



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Muscle Injury Panel 2 (rat) Calibrator Blend

**Catalog Number:** C0180-2

**Lot Number:** A0040252

**Expiration Date:** 31 JAN 2023

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

### Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
TIMP-1	882	ng/mL
CK	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
<b>Review/Approval</b>	Karen Hamilla	Quality		09 NOV 2020