

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Rat NT-proBNP Kit  
**Kit Lot Number:** K0040496  
**Expiration Date:** 31 Dec 2021  
**Kit Catalog Numbers:** K153JKD-Series

**Precision and Accuracy Specifications:**

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

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	NT-proBNP	91600	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Rat NT-proBNP Plate	Z0056717	2-8°C	30 Nov 2022
Anti-rat NT-proBNP Antibody	D0040919	2-8°C	31 Oct 2023
Rat NT-proBNP Calibrator	A0040250	≤ -70°C	30 Apr 2023
Diluent 34	M0340016	≤ -10°C	31 Dec 2021

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		05 MAR 2021