

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Muscle Injury Panel 2 (rat) Kit
Kit Catalog Numbers: K15180C-Series
Kit Lot Number: K0040486
Expiration Date: 30 Sep 2021

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator	
		Recovery	CV
2	TIMP-1	80% - 120%	< 20%
3	CK	80% - 120%	< 20%

Spot	Name	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
2	TIMP-1	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
3	CK	75% - 125%	< 20%	75% - 125%	< 20%	N/A	N/A

Acceptance Data/ Non Specific Binding Specifications:

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

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
2	TIMP-1	20X Calibrator	839	ng/mL
		Muscle Injury Panel 2 (rat) Control 1	2.62	ng/mL
		Muscle Injury Panel 2 (rat) Control 2	0.655	ng/mL
		Muscle Injury Panel 2 (rat) Control 3	0.164	ng/mL
3	CK	20X Calibrator	4930	ng/mL
		Muscle Injury Panel 2 (rat) Control 1	15.4	ng/mL
		Muscle Injury Panel 2 (rat) Control 2	3.85	ng/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

CERTIFICATE OF ANALYSIS

Standard Curve Data:


TIMP-1 - Spot 2			CK - Spot 3		
Conc. ng/mL	ECL Signal	%CV	Conc. ng/mL	ECL Signal	%CV
42.0	1,370,000	2.3	247	94,000	3.7
10.5	575,000	3.7	61.6	15,500	1.7
2.62	97,400	4.9	15.4	3,340	1.7
0.655	13,800	4.8	3.85	814	1.8
0.164	2,380	5.4	0.963	311	3.2
0.0410	613	4.3	0.241	161	6.4
0.0102	239	6.0	0.0602	134	11.0
0.00	123	6.3	0.00	114	4.8

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Muscle Injury Panel 2 (rat) Plate	Z0056623	2-8°C	30 Apr 2022
Anti-rat TIMP-1 Antibody	D0040894	2-8°C	31 Jul 2023
Anti-rat CK Antibody	D0040893	2-8°C	31 Oct 2023
Muscle Injury Panel 2 (rat) Calibrator Blend	A0040233	≤ -70°C	30 Jun 2022
Blocker A	Not Kit Specific	Not Kit Specific	N/A
Diluent 11	M0110044	≤ -10°C	30 Jun 2022
Diluent 5	M0050069	≤ -10°C	31 May 2022
Blocker D-R	R0010656	≤ -10°C	30 Sep 2021
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	Karen Hamilla	Quality		25 JUN 2020