



### CERTIFICATE OF ANALYSIS

**Product Description:** Vascular Injury Panel 2 (human) Calibrator Blend  
**Catalog Number:** C0687-2  
**Lot Number:** A0080367  
**Expiration Date:** 31 DEC 2025  
**Storage Temperature:** ≤ -70°C

**Concentration Assignment:**

Analyte	Stock Concentration	Unit of Measure
SAA	4840000	pg/mL
CRP	3920000	pg/mL
VCAM-1	990000	pg/mL
ICAM-1	1006000	pg/mL

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

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
<b>Review/Approval</b>	Karen Hamilla	Quality		04 AUG 2023