



Extension Memo

Product Description: Vascular Injury Panel 2 (human) Calibrator Blend
Lot Number: A0080367
New Expiration Date: 31 DEC 2026
Catalog Number(s): C0687-2

Based on additional testing, the original date of expiration for this calibrator has been extended by 12 months.

Function	Name	Signature	Date
Quality	Karen Hamilla		12 Nov 2025



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

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
Review/Approval	Karen Hamilla	Quality		04 AUG 2023