

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Rat NT-proBNP Kit
Kit Catalog Numbers: K153JKD-Series
Kit Lot Number: K0040457
Expiration Date: 30 Sep 2020

Acceptance Data/ Percent Recovery 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 Material Tested :

| Spot | Name | Item | Conc | Units |
|------|-----------|-----------------------------|-------|-------|
| 1 | NT-proBNP | 20X Calibrator | 90200 | pg/mL |
| | | Rat NT-proANP/BNP Control 1 | 7901 | pg/mL |
| | | Rat NT-proANP/BNP Control 2 | 990 | pg/mL |
| | | Rat NT-proANP/BNP Control 3 | 186 | pg/mL |

Controls data was generated utilizing internal reference controls.

Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentrations, please refer to the Standard Curve Data

Standard Curve Data:

| NT-proBNP - Spot 1 | | |
|--------------------|---------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 4510 | 355,000 | 2.7 |
| 1130 | 70,500 | 4.0 |
| 282 | 12,000 | 1.6 |
| 70.5 | 2,100 | 2.0 |
| 17.6 | 559 | 2.5 |
| 4.40 | 212 | 3.5 |
| 1.10 | 138 | 2.9 |
| 0.00 | 109 | 5.3 |

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|-----------------------------|------------------|---------------------|-----------------|
| Rat NT-proBNP Plate | Z0056437 | 2-8°C | 30 Jun 2021 |
| Anti-rat NT-proBNP Antibody | D0040820 | 2-8°C | 30 Sep 2022 |
| Rat NT-proBNP Calibrator | A0040207 | ≤ -70°C | 30 Sep 2020 |
| Diluent 34 | M0340015 | ≤ -10°C | 31 Jul 2021 |
| Protease Inhibitor Solution | Not Kit Specific | Not Kit Specific | N/A |
| MSD Read Buffer T | Not Kit Specific | Not Kit Specific | N/A |

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|------------------------|--------------|----------|-------------------------------------------------------------------------------------|-------------|
| Review/Approval | Dawn Stollar | Quality |  | 08 MAY 2019 |