



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Rat NT-proBNP Calibrator

Catalog Number: C03JK-2

Lot Number: A0040207

Expiration Date: 30 SEP 2020

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

Concentration of Materials Tested:

	Assay
	Rat NT-proBNP
Calibrator (pg/mL)	90 200

Concentration: 20X

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	04 SEP 2018