

# MSD<sup>®</sup> MULTI-SPOT Assay System

## Muscle Injury Panel Kits

**cTnl, cTnT, FABP3, Myl3, and sTnl**

|                                  | <b>Catalog No.</b> |
|----------------------------------|--------------------|
| Muscle Injury Panel 1 (rat) Kit  | K15181C            |
| Cardiac Injury Panel 2 (rat) Kit | K15155C            |
| Cardiac Injury Panel 3 (rat) Kit | K15161C            |



# MSD Toxicology Assays

## Muscle Injury Panel 1 (rat) Kit

cTnl, cTnT, FABP3, Myl3, sTnl

## Cardiac Injury Panel 2 (rat) Kit

cTnl, cTnT, FABP3

## Cardiac Injury Panel 3 (rat) Kit

cTnl, cTnT, FABP3, Myl3

*This package insert must be read in its entirety before using this product.*

**FOR RESEARCH USE ONLY.**

**NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

## MESO SCALE DISCOVERY<sup>®</sup>

A division of Meso Scale Diagnostics, LLC.

1601 Research Blvd.

Rockville, MD 20850 USA

[www.mesoscale.com](http://www.mesoscale.com)

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, mesoscale.com, www.mesoscale.com, methodicalmind.com, www.methodicalmind.com, DISCOVERY WORKBENCH, InstrumentLink, MESO, MesoAccess, MesoAdvantage, MesoSphere, Methodical Mind, Methodical Mind Enterprise, MSD GOLD, MULTI-ARRAY, MULTI-SPOT, QuickPlex, Parsec, ProductLink, SECTOR, SECTOR PR, SECTOR HTS, SULFO-TAG, TeamLink, TrueSensitivity, TURBO-BOOST, TURBO-TAG, N-PLEX, R-PLEX, S-PLEX, T-PLEX, U-PLEX, V-PLEX, MSD (design), MSD (luminous design), Methodical Mind (design), Parsec (design), 96 WELL SMALL-SPOT (design), 96 WELL 1-, 4-, 7-, 9-, & 10-SPOT (designs), 384 WELL 1- & 4-SPOT (designs), N-PLEX (design), R-PLEX (design), S-PLEX (design), T-PLEX (design), U-PLEX (design), V-PLEX (design), It's All About U, SPOT THE DIFFERENCE, The Biomarker Company, and The Methodical Mind Experience are trademarks and/or service marks owned by or licensed to Meso Scale Diagnostics, LLC. All other trademarks and service marks are the property of their respective owners.  
©2013, 2018, 2021 Meso Scale Diagnostics, LLC. All rights reserved.

# Table of Contents

|                                                       |    |
|-------------------------------------------------------|----|
| Introduction .....                                    | 4  |
| Principle of the Assay .....                          | 5  |
| Kit Components.....                                   | 6  |
| Additional Materials and Equipment—not supplied ..... | 7  |
| Safety .....                                          | 7  |
| Best Practices.....                                   | 8  |
| Reagent Preparation .....                             | 9  |
| Protocol .....                                        | 12 |
| Analysis of Results .....                             | 12 |
| Assay Qualification and Verification.....             | 12 |
| Typical Data .....                                    | 13 |
| Sensitivity.....                                      | 14 |
| Precision.....                                        | 16 |
| Dilution Linearity .....                              | 16 |
| Spike Recovery .....                                  | 20 |
| Specificity .....                                     | 23 |
| Assay Components .....                                | 25 |
| References.....                                       | 26 |
| Summary Protocol .....                                | 27 |
| Plate Diagrams .....                                  | 27 |

## Contact Information

### MSD Customer Service

Phone: 1-240-314-2795

Fax: 1-301-990-2776

Email: [CustomerService@mesoscale.com](mailto:CustomerService@mesoscale.com)

### MSD Scientific Support

Phone: 1-240-314-2798

Fax: 1-240-632-2219 attn: Scientific Support

Email: [ScientificSupport@mesoscale.com](mailto:ScientificSupport@mesoscale.com)

# Introduction

**Troponin** is a heterotrimer that regulates muscle contraction in skeletal and cardiac muscle (but not in smooth muscle). Troponin acts with intracellular calcium to control the interaction of actin and myosin filaments in striated muscle fibers. Though they perform similar functions, cardiac and skeletal troponins differ in sequence and can be differentiated in immunoassays.

The three subunits of troponin are:

- **Troponin T** is the subunit that interacts with tropomyosin to form the troponin-tropomyosin complex.
- **Troponin I** is an inhibitory subunit that prevents muscle contraction in the absence of calcium. It is responsible for the binding of the troponin-tropomyosin complex to actin. Troponin I exists in three isoforms: slow-twitch (striated) skeletal muscle, fast-twitch (striated) skeletal muscle, and cardiac muscle.
- **Troponin C** binds calcium, producing a conformational change in troponin I and activating the troponin-tropomyosin complex.

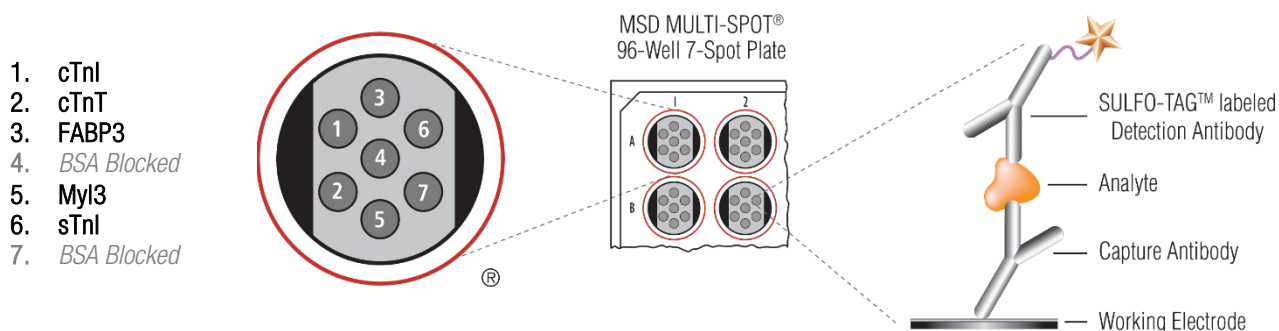
When muscle tissue is damaged, the troponin-tropomyosin complex breaks down and troponin I and troponin T are released into the blood. Cardiac troponin T (cTnT) and cardiac troponin I (cTnI) can be readily distinguished from their skeletal muscle analogs (sTnT and sTnI), allowing confirmation of cardiac muscle tissue damage over skeletal muscle tissue damage. Troponins are excellent biomarkers for research on myocardial injury in cardiotoxicity because of the demonstrated tissue-specificity of cardiac and skeletal troponins.<sup>2</sup>

**Myosin light chain 3 (MyI3)** is an essential light chain of the myosin molecule found in cardiac and slow-twitch skeletal muscle. Myosin is a hexamer ATPase motor protein that is a major constituent of thick muscle filament. The myosin molecule consists of a head domain that “walks” along the actin chain to contract the muscle and a tail domain that is responsible for binding myosin to its cargo. Two heavy chain subunits intertwine to form the head and tail domains. Four light chain subunits—two regulatory light chains with phosphorylation sites (encoded by the MYL2 genes), and two essential light chains (encoded by the MYL3 genes)—bind the heavy chains together in the neck region between the head and tail domains. After damage to muscle tissue, myosin breaks down and MyI3 becomes elevated in the blood. MyI3 can be used during research in conjunction with other toxicity biomarkers to confirm cardiac and slow-twitch skeletal muscle injury.

**Fatty acid binding protein 3 (FABP3)** is a monomeric protein that modulates the uptake of fatty acids in cells. Heart-type fatty acid binding protein is released into circulation after myocardial ischemia and necrosis. FABP3 is mostly present in heart and skeletal muscle but can also be found in brain, liver, and small intestine.

# Principle of the Assay

MSD toxicology assays provide rapid and convenient methods for measuring the levels of protein targets within a single, small-volume sample. The assays in the Muscle Injury Panel 1 (rat) Kit are sandwich immunoassays (Figure 1). MSD provides a plate pre-coated with capture antibodies. The user adds the sample and a solution containing detection antibodies conjugated with electrochemiluminescent MSD SULFO-TAG™ labels throughout one or more incubation periods. Analytes in the sample capture antibodies immobilized on the working electrode surface; recruitment of the detection antibodies by the bound analytes completes the sandwich. The user adds an MSD buffer that provides the appropriate chemical environment for electrochemiluminescence and loads the plate into an MSD® instrument where a voltage applied to the plate electrodes causes the captured labels to emit light. The instrument measures the intensity of emitted light to provide a quantitative measure of analytes in the sample.



