



### CERTIFICATE OF ANALYSIS

**Product Description:** Muscle Injury Panel 2 (rat) Calibrator  
**Catalog Number:** C0180-2  
**Lot Number:** A0040291  
**Expiration Date:** 31 MAR 2027  
**Storage Temperature:** ≤ -70°C

**Concentration Assignment:**

Analyte	Stock Concentration	Unit of Measure
TIMP-1	772	ng/mL
CK	5500	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
<b>Review/Approval</b>	Karen Hamilla	Quality		21 NOV 2024