

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human CRP Kit

**Kit Lot Number:** K0082388

**Expiration Date:** 28 Feb 2025

**Kit Catalog Numbers:** K151ST-Series

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human CRP Plate	Z0057419	2-8°C	30 Apr 2026
Anti-hu CRP Antibody	D0082158	2-8°C	31 Mar 2027
Vascular Injury Panel 2 (human) Calibrator Blend	A0080356	≤ -70°C	31 Jul 2025
Diluent 101	M1010047	≤ -10°C	28 Feb 2025

Additional components may be provided to execute the assay. Additional components are not lot-specific.

### Precision and Accuracy 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	3860000	pg/mL

Dilution to Top of Curve = 20X

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		01 Dec 2023