



### CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C03MB-2

**Lot Number:** A0040191

**Expiration Date:** 31 AUG 2019

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

**Concentration of Materials Tested:**

	Assay
	<b>Rat NT-proANP</b>
Calibrator (ng/mL)	1942

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	06 APR 2017