

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Muscle Injury Panel 2 (rat) Kit
Kit Lot Number: K0040516
Expiration Date: 31 Jan 2023
Kit Catalog Numbers: K15180C-Series

Precision and Accuracy Specifications:

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

Non-Specific Binding Specifications:

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

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
2	TIMP-1	934	ng/mL
3	CK	3880	ng/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Muscle Injury Panel 2 (rat) Plate	Z0056931	2-8°C	31 Aug 2023
Anti-rat TIMP-1 Antibody	D0040951	2-8°C	31 Dec 2024
Anti-rat CK Antibody	D0040950	2-8°C	31 Jan 2025
Muscular Injury Panel 2 (rat) Calibrator Blend	A0040259	≤ -70°C	30 Sep 2023
Diluent 11	M0110050	≤ -10°C	30 Sep 2023
Diluent 5	M0050070	≤ -10°C	31 Jan 2023

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		23 SEP 2021