

**Extension Memo**

Product Description: V-PLEX® and V-PLEX PLUS Vascular Injury Panel 2 (human) Kit
Lot Number: K0082800
New Expiration Date: 31 DEC 2026
Catalog Number(s): K15198-Series, K151SS-Series, K151SR-Series, K151SU-Series

Based on additional testing, the original date of expiration for this kit has been extended by 12 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Vascular Injury Panel 2 (human) Calibrator Blend	A0080367	31 DEC 2025	31 DEC 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		04 Nov 2025

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX PLUS Vascular Injury Panel 2 (human) Kit**Lot Number:** K0082800**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	D0082363	2-8°C	31 AUG 2028
Anti-hu ICAM-1 Antibody	D0082278	2-8°C	30 NOV 2027
Anti-hu SAA Antibody	D0082362	2-8°C	31 MAR 2028
Anti-hu VCAM-1 Antibody	D0082364	2-8°C	31 MAR 2028
Vascular Injury Panel 2 (hu) 96-Well 4-Spot SECTOR Plate	Z0057517	2-8°C	31 MAR 2027
Diluent 101	M1010056	≤ -10°C	30 NOV 2027
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:

		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	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%



CERTIFICATE OF ANALYSIS

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%

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	Karen Hamilla	Quality		03 NOV 2025