

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Rat NT-proBNP Kit
Kit Lot Number: K0040541
Expiration Date: 31 Dec 2024
Kit Catalog Numbers: K153JK-Series

Precision and Accuracy Specifications:

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

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	NT-proBNP	96200	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat NT-proBNP Plate	Z0057265	2-8°C	28 Feb 2025
Anti-rat NT-proBNP Antibody	D0041009	2-8°C	28 Feb 2026
Rat NT-proBNP Calibrator	A0040270	≤ -70°C	31 Dec 2024

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		21 JUN 2023