



<b>Title: Internal Certificate of Analysis</b>			
<b>Revision Code: A</b>	<b>Level: 3</b>	<b>Effective Date: 15MAR2017</b>	<b>Pages: 1 of 1</b>

**Product Description:** Inflammation Panel 1 (rat) Control Pack  
**Catalog Number:** N/A  
**Lot Number:** CP000061  
**Expiration Date:** 30 JUN 2022  
**Storage Temperature:** ≤ -70°C

**Concentration of Materials Tested:**

Units = ng/mL	Assay			
	Lipocalin-2	TSP-1	TIMP-1	MCP-1
Control 1 (A00C0470)	225	9150	66.9	88.1
Control 2 (A00C0471)	20.4	5500	17.6	18.5
Control 3 (A00C0472)	2.12	3290	3.65	3.63

**Concentration: 20X**

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Inflammation Panel 1 (rat) Control 1	A00C0470	≤ -70°C	30 JUN 2022
Inflammation Panel 1 (rat) Control 2	A00C0471	≤ -70°C	30 JUN 2022
Inflammation Panel 1 (rat) Control 3	A00C0472	≤ -70°C	30 JUN 2022

**The above product is intended for Internal Testing Use Only. Not for use by Customers.**

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	26 NOV 2018