



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Rat Skeletal Troponin I Calibrator

Catalog Number: C03IM-2

Lot Number: A0040215

Expiration Date: 31 DEC 2020

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

Concentration of Materials Tested:

	Assay
	rt sTnI
Calibrator (pg/mL)	4 140 000

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>	28 AUG 2018