

## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Kit Lot Number:** K0081801

**Expiration Date:** 30 Nov 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                           | Z0057011   | 2-8°C               | 31 Dec 2023     |
| Anti-hu CRP Antibody                      | D0081726   | 2-8°C               | 28 Feb 2025     |
| Vascular Panel 2 (human) Calibrator Blend | A0080297   | ≤ -70°C             | 30 Nov 2023     |
| Diluent 101                               | M1010043   | ≤ -10°C             | 31 Jan 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  |  | 22 NOV 2021 |