



CERTIFICATE OF ANALYSIS

Control Pack Description: Vascular Injury Control Pack 1
Catalog Number: C4687-1
Lot Number: CP000159
Expiration Date: 31 OCT 2026

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Vascular Injury Control 1	A00C0936	≤ -70°C	31 OCT 2026
Vascular Injury Control 2	A00C0937	≤ -70°C	31 OCT 2026
Vascular Injury Control 3	A00C0938	≤ -70°C	31 OCT 2026

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
SAA	37000	3780	511	pg/mL
CRP	24400	4770	638	pg/mL
VCAM-1	12000	1160	158	pg/mL
ICAM-1	11500	1270	150	pg/mL

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		22 MAY 2024