



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Vascular Injury Panel 2 (human) Calibrator Blend
Catalog Number: C0198-2
Lot Number: A0080212
Expiration Date: 30 NOV 2021
Storage Temperature: ≤ -70°C

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
SAA	4960000	pg/mL
CRP	3920000	pg/mL
VCAM-1	958000	pg/mL
ICAM-1	1002000	pg/mL

Dilution to Top of Curve: 20X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		15 JUL 2019