*Figure 1. Spot diagram showing the placement of analyte capture antibodies. The numbering convention for the different spots is maintained in the software visualization tools, on the plate packaging, and in the data files.*

# Kit Components

*Table 1. Reagents supplied with all kits*

| Product Description                                   | Storage | Catalog No. | Size   | Quantity per Kit |          |           |
|-------------------------------------------------------|---------|-------------|--------|------------------|----------|-----------|
|                                                       |         |             |        | 1 Plate          | 5 Plates | 25 Plates |
| MULTI-SPOT® 96-Well Muscle Injury Panel 1 (rat) Plate | 2–8 °C  | N75181A-1   | 7-spot | 1 plate          | 5 plates | 25 plates |
| Muscle Injury Panel 1 (rat) Calibrator Blend (20X)    | ≤-70 °C | C0181-2     | 20 µL  | 1 vial           | 5 vials  | 25 vials  |
| Diluent 7                                             | ≤-10 °C | R54BB-4     | 5 mL   | 2 bottles        | —        | —         |
|                                                       |         | R54BB-3     | 50 mL  | —                | 1 bottle | 5 bottles |
| Diluent 30                                            | ≤-10 °C | R50AB-4     | 25 mL  | 1 bottle         | 1 bottle | 5 bottles |
| 25 mM DTT                                             | ≤-10 °C | —           | 1 mL   | 1 vial           | 1 vial   | 5 vials   |
| 0.5 M EDTA pH 8.0                                     | RT      | —           | 4 mL   | 1 bottle         | 1 bottle | 5 bottles |
| Read Buffer T (4X)                                    | RT      | R92TC-3     | 50 mL  | 1 bottle         | 1 bottle | 5 bottles |

RT = room temperature

Dash (—) = not applicable

*Table 2. Components that are supplied with specific kits*

| Product Description                     | Storage | Catalog No. | Size   | Quantity per Kit |          |           |
|-----------------------------------------|---------|-------------|--------|------------------|----------|-----------|
|                                         |         |             |        | 1 Plate          | 5 Plates | 25 Plates |
| SULFO-TAG Anti-rat cTnI Antibody (50X)  | 2–8 °C  | D23HR-2     | 75 µL  | 1 vial           | —        | —         |
|                                         |         | D23HR-3     | 375 µL | —                | 1 vial   | 5 vials   |
| SULFO-TAG Anti-rat cTnT Antibody (50X)  | 2–8 °C  | D23EF-2     | 75 µL  | 1 vial           | —        | —         |
|                                         |         | D23EF-3     | 375 µL | —                | 1 vial   | 5 vials   |
| SULFO-TAG Anti-rat FABP3 Antibody (50X) | 2–8 °C  | D23HT-2     | 75 µL  | 1 vial           | —        | —         |
|                                         |         | D23HT-3     | 375 µL | —                | 1 vial   | 5 vials   |
| SULFO-TAG Anti-rat Myl3 Antibody (50X)  | 2–8 °C  | D23JI-2     | 75 µL  | 1 vial           | —        | —         |
|                                         |         | D23JI-3     | 375 µL | —                | 1 vial   | 5 vials   |
| SULFO-TAG Anti-rat sTnI Antibody (50X)  | 2–8 °C  | D23IM-2     | 75 µL  | 1 vial           | —        | —         |
|                                         |         | D23IM-3     | 375 µL | —                | 1 vial   | 5 vials   |

Dash (—) = not applicable

# Additional Materials and Equipment—not supplied

- Appropriately sized tubes for reagent preparation
- Microcentrifuge tubes for preparing serial dilutions
- Phosphate buffered saline plus 0.05% Tween-20 (PBS-T) for plate washing
- MSD Wash Buffer (Catalog No. R61AA-1)
- Liquid handling equipment for the desired throughput, capable of dispensing 10 to 150  $\mu\text{L}$ /well into a 96-well microtiter plate
- Plate washing equipment: automated plate washer or multichannel pipette
- Adhesive plate seals
- Microtiter plate shaker (rotary) capable of shaking at 500–1,000 rpm
- Deionized water
- Vortex mixer

## Safety

Use safe laboratory practices: wear gloves, safety glasses, and lab coats when handling assay components. Handle and dispose of all hazardous samples properly in accordance with local, state, and federal guidelines.

Additional product-specific safety information is available in the applicable safety data sheet(s) (SDS), which can be obtained from MSD Customer Service or at the [www.mesoscale.com](http://www.mesoscale.com)<sup>®</sup> website.

# Best Practices

- Bring frozen diluent to room temperature in a 22-25 °C water bath. Thaw frozen calibrator (when applicable) on wet ice.
- Prepare Calibrator Standards and samples in polypropylene microcentrifuge tubes. Use a fresh pipette tip for each dilution and mix by vortexing after each dilution.
- Avoid prolonged exposure of the detection antibody (stock or diluted) to light. During the antibody incubation step, plates do not need to be shielded from light (except for direct sunlight).
- Avoid bubbles in wells during all pipetting steps as they may lead to variable results. Bubbles introduced when adding Read Buffer may interfere with signal detection.
- Use reverse pipetting when necessary to avoid the introduction of bubbles. For empty wells, pipette gently to the bottom corner. Do not touch the pipette tip on the bottom of the wells when pipetting into the MSD plate.
- Plate shaking should be vigorous, with a rotary motion between 500 and 1,000 rpm. Binding reactions may reach equilibrium sooner if you use shaking at the middle of this range (~700 rpm) or above.
- When using an automated plate washer, rotate the plate 180 degrees between wash steps to improve assay precision.
- Gently tap the plate on a paper towel to remove residual fluid after washing.
- If an incubation step needs to be extended, leave the sample or detection antibody solution in the plate to keep the plate from drying out.
- Remove the plate seal before reading the plate.
- Make sure that the read buffer is at room temperature when added to a plate.
- Do not shake the plate after adding the read buffer.
- To improve interplate precision, keep time intervals consistent between adding the read buffer and reading the plate. Unless otherwise directed, read the plate as soon as possible after adding the read buffer.
- If the sample results are above the top of the calibration curve, dilute the samples and repeat the assay.
- When running a partial plate, seal the unused sectors to avoid contaminating unused wells. Remove all seals before reading. The uncoated wells of a partially used plate may be sealed and stored up to 30 days at 2–8 °C in the original foil pouch with desiccant. You may adjust volumes proportionally when preparing reagents.



# Reagent Preparation

Bring all reagents to room temperature. Thaw the stock calibrator on ice.

**Important:** Upon the first thaw, separate Diluent 7 and Diluent 30 into aliquots appropriate to the size of your needs before refreezing.

## Prepare Diluent 7 + Additives

Samples and calibrators are diluted in Diluent 7 that contains added EDTA and DTT. These two additives must be added into the diluent by the user before each assay is carried out. EDTA and DTT additive stocks are provided at the concentrations in the table below.

*Table 3. EDTA and DDT additives*

| Additive | Stock Conc.    | Final Conc.  |
|----------|----------------|--------------|
| EDTA     | 500 mM (16.7X) | 30 mM (1X)   |
| DTT      | 25 mM (100X)   | 0.25 mM (1X) |

For one plate, combine:

- 540  $\mu$ L of EDTA stock solution
- 90  $\mu$ L of DTT stock solution
- 8,370  $\mu$ L of Diluent 7

## Prepare Calibrator Solutions

The Calibrator blend for the Muscle Injury Panels is supplied at a 20-fold higher concentration than the recommended highest Calibrator. For each assay, a 7-point standard curve is recommended with 3-fold serial dilution steps and a zero Calibrator blank. Signals from the blank should be excluded when generating the curve. Thaw the stock Calibrator and keep it on ice, then add to diluent at room temperature to make the standard curve solutions. For the actual concentration of each Calibrator in the blend, refer to the certificate of analysis (COA) supplied with the kit or available at [www.mesoscale.com](http://www.mesoscale.com).

