



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C0180-2

Lot Number: A0040266

Expiration Date: 31 DEC 2024

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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
TIMP-1	870	ng/mL
CK	4040	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		01 AUG 2022