

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Vascular Injury Control Pack 1
Catalog Number: C4687-1
Lot Number: CP000150
Expiration Date: 31 JAN 2026

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Vascular Injury Control 1	A00C0915	≤ -70°C	31 JAN 2026
Vascular Injury Control 2	A00C0916	≤ -70°C	31 JAN 2026
Vascular Injury Control 3	A00C0917	≤ -70°C	31 JAN 2026

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
SAA	72900	8280	818	pg/mL
CRP	25300	5530	683	pg/mL
VCAM-1	13400	1410	167	pg/mL
ICAM-1	12900	1460	177	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		28 SEP 2023