To prepare 7 standard solutions plus a zero Calibrator blank for up to 5 replicates:

- Prepare the highest standard by adding 12  $\mu$ L of stock Calibrator to 228  $\mu$ L of Diluent 7 + Additives. Mix well.
- Prepare the next standard by transferring 80  $\mu$ L of the highest standard to 160  $\mu$ L of Diluent 7 + Additives. Mix well. Repeat 3-fold serial dilutions 5 additional times to generate 7 standards.
- Use Diluent 7 + Additives as the blank.

After preparation of the Calibrators at the concentrations above, incubate the Calibrator solutions without shaking for 30 minutes at room temperature before addition to the plate.

## Dilute Samples

Rat serum and plasma samples should be run at 4-fold dilution. Diluent 7 + Additives should be used to dilute the samples. To perform sample dilution, add 25  $\mu$ L of sample to 75  $\mu$ L of Diluent 7 + Additives. Depending on the sample set under investigation, higher or lower dilution factors may be necessary. If needed, additional Diluent 7 can be purchased at [www.mesoscale.com](http://www.mesoscale.com).

Diluted samples should be incubated at room temperature without shaking for 30 minutes before adding to the plate.

## Prepare Detection Antibody Solution

MSD provides each detection antibody as a 50X stock solution. The working detection antibody solution is 1X.

For one plate of Muscle Injury Panel 1 (rat), combine:

- 60  $\mu$ L of 50X SULFO-TAG Anti-rat cTnl Antibody
- 60  $\mu$ L of 50X SULFO-TAG Anti-rat cTnT Antibody
- 60  $\mu$ L of 50X SULFO-TAG Anti-rat FABP3 Antibody
- 60  $\mu$ L of 50X SULFO-TAG Anti-rat Myl3 Antibody
- 60  $\mu$ L of 50X SULFO-TAG Anti-rat sTnl Antibody
- 2,700  $\mu$ L of Diluent 30

For one plate of Cardiac Injury Panel 3 (rat), combine:

- 60  $\mu$ L of 50X SULFO-TAG Anti-rat cTnl Antibody
- 60  $\mu$ L of 50X SULFO-TAG Anti-rat cTnT Antibody
- 60  $\mu$ L of 50X SULFO-TAG Anti-rat FABP3 Antibody
- 60  $\mu$ L of 50X SULFO-TAG Anti-rat Myl3 Antibody
- 2,760  $\mu$ L of Diluent 30

For one plate of Cardiac Injury Panel 2 (rat), combine:

- 60  $\mu$ L of 50X SULFO-TAG Anti-rat cTnl Antibody
- 60  $\mu$ L of 50X SULFO-TAG Anti-rat cTnT Antibody
- 60  $\mu$ L of 50X SULFO-TAG Anti-rat FABP3 Antibody
- 2,820  $\mu$ L of Diluent 30

## Prepare Wash Buffer

MSD provides Wash Buffer (Catalog No. R61AA-1) as a 20X stock solution. The working solution is 1X. PBS + 0.05% Tween-20 can be used instead.

For one plate, combine:

- 15 mL of MSD Wash Buffer (20X)
- 285 mL of deionized water
- 1X MSD Wash Buffer can be stored at room temperature for up to two weeks.

## Prepare Read Buffer

MSD provides Read Buffer T as a 4X stock solution. The working solution is 1X.

For one plate, combine:

- 5 mL of Read Buffer T (4X)
- 15 mL of deionized water

You may prepare diluted read buffer in advance and store it at room temperature in a tightly sealed container.

## Prepare MSD Plate

MSD plates are precoated with capture antibodies (Figure 1) and exposed to a proprietary stabilizing treatment to ensure the integrity and stability of the immobilized antibodies. Plates can be used as delivered; no additional preparation (e.g., prewetting) is required.

# Protocol

**Note:** Complete the steps described in **Reagent Preparation** (above) before beginning this assay protocol.

## Step 1: Add Diluent 7 + Additives

- ❑ Add 25  $\mu\text{L}$  of Diluent 7 + Additives solution to each well. Seal the plate with an adhesive plate seal and incubate for 30 minutes with vigorous shaking (500–1,000 rpm) at room temperature.

## STEP 1: Step 2: Add Sample or Calibrator

- ❑ Add 25  $\mu\text{L}$  of sample or Calibrator (which has been preincubated for 30 min following dilution with Diluent 7 + Additives) per well. Seal the plate with an adhesive plate seal and incubate for 2 hours with vigorous shaking (500–1,000 rpm) at room temperature

**Note:** You may prepare detection antibody solution during incubation.

## Step 3: Wash and Add Detection Antibody Solution

- ❑ Wash the plate 3 times with at least 300  $\mu\text{L}$ /well of 1X MSD Wash Buffer or PBS-T.
- ❑ Add 25  $\mu\text{L}$  of 1X detection antibody solution to each well. Seal the plate with an adhesive plate seal and incubate for 2 hours with vigorous shaking (500–1,000 rpm) at room temperature.

**Note:** You may prepare diluted read buffer during incubation.

## Step 4: Wash and Read

- ❑ Wash the plate 3 times with at least 300  $\mu\text{L}$ /well of 1X MSD Wash Buffer or PBS-T.
- ❑ Add 150  $\mu\text{L}$  of 1X Read Buffer T to each well. Analyze the plate on an MSD instrument. Incubation in read buffer is not required before reading the plate.

# Analysis of Results

Run at least one set of calibrators in duplicate to generate the standard curve. The standard curve is modeled using least squares fitting algorithms so that signals from the calibrators can be used to calculate the concentration of analytes in the samples. The assays have a wide dynamic range (3–4 logs) which allows accurate quantification without the need for dilution in many cases. The MSD DISCOVERY WORKBENCH<sup>®</sup> analysis software uses a 4-parameter logistic model (or sigmoidal dose-response) and includes a  $1/Y^2$  weighting function. The weighting function is important because it provides a better fit of data over a wide dynamic range, particularly at the low end of the standard curve.

# Assay Qualification and Verification

The performance of this Kit meets levels of consistency and robustness as determined by methods based on the principles outlined in “Fit-for-Purpose Method Development and Validation for Successful Biomarker Measurement” by Lee, et al.<sup>1</sup>

Bioanalytical and functional characterizations of calibrators, antibodies, and assay components are completed to allow for bridging of reagents between lots. This includes plate coating uniformity and reagent and component specificity testing for individual kit lots.

Control samples for specific matrices are designed and tested to meet the accuracy, precision, and sensitivity criteria for a Kit that has completed the qualification process. Spike recovery and dilution linearity of endogenous samples, pooled and individual matrices are tested across the assay range.

## Sensitivity, Range, and Curve Fitting

- Sample range and assay sensitivity are established from 4-PL fitted calibration curves with  $1/Y^2$  weighting. Percent recovery of calibrators and controls between the upper limit of quantification (ULOQ) and lower limit of quantification (LLOQ) must have a calculated concentration %CV of less than 20% and accuracy within 20% of the expected concentration.
- The limits of quantification defined in the product insert are verified for each lot as part of the lot verification and quality control release.

## Accuracy and Precision

High, mid, and low controls made in the matrix (need to be defined on a kit-by-kit basis) are run to measure accuracy and precision.

- Qualification – Testing on multiple days (>6 days) and multiple runs per day for a total of 15–20 runs of complete kits. Precision is measured for the standard curve for intra- and interday %CV of less than 20%. %CV and accuracy of the controls are measured on all runs and must meet the kit specification as defined in the certificate of analysis (COA). The typical calculated concentration %CV specification is less than 20% and accuracy within 20% of expected concentration and a total error of less than 30%.
- Verification – A multiday (2–3 days), multiple runs per day for a total of 6–12 plates is performed as part of the release testing for each lot. The specifications for release are provided in the COA.

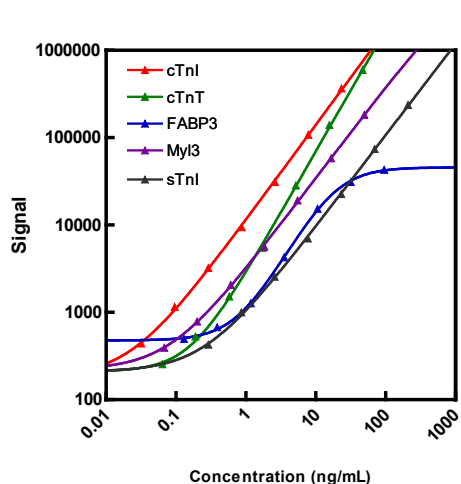
## Robustness and Stability

- Freeze-thaw testing and accelerated stability studies performed during assay development (calibrators, antibodies, controls) are augmented with real-time stability studies on complete kits out to 18 months from the date of manufacture.

