

CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Kit Lot Number: K0081933

Expiration Date: 29 Feb 2024

Kit Catalog Numbers: K151ST-Series

Precision and Accuracy Specifications:

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

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	3860000	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human CRP Plate	Z0057087	2-8°C	30 Jun 2024
Anti-hu CRP Antibody	D0081724	2-8°C	31 Mar 2025
Vascular Panel 2 (human) Calibrator Blend	A0080302	≤ -70°C	29 Feb 2024
Diluent 101	M1010045	≤ -10°C	30 Jun 2024

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		18 APR 2022