

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX Plus Vascular Injury Panel 2 (human) Kit
Lot Number: K0082516
Expiration Date: 31 DEC 2025
Catalog Number: K15198-Series, K151SS-Series, K151SR-Series, K151SU-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu CRP Antibody	D0082298	2-8°C	31 DEC 2027
Anti-hu ICAM-1 Antibody	D0082300	2-8°C	31 JAN 2028
Anti-hu SAA Antibody	D0082297	2-8°C	31 DEC 2027
Anti-hu VCAM-1 Antibody	D0082299	2-8°C	29 FEB 2028
Vascular Injury Panel 2 (human) Plate	Z0057481	2-8°C	31 DEC 2026
Diluent 101	M1010052	≤ -10°C	31 MAY 2026
Vascular Injury Panel 2 (human) Calibrator Blend	A0080367	≤ -70°C	31 DEC 2025

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

Precision, Accuracy, and Sensitivity 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	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%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	SAA	54	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	80% - 120%	< 20%	32700	80% - 120%	< 20%

Non-Specific Binding Specifications:

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




CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	SAA	4840000	pg/mL
2	CRP	3920000	pg/mL
3	VCAM-1	990000	pg/mL
4	ICAM-1	1006000	pg/mL

Dilution to top of Curve: 20X

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

	Name	Function	Signature	Date
Review/Approval	Hamilla, Karen	Quality		27 SEP 2024