

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Rat NT-proANP Kit
Kit Catalog Numbers: K153MBD-Series
Kit Lot Number: K0040411
Expiration Date: 31 Dec 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-proANP	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NT-proANP	20X Calibrator	2020	ng/mL
		Rat NT-proANP/BNP Control 1	173	ng/mL
		Rat NT-proANP/BNP Control 2	44.1	ng/mL
		Rat NT-proANP/BNP Control 3	12.7	ng/mL

Control data was generated utilizing internal reference controls.

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-proANP - Spot 1		
Conc. ng/mL	ECL Signal	%CV
101	271,000	5.2
33.7	144,000	5.7
11.2	41,300	4.2
3.74	7,730	4.3
1.25	1,450	5.4
0.416	334	3.8
0.139	131	3.5
0.00	66	7.5

CERTIFICATE OF ANALYSIS


Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat NT-proANP Plate	Z0056052	2-8°C	28 Feb 2020
Anti-rat NT-proANP Antibody	D0040741	2-8°C	29 Feb 2020
Rat NT-ProANP Calibrator	A0040203	≤ -70°C	28 Feb 2020
Diluent 30	M0300020	≤ -10°C	31 Dec 2019
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		06 MAR 2019