

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Vascular Injury Panel 2 (human) Kit  
**Kit Lot Number:** K0082015  
**Expiration Date:** 29 Feb 2024  
**Kit Catalog Numbers:** K15198-Series; K151SS-Series; K151SR-Series; K151SU-Series

**Precision and Accuracy Specifications:**

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

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	SAA	54.0	80% - 120%	< 20%	138000	80% - 120%	< 20%
2	CRP	27.6	80% - 120%	< 20%	49600	80% - 120%	< 20%
3	VCAM-1	37.6	80% - 120%	< 20%	32000	80% - 120%	< 20%
4	ICAM-1	15.0	80% - 120%	< 20%	32700	80% - 120%	< 20%

**Non-Specific Binding Specifications:**

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

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	SAA	4260000	pg/mL
2	CRP	3860000	pg/mL
3	VCAM-1	896000	pg/mL
4	ICAM-1	928000	pg/mL


**CERTIFICATE OF ANALYSIS**

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Vascular Injury Panel 2 Plate	Z0057076	2-8°C	31 May 2024
Anti-hu SAA Antibody	D0081725	2-8°C	31 Mar 2025
Anti-hu CRP Antibody	D0081724	2-8°C	31 Mar 2025
Anti-hu VCAM-1 Antibody	D0081727	2-8°C	31 Jan 2025
Anti-hu ICAM-1 Antibody	D0081728	2-8°C	28 Feb 2025
Vascular Injury Panel 2 (human) Calibrator Blend	A0080302	≤ -70°C	29 Feb 2024
Diluent 101	M1010045	≤ -10°C	30 Jun 2024

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		21 JUN 2022