



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Rat Skeletal Troponin I Calibrator

**Catalog Number:** C03IM-2

**Lot Number:** A0040267

**Expiration Date:** 30 JUN 2024


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

### Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
sTnI	3960000	pg/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		10 FEB 2022