



CERTIFICATE OF ANALYSIS

Product Description: Rat NT-proBNP Kit
Lot Number: K0040584
Expiration Date: 31 JAN 2027
Catalog Number: K153JKD-Series

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|---|------------|---------------------|-----------------|
| Anti-rat NT-proBNP Antibody | D0041081 | 2-8°C | 31 AUG 2028 |
| Rat NT-proBNP 96-Well Small Spot SECTOR Plate | Z0057561 | 2-8°C | 31 AUG 2027 |
| Diluent 34 | M0340024 | ≤ -10°C | 30 SEP 2027 |
| Rat NT-proBNP Calibrator | A0040288 | ≤ -70°C | 31 JAN 2027 |

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

Precision, Accuracy, and Sensitivity Specifications:


| | | Calibrator | | Rat NT-proBNP Control 1 | | Rat NT-proBNP Control 2 | | Rat NT-proBNP Control 3 | |
|------|-----------|------------|-------|-------------------------|-------|-------------------------|-------|-------------------------|-------|
| Spot | Name | 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 |  | 13 JUN 2025 |