



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Vascular Injury Panel 2 (human) Calibrator Blend

**Catalog Number:** C0198-2

**Lot Number:** A0080234

**Expiration Date:** 31 MAY 2022

**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

### Concentration Assignment:

| Analyte | Stock Concentration | Unit of Measure |
|---------|---------------------|-----------------|
| SAA     | 4860000             | pg/mL           |
| CRP     | 4000000             | pg/mL           |
| VCAM-1  | 988000              | pg/mL           |
| ICAM-1  | 998000              | pg/mL           |

**Dilution to Top of Curve:** 20X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

|                        | Name          | Function | Signature | Date        |
|------------------------|---------------|----------|-----------|-------------|
| <b>Review/Approval</b> | Karen Hamilla | Quality  |           | 02 JAN 2020 |