



RFE MEMO

Page 1 of 1

Product Description(s): Rat NT-proANP Kit
Catalog Number(s) / Size(s): K153MBD-Series
Lot Number: K0040498

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0040498 has been extended to 31 JUL 2022.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Anti-rat NT-proANP Antibody	D0040761	31 DEC 2021	31 JUL 2022
Rat NT-proANP Calibrator	A0040227	31 AUG 2021	31 JUL 2022
Diluent 30	M0300025	31 JUL 2021	31 JUL 2022

Function	Name	Signature	Date
Quality	Karen Hamilla		02 AUG 2021

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

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Rat NT-proANP Kit
Kit Lot Number: K0040498
Expiration Date: 31 Jul 2021
Kit Catalog Numbers: K153MBD-Series

Precision and Accuracy 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%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	NT-proANP	1794	ng/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat NT-proANP Plate	Z0056724	2-8°C	30 Nov 2022
Anti-rat NT-proANP Antibody	D0040761	2-8°C	31 Dec 2021
Rat NT-proANP Calibrator	A0040227	≤ -70°C	31 Aug 2021
Diluent 30	M0300025	≤ -10°C	31 Jul 2021

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		23 NOV 2020