

**CERTIFICATE OF ANALYSIS**

Page 1 of 1

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

**Kit Lot Number:** K0040549

**Expiration Date:** 31 Jul 2025

**Kit Catalog Numbers:** K153MBD-Series

**Precision and Accuracy Specifications:**

		Calibrator		Rat NT-proANP Control 1		Rat NT-proANP Control 2		Rat NT-proANP Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	NT-proANP	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	NT-proANP	2240	ng/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Rat NT-proANP Plate	Z0057348	2-8°C	31 Aug 2025
Anti-rat NT-proANP Antibody	D0041019	2-8°C	30 Nov 2026
Rat NT-proANP Calibrator	A0040281	≤ -70°C	31 Oct 2025
Diluent 30	M0300038	≤ -10°C	31 Jul 2025

Additional components may be provided to execute the assay. Additional components are not lot-specific.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		14 JUL 2023