All acceptance criteria and verification conformance are defined in the COA for all kit lots. Presented below are representative data from the assay qualification for this assay that meets the criteria described above. The actual kit-specific standard curve and measured limits of quantification can be found in the COA enclosed with the kit.

# Typical Data

The following standard curve and data illustrate the dynamic range of the assay. Actual signals will vary. The best quantification of unknown samples will be achieved by generating a standard curve for each plate using a minimum of two replicates of standards. For each kit lot, refer to the COA for the actual concentration of the Calibrator.



| cTnI          |                |      |
|---------------|----------------|------|
| Conc. (ng/mL) | Average Signal | %CV  |
| 0             | 107            | 12.4 |
| 0.0321        | 440            | 6.2  |
| 0.0963        | 1,155          | 7.0  |
| 0.289         | 3,214          | 6.4  |
| 0.867         | 9,433          | 5.5  |
| 2.60          | 31,018         | 4.5  |
| 7.80          | 107,673        | 4.0  |
| 23.4          | 365,820        | 7.0  |

| cTnT          |                |     |
|---------------|----------------|-----|
| Conc. (ng/mL) | Average Signal | %CV |
| 0             | 120            | 8.5 |
| 0.0645        | 254            | 6.8 |
| 0.193         | 529            | 6.7 |
| 0.580         | 1,519          | 5.6 |
| 1.74          | 5,468          | 5.1 |
| 5.22          | 28,325         | 4.1 |
| 15.7          | 139,526        | 5.9 |
| 47.0          | 594,838        | 5.6 |

| FABP3         |                |     |
|---------------|----------------|-----|
| Conc. (ng/mL) | Average Signal | %CV |
| 0             | 421            | 6.8 |
| 0.131         | 496            | 7.4 |
| 0.392         | 675            | 4.9 |
| 1.18          | 1,270          | 7.1 |
| 3.53          | 4,257          | 5.2 |
| 10.6          | 15,362         | 5.7 |
| 31.7          | 31,031         | 6.0 |
| 95.2          | 42,881         | 6.4 |

| MyI3          |                |      |
|---------------|----------------|------|
| Conc. (ng/mL) | Average Signal | %CV  |
| 0             | 167            | 10.8 |
| 0.0682        | 394            | 4.8  |
| 0.205         | 785            | 6.6  |
| 0.614         | 2,066          | 5.1  |
| 1.84          | 5,645          | 6.4  |
| 5.52          | 19,059         | 3.9  |
| 16.6          | 58,134         | 3.0  |
| 49.7          | 182,582        | 3.5  |

| sTnI          |                |     |
|---------------|----------------|-----|
| Conc. (ng/mL) | Average Signal | %CV |
| 0             | 165            | 9.6 |
| 0.286         | 427            | 5.8 |
| 0.858         | 996            | 3.7 |
| 2.58          | 2,556          | 3.0 |
| 7.74          | 7,019          | 3.8 |
| 23.2          | 22,614         | 3.5 |
| 69.7          | 74,092         | 4.5 |
| 209           | 237,149        | 6.1 |

# Sensitivity

The lower limit of detection (LLOD) is a calculated concentration based on a signal 2.5 standard deviations above the background (zero calibrator).

A multiplate, multiday study was performed to measure the reproducibility of the assay. The lower limit of quantification (LLOQ) and upper limit of quantification (ULOQ) were established from the multiple plate run.

The LLOQ is determined as the lowest concentration where the %CV of the calculated concentration is less than 20% and the percent recovery of the standard is between 80% and 120%. For FABP3 and Myl3, the specification on percent recovery was widened to 75% to 125%.

The ULOQ is determined as the highest concentration where the %CV of the calculated concentration is less than 20% and the percent recovery of the standard is between 80% and 120%. For FABP3, the specification on percent recovery was widened to 75% to 125%.

**Table 4. LLOD, LLOQ, and ULOQ for each analyte in the Muscle Injury Panel (rat) Kits**

|      | cTnl    | cTnT    | FABP3   | Myl3    | sTnl    |
|------|---------|---------|---------|---------|---------|
|      | (ng/mL) | (ng/mL) | (ng/mL) | (ng/mL) | (ng/mL) |
| LLOD | 0.00400 | 0.0171  | 0.155   | 0.0865  | 0.0472  |
| LLOQ | 0.0977  | 0.488   | 0.391   | 0.170   | 0.781   |
| ULOQ | 20.0    | 35.0    | 15.0    | 44.0    | 160     |

# Precision

Control samples of high, mid, and low levels of each analyte were measured on each plate. Normal rat serum, rat soleus homogenate, and assay Calibrators are used to make control samples. The high control contains 25% normal rat serum and Calibrators. The mid control contains rat soleus homogenate and Calibrators. The low control contains only the assay Calibrators.

The controls were run in triplicate or quadruplicate on each of 9 plates run across 3 days.

Average Intraplate %CV is the average %CV of the control replicates within an individual plate.

Interplate %CV is the variability of controls across 9 plates over 3 days.

**Table 5. Intraplate and Interplate %CVs for each analyte in the Muscle Injury Panel (rat) Kits**

|       | Control | Plates | Average Conc. (ng/mL) | Average Intraplate %CV | Interplate %CV |
|-------|---------|--------|-----------------------|------------------------|----------------|
| cTnl  | High    | 9      | 9.36                  | 3.8                    | 5.9            |
|       | Mid     | 9      | 1.31                  | 3.9                    | 6.0            |
|       | Low     | 9      | 0.278                 | 4.0                    | 5.1            |
| cTnT  | High    | 9      | 35.4                  | 3.9                    | 5.0            |
|       | Mid     | 9      | 6.46                  | 2.5                    | 4.0            |
|       | Low     | 9      | 1.06                  | 3.3                    | 4.6            |
| FAPB3 | High    | 9      | 13.9                  | 5.9                    | 7.5            |
|       | Mid     | 9      | 8.44                  | 5.5                    | 7.6            |
|       | Low     | 9      | 2.85                  | 2.8                    | 4.8            |
| MyI3  | High    | 9      | 30.3                  | 5.9                    | 7.0            |
|       | Mid     | 9      | 3.01                  | 3.8                    | 6.9            |
|       | Low     | 9      | 0.319                 | 6.0                    | 7.1            |
| sTnl  | High    | 9      | 108                   | 4.3                    | 5.8            |
|       | Mid     | 9      | 17.4                  | 2.4                    | 3.1            |
|       | Low     | 9      | 2.88                  | 3.5                    | 10.5           |

# Dilution Linearity

To assess linearity, serum, EDTA plasma and heparin plasma samples from Sprague-Dawley rats were diluted 2-fold, 4-fold, 8-fold, and 16-fold before testing. All samples were taken from normal rats. The concentrations shown below have been corrected for dilution (concentration = measured × dilution factor). Measurements that were outside of the quantitative range are shown in italics. Percent recovery is calculated as the measured concentration divided by the concentration measured from the previous dilution (expected).

$$\% \text{ recovery} = \frac{\text{measured concentration}}{\text{expected concentration}} \times 100$$



