



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Muscle Injury Panel 3 (mouse) Calibrator Blend

Catalog Number: C0186-2

Lot Number: A0040196

Expiration Date: 30 NOV 2019

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

Concentration of Materials Tested:

	Assay			
	cTnl	FABP3	Myl3	sTnl
Calibrator (ng/mL)	480	1966	1300	3860

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>	04 DEC 2017