



CERTIFICATE OF ANALYSIS

Product Description: Muscle Injury Panel 3 (mouse) Calibrator Blend
Catalog Number: C0186-2
Lot Number: A0040282
Expiration Date: 31 DEC 2025
Storage Temperature: ≤ -70°C

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
cTnI	336	ng/mL
FABP3	678	ng/mL
Myl3	946	ng/mL
Troponin I (fast-twitch)	3860	ng/mL

Dilution to Top of Curve: 20X

This product was manufactured and tested according to MSD documents. This lot meets MSD’s specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		23 AUG 2023