

**CERTIFICATE OF ANALYSIS****Product Description:** Muscle Injury Panel 3 (mouse) Kit**Lot Number:** K0040590**Expiration Date:** 31 JAN 2028**Catalog Number:** K15186C-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-ms FABP3 Antibody	D0041093	2-8°C	28 FEB 2029
Anti-ms Myl3 Antibody	D0041094	2-8°C	28 FEB 2029
Anti-ms Troponin I (cardiac) Antibody	D0041092	2-8°C	31 JAN 2029
Anti-ms Troponin I (fast-twitch) Antibody	D0041095	2-8°C	31 JAN 2029
Muscle Injury Panel 3 (Mouse) 96-Well 7-Spot SECTOR Plate	Z0057601	2-8°C	31 JAN 2028
Diluent 30	M0300041	≤ -10°C	31 MAR 2028
Diluent 33	M0330021	≤ -10°C	30 APR 2028
Muscle Injury Panel 3 Calibrator Blend	A0040290	≤ -70°C	31 JAN 2028

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 3 Control 1		Muscle Injury Panel 3 Control 2		Mouse Injury Panel 3 Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Troponin I (cardiac)	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%*	< 20%	N/A*	N/A*
3	FABP3	80% - 120%	< 20%	75% - 125%*	< 20%	N/A*	N/A*	N/A*	N/A*
5	Myl3	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
6	Troponin I (fast-twitch)	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

*See Additional Comments

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Troponin I (cardiac)	0.070	75% - 125%	< 25%	13.2	80% - 120%*	< 20%
3	FABP3	1.00	75% - 125%	< 25%	26.8	80% - 120%	< 20%*
5	Myl3	0.300	75% - 125%	< 25%	41.4	80% - 120%	< 20%
6	Troponin I (fast-twitch)	0.505	75% - 125%	< 25%	155	80% - 120%	< 20%

*See Additional Comments



CERTIFICATE OF ANALYSIS

Non-Specific Binding Specifications:

Spot	Name	Troponin I (cardiac)	FABP3	Myl3	Troponin I (fast-twitch)
1	Troponin I (cardiac)	NA	< 1.0%	< 1.0%	< 1.0%
3	FABP3	< 1.0%	NA	< 1.0%	< 1.0%
5	Myl3	< 1.0%	< 1.0%	NA	< 1.0%
6	Troponin I (fast-twitch)	< 1.0%	< 1.0%	< 1.0%	NA

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	Troponin I (cardiac)	330	ng/mL
3	FABP3	1412	ng/mL
5	Myl3	1342	ng/mL
6	Troponin I (fast-twitch)	3920	ng/mL

Dilution to top of Curve: 20X

Additional Comments:

Spot	Name	Specification	Comment
1	Troponin I (cardiac)	Control Conc.	The specification for Control 2 % recovery is 75-125%; the actual measured value was 145.2%. Applying the LLOQ at 0.07 ng/mL caused Control 3 to fall outside of the quantifiable range.
3	FABP3	Control Conc.	The specification for Control 1 % recovery is 75-125%; the actual measured value was 134.7%. Applying the LLOQ at 1.00 ng/mL caused Control 2 and 3 to fall outside of the quantifiable range.
1	Troponin I (cardiac)	ULOQ Conc.	Alternate ULOQ @ 13.2ng/mL selected; targeted ULOQ at 20.0 ng/mL not in the dynamic range.
3	FABP3	ULOQ CV	ULOQ inter-plate CV specification (< 20%) at 26.8 ng/mL; Actual measured value was 34.2%.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		05 JAN 2026