



CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-04571

Product Description: Rat NT-proANP Kit

Kit Catalog Numbers: K153MBD-1, K153MBD-2, K153MBD-4

Kit Lot Number: K0040222

Expiration Date: 29 FEB 2016

Acceptance Data:

Metric		Specifications	Assay
			Rat NT-proANP
% Recovery Concentration	Calibrators	80% - 120%	Pass
	Controls	80% - 120%	Pass
% Recovery CV	Calibrators	< 20%	Pass
	Controls	< 20%	Pass

Concentration of Materials Tested:

Units = ng/mL	Assay
Rat NT-proANP	
20X Calibrator	1970
Control 1	39.1
Control 2	3.44

Standard Curve Data from Kit QC:

Rat NT-proANP		
Conc., (ng/mL)	Average Signal	%CV
98.5	818 492	3.6
32.8	431 116	1.6
10.9	136 321	2.1
3.65	28 257	4.4
1.22	5384	5.0
0.405	1000	5.1
0.135	255	3.2
0	60	10.1




CERTIFICATE OF ANALYSIS

Page 2 of 2

COA-04571

Rat NT-proANP Components			
Description	Lot Number	Storage Temperature	Expiration Date
Rat NT-proANP Plate	Z0054760	2-8°C	29 FEB 2016
Anti-rat NT-proANP Antibody	D0040411	2-8°C	30 JUN 2016
Rat NT-proANP Calibrator	A0040081	≤ -70°C	30 NOV 2016
Diluent 30	M0300013	≤ -10°C	31 MAR 2016
MSD Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A

Statement: The above kit lot 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	Dana Short	Quality Control		09 JUL 2014