

**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** Rat NT-proBNP Kit  
**Kit Lot Number:** K0040532  
**Expiration Date:** 30 Nov 2024  
**Kit Catalog Numbers:** K153JKD-Series

**Precision and Accuracy Specifications:**

Spot	Name	Calibrator		Rat NT-proBNP Control 1		Rat NT-proBNP Control 2		Rat NT-proBNP Control 3	
		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	Z0057201	2-8°C	30 Nov 2024
Anti-rat NT-proBNPAntibody	D0040987	2-8°C	30 Nov 2025
Rat NT-proBNP Calibrator	A0040270	≤ -70°C	31 Dec 2024
Diluent 34	M0340021	≤ -10°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		07 SEP 2022