

**CERTIFICATE OF ANALYSIS**

Product Description: Vascular Injury Panel 2 (human) Calibrator Blend
Catalog Number: C0687-2
Lot Number: A0080445
Expiration Date: 31 MAR 2028
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
SAA	3200000	pg/mL
CRP	3800000	pg/mL
VCAM-1	1022000	pg/mL
ICAM-1	1014000	pg/mL

Dilution to Top of Curve: 20 X

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		11 NOV 2025