

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human CRP Kit
Kit Lot Number: K0082048
Expiration Date: 29 Feb 2024
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 | 3860000 | pg/mL |

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|---------------------|-----------------|
| Human CRP Plate | Z0057185 | 2-8°C | 30 Nov 2024 |
| Anti-hu CRP Antibody | D0081870 | 2-8°C | 30 Nov 2025 |
| Vascular Injury Panel 2 (human) Calibrator Blend | A0080302 | ≤ -70°C | 29 Feb 2024 |
| Diluent 101 | M1010046 | ≤ -10°C | 31 Dec 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 |  | 19 AUG 2022 |