



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Vascular Injury Panel 2 (human) Calibrator Blend

Catalog Number: C0198-2

Lot Number: A0080132

Expiration Date: 30 SEP 2019

Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration of Materials Tested:

	Assay			
	hu SAA	hu CRP	hu VCAM-1	hu ICAM-1
Reconstituted Calibrator (pg/mL)	4 800 000	4 040 000	1 038 000	968 000

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	06 JUN 2017