

**CERTIFICATE OF ANALYSIS****Product Description:** Rat NT-proBNP Kit**Lot Number:** K0040591**Expiration Date:** 29 FEB 2028**Catalog Number:** K153JKD-Series**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Anti-rat NT-proBNP Antibody	D0041096	2-8°C	31 MAR 2029
Rat NT-proBNP 96-Well Small Spot SECTOR Plate	Z0057618	2-8°C	31 MAR 2028
Diluent 34	M0340025	≤ -10°C	29 FEB 2028
Rat NT-proBNP Calibrator	A0040299	≤ -70°C	30 APR 2028

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	94400	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
<b>Review/Approval</b>	Karen Hamilla	Quality		01 APR 2026