



**RFE MEMO**

Page 1 of 1

**Product Description:** Vascular Injury Control Pack 1

**Catalog Number:** C4198-1

**Lot Number:** CP000075

Based on additional testing, the original date of expiration for this control pack has been extended by 22 months. The expiration date of CP000075 has been extended to 30 SEP 2022.

The expiration dates for the following controls have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Vascular Injury Control 1	A00C0601	31 OCT 2020	30 SEP 2022
Vascular Injury Control 2	A00C0602	31 OCT 2020	30 SEP 2022
Vascular Injury Control 3	A00C0603	31 OCT 2020	30 SEP 2022

Function	Name	Signature	Date
Quality	Karen Hamilla		29 SEP 2021

This memo supersedes all other statements regarding date of expiration for this product.



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Vascular Injury Control Pack 1

**Catalog Number:** C4198-1

**Lot Number:** CP000075

**Expiration Date:** 31 OCT 2020

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

### Concentration of Materials Tested:

Units = pg/mL	Assay			
	SAA	CRP	VCAM-1	ICAM-1
Control 1 (A00C0601)	49 800	25 100	13 900	12 900
Control 2 (A00C0602)	4640	5060	1290	1320
Control 3 (A00C0603)	597	639	160	160

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Vascular Injury Control 1	A00C0601	$\leq -70^{\circ}\text{C}$	31 OCT 2020
Vascular Injury Control 2	A00C0602	$\leq -70^{\circ}\text{C}$	31 OCT 2020
Vascular Injury Control 3	A00C0603	$\leq -70^{\circ}\text{C}$	31 OCT 2020

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	30 MAY 2018