

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** Muscle Injury Panel 1 (rat) Kit  
**Kit Lot Number:** K0040508  
**Expiration Date:** 31 Oct 2021  
**Kit Catalog Numbers:** K15181C-Series

**Precision and Accuracy Specifications:**

		Calibrator		Muscle Injury Panel 1 Control 1		Muscle Injury Panel 1 Control 2		Muscle Injury Panel 1 Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	cTnI	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	cTnT	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
3	FABP3	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
5	MyI3	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
6	sTnI	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	cTnI	0.0977	80% - 120%	< 20%	20.0	80% - 120%	< 20%
2	cTnT	0.488	80% - 120%	< 20%	35.0	80% - 120%	< 20%
3	FABP3	0.391	75% - 125%	< 25%	15.0	75% - 125%	< 25%
5	MyI3	0.170	75% - 125%	< 25%	44.0	80% - 120%	< 20%
6	sTnI	0.781	80% - 120%	< 20%	160	80% - 120%	< 20%

**Non-Specific Binding Specifications:**

Spot	Name	cTnI	cTnT	FABP3	MyI3	sTnI
1	cTnI	NA	< 1.0%	< 1.0%	< 1.0%	< 65.0%
2	cTnT	< 1.0%	NA	< 1.0%	< 1.0%	< 15.0%*
3	FABP3	< 1.0%	< 1.0%	NA	< 1.0%	< 1.0%
5	MyI3	< 1.0%	< 1.0%	< 1.0%	NA	< 1.0%
6	sTnI	< 1.0%	< 1.0%	< 1.0%	< 1.0%	NA

\* See Additional Comments.

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	cTnI	576	ng/mL
2	cTnT	1180	ng/mL
3	FABP3	1758	ng/mL
5	MyI3	1132	ng/mL
6	sTnI	3440	ng/mL

## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Muscle Injury Panel 1 (rat) Plate	Z0056739	2-8°C	31 Dec 2022
Anti-rcTnI Antibody	D0040848	2-8°C	31 Jan 2023
Anti-rat cTnT Antibody	D0040921	2-8°C	31 Dec 2023
Anti-rat FABP3 Antibody	D0040899	2-8°C	31 Oct 2023
Anti-rat MyI3 Antibody	D0040759	2-8°C	31 Oct 2021
Anti-rat sTnI Antibody	D0040922	2-8°C	31 Dec 2023
Muscle Injury Panel 1 (rat) Calibrator Blend	A0040256	≤ -70°C	30 Jun 2022
Diluent 7	M0070064	≤ -10°C	30 Nov 2022
Diluent 30	M0300031	≤ -10°C	31 Dec 2022

### Additional Comments:

	Assay	Spot	Specification	Comment
1	cTnI	1	< 15.0%	NSB sTnI on cTnT = 19.0%

Additional components may be provided to execute the assay. Additional components are not lot-specific.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		17 NOV 2020