



CERTIFICATE OF ANALYSIS

Product Description: Vascular Injury Panel 2 (human) Calibrator Blend
Catalog Number: C0687-2
Lot Number: A0080396
Expiration Date: 30 SEP 2026
Storage Temperature: ≤ -70°C

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
SAA	3340000	pg/mL
CRP	3820000	pg/mL
VCAM-1	1016000	pg/mL
ICAM-1	974000	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		16 APR 2024