



CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-05029

Product Description: Rat BNP Kit

Kit Catalog Numbers: K153KFD-1, K153KFD-2, K153KFD-4

Kit Lot Number: K0040243

Expiration Date: 30 NOV 2015

Acceptance Data:

| Metric | | Specifications | Assay Rat BNP |
|--------------------------|-------------|----------------|------------------|
| % Recovery Concentration | Calibrators | 80% - 120% | Pass |
| | Controls | 75% - 125% | Pass |
| % Recovery CV | Calibrators | < 20% | Pass |
| | Controls | < 20% | Pass |

Concentration of Materials Tested:

| Units = pg/mL | Assay Rat BNP |
|-----------------------------|------------------|
| 20X Calibrator ¹ | 32 400 |
| Control 1 | 1196 |
| Control 2 | 252 |
| Control 3 | 75.3 |

¹ Concentration provided in this table refers to the stock concentration of the calibrator. For top of the standard curve concentration, please refer to the Standard Curve Data on page 2.



CERTIFICATE OF ANALYSIS

Page 2 of 2

COA-05029

Standard Curve Data from Kit QC:

| Rat BNP | | |
|-------------------|----------------|-----|
| Conc., (pg/mL) | Average Signal | %CV |
| 1620 | 36 599 | 3.7 |
| 540 | 9947 | 1.7 |
| 180 | 2931 | 2.5 |
| 60.0 | 977 | 2.0 |
| 20.0 | 389 | 2.3 |
| 6.67 | 207 | 3.0 |
| 2.22 | 142 | 2.6 |
| 0 | 115 | 6.5 |

| Rat BNP Kit Components | | | |
|--|------------------|---------------------|-----------------|
| Description | Lot Number | Storage Temperature | Expiration Date |
| MULTI-ARRAY® 96-well Streptavidin Gold Plate | Z0020963 | 2-8°C | 31 MAR 2016 |
| Anti-rat BNP Antibody | D0040436 | 2-8°C | 31 OCT 2016 |
| Anti-rat BNP Biotinylated Capture Antibody | D0040437 | 2-8°C | 31 OCT 2016 |
| Rat BNP Calibrator | A0040089 | ≤-70°C | 30 NOV 2015 |
| Diluent 8 | M0080025 | ≤-10°C | 31 MAY 2016 |
| Protease Inhibitor Solution | Not Kit Specific | 2-8°C | N/A |
| MSD Phosphate Buffer (5X) | Not Kit Specific | Room Temperature | N/A |
| MSD Blocker A | Not Kit Specific | Room Temperature | N/A |
| MSD Read Buffer T (4X) | Not Kit Specific | Room Temperature | N/A |

Statement: The above kit lot was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|-----------------|------------|-----------------|-----------|-------------|
| Review/Approval | Dana Short | Quality Control | | 11 DEC 2014 |