



Extension Memo

Page 1 of 1

Product Description: Muscle Injury Panel 2 (rat) Kit
Catalog Number(s): K15180C-Series
Lot Number: K0040568
New Expiration Date: 31 JAN 2026

Based on additional testing, the original date of expiration for this kit has been extended by 12 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 5	M0050071	31 JAN 2025	31 JAN 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		31 JAN 2025



CERTIFICATE OF ANALYSIS

Product Description: Muscle Injury Panel 2 (rat) Kit
Lot Number: K0040568
Expiration Date: 31 JAN 2025
Catalog Number: K15180C-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-rat CK Antibody	D0041048	2-8°C	30 NOV 2027
Anti-rat TIMP-1 Antibody	D0041049	2-8°C	31 OCT 2027
Muscle Injury Panel 2 (rat) Plate	Z0057463	2-8°C	31 OCT 2026
Diluent 11	M0110065	≤ -10°C	31 JUL 2026
Diluent 5	M0050071	≤ -10°C	31 JAN 2025
Muscle Injury Panel 2 (rat) Calibrator Blend	A0040287	≤ -70°C	31 OCT 2026

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

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator		Muscle Injury Panel 2 Control 1		Muscle Injury Panel 2 Control 2		Muscle Injury Panel 2 Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
2	TIMP-1	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
3	CK	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Non-Specific Binding Specifications:

Spot	Name	TIMP-1	CK
2	TIMP-1	NA	< 1.0%
3	CK	< 1.0%	NA

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
2	TIMP-1	700	ng/mL
3	CK	4060	ng/mL

Dilution to top of Curve: 20X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 JUL 2024