



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-04725

Product Description: Vascular Injury Panel 2 (human) Calibrator Blend
Catalog Number: C0198-2
Lot Number: A0040098
Expiration Date: 31 DEC 2016
Storage Temperature: ≤ -70°C

Concentration of Materials Tested:

	Assay			
	SAA	CRP	VCAM-1	ICAM-1
Calibrator (pg/mL)	4 150 000	3 650 000	983 000	919 000

Statement: The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dana Short	Quality Control		11 SEP 2014