

CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Kit Lot Number: K0081646

Expiration Date: 31 May 2022

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	4000000	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	A0080234	≤ -70°C	31 May 2022
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	Carissa B. Hubbard	Quality		25 JUN 2021