



CERTIFICATE OF ANALYSIS

Page 1 of 1

Control Pack Description: Vascular Injury Control Pack 1

Catalog Number: C4198-1

Lot Number: CP000087

Expiration Date: 31 AUG 2021

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
SAA	48500	4720	607	pg/mL
CRP	24600	4800	610	pg/mL
VCAM-1	16600	1600	200	pg/mL
ICAM-1	13000	1350	166	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Vascular Injury Control 1	A00C0604	≤ -70°C	31 AUG 2021
Vascular Injury Control 2	A00C0605	≤ -70°C	31 AUG 2021
Vascular Injury Control 3	A00C0606	≤ -70°C	31 AUG 2021

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		17 SEP 2020