

**CERTIFICATE OF ANALYSIS**

Product Description: Rat NT-proANP Kit
Lot Number: K0040570
Expiration Date: 31 OCT 2025
Catalog Number: K153MBD-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat NT-proANP Antibody	D0041054	2-8°C	30 NOV 2027
Rat NT-proANP Plate	Z0057468	2-8°C	31 OCT 2026
Diluent 30	M0300039	≤ -10°C	31 JAN 2026
Rat NT-proANP Calibrator	A0040281	≤ -70°C	31 OCT 2025

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

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator		Rat NT-proANP Control 1		Rat NT-proANP Control 2		Rat NT-proANP Control 3	
		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

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		17 JUL 2024