



CERTIFICATE OF ANALYSIS

Product Description: Rat NT-proBNP Calibrator
Catalog Number: C03JK-2
Lot Number: A0040288
Expiration Date: 31 JAN 2027
Storage Temperature: ≤ -70°C

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
NT-proBNP	95400	pg/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		11 SEP 2024