



CERTIFICATE OF ANALYSIS

Product Description: Muscle Injury Panel 2 (rat) Kit
Lot Number: K0040587
Expiration Date: 28 FEB 2027
Catalog Number: K15180C-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-rat CK Antibody	D0041083	2-8°C	31 DEC 2028
Anti-rat TIMP-1 Antibody	D0041084	2-8°C	28 FEB 2029
Muscle Injury Panel 2 (rat) Plate	Z0057582	2-8°C	31 OCT 2027
Diluent 11	M0110072	≤ -10°C	30 NOV 2027
Diluent 5	M0050074	≤ -10°C	28 FEB 2027
Muscle Injury Panel 2 (rat) Calibrator Blend	A0040296	≤ -70°C	31 DEC 2027

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	880	ng/mL
3	CK	4800	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		15 OCT 2025