

CERTIFICATE OF ANALYSIS

COA-5942

Product Description: Rat NT-proBNP Kit
Kit Catalog Numbers: K153JKD-1, K153JKD-2, K153JKD-4
Kit Lot Number: K0040308
Expiration Date: 31 Jan 2017

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Control 1		Control 2	
		Recovery	CV	Recovery	CV	Recovery	CV
1	NT-ProBNP	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NT-ProBNP	20X Calibrator	88200	pg/ml
		Control 1	6735	pg/ml
		Control 2	1100	pg/ml

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
4410	386,000	7.4
1100	88,300	4.7
276	14,400	3.4
68.9	2,310	5.5
17.2	498	4.4
4.31	190	2.7
1.08	117	3.5
0.00	95	4.8

CERTIFICATE OF ANALYSIS
Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat NT-ProBNP Plate	Z0055383	2-8°C	31 Dec 2017
Anti-rat NT-proBNP Antibody	D0040540	2-8°C	31 Dec 2017
Rat NT-ProBNP Calibrator	A0040135	≤ -70°C	31 Dec 2017
Diluent 34	M0340007	≤ -10°C	30 Nov 2017
Protease Inhibitor Solution	R0010529	2-8°C	31 Jan 2017
Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A

Statement:

The above kit 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	Monika Jain	Quality Assurance	<i>Monika Jain</i>	04 SEP 2015