

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Rat NT-proBNP Kit
Kit Catalog Numbers: K153JKD-Series
Kit Lot Number: K0040396
Expiration Date: 30 Apr 2019

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	Rat 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	Rat NT-proBNP	20X Calibrator	96800	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 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:

Rat NT-proBNP - Spot 1		
Conc. pg/mL	ECL Signal	%CV
4840	408,000	2.3
1210	80,500	3.8
303	13,100	5.6
75.6	2,150	5.5
18.9	504	7.0
4.73	186	3.8
1.18	113	3.5
0.00	92	3.5

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat NT-proBNP Plate	Z0055972	2-8°C	30 Nov 2019
Anti-rat NT-proBNP Antibody	D0040704	2-8°C	30 Nov 2019
Rat NT-proBNP Calibrator	A0040184	≤ -70°C	30 Apr 2019
Diluent 34	M0340012	≤ -10°C	31 Dec 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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	05 SEP 2017