

CERTIFICATE OF ANALYSIS

Page 1 of 4

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

Kit Catalog Numbers: K15048-Series; K152QR-Series; K152QS-Series; K152QT-Series; K152QU-Series; K152QV-Series

Kit Lot Number: K0081503

Expiration Date: 31 Aug 2022

Acceptance Data/ Percent Recovery Specifications:

		Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IFN- γ	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
2	IL-1 β	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
3	IL-2	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
4	IL-4	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
5	IL-5	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
6	IL-6	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
7	KC/GRO	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
8	IL-10	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
9	IL-12p70	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%
10	TNF- α	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IFN- γ	0.390	75% - 125%	< 25%	570	75% - 125%	< 25%
2	IL-1 β	0.720	75% - 125%	< 25%	1030	75% - 125%	< 25%
3	IL-2	1.03	75% - 125%	< 25%	1570	75% - 125%	< 25%
4	IL-4	0.818	75% - 125%	< 25%	1060	75% - 125%	< 25%
5	IL-5	0.302	75% - 125%	< 25%	590	75% - 125%	< 25%
6	IL-6	7.61	75% - 125%	< 25%	3140	75% - 125%	< 25%
7	KC/GRO	3.29	75% - 125%	< 25%	1230	75% - 125%	< 25%
8	IL-10	7.26	75% - 125%	< 25%	2030	75% - 125%	< 25%
9	IL-12p70	179	75% - 125%	< 25%	20600	75% - 125%	< 25%
10	TNF- α	0.980	75% - 125%	< 25%	403	75% - 125%	< 25%

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-5	< 1.0%
6	IL-6	< 1.0%
7	KC/GRO	< 1.0%
8	IL-10	< 1.0%
9	IL-12p70	< 1.0%
10	TNF- α	< 1.0%

CERTIFICATE OF ANALYSIS

Page 2 of 4

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IFN- γ	Calibrator	764	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	469	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	32	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	2.12	pg/mL
2	IL-1 β	Calibrator	1510	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	722	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	52	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	4.32	pg/mL
3	IL-2	Calibrator	2430	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	908	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	54.5	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	5.57	pg/mL
4	IL-4	Calibrator	1660	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	479	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	62.1	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	14.5	pg/mL
5	IL-5	Calibrator	945	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	356	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	54.2	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	8.3	pg/mL
6	IL-6	Calibrator	5120	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	990	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	152	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	28.9	pg/mL
7	KC/GRO	Calibrator	2150	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	709	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	119	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	17.3	pg/mL
8	IL-10	Calibrator	2630	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	1100	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	331	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	124	pg/mL
9	IL-12p70	Calibrator	29800	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	14400	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	3710	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	822	pg/mL
10	TNF- α	Calibrator	572	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	158	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	28.1	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	4.46	pg/mL

Controls tested may not be the same lot as shipped.

CERTIFICATE OF ANALYSIS

Page 3 of 4

Standard Curve Data:

IFN- γ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
764	1,140,000	2.2
191	306,000	2.5
47.8	74,800	2.7
11.9	18,500	2.8
2.98	4,760	3.8
0.746	1,400	3.3
0.187	504	11.0
0.00	256	32.0

IL-1 β - Spot 2		
Conc. pg/mL	ECL Signal	%CV
1510	524,000	4.0
378	144,000	4.2
94.4	37,800	2.9
23.6	9,380	4.3
5.90	2,570	4.5
1.47	798	4.7
0.369	342	8.2
0.00	226	14.0

IL-2 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
2430	996,000	3.4
608	249,000	4.0
152	62,300	4.5
38.0	14,100	5.4
9.49	3,820	4.9
2.37	1,050	3.4
0.593	411	7.7
0.00	189	14.0

IL-4 - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1660	1,030,000	2.7
415	292,000	2.2
104	77,600	2.0
25.9	20,000	3.0
6.48	5,370	2.6
1.62	1,710	5.7
0.405	734	12.0
0.00	370	19.0

IL-5 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
945	1,190,000	2.3
236	264,000	2.6
59.1	60,000	2.6
14.8	14,100	2.7
3.69	3,820	2.8
0.923	1,150	3.2
0.231	472	5.8
0.00	267	16.0

IL-6 - Spot 6		
Conc. pg/mL	ECL Signal	%CV
5120	638,000	2.1
1280	143,000	2.3
320	34,400	2.1
80.0	8,420	2.1
20.0	2,350	2.1
5.00	767	3.1
1.25	363	4.7
0.00	236	11.0

KC/GRO - Spot 7		
Conc. pg/mL	ECL Signal	%CV
2150	2,150,000	2.1
538	473,000	3.2
134	100,000	3.6
33.6	21,900	4.6
8.40	5,550	2.7
2.10	1,540	5.1
0.525	537	6.0
0.00	187	19.0

IL-10 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
2630	594,000	4.2
658	132,000	3.4
164	32,700	3.2
41.1	7,960	4.2
10.3	2,350	2.7
2.57	882	4.8
0.642	506	6.9
0.00	387	10.0

IL-12p70 - Spot 9		
Conc. pg/mL	ECL Signal	%CV
29800	1,250,000	2.2
7450	283,000	2.5
1860	51,000	2.1
466	9,570	2.8
116	2,410	4.0
29.1	768	4.3
7.28	373	7.9
0.00	248	9.5

CERTIFICATE OF ANALYSIS

Page 4 of 4

TNF- α - Spot 10		
Conc. pg/mL	ECL Signal	%CV
572	390,000	3.0
143	95,800	3.3
35.8	24,100	2.5
8.94	5,860	3.6
2.23	1,690	1.9
0.559	578	5.0
0.140	299	9.9
0.00	207	8.9

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (mouse) Plate	Z0047349	2-8°C	30 Nov 2022
Anti-ms IFN- γ Antibody	D0081487	2-8°C	31 Oct 2023
Anti-ms IL-1 β Antibody	D0081439	2-8°C	30 Jun 2023
Anti-ms IL-2 Antibody	D0081518	2-8°C	31 Oct 2023
Anti-ms IL-4 Antibody	D0081441	2-8°C	30 Jun 2023
Anti-ms IL-5 Antibody	D0081290	2-8°C	31 Dec 2022
Anti-ms IL-6 Antibody	D0081443	2-8°C	31 Jul 2023
Anti-ms KC/GRO Antibody	D0081444	2-8°C	30 Jun 2023
Anti-ms IL-10 Antibody	D0081291	2-8°C	31 Mar 2023
Anti-ms IL-12p70 Antibody	D0081446	2-8°C	31 Aug 2023
Anti-ms TNF- α Antibody	D0081359	2-8°C	30 Apr 2023
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080250	2-8°C	28 Feb 2025
Diluent 41	M0410032	$\leq -10^{\circ}\text{C}$	31 Aug 2022
Diluent 45	M0450033	$\leq -10^{\circ}\text{C}$	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 (mouse) Control 1	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 2	Not Kit Specific
Proinflammatory Panel 1 (mouse) 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		04 AUG 2020