



Extension Memo

Product Description: V-PLEX and V-PLEX Plus Vascular Injury Panel 2 (human) Kit
Lot Number: K0082671
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 7 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	A0080396	30 SEP 2026	31 DEC 2026
Diluent 101	M1010052	31 MAY 2026	31 DEC 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		18 Mar 2026



CERTIFICATE OF ANALYSIS

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

Lot Number: K0082671

Expiration Date: 31 MAY 2026

Catalog Number: K15198-Series, K151SS-Series, K151SR-Series, K151SU-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu CRP Antibody	D0082436	2-8°C	31 AUG 2028
Anti-hu ICAM-1 Antibody	D0082438	2-8°C	31 AUG 2028
Anti-hu SAA Antibody	D0082435	2-8°C	31 OCT 2028
Anti-hu VCAM-1 Antibody	D0082437	2-8°C	31 AUG 2028
Vascular Injury Panel 2 (hu) 96-Well 4-Spot SECTOR Plate	Z0057560	2-8°C	31 AUG 2027
Diluent 101	M1010052	≤ -10°C	31 MAY 2026
Vascular Injury Panel 2 (human) Calibrator Blend	A0080396	≤ -70°C	30 SEP 2026

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%



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	3340000	pg/mL
2	CRP	3820000	pg/mL
3	VCAM-1	1016000	pg/mL
4	ICAM-1	974000	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		24 JUL 2025