


**Extension Memo**

**Product Description:** V-PLEX® and V-PLEX Plus Human CRP Kit  
**Catalog Number(s):** K151ST-Series  
**Lot Number:** K0082564  
**New Expiration Date:** 31 DEC 2026

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 Calibrator Blend	A0080367	31 DEC 2025	31 DEC 2026
Diluent 101	M1010052	31 MAY 2026	31 DEC 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		23 Sep 2025



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX Plus Human CRP Kit  
**Lot Number:** K0082564  
**Expiration Date:** 31 DEC 2025  
**Catalog Number:** K151ST-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu CRP Antibody	D0082298	2-8°C	31 DEC 2027
Human CRP 96-Well 4-Spot SECTOR Plate	Z0057511	2-8°C	31 MAR 2027
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:**

		Calibrator		Vascular Injury Control 1		Vascular Injury Control 2		Vascular Injury Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	CRP	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%


		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	CRP	27.6	80% - 120%	< 20%	49600	80% - 120%	< 20%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	CRP	3920000	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	Michelle Likiam	Quality		03 Dec 2024