



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C0180-2

Lot Number: A0040211

Expiration Date: 31 DEC 2020

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

Concentration of Materials Tested:

	Assay	
	TIMP-1	CK
Calibrator (ng/mL)	782	4600

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>	27 JUL 2018