

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Rat BNP Kit
Kit Lot Number: K0040544
Expiration Date: 28 Feb 2025
Kit Catalog Numbers: K153KFD-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Rat BNPControl 1		Rat BNPControl 2		Rat BNPControl 3	
		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

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
MSD GOLD™ 96-well Streptavidin Plate	Z0021960	2-8°C	28 Feb 2025
Anti-rat BNP Antibody	D0041013	2-8°C	30 Nov 2026
Rat BNP Calibrator	A0040276	≤ -70°C	30 Apr 2025
Anti-rat BNP Biotinylated Capture Antibody	D0041012	2-8°C	31 Mar 2026

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		19 JUN 2023