Table 6: Dilution Linearity: cTnI and cTnT

| Sample               | Fold Dilution | cTnI          |           |            | cTnT          |           |            |
|----------------------|---------------|---------------|-----------|------------|---------------|-----------|------------|
|                      |               | Conc. (ng/mL) | Conc. %CV | % Recovery | Conc. (ng/mL) | Conc. %CV | % Recovery |
| Serum No. 1          | 2             | 19.4          | 0.2       | —          | 13.2          | 2.9       | —          |
|                      | 4             | 17.3          | 3.8       | 89         | 12.5          | 1.4       | 95         |
|                      | 8             | 16.8          | 7.1       | 97         | 12.4          | 2.2       | 99         |
|                      | 16            | 15.7          | 6.2       | 94         | 11.4          | 5.1       | 93         |
| Serum No. 2          | 2             | 5.82          | 7.6       | —          | 3.7           | 11.0      | —          |
|                      | 4             | 5.99          | 4.2       | 103        | 3.3           | 1.8       | 89         |
|                      | 8             | 6.32          | 1.4       | 105        | < LLOQ        | —         | —          |
|                      | 16            | 5.72          | 2.4       | 91         | < LLOQ        | —         | —          |
| Serum No. 3          | 2             | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
|                      | 4             | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
|                      | 8             | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
|                      | 16            | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
| EDTA Plasma No. 1    | 2             | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
|                      | 4             | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
|                      | 8             | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
|                      | 16            | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
| EDTA Plasma No. 2    | 2             | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
|                      | 4             | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
|                      | 8             | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
|                      | 16            | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
| EDTA Plasma No. 3    | 2             | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
|                      | 4             | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
|                      | 8             | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
|                      | 16            | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
| Heparin Plasma No. 1 | 2             | 1.62          | 2.0       | —          | < LLOQ        | —         | —          |
|                      | 4             | 1.57          | 2.8       | 97         | < LLOQ        | —         | —          |
|                      | 8             | 1.52          | 2.6       | 97         | < LLOQ        | —         | —          |
|                      | 16            | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
| Heparin Plasma No. 2 | 2             | 1.30          | 4.2       | —          | < LLOQ        | —         | —          |
|                      | 4             | 1.25          | 5.1       | 96         | < LLOQ        | —         | —          |
|                      | 8             | 1.29          | 1.2       | 103        | < LLOQ        | —         | —          |
|                      | 16            | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
| Heparin Plasma No. 3 | 2             | 1.30          | 5.0       | —          | < LLOQ        | —         | —          |
|                      | 4             | 1.27          | 5.7       | 98         | < LLOQ        | —         | —          |
|                      | 8             | 1.24          | 2.7       | 97         | < LLOQ        | —         | —          |
|                      | 16            | < LLOQ        | —         | —          | < LLOQ        | —         | —          |

LLOQ = lower limit of quantification

Dash (-) = not applicable

Table 7. Dilution Linearity: FABP3 and Myl3

| Sample               | Fold Dilution | FABP3         |           |            | Myl3          |           |            |
|----------------------|---------------|---------------|-----------|------------|---------------|-----------|------------|
|                      |               | Conc. (ng/mL) | Conc. %CV | % Recovery | Conc. (ng/mL) | Conc. %CV | % Recovery |
| Serum No. 1          | 2             | > ULOQ        | —         | —          | 15.5          | 5.5       | —          |
|                      | 4             | > ULOQ        | —         | —          | 13.4          | 0.5       | 86         |
|                      | 8             | > ULOQ        | —         | —          | 13.8          | 4.6       | 103        |
|                      | 16            | > ULOQ        | —         | —          | 12.9          | 2.2       | 93         |
| Serum No. 2          | 2             | > ULOQ        | —         | —          | 8.32          | 11.3      | —          |
|                      | 4             | > ULOQ        | —         | —          | 7.53          | 2.5       | 90         |
|                      | 8             | > ULOQ        | —         | —          | 7.23          | 3.5       | 96         |
|                      | 16            | 124           | 10.9      | —          | 6.57          | 4.7       | 91         |
| Serum No. 3          | 2             | 3.24          | 6.0       | —          | 0.348         | 7.0       | —          |
|                      | 4             | 3.05          | 3.7       | 94         | 0.306         | 5.1       | 88         |
|                      | 8             | 3.28          | 6.3       | 108        | < LLOQ        | —         | —          |
|                      | 16            | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
| EDTA Plasma No. 1    | 2             | > ULOQ        | —         | —          | 1.93          | 1.5       | —          |
|                      | 4             | 54.7          | 10.4      | 111        | 1.88          | 3.5       | 98         |
|                      | 8             | 55.6          | 3.7       | 102        | 1.87          | 3.9       | 99         |
|                      | 16            | 52.1          | 6.5       | 94         | 1.47          | 5.1       | 79         |
| EDTA Plasma No. 2    | 2             | > ULOQ        | —         | —          | 2.94          | 7.2       | —          |
|                      | 4             | 40.8          | 13.8      | 100        | 2.67          | 3.5       | 91         |
|                      | 8             | 38.1          | 10.3      | 93         | 2.56          | 6.2       | 96         |
|                      | 16            | 37.7          | 4.4       | 99         | 1.91          | 8.9       | 74         |
| EDTA Plasma No. 3    | 2             | 6.00          | 2.8       | —          | 0.538         | 8.5       | —          |
|                      | 4             | 5.47          | 5.1       | 91         | 0.478         | 8.5       | 89         |
|                      | 8             | 5.29          | 6.5       | 97         | 0.507         | 12.4      | 106        |
|                      | 16            | < LLOQ        | —         | —          | < LLOQ        | —         | —          |
| Heparin Plasma No. 1 | 2             | > ULOQ        | —         | —          | 3.08          | 3.7       | —          |
|                      | 4             | > ULOQ        | —         | —          | 2.84          | 3.3       | 92         |
|                      | 8             | 66.8          | 6.3       | —          | 2.58          | 8.9       | 91         |
|                      | 16            | 67.2          | 7.2       | 101        | 2.10          | 8.3       | 82         |
| Heparin Plasma No. 2 | 2             | > ULOQ        | —         | —          | 3.43          | 3.1       | —          |
|                      | 4             | 56.1          | 5.4       | —          | 3.00          | 4.4       | 87         |
|                      | 8             | 55.7          | 1.0       | 99         | 2.90          | 5.9       | 97         |
|                      | 16            | 50.6          | 6.5       | 91         | 2.48          | 8.8       | 86         |
| Heparin Plasma No. 3 | 2             | > ULOQ        | —         | —          | 3.48          | 5.0       | —          |
|                      | 4             | 59.4          | 3.2       | —          | 3.12          | 3.1       | 90         |
|                      | 8             | 57.9          | 2.9       | 98         | 3.10          | 3.4       | 100        |
|                      | 16            | 51.8          | 1.6       | 89         | 2.60          | 0.4       | 84         |

ULOQ = upper limit of quantification

LLOQ = lower limit of quantification

Dash (—) = not applicable

Table 8. Dilution Linearity: sTnI

| Sample               | Fold Dilution | sTnI                  |                   |                  |
|----------------------|---------------|-----------------------|-------------------|------------------|
|                      |               | Concentration (ng/mL) | Concentration %CV | Percent Recovery |
| Serum No. 1          | 2             | 23.1                  | 3.4               | —                |
|                      | 4             | 22.4                  | 5.4               | 97               |
|                      | 8             | 22.0                  | 3.9               | 98               |
|                      | 16            | 21.5                  | 3.3               | 98               |
| Serum No. 2          | 2             | 14.9                  | 7.8               | —                |
|                      | 4             | 14.2                  | 3.7               | 95               |
|                      | 8             | 13.9                  | 6.1               | 98               |
|                      | 16            | 13.1                  | 1.1               | 95               |
| Serum No. 3          | 2             | < LLOQ                | —                 | —                |
|                      | 4             | < LLOQ                | —                 | —                |
|                      | 8             | < LLOQ                | —                 | —                |
|                      | 16            | < LLOQ                | —                 | —                |
| EDTA Plasma No. 1    | 2             | 20.6                  | 6.1               | —                |
|                      | 4             | 20.4                  | 2.8               | 99               |
|                      | 8             | 19.5                  | 6.8               | 95               |
|                      | 16            | 19.3                  | 9.0               | 99               |
| EDTA Plasma No. 2    | 2             | 13.0                  | 11.7              | —                |
|                      | 4             | 13.3                  | 6.5               | 102              |
|                      | 8             | 13.1                  | 7.5               | 98               |
|                      | 16            | < LLOQ                | —                 | —                |
| EDTA Plasma No. 3    | 2             | < LLOQ                | —                 | —                |
|                      | 4             | < LLOQ                | —                 | —                |
|                      | 8             | < LLOQ                | —                 | —                |
|                      | 16            | < LLOQ                | —                 | —                |
| Heparin Plasma No. 1 | 2             | 16.4                  | 2.7               | —                |
|                      | 4             | 16.0                  | 2.6               | 97               |
|                      | 8             | 15.3                  | 2.4               | 96               |
|                      | 16            | 15.3                  | 6.4               | 100              |
| Heparin Plasma No. 2 | 2             | 14.1                  | 8.1               | —                |
|                      | 4             | 13.8                  | 2.3               | 98               |
|                      | 8             | 13.3                  | 2.3               | 97               |
|                      | 16            | < LLOQ                | —                 | —                |
| Heparin Plasma No. 3 | 2             | 13.9                  | 3.9               | —                |
|                      | 4             | 14.0                  | 5.8               | 101              |
|                      | 8             | 13.7                  | 2.0               | 98               |
|                      | 16            | 12.7                  | 4.9               | 93               |

