



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Muscle Injury Panel 1 (rat) Calibrator Blend

Catalog Number: C0181-2

Lot Number: A0040216

Expiration Date: 31 MAR 2021

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

Concentration of Materials Tested:

	Assay				
	cTnI	cTnT	FABP3	Myl3	sTnI
Calibrator (ng/mL)	522	1010	1916	990	3900

Concentration: 20X

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	28 JAN 2019