



September 2022

Re: Vascular Injury Control Pack 1 (Catalog No. C4198-1) Lot No. CP000123, Revised Concentrations

Dear Valued Customer,

As part of our process for continuous product improvement, we periodically review the performance specifications of V-PLEX® products. As we accumulate data from additional lots, if the larger data set indicates a change to a more accurate value, we update the product accordingly.

We recently reviewed the data for the Controls in our Vascular Injury Control Pack 1. Based on the larger data set, we have decided to reassign the values of these controls for all of the analytes in this panel. We do not think that the controls have changed. Rather, the larger data set provides more accurate assignments. The table below shows the revised and original quantitation.

Units = pg/mL	Assay			
	hu SAA	hu CRP	hu VCAM-1	hu ICAM-1
Control 1 (A00C0789)	48 300	23 500	13 100	12 200
Control 2 (A00C0790)	4920	4530	1160	1300
Control 3 (A00C0791)	856	550	152	159

Table 1. Revised concentrations

Units = pg/mL	Assay			
	hu SAA	hu CRP	hu VCAM-1	hu ICAM-1
Control 1 (A00C0789)	49 600	23 700	13 400	12 200
Control 2 (A00C0790)	5310	4680	1240	1370
Control 3 (A00C0791)	939	575	158	175

Table 2. Original concentrations

If you have any questions or concerns, please contact MSD Scientific Support at +1 240-314-2798 or scientificsupport@mesoscale.com.

Sincerely,

MSD Quality Assurance



CERTIFICATE OF ANALYSIS

Page 1 of 1

Control Pack Description: Vascular Injury Control Pack 1

Catalog Number: C4198-1

Lot Number: CP000123

Expiration Date: 29 FEB 2024

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
SAA	48300	4920	856	pg/mL
CRP	23500	4530	550	pg/mL
VCAM-1	13100	1160	152	pg/mL
ICAM-1	12200	1300	159	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Vascular Injury Control 1	A00C0789	≤ -70°C	29 FEB 2024
Vascular Injury Control 2	A00C0790	≤ -70°C	29 FEB 2024
Vascular Injury Control 3	A00C0791	≤ -70°C	29 FEB 2024

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		27 SEP 2022