



RFE MEMO

Page 1 of 1

Product Description(s): Rat NT-proANP Kit
Catalog Number(s) / Size(s): K153MBD-Series
Lot Number: K0040424

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0040424 has been extended to 28 Feb 2021.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Rat NT-proANP Plate	Z0056183	31 Aug 2020	28 Feb 2021
Rat NT-proANP Calibrator	A0040203	28 Feb 2020	28 Feb 2021
Diluent 30	M0300022	30 Nov 2020	28 Feb 2021

Function	Name	Signature	Date
Quality	Karen Hamilla		14 FEB 2020

This memo supersedes all other statements regarding date of expiration for this product.

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Rat NT-proANP Kit
Kit Catalog Numbers: K153MBD-Series
Kit Lot Number: K0040424
Expiration Date: 28 Feb 2020

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	rt 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	rt 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:

rt NT-proANP - Spot 1		
Conc. ng/mL	ECL Signal	%CV
101	364,000	3.9
33.7	189,000	4.9
11.2	56,200	3.6
3.74	10,400	3.4
1.25	1,900	4.5
0.416	396	3.5
0.139	139	3.6
0.00	68	7.9

CERTIFICATE OF ANALYSIS


Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat NT-proANP Plate	Z0056183	2-8°C	31 Aug 2020
Anti-rat NT-proANP Antibody	D0040761	2-8°C	31 Dec 2021
Rat NT-proANP Calibrator	A0040203	2-8°C	28 Feb 2020
Diluent 30	M0300022	≤ -10°C	30 Nov 2020
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	Karen Hamilla	Quality		14 FEB 2020