



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C0198-2

Lot Number: A0080341

Expiration Date: 28 FEB 2025

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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
SAA	4380000	pg/mL
CRP	3780000	pg/mL
VCAM-1	1002000	pg/mL
ICAM-1	936000	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
Review/Approval	Karen Hamilla	Quality		11 OCT 2022