



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C0181-2

Lot Number: A0040253

Expiration Date: 30 APR 2022

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

Concentration Assignment:

| Analyte | Stock Concentration | Unit of Measure |
|---------|---------------------|-----------------|
| cTnI | 386 | ng/mL |
| cTnT | 1118 | ng/mL |
| FABP3 | 1830 | ng/mL |
| MyI3 | 1070 | ng/mL |
| sTnI | 3820 | 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 | | 03 SEP 2020 |