control 3	44.2	pg/mL
2	IL-1 β	Calibrator	583	pg/mL
		Proinflammatory Panel 1 (human) Control 1	235	pg/mL
		Proinflammatory Panel 1 (human) Control 2	55.9	pg/mL
		Proinflammatory Panel 1 (human) Control 3	11.9	pg/mL
3	IL-2	Calibrator	1250	pg/mL
		Proinflammatory Panel 1 (human) Control 1	390	pg/mL
		Proinflammatory Panel 1 (human) Control 2	39.8	pg/mL
		Proinflammatory Panel 1 (human) Control 3	3.46	pg/mL
4	IL-4	Calibrator	213	pg/mL
		Proinflammatory Panel 1 (human) Control 1	94.3	pg/mL
		Proinflammatory Panel 1 (human) Control 2	20	pg/mL
		Proinflammatory Panel 1 (human) Control 3	2.8	pg/mL
5	IL-6	Calibrator	663	pg/mL
		Proinflammatory Panel 1 (human) Control 1	255	pg/mL
		Proinflammatory Panel 1 (human) Control 2	47.2	pg/mL
		Proinflammatory Panel 1 (human) Control 3	7.83	pg/mL
6	IL-8	Calibrator	575	pg/mL
		Proinflammatory Panel 1 (human) Control 1	233	pg/mL
		Proinflammatory Panel 1 (human) Control 2	36.8	pg/mL
		Proinflammatory Panel 1 (human) Control 3	6.29	pg/mL
7	IL-10	Calibrator	346	pg/mL
		Proinflammatory Panel 1 (human) Control 1	134	pg/mL
		Proinflammatory Panel 1 (human) Control 2	25.5	pg/mL
		Proinflammatory Panel 1 (human) Control 3	3.93	pg/mL
8	IL-12p70	Calibrator	447	pg/mL
		Proinflammatory Panel 1 (human) Control 1	177	pg/mL
		Proinflammatory Panel 1 (human) Control 2	31.9	pg/mL
		Proinflammatory Panel 1 (human) Control 3	6.39	pg/mL
9	IL-13	Calibrator	466	pg/mL
		Proinflammatory Panel 1 (human) Control 1	164	pg/mL
		Proinflammatory Panel 1 (human) Control 2	45.2	pg/mL
		Proinflammatory Panel 1 (human) Control 3	12.1	pg/mL
10	TNF- α	Calibrator	338	pg/mL
		Proinflammatory Panel 1 (human) Control 1	96.2	pg/mL
		Proinflammatory Panel 1 (human) Control 2	16.9	pg/mL
		Proinflammatory Panel 1 (human) Control 3	2.32	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
1560	408,000	1.7
390	112,000	2.9
97.5	28,600	2.2
24.4	7,230	3.2
6.09	2,030	3.2
1.52	711	6.3
0.381	371	8.4
0.00	280	24.0

IL-1 β - Spot 2		
Conc. pg/mL	ECL Signal	%CV
583	1,640,000	1.7
146	460,000	1.9
36.4	117,000	1.1
9.11	28,800	1.9
2.28	8,200	2.1
0.569	2,820	3.0
0.142	1,430	4.2
0.00	1,080	16.0

IL-2 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
1250	1,100,000	2.6
313	318,000	1.8
78.1	85,600	3.1
19.5	21,500	2.9
4.88	5,720	2.4
1.22	1,640	2.9
0.305	637	5.7
0.00	292	11.0

IL-4 - Spot 4		
Conc. pg/mL	ECL Signal	%CV
213	869,000	2.7
53.3	265,000	3.1
13.3	68,400	3.6
3.33	17,800	3.2
0.832	4,530	6.2
0.208	1,400	12.0
0.0520	490	15.0
0.00	188	16.0

IL-6 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
663	980,000	2.4
166	226,000	2.4
41.4	47,800	2.2
10.4	11,200	1.7
2.59	2,880	2.7
0.647	865	5.0
0.162	347	5.3
0.00	174	11.0

IL-8 - Spot 6		
Conc. pg/mL	ECL Signal	%CV
575	1,020,000	1.9
144	259,000	2.0
35.9	64,100	1.6
8.98	15,400	1.5
2.25	4,200	1.6
0.562	1,220	2.5
0.140	478	6.6
0.00	226	12.0

IL-10 - Spot 7		
Conc. pg/mL	ECL Signal	%CV
346	692,000	1.7
86.5	207,000	1.4
21.6	54,100	2.2
5.41	13,200	1.9
1.35	3,540	1.6
0.338	1,030	4.1
0.0845	391	7.7
0.00	176	15.0

IL-12p70 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
447	344,000	5.0
112	105,000	2.9
27.9	27,300	3.9
6.98	6,940	4.6
1.75	1,990	4.5
0.437	732	10.0
0.109	365	10.0
0.00	294	13.0

IL-13 - Spot 9		
Conc. pg/mL	ECL Signal	%CV
466	856,000	2.7
117	152,000	1.5
29.1	20,300	3.4
7.28	2,980	2.6
1.82	872	3.6
0.455	343	4.3
0.114	216	10.0
0.00	180	11.0

CERTIFICATE OF ANALYSIS

Page 4 of 4

TNF- α - Spot 10		
Conc. pg/mL	ECL Signal	%CV
338	409,000	2.6
84.5	104,000	3.9
21.1	25,500	3.3
5.28	6,340	4.7
1.32	1,800	2.4
0.330	601	5.1
0.0825	298	6.2
0.00	202	11.0

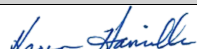
Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Plate	Z0047334	2-8°C	30 Nov 2022
Anti-hu IFN- γ Antibody	D0081474	2-8°C	30 Nov 2023
Anti-hu IL-1 β Antibody	D0081380	2-8°C	31 Mar 2023
Anti-hu IL-2 Antibody	D0081510	2-8°C	30 Sep 2023
Anti-hu IL-4 Antibody	D0081342	2-8°C	31 May 2023
Anti-hu IL-6 Antibody	D0081420	2-8°C	30 Apr 2023
Anti-hu IL-8 Antibody	D0081511	2-8°C	30 Sep 2023
Anti-hu IL-10 Antibody	D0081496	2-8°C	31 Oct 2023
Anti-hu 12p70 Antibody	D0081339	2-8°C	31 Jan 2023
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	A0080262	2-8°C	30 Jun 2025
Diluent 2	M0020113	\leq -10°C	30 Nov 2022
Diluent 3	M0030100	\leq -10°C	31 Oct 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		29 JUL 2020