LLOQ = lower limit of quantification  
 Dash (—) = not applicable

# Spike Recovery

Normal serum, heparin plasma, and EDTA plasma were spiked with the Calibrators at multiple values throughout the range of the assay. Spikes were made into neat samples. Values in italics are outside of the range of quantitation.

$$\% \text{ recovery} = \frac{\text{measured concentration}}{\text{expected concentration}} \times 100$$

Table 9. Spike recovery: cTnI and cTnT

| Sample            | cTnI                |                        |                    |            | cTnT                |                        |                    |            |
|-------------------|---------------------|------------------------|--------------------|------------|---------------------|------------------------|--------------------|------------|
|                   | Spike Conc. (ng/mL) | Measured Conc. (ng/mL) | Measured Conc. %CV | % Recovery | Spike Conc. (ng/mL) | Measured Conc. (ng/mL) | Measured Conc. %CV | % Recovery |
| Serum #1          | 0                   | 1.50                   | 2.6                | —          | 0                   | 0.986                  | 5.2                | —          |
|                   | 0.260               | 1.80                   | 5.2                | 102        | 0.522               | 1.55                   | 2.4                | 103        |
|                   | 0.780               | 2.37                   | 3.8                | 104        | 1.57                | 2.53                   | 5.4                | 99         |
|                   | 2.34                | 3.90                   | 8.3                | 102        | 4.70                | 5.84                   | 3.8                | 103        |
| Serum #2          | 0                   | 0                      | —                  | —          | 0                   | 0                      | —                  | —          |
|                   | 0.260               | 0.323                  | 2.5                | 124        | 0.522               | 0.396                  | 4.3                | 76         |
|                   | 0.780               | 0.925                  | 3.9                | 119        | 1.57                | 1.27                   | 3.8                | 81         |
|                   | 2.34                | 2.67                   | 1.9                | 114        | 4.70                | 4.20                   | 2.1                | 89         |
| Serum #3          | 0                   | 0                      | —                  | —          | 0                   | 0                      | —                  | —          |
|                   | 0.260               | 0.332                  | 7.0                | 128        | 0.522               | 0.443                  | 6.0                | 85         |
|                   | 0.780               | 0.982                  | 2.3                | 126        | 1.57                | 1.37                   | 0.9                | 88         |
|                   | 2.34                | 2.80                   | 5.6                | 120        | 4.70                | 4.31                   | 4.0                | 92         |
| EDTA Plasma #1    | 0                   | 0.0388                 | 1.5                | —          | 0                   | 0                      | —                  | —          |
|                   | 0.260               | 0.374                  | 4.0                | 125        | 0.522               | 0.459                  | 3.7                | 88         |
|                   | 0.780               | 1.04                   | 2.6                | 127        | 1.57                | 1.44                   | 1.2                | 92         |
|                   | 2.34                | 2.74                   | 3.3                | 115        | 4.70                | 4.56                   | 1.9                | 97         |
| EDTA Plasma #2    | 0                   | 0.0262                 | 5.1                | —          | 0                   | 0                      | —                  | —          |
|                   | 0.260               | 0.373                  | 5.5                | 130        | 0.522               | 0.457                  | 8.8                | 87         |
|                   | 0.780               | 0.975                  | 4.1                | 121        | 1.57                | 1.35                   | 1.0                | 86         |
|                   | 2.34                | 2.74                   | 3.7                | 116        | 4.70                | 4.42                   | 3.1                | 94         |
| EDTA Plasma #3    | 0                   | 0                      | —                  | —          | 0                   | 0                      | —                  | —          |
|                   | 0.260               | 0.350                  | 1.1                | 135        | 0.522               | 0.443                  | 5.7                | 85         |
|                   | 0.780               | 1.00                   | 2.0                | 128        | 1.57                | 1.34                   | 1.1                | 86         |
|                   | 2.34                | 2.77                   | 2.4                | 118        | 4.70                | 4.32                   | 0.7                | 92         |
| Heparin Plasma #1 | 0                   | 0.269                  | 1.8                | —          | 0                   | 0.0485                 | 35.0               | —          |
|                   | 0.260               | 0.590                  | 5.9                | 111        | 0.522               | 0.577                  | 3.6                | 101        |
|                   | 0.780               | 1.20                   | 3.9                | 114        | 1.57                | 1.54                   | 3.1                | 95         |
|                   | 2.34                | 2.99                   | 3.6                | 115        | 4.70                | 4.70                   | 3.3                | 99         |
| Heparin Plasma #2 | 0                   | 0.276                  | 6.4                | —          | 0                   | 0.0576                 | 86.7               | —          |
|                   | 0.260               | 0.611                  | 3.2                | 114        | 0.522               | 0.588                  | 3.7                | 101        |
|                   | 0.780               | 1.27                   | 0.2                | 120        | 1.57                | 1.61                   | 0.9                | 99         |
|                   | 2.34                | 2.82                   | 8.4                | 108        | 4.70                | 4.40                   | 9.4                | 92         |
| Heparin Plasma #3 | 0                   | 0.325                  | 3.1                | —          | 0                   | 0.133                  | 10.4               | —          |
|                   | 0.260               | 0.612                  | 2.2                | 105        | 0.522               | 0.665                  | 2.2                | 102        |
|                   | 0.780               | 1.10                   | 1.4                | 100        | 1.57                | 1.58                   | 3.9                | 93         |
|                   | 2.34                | 2.71                   | 2.2                | 102        | 4.70                | 4.84                   | 3.7                | 100        |

