

**CERTIFICATE OF ANALYSIS**

Page 1 of 3

**Product Description:** V-PLEX® and V-PLEX Plus Vascular Injury Panel 2 (human) Kit

**Kit Catalog Numbers:** K15198-Series

**Kit Lot Number:** K0080911

**Expiration Date:** 30 Sep 2019

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		Vascular Injury Control 1		Vascular Injury Control 2		Vascular Injury Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu SAA	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	hu CRP	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
3	hu VCAM-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	hu ICAM-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu SAA	54.0	80% - 120%	< 20%	138000	80% - 120%	< 20%
2	hu CRP	27.6	80% - 120%	< 20%	49600	80% - 120%	< 20%
3	hu VCAM-1	37.6	80% - 120%	< 20%	32000	80% - 120%	< 20%
4	hu ICAM-1	6.40	80% - 120%	< 20%	32700	80% - 120%	< 20%

**Acceptance Data/ Non Specific Binding Specifications:**

Spot	Name	NSB Spec
1	hu SAA	< 1.0%
2	hu CRP	< 1.0%
3	hu VCAM-1	< 1.0%
4	hu ICAM-1	< 1.0%

**CERTIFICATE OF ANALYSIS**

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	hu SAA	20X Calibrator	4800000	pg/mL
		Vascular Injury Control 1	70500	pg/mL
		Vascular Injury Control 2	6620	pg/mL
		Vascular Injury Control 3	833	pg/mL
2	hu CRP	20X Calibrator	4040000	pg/mL
		Vascular Injury Control 1	24300	pg/mL
		Vascular Injury Control 2	4900	pg/mL
		Vascular Injury Control 3	604	pg/mL
3	hu VCAM-1	20X Calibrator	1038000	pg/mL
		Vascular Injury Control 1	13900	pg/mL
		Vascular Injury Control 2	1331	pg/mL
		Vascular Injury Control 3	165	pg/mL
4	hu ICAM-1	20X Calibrator	968000	pg/mL
		Vascular Injury Control 1	16300	pg/mL
		Vascular Injury Control 2	1720	pg/mL
		Vascular Injury Control 3	209	pg/mL

Controls tested may not be the same lot as shipped.

Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentrations, please refer to the Standard Curve Data

**Standard Curve Data:**

hu SAA - Spot 1			hu CRP - Spot 2			hu VCAM-1 - Spot 3		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
240000	798,000	3.9	202000	1,120,000	2.3	51900	486,000	2.9
48000	161,000	3.8	40400	445,000	2.3	10400	121,000	2.3
9600	29,000	3.4	8080	104,000	2.8	2080	20,600	3.1
1920	5,670	3.8	1620	21,900	3.1	415	3,220	3.5
384	1,200	3.3	323	4,470	3.6	83.0	673	1.9
76.8	296	4.3	64.6	976	3.0	16.6	184	4.0
15.4	106	5.0	12.9	251	3.8	3.32	87	6.0
0.00	60	8.8	0.00	69	12.0	0.00	64	8.4

**CERTIFICATE OF ANALYSIS**

Page 3 of 3

hu ICAM-1 - Spot 4		
Conc. pg/mL	ECL Signal	%CV
48400	612,000	3.5
9680	132,000	4.2
1940	27,600	4.0
387	5,560	6.1
77.4	1,230	5.6
15.5	295	5.5
3.10	125	12.0
0.00	73	19.0


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Vascular Injury Panel 2 Plate	Z0056102	2-8°C	30 Apr 2020
Anti-hu SAA Antibody	D0081042	2-8°C	30 Apr 2020
Anti-hu CRP Antibody	D0081021	2-8°C	31 Jan 2020
Anti-hu VCAM-1 Antibody	D0080980	2-8°C	30 Sep 2019
Anti-hu ICAM-1 Antibody	D0081043	2-8°C	30 Apr 2020
Vascular Injury Panel 2 (human) Calibrator Blend	A0080132	≤ -70°C	30 Sep 2019
Diluent 101	M1010029	≤ -10°C	31 Mar 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)	Not Kit Specific
Vascular Injury Control 1	Not Kit Specific
Vascular Injury Control 2	Not Kit Specific
Vascular Injury 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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		16 JAN 2018