



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Rat NT-proBNP Calibrator

**Catalog Number:** C03JK-2

**Lot Number:** A0040270

**Expiration Date:** 31 DEC 2024


**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

### Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
NT-proBNP	96200	pg/ml

**Dilution to Top of Curve:** 20X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		09 AUG 2022