



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0186-2

**Lot Number:** A0040244

**Expiration Date:** 30 JUN 2022


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

### Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
cTnI	618	ng/mL
FABP3	1550	ng/mL
MyI3	866	ng/mL
sTnI	4680	ng/mL

**Dilution to Top of Curve:** 20X

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		16 JAN 2020