

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** Rat BNP Kit  
**Kit Catalog Numbers:** K153KFD-Series  
**Kit Lot Number:** K0040456  
**Expiration Date:** 30 Nov 2020

**Acceptance Data/ Percent Recovery 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 Material Tested :**

Spot	Name	Item	Conc	Units
1	BNP	20X Calibrator	41600	pg/mL
		Rat BNP Control 1	6700	pg/mL
		Rat BNP Control 2	3840	pg/mL
		Rat BNP Control 3	1000	pg/mL

Control data was generated utilizing internal 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:**

BNP - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2080	57,700	1.8
693	16,800	1.1
231	5,490	1.1
77.0	2,100	1.0
25.7	1,080	1.7
8.56	749	2.7
2.85	645	2.3
0.00	589	2.1

## CERTIFICATE OF ANALYSIS


Page 2 of 2

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
MSD Gold Plate	Z0021440	2-8°C	30 Nov 2020
Anti-rat BNP Antibody	D0040818	2-8°C	31 Aug 2022
Rat BNP Calibrator	A0040226	≤ -70°C	30 Jun 2021
Blocker A	Not Kit Specific	Not Kit Specific	N/A
Diluent 8	M0080034	≤ -10°C	31 Jul 2021
Anti-rat BNP Biotinylated Capture Antibody	D0040819	2-8°C	31 Aug 2022
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
<b>Review/Approval</b>	Dawn Stollar	Quality		12 APR 2019