



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C0181-2

Lot Number: A0040206

Expiration Date: 31 OCT 2020

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

Concentration of Materials Tested:

| | Assay | | | | |
|-----------------------|-------|------|-------|------|------|
| | cTnI | cTnT | FABP3 | Myl3 | sTnI |
| Calibrator (ng/mL) | 608 | 1328 | 1698 | 992 | 4120 |

Concentration: 20X

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

| | Name | Function | Signature | Date |
|------------------------|--------------|-------------------|---------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 22 JUN 2018 |