



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Control Pack Description:** Vascular Injury Control Pack 1

**Catalog Number:** C4198-1

**Lot Number:** CP000111

**Expiration Date:** 30 JUN 2023

**Concentration Assignment:**

Analyte	Control 1	Control 2	Control 3	Unit of Measure
SAA	48900	4460	548	pg/mL
CRP	26100	4890	630	pg/mL
VCAM-1	13500	1150	141	pg/mL
ICAM-1	12700	1260	144	pg/mL

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Vascular Injury Control 1	A00C0749	≤ -70°C	30 JUN 2023
Vascular Injury Control 2	A00C0750	≤ -70°C	30 JUN 2023
Vascular Injury Control 3	A00C0751	≤ -70°C	30 JUN 2023

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		13 JAN 2021