



CERTIFICATE OF ANALYSIS

Product Description: Rat NT-proANP Kit
Lot Number: K0040583
Expiration Date: 31 DEC 2026
Catalog Number: K153MBD-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-rat NT-proANP Antibody	D0041080	2-8°C	31 AUG 2028
Rat NT-proANP Plate	Z0057554	2-8°C	31 JUL 2027
Diluent 30	M0300040	≤ -10°C	31 DEC 2026
Rat NT-proANP Calibrator	A0040294	≤ -70°C	30 SEP 2027

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

Precision, Accuracy, and Sensitivity 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

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		09 SEP 2025