

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Vascular Injury Panel 2 (human) Kit**Lot Number:** K0082862**Expiration Date:** 30 NOV 2027**Catalog Number:** K15198-Series, K151SS-Series, K151SR-Series, K151SU-Series**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu CRP Antibody	D0082525	2-8°C	31 MAR 2029
Anti-hu ICAM-1 Antibody	D0082527	2-8°C	31 MAR 2029
Anti-hu SAA Antibody	D0082275	2-8°C	30 NOV 2027
Anti-hu VCAM-1 Antibody	D0082526	2-8°C	30 APR 2029
Vascular Injury Panel 2 (hu) 96-Well 4-Spot SECTOR Plate	Z0057611	2-8°C	29 FEB 2028
Diluent 101	M1010057	≤ -10°C	31 MAR 2028
Vascular Injury Panel 2 (human) Calibrator Blend	A0080445	≤ -70°C	31 MAR 2028

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	54.0	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%

*See Additional Comments

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	3200000	pg/mL
2	CRP	3800000	pg/mL
3	VCAM-1	1022000	pg/mL
4	ICAM-1	1014000	pg/mL

Dilution to top of Curve: 20X

Additional Comments:

Spot	Name	Specification	Comment
1	SAA	LLOQ Conc.	% LLOQ measured 140% (specification 80-120%) at target 54.0pg/mL, was fully conforming at 366pg/mL.
3	VCAM-1	LLOQ Conc.	% LLOQ measured 131% (specification 80-120%) at target 37.6pg/mL, was fully conforming at 39.3pg/mL.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		06 MAR 2026