



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C0186-2

Lot Number: A0040249

Expiration Date: 30 SEP 2022

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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
cTnI	662	ng/mL
FABP3	1706	ng/mL
MyI3	956	ng/mL
sTnI	4820	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 JUL 2020