

# CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** Muscle Injury Panel 2 (rat) Kit

**Kit Catalog Numbers:** K15180C-Series

**Kit Lot Number:** K0040426

**Expiration Date:** 31 Mar 2019

## Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Muscle Injury Panel 2 (rat) Control-1		Muscle Injury Panel 2 (rat) Control-2		Muscle Injury Panel 2 (rat) Control-3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
2	TIMP-1	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	N/A	N/A
3	CK	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

## Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
2	Rat TIMP-1	< 1.0%
3	Rat CK	< 1.0%

## Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
2	TIMP-1	20X Calibrator	782	ng/mL
		Muscle Injury Panel 2 (rat) Control-1	203	ng/mL
		Muscle Injury Panel 2 (rat) Control-2	8.37	ng/mL
3	CK	20X Calibrator	4600	ng/mL
		Muscle Injury Panel 2 (rat) Control-1	915	ng/mL
		Muscle Injury Panel 2 (rat) Control-2	154	ng/mL
		Muscle Injury Panel 2 (rat) Control-3	53.3	ng/mL

Controls tested may not be the same lot as shipped.

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

## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Standard Curve Data:


TIMP-1 - Spot 2			CK - Spot 3		
Conc. ng/mL	ECL Signal	%CV	Conc. ng/mL	ECL Signal	%CV
39.1	1,370,000	1.2	230	268,000	3.9
9.78	783,000	6.0	57.5	35,500	3.1
2.44	166,000	5.3	14.4	5,610	3.4
0.611	22,900	6.1	3.59	1,330	3.8
0.153	3,380	4.7	0.898	495	3.3
0.0382	816	4.7	0.225	268	4.0
0.00955	352	3.6	0.0562	224	5.1
0.00	208	8.5	0.00	202	9.9

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Muscle Injury Panel 2 (rat) Plate	Z0056233	2-8°C	31 Oct 2020
Anti-rat TIMP-1 Antibody	D0040765	2-8°C	30 Nov 2021
Anti-rat CK Antibody	D0040764	2-8°C	30 Sep 2021
Muscle Injury Panel 2 (rat) Calibrator Blend	A0040211	≤ -70°C	31 Dec 2020
Blocker A Kit	Not Kit Specific	Not Kit Specific	N/A
Diluent 11	M0110030	≤ -10°C	31 Mar 2019
Diluent 5	M0050063	≤ -10°C	30 Nov 2020
Blocker D-R	R0060022	≤ -10°C	31 Oct 2019
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
Review/Approval	Dawn Stollar	Quality Assurance		27 JUL 2018