



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0198-2

**Lot Number:** A0080284

**Expiration Date:** 31 OCT 2023

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

### Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
SAA	3880000	pg/mL
CRP	3720000	pg/mL
VCAM-1	964000	pg/mL
ICAM-1	914000	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		19 MAY 2021