



### 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:

Spot	Name	Calibrator		Rat BNP Control 1		Rat BNP Control 2		Rat BNP Control 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

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