



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C0181-2

Lot Number: A0040258

Expiration Date: 30 SEP 2022

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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
cTnI	354	ng/mL
cTnT	1036	ng/mL
FABP3	2060	ng/mL
Myl3	1402	ng/mL
sTnI	4300	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		25 FEB 2021