



CERTIFICATE OF ANALYSIS

Product Description: Muscle Injury Panel 1 (rat) Calibrator Blend
Catalog Number: C0181-2
Lot Number: A0040277
Expiration Date: 30 JUN 2024
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
cTnl	654	ng/mL
cTnT	942	ng/mL
FABP3	599	ng/mL
Myl3	808	ng/mL
sTnl	3600	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		10 JAN 2023