Dash (—) = not applicable

Table 10. Spike recovery: FABP3 and Myl3

| Sample            | FABP3               |                        |                    |            | Myl3                |                        |                    |            |
|-------------------|---------------------|------------------------|--------------------|------------|---------------------|------------------------|--------------------|------------|
|                   | Spike Conc. (ng/mL) | Measured Conc. (ng/mL) | Measured Conc. %CV | % Recovery | Spike Conc. (ng/mL) | Measured Conc. (ng/mL) | Measured Conc. %CV | % Recovery |
| Serum #1          | 0                   | 28.4                   | 10.1               | —          | 0                   | 1.51                   | 4.2                | —          |
|                   | 1.06                | 26.9                   | 7.6                | 91         | 0.552               | 2.28                   | 3.5                | 111        |
|                   | 3.17                | 30.6                   | 8.8                | 97         | 1.66                | 3.47                   | 4.2                | 110        |
|                   | 9.52                | 32.0                   | 21.7               | 84         | 4.97                | 6.81                   | 6.1                | 105        |
| Serum #2          | 0                   | 0.643                  | 4.8                | —          | 0                   | 0.0492                 | 40.1               | —          |
|                   | 1.06                | 1.84                   | 2.2                | 108        | 0.552               | 0.692                  | 6.9                | 115        |
|                   | 3.17                | 4.51                   | 4.6                | 118        | 1.66                | 2.05                   | 5.2                | 120        |
|                   | 9.52                | 12.1                   | 10.8               | 119        | 4.97                | 6.14                   | 6.5                | 122        |
| Serum #3          | 0                   | 1.05                   | 10.2               | —          | 0                   | 0.0557                 | 8.5                | —          |
|                   | 1.06                | 2.19                   | 4.8                | 104        | 0.552               | 0.734                  | 4.4                | 121        |
|                   | 3.17                | 4.86                   | 6.1                | 115        | 1.66                | 2.06                   | 4.3                | 120        |
|                   | 9.52                | 12.6                   | 2.3                | 119        | 4.97                | 6.14                   | 4.4                | 122        |
| EDTA Plasma #1    | 0                   | 11.3                   | 5.9                | —          | 0                   | 0.395                  | 2.9                | —          |
|                   | 1.06                | 12.5                   | 11.4               | 101        | 0.552               | 1.08                   | 6.4                | 114        |
|                   | 3.17                | 14.2                   | 1.2                | 98         | 1.66                | 2.39                   | 0.6                | 116        |
|                   | 9.52                | 19.0                   | 2.5                | 91         | 4.97                | 6.29                   | 2.5                | 117        |
| EDTA Plasma #2    | 0                   | 8.66                   | 4.6                | —          | 0                   | 0.493                  | 6.2                | —          |
|                   | 1.06                | 10.1                   | 6.6                | 104        | 0.552               | 1.15                   | 8.1                | 110        |
|                   | 3.17                | 12.7                   | 4.1                | 107        | 1.66                | 2.46                   | 3.2                | 114        |
|                   | 9.52                | 17.4                   | 5.6                | 96         | 4.97                | 6.21                   | 3.6                | 114        |
| EDTA Plasma #3    | 0                   | 1.33                   | 8.3                | —          | 0                   | 0.0813                 | 19.2               | —          |
|                   | 1.06                | 2.51                   | 4.0                | 105        | 0.552               | 0.772                  | 5.6                | 122        |
|                   | 3.17                | 5.15                   | 1.1                | 114        | 1.66                | 2.16                   | 2.6                | 124        |
|                   | 9.52                | 12.1                   | 6.5                | 111        | 4.97                | 6.28                   | 3.6                | 124        |
| Heparin Plasma #1 | 0                   | 12.1                   | 7.4                | —          | 1.92                | 0.516                  | 11.4               | —          |
|                   | 1.06                | 13.8                   | 4.6                | 105        | 4.50                | 1.24                   | 11.3               | 116        |
|                   | 3.17                | 16.9                   | 12.0               | 110        | 9.22                | 2.53                   | 7.9                | 116        |
|                   | 9.52                | 22.4                   | 8.0                | 103        | 23.5                | 6.28                   | 3.7                | 115        |
| Heparin Plasma #2 | 0                   | 13.1                   | 30.7               | —          | 1.98                | 0.535                  | 5.5                | —          |
|                   | 1.06                | 12.4                   | 4.1                | 88         | 4.48                | 1.24                   | 1.5                | 114        |
|                   | 3.17                | 14.8                   | 6.9                | 91         | 9.42                | 2.58                   | 3.6                | 118        |
|                   | 9.52                | 20.6                   | 8.3                | 91         | 24.3                | 6.48                   | 2.8                | 118        |
| Heparin Plasma #3 | 0                   | 12.8                   | 0.9                | —          | 0                   | 0.671                  | 3.4                | —          |
|                   | 1.06                | 14.6                   | 8.4                | 106        | 0.552               | 1.43                   | 3.2                | 117        |
|                   | 3.17                | 16.4                   | 1.2                | 103        | 1.66                | 2.66                   | 4.3                | 114        |
|                   | 9.52                | 23.8                   | 9.2                | 107        | 4.97                | 6.69                   | 0.5                | 119        |

Dash (—) = not applicable

Table 11. Spike recovery: sTnI

| Sample            | sTnI                |                        |                    |            |
|-------------------|---------------------|------------------------|--------------------|------------|
|                   | Spike Conc. (ng/mL) | Measured Conc. (ng/mL) | Measured Conc. %CV | % Recovery |
| Serum #1          | 0                   | 3.50                   | 1.6                | —          |
|                   | 2.32                | 6.14                   | 2.8                | 106        |
|                   | 6.95                | 11.0                   | 3.2                | 105        |
|                   | 20.9                | 26.1                   | 3.2                | 107        |
| Serum #2          | 0                   | 0.0288                 | 64.0               | —          |
|                   | 2.32                | 2.48                   | 5.3                | 106        |
|                   | 6.95                | 7.20                   | 5.9                | 103        |
|                   | 20.9                | 20.6                   | 4.3                | 98         |
| Serum #3          | 0                   | 0.0136                 | 71.4               | —          |
|                   | 2.32                | 2.44                   | 1.5                | 105        |
|                   | 6.95                | 7.25                   | 2.4                | 104        |
|                   | 20.9                | 21.7                   | 3.6                | 104        |
| EDTA Plasma #1    | 0                   | 4.40                   | 2.7                | —          |
|                   | 2.32                | 6.92                   | 6.1                | 103        |
|                   | 6.95                | 12.0                   | 5.1                | 106        |
|                   | 20.9                | 25.4                   | 2.5                | 101        |
| EDTA Plasma #2    | 0                   | 3.19                   | 5.6                | —          |
|                   | 2.32                | 5.79                   | 4.1                | 105        |
|                   | 6.95                | 11.1                   | 1.7                | 109        |
|                   | 20.9                | 24.3                   | 3.3                | 101        |
| EDTA Plasma #3    | 0                   | 0.0838                 | 63.1               | —          |
|                   | 2.32                | 2.59                   | 3.0                | 108        |
|                   | 6.95                | 7.26                   | 1.5                | 103        |
|                   | 20.9                | 20.2                   | 3.1                | 96         |
| Heparin Plasma #1 | 0                   | 3.23                   | 8.2                | —          |
|                   | 2.32                | 5.70                   | 7.6                | 103        |
|                   | 6.95                | 10.8                   | 5.4                | 106        |
|                   | 20.9                | 24.3                   | 6.2                | 101        |
| Heparin Plasma #2 | 0                   | 3.00                   | 3.3                | —          |
|                   | 2.32                | 5.72                   | 3.0                | 108        |
|                   | 6.95                | 11.3                   | 6.2                | 114        |
|                   | 20.9                | 26.0                   | 5.4                | 109        |
| Heparin Plasma #3 | 0                   | 3.28                   | 0.4                | —          |
|                   | 2.32                | 5.79                   | 3.5                | 103        |
|                   | 6.95                | 10.2                   | 3.3                | 100        |
|                   | 20.9                | 24.3                   | 2.4                | 101        |

Dash (—) = not applicable

# Specificity

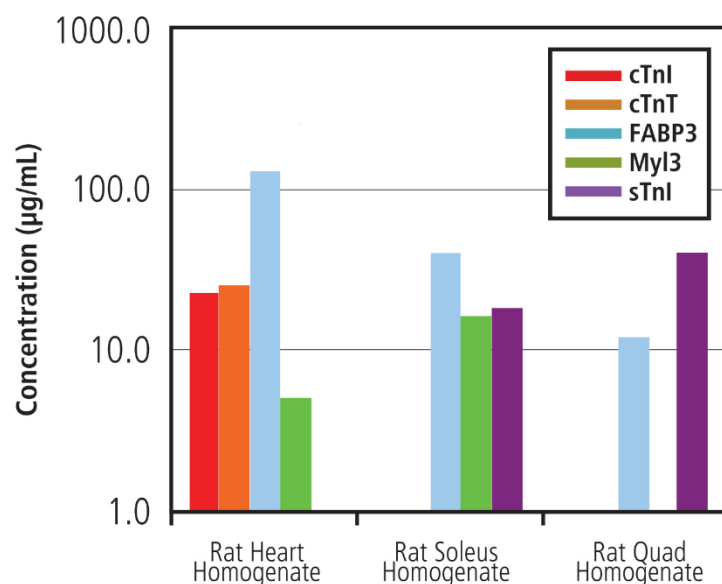
The assays on the Muscle Injury Panel 1 (rat) are all highly specific for their respective analytes. Specificity was demonstrated by testing rat muscle homogenates, and by running the assay with single Calibrators and single detection antibodies. Results are detailed in the sections below.

## Specificity of the assays for muscle homogenates

Tissue homogenates from heart, fast-twitch, and slow-twitch muscle were tested at 100-fold, 1000-fold, and 10,000-fold sample dilution. The assays for cardiac troponins were positive for cardiac homogenates and negative for other muscle homogenates, demonstrating specificity for cardiac tissue. The assay for skeletal Troponin I was specific for fast and slow-twitch skeletal muscle. The assay measured FABP3 in cardiac muscle and skeletal muscle. The slow-twitch muscle was positive for Myl3, and approximately 200-fold less Myl3 was measured in fast-twitch.

