

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Vascular Injury Panel 2 (human) Kit
Kit Lot Number: K0081767
Expiration Date: 31 Aug 2023
Kit Catalog Numbers: K15198-Series; K151SS-Series, K151SR-Series; K151SU-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 | SAA | 80% - 120% | < 20% | 75% - 125% | < 20% | 75% - 125% | < 20% | 75% - 125% | < 20% |
| 2 | CRP | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |
| 3 | VCAM-1 | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |
| 4 | ICAM-1 | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |

| | | LLOQ | | | ULOQ | | |
|------|--------|-------|------------|-------|--------|------------|-------|
| Spot | Name | Value | Recovery | CV | Value | Recovery | CV |
| 1 | SAA | 54.0 | 80% - 120% | < 20% | 138000 | 80% - 120% | < 20% |
| 2 | CRP | 27.6 | 80% - 120% | < 20% | 49600 | 80% - 120% | < 20% |
| 3 | VCAM-1 | 37.6 | 80% - 120% | < 20% | 32000 | 80% - 120% | < 20% |
| 4 | ICAM-1 | 15.0 | 80% - 120% | < 20% | 32700 | 80% - 120% | < 20% |

Non-Specific Binding Specifications:

| Spot | Name | NSB Spec |
|------|--------|----------|
| 1 | SAA | < 1.0% |
| 2 | CRP | < 1.0% |
| 3 | VCAM-1 | < 1.0% |
| 4 | ICAM-1 | < 1.0% |

Stock Concentration of Calibrator:

| Spot | Name | Conc. | Units |
|------|--------|---------|-------|
| 1 | SAA | 3880000 | pg/mL |
| 2 | CRP | 3720000 | pg/mL |
| 3 | VCAM-1 | 964000 | pg/mL |
| 4 | ICAM-1 | 914000 | pg/mL |


CERTIFICATE OF ANALYSIS

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|---------------------|-----------------|
| Vascular Injury Panel 2 Plate | Z0056928 | 2-8°C | 31 Aug 2023 |
| Anti-hu SAA Antibody | D0081657 | 2-8°C | 31 Aug 2024 |
| Anti-hu CRP Antibody | D0081620 | 2-8°C | 31 Jul 2024 |
| Anti-hu VCAM-1 Antibody | D0081658 | 2-8°C | 31 Oct 2024 |
| Anti-hu ICAM-1 Antibody | D0081659 | 2-8°C | 30 Sep 2024 |
| 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 |  | 23 AUG 2021 |