

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Rat NT-proBNP Kit
Kit Catalog Numbers: K153JKD-Series
Kit Lot Number: K0040389
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	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
		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:

NT-ProBNP - Spot 1		
Conc. pg/ml	ECL Signal	%CV
4840	440,000	2.1
1210	88,300	2.4
303	15,300	2.9
75.6	2,430	3.0
18.9	623	3.0
4.73	224	19.0
1.18	120	3.8
0.00	90	3.5

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat NT-proBNP	Z0055917	2-8°C	31 Oct 2019
Anti-rat NT-proBNP Antibody	D0040684	2-8°C	31 Jul 2019
Rat NT-proBNP Calibrator	A0040184	≤ -70°C	30 Apr 2019
Diluent 34	M0340011	≤ -10°C	31 Oct 2019
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 Assurance	<i>Dawn Stollar</i>	05 JUN 2017