



## CERTIFICATE OF ANALYSIS

**Product Description:** Muscle Injury Panel 2 (rat) Kit  
**Lot Number:** K0040586  
**Expiration Date:** 31 MAR 2027  
**Catalog Number:** K15180C-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-rat CK Antibody	D0041071	2-8°C	31 MAR 2028
Anti-rat TIMP-1 Antibody	D0041072	2-8°C	31 MAR 2028
Muscle Injury Panel 2 (rat) Plate	Z0057509	2-8°C	31 MAR 2027
Diluent 11	M0110071	≤ -10°C	31 JUL 2027
Diluent 5	M0050075	≤ -10°C	30 JUN 2027
Muscle Injury Panel 2 (rat) Calibrator Blend	A0040291	≤ -70°C	31 MAR 2027

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

**Precision, Accuracy, and Sensitivity 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	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	772	ng/mL
3	CK	5500	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		17 APR 2025