

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Muscle Injury Panel 2 (rat) Kit
Kit Catalog Numbers: K15180C-Series
Kit Lot Number: K0040467
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

* See Additional Comments.

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	2.44	ng/mL
		Muscle Injury Panel 2 (rat) Control 2	0.611	ng/mL
		Muscle Injury Panel 2 (rat) Control 3	0.153	ng/mL
3	CK	20X Calibrator	4930	ng/mL
		Muscle Injury Panel 2 (rat) Control 1	14.4	ng/mL
		Muscle Injury Panel 2 (rat) Control 2	3.59	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

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
42.0	1,660,000	1.4	247	200,000	6.9
10.5	1,220,000	3.2	61.6	30,200	4.2
2.62	256,000	6.1	15.4	5,610	2.5
0.655	30,900	4.7	3.85	1,270	5.0
0.164	4,090	4.9	0.963	432	3.2
0.0410	857	4.0	0.241	178	3.7
0.0102	303	7.7	0.0602	120	4.7
0.00	147	18.0	0.00	97	5.1

Components:


Description	Lot Number	Storage Temperature	Expiration Date
Muscle Injury Panel 2 (rat) Plate	Z0056499	2-8°C	30 Sep 2021
Anti-rat TIMP-1 Antibody	D0040839	2-8°C	30 Sep 2022
Anti-rat CK Antibody	D0040838	2-8°C	31 Mar 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	M0050066	≤ -10°C	31 Oct 2021
Blocker D-R	R0010656	≤ -10°C	30 Sep 2021
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Comments:

Assay	Spot	Specification	Comment
TIMP-1	2	Calibrator Conc.	TIMP-1 Cal-2 recovery 123.3% (spec 80-120%)

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		17 APR 2020