*Table 12. Muscle homogenate samples tested for assay specificity*

| Sample Group                        | cTnI            |               | cTnT            |               | FABP3           |               | Myl3            |               | sTnI            |               |
|-------------------------------------|-----------------|---------------|-----------------|---------------|-----------------|---------------|-----------------|---------------|-----------------|---------------|
|                                     | Sample Dilution | Conc. (µg/mL) | Sample Dilution | Conc. (µg/mL) | Sample Dilution | Conc. (µg/mL) | Sample Dilution | Conc. (µg/mL) | Sample Dilution | Conc. (µg/mL) |
| Rat Heart Homogenate                | 1,000           | 22.6          | 1,000           | 25.1          | 10,000          | 125.2         | 1,000           | 5.00          | 100             | < LLOD        |
| Rat Soleus Homogenate (slow-twitch) | 100             | < LLOD        | 100             | < LLOD        | 10,000          | 38.8          | 1,000           | 16.4          | 1,000           | 18.1          |
| Rat Quad Homogenate (fast-twitch)   | 100             | < LLOD        | 100             | < LLOD        | 1000            | 12.2          | 100             | 0.0837        | 1,000           | 40.9          |



*Figure 2. Muscle homogenate samples tested for assay specificity.*

## Specificity of the Assays for Individual Calibrators:

The Muscle Injury Panel 1 assay was run with each single Calibrator at a high level to assess specificity. The table below shows the % cross-reactivity of each assay for each Calibrator. The % cross-reactivity is calculated as the ratio of the nonspecific signal to the specific signal.

*Table 13. Assay specificity for individual calibrators*

| Blended Detection Antibodies and Single Calibrator % Cross-Reactivity |       |       |       |       |       |
|-----------------------------------------------------------------------|-------|-------|-------|-------|-------|
| Spot                                                                  | cTnl  | cTnT  | FABP3 | Myl3  | sTnl  |
| cTnl                                                                  | 100   | 0.55  | < 0.1 | < 0.1 | < 0.1 |
| cTnT                                                                  | 0.19  | 100   | < 0.1 | < 0.1 | 1.93  |
| FABP3                                                                 | < 0.1 | < 0.1 | 100   | < 0.1 | < 0.1 |
| Myl3                                                                  | < 0.1 | < 0.1 | < 0.1 | 100   | < 0.1 |
| sTnl                                                                  | < 0.1 | < 0.1 | < 0.1 | < 0.1 | 100   |

## Specificity of the Detection Antibodies:

To assess specificity of the detection antibodies, the Muscle Injury Panel 1 panel was run with blended Calibrator diluted to STD-02, and single detection antibodies. The table below shows the % cross-reactivity for each individual detection antibody. The data shows that the presence of cTnl Calibrator on the cTnl spot is reported out by the cTnl, cTnT, and sTnl detection antibodies.

*Table 14. Detection antibody specificity*

| Blended Calibrator and Single Detection Antibody % Cross-Reactivity |       |       |       |       |      |
|---------------------------------------------------------------------|-------|-------|-------|-------|------|
| Spot                                                                | cTnl  | cTnT  | FABP3 | Myl3  | sTnl |
| cTnl                                                                | 100   | < 0.1 | < 0.1 | < 0.1 | 50.9 |
| cTnT                                                                | 0.44  | 100   | < 0.1 | 0.17  | 7.89 |
| FABP3                                                               | < 0.1 | < 0.1 | 100   | < 0.1 | 0.11 |
| Myl3                                                                | < 0.1 | < 0.1 | < 0.1 | 100   | 0.11 |
| sTnl                                                                | 0.24  | < 0.1 | < 0.1 | < 0.1 | 100  |

While these detection antibodies are not highly specific, the assays themselves are highly specific due to the specificity of the capture antibodies.



## Tested Samples

Serum, EDTA plasma, and heparin plasma samples collected from normal Sprague-Dawley rats were tested at 2-fold dilution on the Muscle Injury Panel 1 (rat). Shown below are the median and range of concentrations for each sample set. Skeletal Troponin I was below the quantitative range for all samples.

*Table 15. Samples tested in the Muscle Injury Panel (rat) Kits*

| Sample         | Statistic      | cTnI       | cTnT         | FABP3       | MyI3       | sTnI   |
|----------------|----------------|------------|--------------|-------------|------------|--------|
| Serum          | Median (ng/mL) | 1.59       | 0.83         | 19.84       | 0.61       | < 1.56 |
|                | Range (ng/mL)  | 0.436–3.53 | < 0.976–2.04 | 5.28–> 30.0 | 0.25–1.05  | < 1.56 |
|                | N              | 10         | 10           | 10          | 10         | 10     |
| EDTA Plasma    | Median (ng/mL) | 2.31       | 1.28         | 35.63       | 0.81       | < 1.56 |
|                | Range (ng/mL)  | 1.45–4.05  | < 0.976–2.80 | 19.6–> 30.0 | 0.432–1.17 | < 1.56 |
|                | N              | 10         | 10           | 10          | 10         | 10     |
| Heparin Plasma | Median (ng/mL) | 1.81       | 1.06         | 28.35       | 0.55       | < 1.56 |
|                | Range (ng/mL)  | 0.372–4.10 | < 0.976–2.79 | 4.12–> 30.0 | 0.162–1.16 | < 1.56 |
|                | N              | 10         | 10           | 10          | 10         | 10     |

# Assay Components

## Calibrators

Rat cardiac troponin I (cTnI), rat cardiac troponin T (cTnT), and rat fatty acid binding protein 3 (FABP3) were purified from rat heart tissue. Rat skeletal troponin I was purified from rat skeletal muscle. Full-length recombinant rat myosin light chain 3 (MyI3) with an N-terminal 10XHis-tag was expressed in *E. coli*. These analytes were calibrated against internal controls, diluted, and pooled to make the Muscle Injury Panel 1 (rat) Calibrator Blend.

## Antibodies

Table 16. Antibody source species

|         | Source Species       |                        |
|---------|----------------------|------------------------|
| Analyte | MSD Capture Antibody | MSD Detection Antibody |
| cTnI    | Mouse Monoclonal     | Mouse Monoclonal       |
| cTnT    | Mouse Monoclonal     | Mouse Monoclonal       |
| FABP3   | Chicken Polyclonal   | Mouse Monoclonal       |
| MyI3    | Mouse Monoclonal     | Mouse Monoclonal       |
| sTnI    | Mouse Monoclonal     | Mouse Monoclonal       |

## References

1. Lee JW, et al. Fit-for-purpose method development and validation for successful biomarker measurement. *Pharm Res.* 2006 Feb;23(2):312-28. doi: 10.1007/s11095-005-9045-3.
2. Babuin L, Jaffe AS. Troponin: the biomarker of choice for the detection of cardiac injury. *CMAJ.* 2005 Nov 8;173(10):1191-202. doi: 10.1503/cmaj/051291.

# Summary Protocol

## Muscle Injury Panel Kits

MSD provides this summary protocol for your convenience.

Please read the entire detailed protocol before performing the Muscle Injury Panel assays.

## Sample and Reagent Preparation

- Bring all reagents to room temperature and thaw the Calibrator on ice.
- Prepare Diluent 7 + Additives by diluting the provided DTT (100X) and EDTA (16.7X) stock solutions to 1X concentrations in Diluent 7.
- Prepare 8 calibration solutions using the supplied Calibrator:
  - Dilute the stock Calibrator 20-fold in Diluent 7 + Additives.
  - Perform a series of 3-fold dilution steps and prepare a zero Calibrator.
- Dilute serum and plasma samples 4-fold in Diluent 7 + Additives before adding to the plate.
- Incubate the diluted Calibrators and diluted samples at room temperature for 30 minutes before adding them to the plate.
- Prepare combined detection antibody solution by diluting each stock detection antibody 50-fold in Diluent 30.
- Prepare 20 mL of 1X Read Buffer T by diluting 4X Read Buffer T 4-fold with deionized water.

### STEP 1: Add Diluent 7 + Additives

- Add 25  $\mu$ L/well of Diluent 7 + Additives.
- Incubate at room temperature with vigorous shaking (500–1,000 rpm) for 30 minutes.

### STEP 2: Add Sample or Calibrator

- Add 25  $\mu$ L/well of sample or Calibrator.
- Incubate at room temperature with vigorous shaking (500–1,000 rpm) for 2 hours.

### STEP 3: Wash and Add Detection Antibody Solution

- Wash plate 3 times with at least 150  $\mu$ L/well of 1X MSD Wash Buffer or PBS-T.
- Add 25  $\mu$ L/well of 1X detection antibody solution.
- Incubate at room temperature with vigorous shaking (500–1,000 rpm) for 2 hours.

### STEP 4: Wash and Read Plate

- Wash plate 3 times with at least 150  $\mu$ L/well of 1X MSD Wash Buffer or PBS-T.
- Add 150  $\mu$ L/well of 1X Read Buffer T.
- Analyze plate on an MSD instrument.

# Plate Diagrams

