

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human CRP Kit  
**Kit Lot Number:** K0081852  
**Expiration Date:** 31 May 2023  
**Kit Catalog Numbers:** K151ST-Series

**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	3720000	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human CRP Plate	Z0056868	2-8°C	31 Jul 2023
Anti-hu CRP Antibody	D0081404	2-8°C	31 May 2023
Vascular Injury Panel 2 (human) Calibrator Blend	A0080284	≤ -70°C	31 Oct 2023
Diluent 101	M1010042	≤ -10°C	30 Nov 2023

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		07 OCT 2021