



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C0186-2

Lot Number: A0040254

Expiration Date: 28 FEB 2023

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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
cTnI	928	ng/mL
FABP3	1776	ng/mL
MyI3	928	ng/mL
sTnI	5720	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
Review/Approval	Karen Hamilla	Quality		16 DEC 2020