



**RFE MEMO**

Page 1 of 1

**Product Description(s):** Rat NT-proANP Kit  
**Catalog Number(s) / Size(s):** K153MBD- Series  
**Lot Number:** K0040510

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0040510 has been extended to 31 May 2024.

The expiration date for the following kit component has been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Rat NT-proANP Plate	Z0056837	31 May 2023	31 May 2024
Rat NT-proANP Calibrator	A0040262	30 Nov 2023	31 May 2024
Diluent 30	M0300032	30 Sep 2023	31 May 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		04 APR 2023

This memo supersedes all other statements regarding date of expiration for this product.

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Rat NT-proANP Kit  
**Kit Lot Number:** K0040510  
**Expiration Date:** 31 May 2023  
**Kit Catalog Numbers:** K153MBD-Series

**Precision and Accuracy Specifications:**

		Calibrator		Control 1		Control 2		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	2000	ng/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Rat NT-proANP Plate	Z0056837	2-8°C	31 May 2023
Anti-rat NT-proANP Antibody	D0040940	2-8°C	31 May 2024
Rat NT-proANP Calibrator	A0040262	≤ -70°C	30 Nov 2023
Diluent 30	M0300032	≤ -10°C	30 Sep 2023

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	Xiomara Talavera	Quality		27 MAY 2022