

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Human CRP Kit**Lot Number:** K0082940**Expiration Date:** 30 NOV 2027**Catalog Number:** K151ST-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	D0082363	2-8°C	31 AUG 2028
Human CRP 96-Well 4-Spot SECTOR Plate	Z0057690	2-8°C	30 SEP 2028
Diluent 101	M1010056	≤ -10°C	30 NOV 2027
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	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	3800000	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
<b>Review/Approval</b>	Karen Hamilla	Quality		15 MAY 2026