



CERTIFICATE OF ANALYSIS

Product Description: Rat NT-proBNP Kit
Lot Number: K0040573
Expiration Date: 31 JUL 2026
Catalog Number: K153JKD-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-rat NT-proBNP Antibody	D0041061	2-8°C	31 JAN 2028
Rat NT-proBNP Plate	Z0057483	2-8°C	31 DEC 2026
Rat NT-proBNP Calibrator	A0040288	≤ -70°C	31 JAN 2027
Diluent 34	M0340023	≤ -10°C	31 JUL 2026

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

Precision, Accuracy, and Sensitivity 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	95400	pg/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		27 SEP 2024