



**CERTIFICATE OF ANALYSIS**

COA-04515

**Product Description:** Muscle Injury Panel 1 (rat) Kit  
**Kit Catalog Numbers:** K15181C-1, K15181C-2, K15181C-4  
**Kit Lot Number:** K0040261  
**Expiration Date:** 31 JAN 2015

**Acceptance Data:**

Metric		Assay					
		cTnI		cTnT		FABP3	
		Specification	Pass/ Fail	Specification	Pass/ Fail	Specification	Pass/ Fail
% Recovery Concentration	Calibrators	80% -120%	Pass	80% -120%	Pass	80% -120%	Pass
	Controls	75% - 125%	Pass	75% - 125%	Pass	75% - 125%	Pass
	LLOQ	80% -120%	Pass	80% -120%	Pass	75% - 125%	Pass
	ULOQ	80% -120%	Pass	80% -120%	Pass	75% - 125%	Pass
% Recovery CV	Calibrators	<20%	Pass	<20%	Pass	<20%	Pass
	Controls	<20%	Pass	<20%	Pass	<20%	Pass
	LLOQ	<20%	Pass	<20%	Pass	<25%	Pass
	ULOQ	<20%	Pass	<20%	Pass	<25%	Pass

Metric		Assay			
		Myl3		sTnI	
		Specification	Pass/ Fail	Specification	Pass/ Fail
% Recovery Concentration	Calibrators	80% -120%	Pass	80% -120%	Pass
	Controls	75% - 125%	Pass	75% - 125%	Pass
	LLOQ	75% - 125%	Pass	80% -120%	Pass
	ULOQ	80% -120%	Pass	80% -120%	Pass
% Recovery CV	Calibrators	<20%	Pass	<20%	Pass
	Controls	<20%	Pass	<20%	Pass
	LLOQ	<25%	Pass	<20%	Pass
	ULOQ	<20%	Pass	<20%	Pass

**Concentration of Materials Tested:**

Units = ng/mL	Assay				
	cTnI	cTnT	FABP3	Myl3	sTnI
20X Calibrator	440	911	1910	822	4000
Control 1	23.7	84.7	84.6	15.0	92.3
Control 2	2.57	11.7	22.0	1.69	9.40
ULOQ	20.0	41.4	15.9	45.0	182
LLOQ	0.0900	0.485	0.390	0.0530	0.781



**CERTIFICATE OF ANALYSIS**

COA-04515

**Standard Curve Data from Kit QC:**

cTnI		
Conc., (ng/mL)	Average Signal	%CV
22.0	380 405	5.5
7.33	101 978	5.9
2.44	30 508	5.1
0.815	9406	5.2
0.272	3173	6.7
0.0905	1136	6.0
0.0302	445	4.4
0	102	9.6

cTnT		
Conc., (ng/mL)	Average Signal	%CV
45.6	331 242	6.4
15.2	55 725	6.4
5.06	9483	4.1
1.69	1931	5.5
0.562	649	5.8
0.187	299	6.6
0.0625	199	8.0
0	161	7.8

FABP3		
Conc., (ng/mL)	Average Signal	%CV
95.5	49 035	4.4
31.8	37 355	5.1
10.6	22 535	4.3
3.54	8528	5.0
1.18	2742	4.1
0.393	1359	4.5
0.131	1011	3.2
0	865	4.4

MyI3		
Conc., (ng/mL)	Average Signal	%CV
41.1	153 619	8.2
13.7	45 836	7.8
4.57	14 791	6.1
1.52	4510	7.4
0.507	1689	8.1
0.169	658	6.8
0.0564	331	8.4
0	174	15.8

sTnI		
Conc., (ng/mL)	Average Signal	%CV
200	280 865	3.4
66.7	78 539	3.4
22.2	23 212	3.0
7.41	7143	3.2
2.47	2732	3.3
0.823	1068	4.3
0.274	497	5.1
0	220	6.6


Muscle Injury Panel 1 (rat) Components			
Description	Lot Number	Storage Temperature	Expiration Date
Muscle Injury Panel Plate	Z0054900	2-8°C	31 JUL 2016
Anti-rat cTnI Antibody	D0040396	2-8°C	31 MAR 2016
Anti-rat cTnT Antibody	D0040397	2-8°C	31 MAR 2016
Anti-rat FABP3 Antibody	D0040398	2-8°C	31 MAR 2016
Anti-rat MyI3 Antibody	D0040399	2-8°C	29 FEB 2016
Anti-rat sTnI Antibody	D0040400	2-8°C	30 APR 2016
Muscle Injury Panel 1 (rat) Calibrator Blend	A0040103	≤ -70°C	31 JAN 2015
Diluent 7	M0070042	≤ -10°C	31 MAY 2016
Diluent 30	M0300013	≤ -10°C	31 MAR 2016
EDTA pH 8.0	R0010433	Room Temperature	31 MAR 2016
DTT	R0010432	≤ -10°C	31 MAR 2016
MSD Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A



**CERTIFICATE OF ANALYSIS**

COA-04515

**Statement:** The above kit lot was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Dana Short	Quality Control		27 JUN 2014