



## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** Rat NT-proBNP Kit  
**Kit Catalog Numbers:** K153JKD-Series  
**Kit Lot Number:** K0040374  
**Expiration Date:** 30 Apr 2019

### Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Control 1		Control 2		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	96800	pg/mL
		Control 1	7901	pg/mL
		Control 2	990	pg/mL
		Control 3	186	pg/mL

Controls tested may not be the same lot as shipped.

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
4840	477,000	1.8
1210	105,000	2.3
303	17,400	2.4
75.6	2,850	2.4
18.9	634	4.2
4.73	210	2.7
1.18	122	4.7
0.00	92	6.7

**CERTIFICATE OF ANALYSIS**

Page 2 of 2

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Rat NT-proBNP Plate	Z0055746	2-8°C	30 Apr 2019
Anti-Rat NT-proBNP Antibody	D0040655	2-8°C	31 May 2019
Rat NT-proBNP Calibrator	A0040184	≤ -70°C	30 Apr 2019
Diluent 34	M0340010	≤ -10°C	31 May 2019
Protease Inhibitor	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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	19 JAN 2017