

**CERTIFICATE OF ANALYSIS**


Product Description: Rat NT-proANP Calibrator
Catalog Number: C03MB-2
Lot Number: A0040294
Expiration Date: 30 SEP 2027
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
NT-proANP	2240	ng/mL

Dilution to Top of Curv 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		04 JUN 2025