

**CERTIFICATE OF ANALYSIS**


Product Description: Muscle Injury Panel 2 (rat) Calibrator
Catalog Number: C0180-2
Lot Number: A0040296
Expiration Date: 31 DEC 2027
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
TIMP-1	880	ng/mL
CK	4800	ng/mL

Dilution to Top of Curve: 20 X

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		08 AUG 2025