



Meso Scale Diagnostics, LLC.

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Rat Skeletal Troponin I Kit
Kit Lot Number: K0040537
Expiration Date: 30 Sep 2023
Kit Catalog Numbers: K153IMC-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Muscle Injury Panel 1 Control 1		Muscle Injury Panel 1 Control 2		Muscle Injury Panel 1 Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
6	sTnI	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
6	sTnI	781	80% - 120%	< 20%	160000	80% - 120%	< 25%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
6	sTnI	3960000	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Muscle Injury Panel 1 (rat) Plate	Z0056967	2-8°C	30 Nov 2023
Anti-rat s TnI Antibody	D0040960	2-8°C	30 Nov 2025
Rat Skeletal Troponin I Calibrator	A0040267	≤ -70°C	30 Jun 2024
Diluent 30	M0300035	≤ -10°C	30 Jun 2024
Diluent 7	M0070067	≤ -10°C	30 Sep 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		18 AUG 2022