

**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:**

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

Spot	Name	LLOQ			ULOQ		
		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