



## Extension Memo


Page 1 of 1

**Product Description:** Rat BNP Kit  
**Catalog Number(s):** K153KFD-Series  
**Lot Number:** K0040571  
**New Expiration Date:** 31 May 2026

Based on additional testing, the original date of expiration for this kit has been extended by 16 months.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Rat BNP Calibrator	A0040276	30 APR 2025	31 May 2026
Diluent 8	M0080043	31 JAN 2025	31 May 2026
Anti-rat BNP Biotinylated Capture Antibody	D0041012	31 MAR 2026	31 May 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		03 Jun 2025



## CERTIFICATE OF ANALYSIS

**Product Description:** Rat BNP Kit  
**Lot Number:** K0040571  
**Expiration Date:** 31 JAN 2025  
**Catalog Number:** K153KFD-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-rat BNP Antibody	D0041013	2-8°C	30 NOV 2026
Anti-rat BNP Biotinylated Capture Antibody	D0041012	2-8°C	31 MAR 2026
MSD GOLD™ 96-well Streptavidin SECTOR Plate	Z0022275	2-8°C	30 SEP 2026
Diluent 8	M0080043	≤ -10°C	31 JAN 2025
Rat BNP Calibrator	A0040276	≤ -70°C	30 APR 2025

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

**Precision, Accuracy, and Sensitivity Specifications:**


		Calibrator		Rat BNP Control 1		Rat BNP Control 2		Rat BNP Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	BNP	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	BNP	40400	pg/mL

Dilution to top of Curve: 20X

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		27 JUN 2024