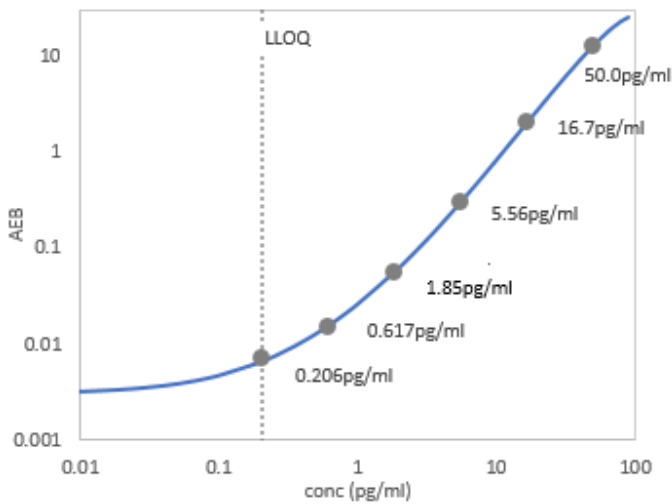


Description

Natriuretic peptides are produced primarily within the heart and released into the circulation in response to increased wall tension, reflecting increased volume or pressure overload.¹ Under pathologic stimuli, the prohormone of BNP is synthesized, cleaved to BNP, the biologically-active peptide, and inactive fragment NT-proBNP, a 76aa amino terminal peptide.² Both BNP and NT-proBNP are released predominantly by the ventricles in response to stretch, and are used for the diagnosis of systolic heart failure.³ NT-proBNP circulates at higher plasma concentrations and has a longer half-life when compared with BNP.⁴ Both peptides have proved equally useful for the diagnosis of ambulatory patients with heart failure and left ventricular dysfunction.⁵

Calibration Curve: Calibrator concentrations and Lower Limit of Quantification depicted (Cubic fitting).



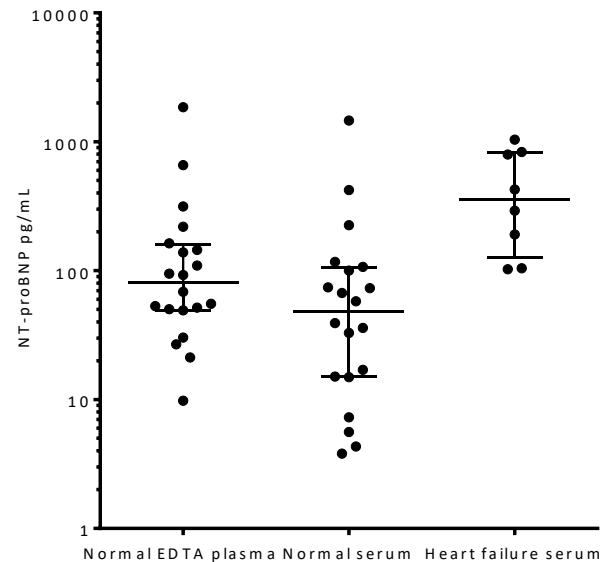
Lower Limit of Quantification (LLOQ): Triplicate measurements of serially diluted calibrator were read back on the calibration curve for 5 runs using 1 reagent lot on 3 instruments. The LLOQ is determined as the lowest dilution with a pooled CV ≤ 20% and sample concentration recovery between 80-120% of the expected.

Limit of Detection (LOD): Calculated as 2.5 standard deviations from the mean of background signal read back on each calibration over 5 runs for 1 reagent lot on 3 instruments.

| | |
|-------------------------------|---|
| LLOQ | 0.206pg/mL pooled CV 15.8%, mean recovery 101.9% |
| LOD | 0.0433 pg/mL range 0.0071-0.1190 pg/mL |
| Sample range | 0–500 pg/mL |
| Diluted Sample volume* | 100 µL per measurement |
| Tests per kit | 96 |

***See Kit Instruction for details**

Endogenous Sample Reading: Healthy donor matched EDTA plasma (n=20) and serum (n=20), and Heart failure samples (n=8) were measured. Bars depict median with interquartile range.



| Human samples | Mean pg/mL | Median pg/mL | % Above LOD |
|---------------------|------------|--------------|-------------|
| EDTA plasma | 265.2 | 93.7 | 100% |
| Serum | 200.9 | 70.6 | 100% |
| Heart failure serum | 473.4 | 359.2 | 100% |

Precision: Measurements of 3 serum or plasma based panels. Triplicate measurements were made for 5 runs using 1 reagent lot and 3 instruments (5 runs total, 15 measurements).

| Sample | Mean (pg/mL) | Within run CV | Between run CV |
|---------|--------------|---------------|----------------|
| Panel 1 | 7.347 | 6.0% | 7.7% |
| Panel 2 | 30.67 | 2.7% | 8.7% |
| Panel 3 | 172.6 | 2.7% | 7.5% |

Spike and Recovery: 4 serum samples were spiked at high and low concentrations within the range of the assay.

Dilution Linearity: 1 endogenous plasma sample was diluted 2x serially from MRD (10x) to 320x with Sample Diluent.

| | |
|--|------------------------------------|
| Spike and Recovery | 96.4% Range 87.8%-109.5% |
| Dilution Linearity (Plasma, 320x) | 84.3% Range: 72.0%-98.4% |

Specificity: Normal serum (n=3) were directly incubated with 20X capture beads and run at MRD. Average knock-down was **93.3%** with a range of **92.2% -94%**.

References:

1. Kinnunen P, Vuolteenaho O, Ruskoaho H. Mechanisms of atrial and brain natriuretic peptide release from rat ventricular myocardium: effect of stretching. *Endocrinology*1993;**132**:1961–70
2. Agency for Health Care Policy and Research *Heart failure evaluation and care of patients with left ventricular systolic dysfunction*1994Rockville, MD Agency for Health Care Policy and Research.
3. Mair J, Hammerer-Lercher A, Puschendorf B. The impact of cardiac natriuretic peptide determination on the diagnosis and management of heart failure. *Clin Chem Lab Med* 20001; 39:571-588
4. Downie PF, Talwar S, Squire IB, Davies JE, Barnett DB, Ng LL. Assessment of the stability of N-terminal pro-brain natriuretic peptide in vitro: implications for assessment of left ventricular dysfunction. *Clin Sci (LOND)* 1999; 97:255-258
5. Masson S, Vago T, Baldi G. Comparative measurement of N-terminal pro-brain natriuretic peptide and brain natriuretic peptide in ambulatory patients with heart failure *Clin Chem Lab Med*2002**40**:761-763