



RFE MEMO

Page 1 of 1

Product Description(s): Rat NT-proBNP Kit
Catalog Number(s) / Size(s): K153JKD- series
Lot Number: K0040422

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0040422 has been extended to 30 APR 2020.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Rat NT-proBNP Calibrator	A0040184	30 APR 2019	30 APR 2020

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar		27 MAR 2019

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

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Rat NT-proBNP Kit
Kit Catalog Numbers: K153JKD-Series
Kit Lot Number: K0040422
Expiration Date: 30 Apr 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Rat NTproANP/BNP Control 1		Rat NTproANP/BNP Control 2		Rat NTproANP/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 NTproANP/BNP Control 1	7901	pg/mL
		Rat NTproANP/BNP Control 2	990	pg/mL
		Rat NTproANP/BNP Control 3	186	pg/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:

Rat NT-ProBNP - Spot 1		
Conc. pg/mL	ECL Signal	%CV
4840	393,000	2.0
1210	75,000	3.0
303	11,500	3.4
75.6	1,920	5.9
18.9	472	3.7
4.73	185	3.3
1.18	121	6.0
0.00	100	4.3

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat NT-proBNP Plate	Z0056184	2-8°C	31 Aug 2020
Anti-rat NT-proBNP Antibody	D0040754	2-8°C	31 Jul 2020
Rat NT-proBNP Calibrator	A0040184	≤ -70°C	30 Apr 2019
Diluent 34	M0340014	≤ -10°C	31 Aug 2020
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>	28 FEB 2019