

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** Rat Skeletal Troponin I Kit  
**Kit Catalog Numbers:** K153IMC-series  
**Kit Lot Number:** K0040430  
**Expiration Date:** 31 Jul 2019

### Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Muscle/Cardiac Panel Control 1 (20X)		Muscle/Cardiac Panel Control 2 (20X)		Muscle/Cardiac Panel Control 3 (20X)	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	rat sTnI	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	rat sTnI	781	80% - 120%	< 20%	16000	80% - 120%	< 25%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	rat sTnI	20X Calibrator	4140000	pg/mL
		Muscle/Cardiac Panel Control 1 (20X)	2000000	pg/mL
		Muscle/Cardiac Panel Control 2 (20X)	297000	pg/mL
		Muscle/Cardiac Panel Control 3 (20X)	32700	pg/mL

Control data was generated utilizing internal reference controls

Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentrations, please refer to the Standard Curve Data

### Standard Curve Data:

rat sTnI - Spot 1		
Conc. pg/mL	ECL Signal	%CV
207000	327,000	3.4
69000	91,500	6.3
23000	28,000	3.6
7670	7,920	5.3
2560	2,990	6.4
852	1,060	4.2
284	474	3.4
0.00	166	4.3


**CERTIFICATE OF ANALYSIS**

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Rat Skeletal Troponin I Plate	Z0056285	2-8°C	30 Nov 2020
Anti-rat sTnI Antibody	D0040772	2-8°C	30 Nov 2021
Rat Skeletal Troponin I Calibrator	A0040215	≤ -70°C	31 Dec 2020
Diluent 7	M0070055	≤ -10°C	31 Dec 2020
Diluent 30	M0300022	≤ -10°C	30 Nov 2020
DTT	R0010583	≤ -10°C	31 Jul 2019
EDTA, pH 8.0	R0010621	Room Temperature	31 Aug 2020
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

**Statement:**

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		25 SEP 2018