



### CERTIFICATE OF ANALYSIS

**Product Description:** Rat NT-proANP Calibrator  
**Catalog Number:** C03MB-2  
**Lot Number:** A0040281  
**Expiration Date:** 31 OCT 2025  
**Storage Temperature:** ≤ -70°C

**Concentration Assignment:**

Analyte	Stock Concentration	Unit of Measure
NT-proANP	2240	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
<b>Review/Approval</b>	Karen Hamilla	Quality	<i>Karen Hamilla</i>	12 JUN 2023