



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

Product Description: Inflammation Panel 1 (rat) Calibrator Blend
Catalog Number: C0179-2
Lot Number: A0040189
Expiration Date: 31 DEC 2019
Storage Temperature: ≤ -70°C

Concentration of Materials Tested:

	Assay			
	Lipocalin-2	TSP-1	TIMP-1	MCP-1
Calibrator (ng/mL)	1032	85900	193	292

Concentration: 20X

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

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