



Document #: **QA-FM-189**

<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:** Cardiac Injury Panel 3 (rat) Calibrator Blend  
**Catalog Number:** C0161-2  
**Lot Number:** A0040218  
**Expiration Date:** 31 MAR 2021  
**Storage Temperature:** ≤ -70°C

**Concentration of Materials Tested:**

	Assay			
	cTnl	cTnT	FABP3	MyI3
Calibrator (ng/mL)	722	1210	1696	1154

**Concentration:** 20X

**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>	24 JAN 2019