



**CERTIFICATE OF ANALYSIS**

**Product Description:** Muscle Injury Panel 2 (rat) Calibrator  
**Catalog Number:** C0180-2  
**Lot Number:** A0040287  
**Expiration Date:** 31 OCT 2026  
**Storage Temperature:** ≤ -70°C

**Concentration Assignment:**

Analyte	Stock Concentration	Unit of Measure
TIMP-1	700	ng/mL
CK	4060	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	<i>Karen Hamilla</i>	05 JUN 2024