

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Rat NT-proBNP Kit
Kit Lot Number: K0040531
Expiration Date: 30 Apr 2023
Kit Catalog Numbers: K153JKD-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Rat NT-proANP/BNP Control 1		Rat NT-proANP/BNP Control 2		Rat NT-proANP/BNP 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	91600	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat NT-proBNP Plate	Z0057157	2-8°C	30 Sep 2024
Anti-rat NT-proBNP Antibody	D0040949	2-8°C	31 Mar 2025
Rat NT-proBNP Calibrator	A0040250	≤ -70°C	30 Apr 2023
Diluent 34	M0340020	≤ -10°C	30 Nov 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		24 JUN 2022